

Meta Imaging Series[®] Software MetaMorph[®] Software

Visual Basic Reference Guide

Version 7.6.5 for Microsoft Windows XP and Microsoft Windows Vista

1020 2201-02

This document is provided to customers who have purchased Molecular Devices, Inc. ("Molecular Devices") equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Equipment, software, reagents, and consumables that may be described in this document are protected under one or more patents filed in the United States, Canada, and other countries. Additional patents are pending.

Software that may be described in this document is furnished under a license agreement. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Molecular Devices makes no warranties or representations as to the fitness of this equipment for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom.

For research use only. Not for use in diagnostic procedures.

META IMAGING SERIES, METAMORPH, METAFLUOR and METAVUE are registered trademarks of Molecular Devices, Inc. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, Inc.

Equipment built by Molecular Devices, Inc. 1311 Orleans Drive, Sunnyvale, California, United States of America 94089. Molecular Devices, Inc. is ISO 9001 registered. © 2010 Molecular Devices, Inc. All rights reserved. Printed in the USA.

Contents

Chapter 1: Introduction	
Overview	
Conventions Used in This Manual	4
The Run User Program Command	
Visual Basic and User Programs	
Creating a User Program	
Data Types and Arrays	
Chapter 2: Performing Serial and Digital I/O Communication	
Overview	
Performing Serial Data Transmission	
Communicating with a Digital I/O Device	
Chapter 3: Executing Commands and Journals	
Overview	
Executing Commands and Journals	
-	
Chapter 4: Reading and Manipulating Images and Image Windows	
Overview	
Loading, Creating, Copying, and Closing Images	
Finding Loaded Images	
Manipulating Image Windows	
Reading and Using Image Properties	
Chapter 5: Adjusting Image Display	
Overview	60
Updating the Image After Changing the Display	61
Adjusting Brightness and Contrast	
Autoscaling 16-Bit Images	
Working with Look-up Tables and Palettes	71
Chapter 6: Reading and Using Image Pixel Data	
Overview	77
Applying Thresholding	
Reading and Manipulating Image Data	
Chapter 7: Working with Regions of Interest	
Overview	
Creating and Removing Regions	
Finding Regions	
Reading and Manipulating Region Properties	
Reading Image Data from Regions	
Chapter 8: Performing Morphometry	
Overview	
Configuring Measurement Preferences	
Configuring Object Measurements	
Configuring Classifier Filters	
Measuring All Objects in an Image	
Measuring Single Objects	

Figures

Introducti	on	
1.1	The Run User Program Dialog Box	6
1.2	Class Module UserMethods Section Code	14
Reading a	nd Manipulating Images and Image Windows	
4.1	Image "Get" Function Programming Example	56
Working v	with Regions of Interest	
7.1	Region Property "Get" Function Programming Example	
7.2	Region Data "Get" Function Programming Example	
Performin	g Morphometry	
8.1	Single Object Data "Get" Function Programming Example	

Tables

Introduction

1.1	Creating a User Program with Visual Basic .NET 2005/2008	.11
1.2	Creating a User Program with Visual Basic .NET 2002/2003	.12
1.3	Creating a User Program with Visual Basic Version 5 or 6	15
1.4	Data Types and Arrays	.18
Performing Se	rial and Digital I/O Communication	
2.1	Syntax Rules for Serial Data Transmission	.23
2.2	ASCII Control Codes	.24

Chapter 1 – Introduction

1.1 Overview

Introduction

Welcome to the Visual Basic Reference Guide for the MetaMorph[®] software for image processing and analysis, and the Meta Imaging Series[®] software system. This manual has been designed to serve as a working reference for the use of the extended set of Visual Basic programming functions by the Run User Program drop-in command in the MetaMorph[®] software. The Run User Program command, located in the File menu, provides you with the ability to run your own Visual Basic programs from within the MetaMorph program. This will allow you to process and analyze images with greater flexibility.

The full power of Microsoft Visual Basic is available to you, with the ability to run conditional "If...Then...Else" routines, to create nested subroutines and loops, and to pass and return values with the imaging system. It is assumed that you already have a working knowledge of Visual Basic and its application. This manual describes the extended set of programming functions that MetaMorph provides, and is intended to serve as a supplement to any Visual Basic references that you may already have.

Note: The following versions of Microsoft Visual Basic can be used to create programs to run in version 7.0 of the MetaMorph software:

- Microsoft Visual Basic .NET 2008 Professional or Enterprise editions
- Microsoft Visual Basic .NET 2005 Professional or Enterprise editions
- Microsoft Visual Basic .NET 2003 Professional or Enterprise editions
- Microsoft Visual Basic .NET 2002 Professional or Enterprise editions
- Microsoft Visual Basic 6.0 Professional or Enterprise editions
- Microsoft Visual Basic 5.0 Professional or Enterprise editions

Note that the Standard and Express editions of the above versions are NOT supported.

This *Reference Guide* starts with some fundamental concepts about Visual Basic user programs and the Run User Program command. The rest of this manual is devoted to a description of the extended set of functions available for use in MetaMorph. Each entry starts with a short description of what the function does. The full syntax of the expression is given, and the involved parameters and returns are explained. For many entries, a "See Also" list of related or otherwise relevant functions is provided. For the most part, these functions will be in the same section of the manual as the function being described. For those located elsewhere in the manual, the pertinent section will be given in parentheses.

The chapters in this guide are roughly arranged in the order that you might expect to use them through the course of an experiment and thereafter:

- Chapter 1 describes how to install the RUNUSER drop-in and create your own user programs with Microsoft Visual Basic.
- Chapter 2 describes functions that are used to communicate with serial or digital devices.
- Chapter 3 covers execution of commands and journals.
- Chapters 4 and 5 deal with the use of image windows and adjusting their display.
- Chapter 6 concerns thresholding and performing basic densitometric procedures.
- Chapter 7 describes how to work with regions of interest.
- And finally, Chapter 8 ends the manual with a description of the morphometric functions that are available.

1.1 Overview, continued

In this chapter This chapter contains the following topics:

Торіс	See Page
Conventions Used in This Manual	4
The Run User Program Command	6
Visual Basic and User Programs	8
Creating a User Program 11	
Data Types and Arrays	18

1.2 Conventions Used in This Manual

Introduction Some definitions	Before we proceed, it may be useful first to define some of the terms used in this manual and to review some of the typographical conventions that are used.			
	<i>Array</i> A multidimensional (usually one- or two-dimensional) "table" that stores values. For example, you can define an array and read into it the X and Y coordinates of an object's outline by using the MorphGetVertexList function. In Visual Basic, you can declare arrays of up to 60 dimensions. (See Section 1.6, <i>Data Types and Arrays.</i>)			
	<i>Boolean</i> In its widest sense, this is a logical operand such as AND, OR, or NOT. When used to describe a variable, a Boolean indicates a logical state, such as TRUE or FALSE. As integers, FALSE is represented by 0 and TRUE is represented by -1 or some other nonzero value.			
	<i>Double</i> A double-precision (64-bit) floating-point number that encodes the value of a variable. Numbers can range from approximately -1.8×10^{308} through -4.9×10^{-324} for negative values and 4.9×10^{-324} through 1.8×10^{308} for positive values.			
	<i>Function</i> A programming "command" that carries out some procedure. For example, the PrintMsg function prompts MetaMorph to display a message box that you have previously configured.			
	<i>Handle</i> An index number used by the program to deal with and keep track of images, image windows, functions, regions of interest, and MetaDevices.			
	<i>Integer</i> In Visual Basic, a 16-bit binary signed (negative or positive) whole number that encodes the value of a variable. Integers can range from -32,767 through 32,767.			
	<i>Long</i> A 32-bit signed whole number that encodes the value of a variable. Numbers can range from -2,147,483,647 through 2,147,483,647.			
	<i>Parameter</i> A value or an identifying element that is passed by the user to the program, such as the handle of an image you want to measure, or a set of coordinates that you want to use to position an image window.			
	<i>Return value</i> A value or state returned by the program in response to a query. For example, the upper and lower limits of a threshold range, expressed as integers, are the return values for the GetThresholdRange function.			
	<i>Single</i> A single-precision (32-bit) floating point number that encodes the value of a variable. Numbers can range from approximately -3.4×10^{38} through -1.4×10^{-45} for negative values and 1.4×10^{-45} through 3.4×10^{38} for positive values.			
	String A segment of alphanumeric text.			
	<i>Variable</i> A placeholder that represents something that is acted upon by a function. For example, the variable <i>nObjectID</i> is used to represent the object in an image that is to be deleted by the MorphDeleteObject function.			
	<i>Variant</i> A variable that is not explicitly assigned to a particular data type. Variants can store different types of data—integers, floating point numbers, or character strings—depending on the need at the time.			
	Continued on next page			

1.2 Conventions Used in This Manual, continued

Typographical
conventionsThe following table lists the typographical conventions used in the Visual Basic
Reference Guide.

This:	Represents
Italics	Variables Examples: <i>hImage, nXPos</i>
Bold	Functions Examples: LoadImage, SetRegionSize
ALL CAPITALS	Returned or passed states or Booleans Examples: EXCLUSIVE, TRUE

1.3 The Run User Program Command

Installing the RUNUSER drop-in	MetaMorph allows you to run Visual Basic functions by using the Run User Program drop-in command. As with all other drop-ins, you will need to use the Configure Drop-ins command in the Meta Imaging Series Administrator to load the RUNUSER drop-in prior to starting MetaMorph. This will place the Run User Program command in your File menu, below the Run Program command (don't confuse the two!).
The Run User Program dialog box	When you choose Run User Program from the File menu, a dialog box will appear which contains a drop-down list box, a text box, a check box, and four command buttons (see Fig. 1.1). This dialog box provides an interface from within MetaMorph for you to pass parameters to the user program which you will have written in Visual Basic.



Run User Program	×
Program name:	
Command Line:	Browse
	Remove
Keep program in memory after execution	Cancel

Run User Program dialog box options

Program Name

Contains a list of all of your user programs that are currently registered with the system. The entries in the list will be descriptions that were entered in Visual Basic when you created your program. If there is no description, the project name entered in Visual Basic will appear instead.

Command Line

This is a text field which will be passed to your user program as the parameter for the **Startup** and **DoCommand** functions (see Section 1.4). You might use this field, for example, to specify the name of an image that your program will then load, convolve with an image filter, threshold, measure, log measurement data from, and close.

1.3 The Run User Program Command, continued

Run User Program dialog box options (continued)

Keep Program in Memory After Execution

This check box determines whether your user program will stay in memory after running, or if it will be unloaded when the routine is completed. If you select the check box, the program will stay in memory, and will therefore run more quickly on subsequent runs. If you clear this check box, the program will be unloaded after running. This may be useful when you are debugging your program, as you can leave both MetaMorph and Visual Basic running at the same time, alternating between running your program and editing it. Visual Basic would not be able to recompile your program if the check box were still selected.

When the *Keep Program in Memory* check box is selected, the **Startup** function will be called the first time your program is run after being loaded. Subsequent runs will call the **DoCommand** function. When this check box is cleared, the **Shutdown** function will be called after the program finishes. If this check box is cleared before you run the program for the first time, the **Startup** and **Shutdown** functions will be called each time you run the program (see Section 1.4 for more about these three functions).

Browse

If the program you want to run does not appear in the *Program Name* list, this command button will allow you to search your system for it. This button opens the Select Start File dialog box, which is a standard file-selection dialog box that has a *Look In* drop-down list, Up One Level icon button, *File Name* text box, and a table that displays the files in the current folder.

When you create a user program, Visual Basic registers it with the system when it is compiled, and it should then be available in the *Program Name* list. However, if you obtain a program that was compiled elsewhere, it will not have been registered on your system. The *Browse* command button will register the program and insert it in the *Program Name* list.

NOTE- this only works for user programs created with Visual Basic 6 and earlier.

Remove

Choosing this button will remove the currently highlighted user program from the *Program Name* list and unregisters it with the system.

NOTE- this only works for user programs created with Visual Basic 6 and earlier.

ок

Loads the program selected in the *Program Name* list, passes the parameter you specify in the *Command Line* text box, runs the user program, and closes the Run User Program dialog box.

Cancel

Cancels any changes made in the Run User Program dialog box and closes it.

1.4 Visual Basic and User Programs

About user programs	A user program is essentially an ActiveX object that communicates with the MetaMorph application using Microsoft's object linking and embedding (OLE) interface. However, it is not necessary to be familiar with the details of this communication when using Microsoft Visual Basic, the language supported by the MetaMorph software to write user programs. Note: The following versions of Microsoft Visual Basic can be used to create programs to run in version 7.0 of the MetaMorph software: Microsoft Visual Basic .NET 2008 Professional or Enterprise editions Microsoft Visual Basic .NET 2005 Professional or Enterprise editions Microsoft Visual Basic .NET 2003 Professional or Enterprise editions Microsoft Visual Basic .NET 2002 Professional or Enterprise editions Microsoft Visual Basic 5.0 Professional or Enterprise editions Nicrosoft Visual Basic 5.0 Professional or Enterprise editions	
	Note that the Standard and Express editions of the above versions are NOT supported.	
Variables and functions required by the MetaMorph software	To be recognized by the MetaMorph software, a Visual Basic user program must contain three variables, <i>MM</i> , <i>gParentWnd</i> and <i>gUserID</i> , and three functions, Startup , DoCommand , and Shutdown . There are several other more technical requirements, but these will be covered in the discussion of the use of specific versions of Visual Basic (Section 1.5). The two variables, <i>gParentWnd</i> and <i>gUserID</i> , are of type Long, and <i>MM</i> is a variant:	
	MM MM is a variant that is used for all of your communication with the MetaMorph software. For example, to put the handle of the current image in the MetaMorph software into the variable <i>sourceImage</i> , you would use the following:	
	Public sourceImage As Long MM.GetCurrentImage sourceImage	
	gParentWnd	
	This will have the handle of the MetaMorph program's main window placed in it. It is primarily useful for languages other than Visual Basic which are not yet supported.	
	gUserID	
	This variable is for future expansion, so you won't be using it immediately, but it must be present nonetheless.	
	Continuea on next page	

1.4 Visual Basic and User Programs, continued

Variables and functions required by the MetaMorph software software (continued)

The three functions are Startup, DoCommand, and Shutdown.

```
Startup(cmdLine As String) As Long (or Integer)
```

This function is called when the program needs to be loaded into memory to run. *cmdLine* will contain the text entered in the *Command Line* text box of the Run User Program dialog box. **Startup** should return a value of 0 on success and a nonzero value on failure. The return value is currently ignored in MetaMorph, but in the future it may be used.

DoCommand(cmdLine As String) As Long (or Integer)

This function is called when the program is run after it has already been loaded into memory. *cmdLine* will contain the text entered in the *Command Line* text box of the Run User Program dialog box. **DoCommand** should return a value of 0 on success and a nonzero value on failure. As with the **Startup** function return, this return value is currently ignored in MetaMorph, but in the future it may be used. If your program behaves the same way whether it is already loaded or not, you may want to have **Startup** simply call **DoCommand**, and put the code that does the work in **DoCommand**.

Shutdown() As Long (or Integer)

This function is called when the user program is unloaded from memory. **Shutdown** should return a value of 0 on success and a nonzero value on failure. The return value is currently ignored in MetaMorph but in the future it may be used.

1.4 Visual Basic and User Programs, continued

Class modules	Note : the Class modules section below applies only to Visual Basic versions 5 and 6. It does NOT apply either version of Visual Studio .NET.		
	All of the above functions and variables must be in a Class Module named UserMethods.		
	To access the <i>MM</i> variable from other modules, you must declare a Public Variant in one of the other modules, and then in Startup and DoCommand , use Set to assign <i>MM</i> to that variable. For example, suppose your program consists of the Class Module UserMethods, and two regular Modules, Mod1 and Mod2:		
	Modl: Public <i>pubMM</i> As Variant		
	UserMethods: Public <i>MM</i> as MMAppLib.UserCall		
	Function Startup (<i>cmdLine</i> As String) As Long DoCommand <i>cmdLine</i> End Function		
	<pre>Function DoCommand(cmdLine As String) As Long Set pubMM = MM</pre>		
	Doit End Function		
	Mod2: Function Doit Dim <i>i</i> As Long		
	pubMM.GetCurrentImage i End Function		
	You see that <i>pubMM</i> is visible to the whole program. Set must be used for the assignment, since <i>MM</i> is a variant. If Set is not used, problems will occur.		

1.5 Creating a User Program

Introduction	Note: The following versions of Microsoft Visual Basic can be used to create programs to run in MetaMorph 7.0:	
	Note that	Microsoft Visual Basic .NET 2008 Professional or Enterprise editions Microsoft Visual Basic .NET 2005 Professional or Enterprise editions Microsoft Visual Basic .NET 2003 Professional or Enterprise editions Microsoft Visual Basic .NET 2002 Professional or Enterprise editions Microsoft Visual Basic 6.0 Professional or Enterprise editions Microsoft Visual Basic 5.0 Professional or Enterprise editions Microsoft Visual Basic 5.0 Professional or Enterprise editions
		the Standard and Express editions of the above versions are 1001 supported.
Note	One limitation of user programs written in Visual Basic versions 5 and 6 must be noted: because of the manner in which the programs run, the user will not be able to use non-modal dialog boxes. That is, you will be able to work in only one command dialog box at a time. Any attempt to run a program with non-modal dialog boxes will return an error message from the Visual Basic runtime controller.	
	Visual Stu has not be Molecular	adio .NET enables the use of non-modal dialog boxes; however, their use een tested with the MetaMorph software and they are not supported by r Devices. Use at your own risk.
Creating a user program with Visual Basic .NET	To create one of the	the framework for a user program using Microsoft Visual Basic .NET, use procedures presented in the following tables.
	Table 1.1	Creating a User Program with Visual Basic .NET 2005 or 2008
	Step	Action
	1	Open Visual Studio and create a new project of type Visual Basic: Class Library.
	2	Select My Project in the Solution Explorer Select the App Tab and Press "Assembly Info" button Check off "MakeAssemblyCom – visible Press OK
	3	Select the References Tab and click "Add" button Select COM Tab. Find the MetaMorph Type Library and select it. Press OK - Adds MetaMorph to the Reference List
	4	In the Solution Explorer, choose the project and right click on it Choose Add > New Item Choose "module" and click Add
	5	Add the code in Figure 1.2 (below) to Module1.vb.
	6	Add any code for your VB application.

7	If deploying this user program to multiple computers with MetaMorph, you must perform the following:	
	1.	In Solution Explorer, highlight the solution, right click on it and choose Add > New Project
	2.	Find and choose "set up project" and press OK
	3.	Right click on the set up project and choose Add > Project Output
	4.	Choose the name of your project from the drop down menu
	5.	Select Primary Out from the list in the window
	6.	Choose Configure > Release
	7.	Press OK
8	From the list of Detected Dependencies, choose "mmapp.exe" and right click on it. Select "Exclude"	
9	Under the Build menu, select Build Solution. Once the program has been compiled without any errors, it is ready to be run from the MetaMorph application	

Note

The default condition for .NET 2005/2008 for Assembly Name and Root Namespace MUST be the same as the project name. DO NOT CHANGE these names or this will not work with the MetaMorph application.

Table 1.2Creating a User Program with Visual Basic .NET 2002 or 2003

Step	Action				
1	Open Visual Studio and create a new project of type Visual Basic: Windows Application.				
2	Add the following lines to the AssemblyInfo.vb file:				
	<assembly: comvisible(true)=""></assembly:>				
	<pre><assembly: classinterface(classinterfacetype.autodual)=""></assembly:></pre>				
3	From the Project menu, choose Properties, and go to the General pane in the window that appears (it should start with the General pane automatically).				
4	Bring up the properties for the project, select Common Properties, General. Under Output Type, select Class Library. Under Root Namespace, type in the name you want to appear in MetaMorph in the Run User Program dialog for your project. Note that if you are running Visual Basic.NET Standard Edition, you will not have a Class Library option. In this case, close Visual Studio, and open the folder that contains your project. Find the file with the .vbproj extension and edit it using Notepad. Change the line that says OutputType = "WinEXE" to OutputType = "Library" Change the line that says StartupObject = "YourApplicationName.YourFormName" to StartupObject = ""				
	continue with these steps.				
5	Bring up the properties for the project, select Configuration Properties, then Build. Check the <i>Register for COM Interop</i> checkbox.				

6	On the Project menu, select Add Reference, select the COM pane, click the Browse button, and find your MMApp.exe file and select it. Then click <i>OK</i> to close the Add Reference window.			
7	Add the code in Figure 1.2 (below) to Module1.vb.			
8	Add any code for your VB application.			
9	Under the Build menu, select Build Solution. Once the program has been compiled without any errors, it is ready to be run from MetaMorph			

```
Option Strict Off
Option Explicit On
Public Interface IUserMethods
   Property mm() As MMAppLib.UserCall
    Property gParentWnd() As Integer
   Property gUserID() As Integer
   Function Startup(ByRef cmdLine As String) As Integer
   Function Docommand(ByRef cmdLin As String) As Integer
   Function Shutdown() As Integer
End Interface
<ComClass(UserMethods.ClassId, UserMethods.InterfaceId)> _
Public Class UserMethods
   Implements IUserMethods
   Private mygParentWnd As Integer
   Private mygUserID As Integer
   Public mymm As MMAppLib.UserCall
    Public Const ClassId As String = "832F34A5-5CF5-403f-B4A8-428C8351FD02"
   Public Const InterfaceId As String = "3D8B5BA4-FB8C-5ff8-8468-
11BF6BD5CF91"
    Property mm() As MMAppLib.UserCall Implements IUserMethods.mm
        Get
           Return mymm
        End Get
        Set(ByVal Value As MMAppLib.UserCall)
           mymm = Value
        End Set
   End Property
    Property gParentWnd() As Integer Implements IUserMethods.gParentWnd
       Get
           Return mygParentWnd
        End Get
        Set(ByVal Value As Integer)
           mygParentWnd = Value
       End Set
    End Property
   Property gUserID() As Integer Implements IUserMethods.gUserID
       Get
           Return mygUserID
        End Get
        Set(ByVal Value As Integer)
           mygUserID = Value
        End Set
   End Property
   Public Function Startup(ByRef cmdLine As String) As Integer Implements
IUserMethods.Startup
       Docommand(cmdLine)
   End Function
   Public Function Docommand(ByRef cmdLine As String) As Integer Implements
IUserMethods.Docommand
   End Function
   Public Function Shutdown() As Integer Implements IUserMethods.Shutdown
   End Function
End Class
```

1.5 Creating a User Program, continued

Creating a user program with Visual Basic 5.0 or 6.0

To create the framework for a user program using Microsoft Visual Basic version 5.0 or 6.0, use the procedure presented in the following table.

Step	Action			
1	Start the Microsoft Visual Basic program and create a new project by choosing New Project from the File menu.			
2	For the type of project to create, select ActiveX DLL.			
3	From the Project menu, choose Properties, and go to the General pane in the window that appears (it should start with the General pane automatically).			
4	From the Startup Object drop-down list, select Sub Main.			
5	In the <i>Project Name</i> text box, type the name you want to appear in MetaMorph when selecting a user program to run.			
6	For <i>Project Description</i> , type the description that you want to appear in MetaMorph when selecting a user program to run.			
	Note: The project description is not currently recognized by MetaMorph when using VB 5.0/6.0. MetaMorph will display the <i>Project Name</i> in the Run User Program dialog box.			
7	Choose OK.			
8	From the View menu, choose Properties Window.			

Table 1.3Creating a User Program with Visual Basic 5.0 or 6.0

1.5 Creating a User Program, continued

Creating a user program with Visual Basic 5.0 or 6.0 or 6.0 (continued)

Step	Action				
9	In the Properties window that appears,				
	 Set Instancing to 5 – MultiUse, and Set (Name) to UserMethods. 				
10	From the Project menu, choose References. Then find <i>MetaMorph Type Library</i> in the list and select its check box. If you don't see <i>MetaMorph Type Library</i> , choose the <i>Browse</i> button, select Files of Type Executable, navigate to your MetaMorph directory, and select MMAPP.EXE.				
11	Once <i>MetaMorph Type Library</i> has been selected, you can choose Object Browser from the View menu and select <i>MMAppLib</i> from the top drop-down list. Then select <i>UserCall</i> from the Classes list and you will get a list of all the available functions in MetaMorph that you can call with the MM variable, along with their parameters. When you start to type the name of the object in the module window, a list will appear on the screen with all of the available functions for that object, along with their parameters.				
12	In the Class Module UserMethods section, insert the code shown in Fig. 1.3, which follows this procedure.				
13	From the Project menu, choose Add Module.				
14	In the module you just inserted, add the following code: Sub Main() End Sub				

Compiling the user program

That's the framework for your program. You can now add your personal code to the program, as described in Section 1.4. When you have finished, you are ready to compile your program. From the File menu, choose Make YourProjectName. You can select a name and location for your compiled program, but note that the name you select in this dialog will not be the name that you will use to reference your program in MetaMorph. In MetaMorph, your program will be referenced by the name or description you entered in the *Project Name* text box (Step 5 in table 1.2). Once the program has been compiled without any errors, it is ready to be run from MetaMorph.

1.5 Creating a User Program, continued

```
Class Module UserMethods Section Code
Figure 1.3
Option
         explicit
Public
         gUserID As Long
Public gParentWnd As Long
Public MM As MMAppLib.UserCall
Public
        Function Startup(cmdLine As String) As Long
End Function
Public
         Function DoCommand(cmdLine As String) As Long
End Function
Public Function Shutdown() As Long
End Function
```

1.6 Data Types and Arrays

Introduction Visual Basic functions in MetaMorph frequently make use of arrays to handle image pixel data, such as edgelist coordinates and intensity or color values. There is a complex interplay between the size of an array, the image bit-depth, and the type of data (Byte, Integer, or Long) being passed. The number of elements in the array and the range of data that can be stored will be determined by both the bit-depth of the image (1, 8, 16, or 24) and by the data type.

Data Types
and ArraysThe following table indicates the number of elements in the array as a function of the
image depth, data depth, and data type.

Image Bit-Depth	Data Bit-Depth	Data Type	# Elements in Array
1	1	Byte	number of pixels / 8
		Integer	number of pixels / 16
		Long	number of pixels / 32
	8	Byte	number of pixels ¹
	16	Integer	number of pixels ²
1	24	Long	number of pixels ³
8	1	same as for 1-bit images ⁴	
	8	Byte	number of pixels
	16	Integer	number of pixels
	24	Long	number of pixels ³
16	1	same as for 1-bit images ⁴	
	8	Byte	number of pixels ⁵
	16	Integer	number of pixels
	24	Long	number of pixels ⁶

Table 1.3Data Types and Arrays

Data Types and Arrays

(continued)

Image Bit-Depth	Data Bit-Depth	Data Type	# Elements in Array
24	1	same as for	1-bit images ⁴
	8	Byte	number of pixels ⁷
	16	Integer	number of pixels ⁷
	24	Long	number of pixels

¹ Pixel values will be converted to 0 or 255.

² Pixel values will be converted to 0 or 65535.

³ Each pixel's values will be packed into three bytes, but each of the 32-bit elements in the array will contain four bytes of data such that element one will contain the three values (red, green, blue) of pixel one and the first value of pixel two, the second element will contain the next two values for pixel two and the first two values of pixel three, and so on. Data from 1-bit (binary) images will be expressed as values of 0,0,0 or 255,255,255. Data from 8-bit images will be expressed as a triplet of the pixel value (Value, Value, Value). Data from 16-bit images will be expressed as a triplet of the low byte value of the pixel (Low,Low,Low).

⁵ Only the low byte of each pixel will be stored.

⁶ Only the low byte from each of the three values (red, green, blue) will be stored.

⁷ Only the intensity value will be stored. This value will be a numeric average of the red, green, and blue intensities.

Chapter 2 – Performing Serial and Digital I/O Communication

2.1 Overview

Introduction	Peripheral devices can be controlled from within MetaMorph by a number of user program functions. This chapter describes two types of functions used for communication with a peripheral device—those that send and receive data streams over a serial port and those that control a digital device.		
In this chapter	This chapter contains the following topics:		
	Торіс	See Page	
	Performing Serial Data Transmission	21	
	Communicating with a Digital I/O Device	26	

Introduction	You can control some devices such as VCRs by sending and receiving sequential
	streams of data over a serial port. A pair of Visual Basic functions, SendSerialData
	and WaitForSerialData, can perform the sending and receiving procedures. You will
	need to install the CUSTOMIO drop-in with the MetaMorph Drop-in Manager and to
	install and configure a Data Stream MetaDevice before you can take advantage of
	these two functions. This section describes the two functions, and provides a list of
	the syntax rules and a table of the serial command codes that are used for serial
	communication.

SendSerialData

Description	Carries out the Custom I/O: Send Serial Data command, which sends a sequential stream of data from the computer to another device via a serial port.		
Syntax	SendSerialData(sData As String, lTimeout As Long, bSendWithEcho As Boolean) As Long		
Remarks	This function is a shortcut to the "Send Serial Data" function of the MetaMorph CUSTOMIO drop-in, which provides control over devices that use Data Stream MetaDevices. SendSerialData runs without displaying a dialog box. The CUSTOMIO drop-in must be loaded for this function to run.		
Parameters	sData Gives the string to be sent to the serial port.		
	<i>lTimeout</i> Specifies a maximum time, in seconds, that MetaMorph should wait if no echo has been returned before continuing.		
	<i>bSendWithEcho</i> Determines whether or not MetaMorph is to wait for an echo from the serial device before sending the next character and to warn you if the character is not received. If <i>bSendWithEcho</i> is set to TRUE, MetaMorph will wait until the device returns an echo, or until <i>lTimeout</i> seconds have transpired, whichever comes first. If <i>bSendWithEcho</i> is set to FALSE, MetaMorph will return immediately.		
Example	' Send the string "TestString" to the serial port with a ' timeout of 5. Do not return until the string has been ' echoed or the timeout has expired. MM.SendSerialData "TestString", 5, TRUE		
See also:	WaitForSerialData		

WaitForSerialData

Description	Carries out the Custom I/O: Wait for Serial Data command, which waits for a sequential stream of data from another device by way of a serial port.			
Syntax	WaitForSerialData(sData As String, lTimeout As Long, bLogData As Boolean, sLogFormat As String) As Long			
Remarks	This function is a shortcut to the "Wait for Serial Data" function of the MetaMorph CUSTOMIO drop-in, which provides control over devices that use Data Stream MetaDevices. It runs without displaying a dialog. The CUSTOMIO drop-in must be loaded for this function to run.			
Parameters	sData Specifies the string from the serial port for which to wait.			
	<i>lTimeout</i> Specifies the number of seconds WaitForSerialData will wait for the string before returning.			
	<i>bLogData</i> Determines whether or not the received string will be written to a log file. If <i>bLogData</i> is set to TRUE and the string that is received matches the one specified in <i>sData</i> , the message that you specify with <i>sLogFormat</i> will be written to the log file (assuming it is open).			
	<i>sLogFormat</i> Specifies a message that will be sent to an open log file if the string that is received matches the one specified in <i>sData</i> and <i>bLogData</i> has been set to TRUE.			
Example	' Wait for 10 seconds for the string "ok" from the serial port. ' Don't log it to the data file. <i>MM.</i> WaitForSerialData "ok", 10, FALSE, ""			
See also:	SendSerialData			

2.2 Performing Serial Data Transmission, continued

Syntax rules for serial data transmission Table 2.1 describes the syntax rules for the command codes used in serial communication. Be sure to consult Table 2.2 for the codes themselves.

Code	Result				
\$	"Escape" character, which is ASCII 027 in decimal.				
^A thru ^Z	"Control" character, which is ASCII 001 for ^A through ASCII 26 for ^Z.				
/c	Sends the character after the slash. In this example, the character "c" would be sent. Useful for sending \land , \backslash , or \$ characters.				
\ddd	Sends ASCII digits in decimal. EXAMPLE: \192.				
\xdd	Sends ASCII digits in hexadecimal. EXAMPLE: \x27.				
(dddd)	Delays for specified number of milliseconds. EXAMPLE: (1000)				

Table 2.1Syntax Rules for Serial Data Transmission

Performing Serial Data Transmission, continued 2.2

ASCII control The following table provides the code strings used in serial communication.

Hex	Dec	Key	Name	Description
00	0	^@	NUL	Null
01	1	^A	SOH	Start of Header
02	2	^B	STX	Start of Text
03	3	^C	ETX	End of Text
04	4	^D	EOT	End of Transmission
05	5	^E	ENQ	Inquiry
06	6	^F	ACK	Acknowledge
07	7	^G	BEL	Bell
08	8	^H	BS	Backspace
09	9	~	HT	Horizontal Tab
0A	10	~7	LF	Line Feed
0B	11	٨K	VT	Vertical Tab
0C	12	۸L	FF	Form Feed
0D	13	^M	CR	Carriage Return
0E	14	^N	SO	Shift Out
0F	15	^O	SI	Shift In
10	16	۸P	DLE	Data Link Escape
11	17	^Q	DC1	Device Control 1
12	18	^R	DC2	Device Control 2
13	19	^S	DC3	Device Control 3
14	20	^T	DC4	Device Control 4
15	21	^U	NAK	Negative Acknowledge
16	22	^V	SYN	Synchronous Idle

Continued on next page

codes

2.2 Performing Serial Data Transmission, continued

ASCII control codes

(continued)

Hex	Dec	Key	Name	Description
17	23	٨W	ЕТВ	End Transmission Block
18	24	^X	CAN	Cancel
19	25	۸Y	EM	End of Medium
1A	26	^Z	SUB	Substitute
1B	27		ESC	Escape
1C	28		FS	File Separator
1D	29		GS	Group Separator
1E	30		RS	Record Separator
1F	31		US	Unit Separator

2.3 Communicating with a Digital I/O Device

Introduction	MetaMorph allows you to control digital devices by sending and receiving TTL-level voltage signals through a parallel port or with the use of a digital I/O board. To do so with a user program, you will need to install the CUSTOMIO drop-in with the MetaMorph Drop-in Manager and to install and configure a Digital I/O MetaDevice.
DICotEirot	
DIGerrist	
Description	Obtains the handle of the first Digital I/O MetaDevice in the current list of Digital I/O MetaDevices.
Syntax	DIGetFirst(hRetDevice As Long) As Long
Return values	<i>hRetDevice</i> Returns the handle of the first Digital I/O MetaDevice in the current list. If there are no Digital I/O MetaDevices, <i>hRetDevice</i> will return a value of -1.
Example	' 'dev' will hold the device handle Dim <i>dev</i> As Long
	' get the handle of the first device and put it in 'dev' MM.DIGetFirst dev
See also:	DIGetNext

DIGetIOStatus

Description	Determines whether a specified I/O line of the given device is an input line or an output line.
Syntax	DIGetIOStatus (<i>hDevice</i> As Long, <i>nLineNumber</i> As Integer, <i>nRetInOut</i> As Integer) As Long
Parameters	<i>hDevice</i> Specifies the handle of the device.
	<i>nLineNumber</i> Specifies the number of the I/O line in question. This number will correspond to the pin number on the line's connector.
Return values	<i>nRetInOut</i> Returns a value corresponding to whether the specified line is an input line or an output line. If it is an input line, a value of 0 will be returned. If it is an output line, a value of 1 will be returned.

2.3 Communicating with a Digital I/O Device, continued

DIGetIOStatus (continued)	
Example	Dim <i>dev</i> As Long Dim <i>inout</i> As Integer
	' Get the IO status of line 2 of a device and place it in ' 'inout'. 'dev' must have been previously obtained using ' DIGetFirst or DIGetNext <i>MM</i> . DIGetIOStatus <i>dev</i> , 2, <i>inout</i>
See also:	DIGetFirst, DIGetLineCount, DIGetNext

DIGetLineCount

Description	Obtains the number of I/O lines for a given device.
Syntax	DIGetLineCount(hDevice As Long, nRetLines As Integer) As Long
Parameters	<i>hDevice</i> Specifies the handle of the device.
Return values	<i>nRetLines</i> Returns the number of I/O lines.
Example	Dim <i>dev</i> As Long Dim <i>numLines</i> As Integer
	' Get the number of lines on a device and place it in ' 'numlines'. 'dev' must have been previously obtained using ' DIGetFirst or DIGetNext <i>MM.</i> DIGetLineCount <i>dev</i> , <i>numLines</i>
See also:	DIGetFirst, DIGetNext

DIGetLineState

Description	Obtains the state of a specified I/O line (Low vs. High).
Syntax	DIGetLineState (<i>hDevice</i> As Long, <i>nLineNumber</i> As Integer, <i>nRetState</i> As Integer) As Long
Parameters	<i>hDevice</i> Specifies the handle of the device. <i>nLineNumber</i> Specifies the number of the I/O line in question. This number will correspond to the pin number on the line's connector.
Return values	<i>nRetState</i> Returns a value corresponding to the state of the I/O line. If the line is Low, a value of 0 will be returned. If the line is High, a value of 1 will be returned.
Example	Dim <i>dev</i> As Long Dim <i>state</i> As Integer
	' Get the line state of line 1 of a device and place it in ' 'state'. 'dev' must have been previously obtained using ' DIGetFirst or DIGetNext . MM. DIGetLineState dev, 1, state
See also:	DIGetFirst, DIGetNext

DIGetName

Description	Obtains the name of the MetaDevice whose handle you pass to it.
Syntax	DIGetName(hDevice As Long, sDeviceName As String) As Long
Parameters	<i>hDevice</i> Specifies the handle of the device.
Return values	<i>sDeviceName</i> Returns the name of the selected MetaDevice. This will be given as a text string.
Example	Dim <i>dev</i> As Long Dim <i>name</i> As String
	' Get the name of a device and place it in 'name'. 'dev' must ' have been previously obtained using DIGetFirst or DIGetNext . MM. DIGetName dev, name
See also:	DIGetFirst, DIGetNext

2.3 Communicating with a Digital I/O Device, continued

DIGetNext

Description	Obtains the handle of the Digital I/O MetaDevice that follows the last one that was read (using either DIGetNext or DIGetFirst).
Syntax	DIGetNext(hRetDevice As Long) As Long
Return values	<i>hRetDevice</i> Returns the handle of the Digital I/O MetaDevice that follows the last one read. If there are no more Digital I/O MetaDevices, <i>hRetDevice</i> will return a value of -1.
Example	Dim dev As Long
	' Get the third device and place it in 'dev'.
	<pre>' Get the first device MM.DIGetFirst dev ' Get the device after the first device (the second device). MM.DIGetNext dev</pre>
	' Get the device after the second device. MM. DIGetNext dev
See also:	DIGetFirst

SetDigitalIO

Description	Carries out the Custom I/O: Set Digital I/O command, which controls a peripheral digital I/O device by sending signals to it from MetaMorph.
Syntax	SetDigitalIO(lBitsToSet As Long, lState As Long) As Long
Remarks	This function is a shortcut to the "Set Digital I/O" function of the MetaMorph CUSTOMIO drop-in. It runs without displaying a dialog box. The CUSTOMIO drop-in must be loaded for this function to run.
Parameters	<i>lBitsToSet</i> Provides a 32-bit binary bit pattern that indicates which digital I/O lines will be active. For example, if bit 3 of <i>lBitsToSet</i> is set to 1, digital I/O line 3 will be enabled to write to the device. If it is set to 0, line 3 will be disabled from writing.
	<i>lState</i> Provides a 32-bit binary bit pattern that sets the state (High vs. Low) of each I/O line identified by the bit pattern in <i>lBitsToSet</i> . To continue the preceding example, if bit 3 of <i>lState</i> is set to 0 and <i>lBitsToSet</i> has assigned a value of 1 to bit 3, this will set the line state for bit 3 to Low. If <i>lState</i> assigns a value of 1 to bit 3, this will set the line state for bit 3 to High. If <i>lBitsToSet</i> has assigned a value of 0 to bit 3, the setting of bit 3 for <i>lState</i> will be ignored, and the line state of bit 3 will be unaffected.

SetDigitalIO (continued)	
Example	' Set bits 3 and 4 of the DIO lines High. 24 in binary is ' 00011000 and 255 is 11111111. Since the <i>lState</i> argument had ' ones in all the first 8 lines, it will set to High whichever ' of those lines were selected by the <i>lBitsToSet</i> argument. <i>MM.SetDigitalIO</i> 24, 255
See also:	WaitForDigitalIO

2.3 Communicating with a Digital I/O Device, continued

WaitForDigitalIO

Description	Carries out the Custom I/O: Wait for Digital I/O command, which configures MetaMorph to wait for specific signals from a peripheral digital I/O device.
Syntax	WaitForDigitalIO(<i>lBitsToRead</i> As Long, <i>lState</i> As Long, <i>lMilliseconds</i> As Long) As Long
Remarks	This function is a shortcut to the "Wait for Digital I/O" function of the MetaMorph CUSTOMIO drop-in. It runs without displaying a dialog. The CUSTOMIO drop-in must be loaded for this function to run.
Parameters	<i>lBitsToRead</i> Provides a 32-bit binary bit pattern that indicates which bits in <i>lState</i> will be compared to data read from the I/O lines. For example, if bit 3 of <i>lBitsToRead</i> is set to 1, WaitForDigitalIO will not return until either digital I/O line 3 matches the state (High vs. Low) that has been set for bit 3 by <i>lState</i> , or until <i>lMilliseconds</i> have elapsed—whichever occurs first.
	<i>lState</i> Provides a 32-bit binary bit pattern against which the state (High vs. Low) of each I/O line identified by the bit pattern in <i>lBitsToRead</i> will be compared. To continue the preceding example, if bit 3 of <i>lState</i> is set to 0 and <i>lBitsToRead</i> has assigned a value of 1 to bit 3, this will set the condition to wait until line 3 switches to Low. If <i>lState</i> assigns a value of 1 to bit 3, this will set the condition to wait for line 3 to switch to High.
	<i>Milliseconds</i> Specifies the number of milliseconds to wait before timing out.
Example	' Wait for bits 3, 4, and 7 to be High, Low, and Low, ' respectively. Timeout if this takes longer than 500 ' milliseconds. 98 is 10011000 in binary, and 128 is 10000000. MM.WaitForDigitalIO 98, 128, 500
See also:	SetDigitalIO

Chapter 3 – Executing Commands and Journals

3.1 Overview

Introduction

Fundamental to the successful deployment of user programs is the ability to set variables and pass parameters to the functions, and then run the functions. Occasionally, you will need to determine a function's handle. You may also want to use message boxes to provide feedback while running or troubleshooting the program. This chapter describes the functions you will need for performing all of these procedures.



The parts of the interface that are accessed through the **RunFunction**, **RunFunctionEx**, and **SetFunctionVariable** functions are subject to change. Care should be taken when considering whether to use them. These functions should be used only when there is no alternative way to accomplish your task. Because of this, and because of the great number of functions and variables, not all of the functions and variables that can be accessed through the **RunFunction** and **SetFunctionVariable** functions have been documented. Information for individual functions and their variables will be provided by e-mail on a case-by-case basis. Send inquiries to: **support.dtn@moldev.com**.

In this chapter This chapter contains the following topics: Topic See Page Executing Commands and Journals 32
GetFunctionHandle

Description	Obtains the handle of a specified function.		
Syntax	GetFunctionHandle(sName As String, lRetFunctionHandle As Long) As Long		
Remarks	You can use the function handles returned by this function when calling RunFunctionEx and SetFunctionVariable .		
Parameters	sName Supplies the name of the function.		
Return values	<i>lRetFunctionHandle</i> Returns the function handle.		
Example	<pre>' Run the MetaMorph Deinterlace function on the image 'src'. ' Put the odd and even images in 'destOdd' and 'destEven'. ' 'src', 'destOdd', and 'destEven' should have been previously ' set to valid existing images. Dim deinterHandle As Long MM.GetFunctionHandle "Deinterlace", deinterHandle MM.SetFunctionVariable deinterHandle, "imSource", src MM.SetFunctionVariable deinterHandle, "imDestOdd", destOdd MM.SetFunctionVariable deinterHandle, "imDestEven", destEven MM.RunFunctionEx deinterHandle, 1</pre>		
	' Run Deinterlace again, this time using the current desktop ' image as the source. This time, use RunFunction to run the ' function instead of RunFunctionEx . MM.GetCurrentImage src MM.RunFunction "Deinterlace", 1		
See also:	RunFunctionEx, SetFunctionVariable		

PrintMsg

Description	Displays a configured message in a diagnostic window in MetaMorph.		
Syntax	PrintMsg(sMessage As String) As Long		
Remarks	This function is useful for the purposes of debugging. You can print diagnostics in a message box without pausing execution. The SetPrintMsgSizeAndPosition function allows you to set properties of the window in which the messages appear.		

3.2 Executing Commands and Journals, continued

PrintMsg (continued)	
Parameters	sMessage Specifies the message that is to appear in the message box.
Example	' Print the string "Hello" in the print message window. Place ' the window in the upper left corner of the screen and ' make it 100 pixels wide and 30 high. MM.PrintMsg "Hello" MM.SetPrintMsgWindowPositionAndSize 1, 1, 100, 30
See also:	SetPrintMsgSizeAndPosition

RunFunction

Description	Carries out the function named by the function name <i>sFunctionName</i> as if you had chosen it from a menu.		
Syntax	RunFunction(sFunctionName As String, nMode As Integer) As Long		
Remarks	This function differs from RunFunctionEx in its use of the function's name. RunFunctionEx uses the function's handle.		
Parameters	<i>sFunctionName</i> Specifies the name of the function to be run. <i>nMode</i> Determines the execution mode. If <i>nMode</i> is 0, the function will run normally. If <i>nMode</i> is 1, the function will run in journal playback mode—for most functions this means no user interface will be displayed.		
Example	' Run the MetaMorph Deinterlace function on the image 'src'. MM.GetCurrentImage src MM.RunFunction "Deinterlace", 1		

RunFunctionEx

Description	Carries out the function named by the function handle <i>lFunctionHandle</i> as if you had chosen it from the menu.		
Syntax	RunFunctionEx(lFunctionHandle As Long, nMode As Integer) As Long		
Remarks	This function differs from RunFunction in its use of the function's handle. RunFunction uses the function's name. Function handles are obtained for use with RunFunctionEx by calling GetFunctionHandle .		
Parameters	<i>lFunctionHandle</i> Specifies the handle of the function to be run.		
	nMode Determines the execution mode. If $nMode$ is 0, the function will run normally. If $nMode$ is 1, the function will run in journal playback mode—for most functions this means no user interface will be displayed.		
Example	<pre>' Run the MetaMorph Deinterlace function on the image 'src'. ' Put the odd and even images in 'destOdd' and 'destEven'. ' 'src', 'destOdd', and 'destEven' should have been previously ' set to valid existing images. Dim deinterHandle As Long MM.GetFunctionHandle "Deinterlace", deinterHandle MM.SetFunctionVariable deinterHandle, "imSource", src MM.SetFunctionVariable deinterHandle, "imDestOdd", destOdd MM.SetFunctionVariable deinterHandle, "imDestEven", destEven MM.RunFunctionEx deinterHandle, 1</pre>		
See also:	GetFunctionHandle		

3.2 Executing Commands and Journals, continued

RunJournal

Description	Runs the specified journal.		
Syntax	RunJournal(sPath As String) As Long		
Remarks	<i>sPath</i> should contain the full path of the journal along with the extension, ".jnl".		
Parameters	<i>sPath</i> Specifies the path and name of the journal to be run.		
Example	' Run a journal named 'closeall.jnl' that resides in the ' directory C:\MM\Journals. MM.RunJournal "c:\mm\journals\closeall.jnl"		

SetFunctionVariable

Description	Sets the value of a variable in a MetaMorph function.			
Syntax	SetFunctionVariable (<i>lFunctionHandle</i> As Long, <i>sVariableName</i> As String, <i>value</i> As Variant) As Long			
Remarks	You can obtain the handle of the function whose variable you want to set by using GetFunctionHandle .			
Parameters	<i>lFunctionHandle</i> Gives the handle of the function whose variable you want to set. <i>sVariableName</i> Gives the name of the function's variable that you want to set. <i>value</i> Specifies the new value of the variable—this must be of the same data type as the variable you are setting.			

3.2 Executing Commands and Journals, continued

SetFunctionVariable

(continued)

Example	<pre>' Run the MetaMorph Deinterlace function on the image 'src'. ' Put the odd and even images in 'destOdd' and 'destEven'. ' 'src', 'destOdd', and 'destEven' should have been previously ' set to valid existing images. Dim deinterHandle As Long MM.GetFunctionHandle "Deinterlace", deinterHandle MM.SetFunctionVariable deinterHandle, "imSource", src MM.SetFunctionVariable deinterHandle, "imDestOdd", destOdd MM.SetFunctionVariable deinterHandle, "imDestEven", destEven MM.RunFunctionEx deinterHandle, 1</pre>
	<pre>' Run Deinterlace again, this time using the current desktop ' image as the source. This time, use RunFunction to run the ' function instead of RunFunctionEx. MM.GetCurrentImage src MM.RunFunction "Deinterlace", 1 CetFunctionHandle</pre>
See also:	GetFunctionHandle

SetPrintMsgWindowPositionAndSize

Description	Sets the position and size of the window used by PrintMsg to display messages.	
Syntax	SetPrintMsgWindowPositionAndSize (<i>nXPos</i> As Integer, <i>nYPos</i> As Integer, <i>nXSize</i> As Integer, <i>nYSize</i> As Integer) As Long	
Parameters	<i>nXPos</i> Specifies the X-coordinate for placement of the upper left corner of the message box.	
	<i>nYPos</i> Specifies the Y-coordinate for placement of the upper left corner of the message box.	
	nXSize Specifies the width of the message box, in pixels.	
	<i>nYSize</i> Specifies the height of the message box, in pixels.	
Example	<pre>' Print the string "Hello" in the print message window. Place ' the window in the upper left corner of the screen and ' make it 100 pixels wide and 30 high. MM.PrintMsg "Hello" MM.SetPrintMsgWindowPositionAndSize 1, 1, 100, 30</pre>	
See also:	PrintMsg	

SetMMVariable

Description	Sets the value of a MetaMorph custom variable or certain writable built in variables.
Syntax	SetMMVariable VariableName, value
Parameters	VariableName can be any existing or undefined custom variable, or a built in variable name, such as <i>\$Camera.Digital.Exposure\$</i> . If it is an undefined custom variable, that variable will be created. <i>value</i> can be a string or a number. If it is not the same type as an existing variable, it will not be set.
Examples	<pre>mm.SetMMVariable "MyComment", "Experiment number 7" mm.SetMMVariable "\$Camera.Digital.Exposure\$", 1250</pre>
See also:	GetMMVariable

GetMMVariable

Description	Gets the value of an existing MetaMorph custom or built in variable.	
Syntax	GetMMVariable VariableName, var	
Parameters	<i>VariableName</i> can be any existing custom variable or a built in variable. <i>var</i> is a VB variable. If the type of <i>var</i> does not match the type of the MetaMorph variable, it will try to convert the type before it assigns it. So, for example, the number 123 assigned to a string will end up as "123".	
Examples	<pre>Dim planeNum as Integer mm.GetMMVariable "\$Image.ActivePlane\$", planeNum Dim comment as String mm.GetMMVariable "MyComment", comment</pre>	
See also:	SetMMVariable	

Chapter 4 – Reading and Manipulating Images and Image Windows

4.1 Overview

Introduction	This chapter deals with the functions that you need for managing image windows and their properties. These functions will be central to any user program that you create.		
In this chapter	This chapter contains the following topics:		
	Торіс	See Page	
	Loading, Creating, Copying, and Closing Images	39	
	Finding Loaded Images	46	
	Manipulating Image Windows	49	
	Reading and Using Image Properties	53	

4.2 Loading, Creating, Copying, and Closing Images

Introduction	Image loading, copying, saving, and closing are vital steps in any user program that
	you will run in MetaMorph. This section describes the functions that are used for
	these procedures. In particular, you will make frequent use of the LoadImage,
	SaveImage, and CloseImage functions.

CloneImage

Description	Creates a copy of a specified image.		
Syntax	CloneImage(hImage As Long, hRetNewImage As Long) As Long		
Remarks	This function differs from CopyImage in that CopyImage requires the existence of a destination image that will be overwritten. The information that will be copied by CloneImage includes the image data, the annotation, zoom factor, calibration setting, thresholding, position, wavelength, scaling, and image times.		
Parameters	<i>hImage</i> Specifies the handle of the image to be copied.		
Return values	hRetNewImage Returns the handle for the new image.		
Example	' Create a copy of the current desktop image and place it in ' 'dest' Dim <i>src</i> As Long, <i>dest</i> As Long <i>MM</i> .GetCurrentImage <i>src</i> <i>MM</i> .CloneImage <i>src</i> , <i>dest</i>		
See also:	CopyImage, GetCurrentImage (Section 4.3), GetImage (Section 4.3)		

4.2 Loading, Creating, Copying, and Closing Images, continued

CloseImage

Description	Closes a specified image and removes it from the desktop.		
Syntax	CloseImage(hImage As Long) As Long		
Remarks	If the image contains unsaved information, you will be prompted before the image is closed. In this regard, CloseImage differs from ForceCloseImage , which will close a modified image even if it has not been saved first. This function does not affect any disk-based copies of the image.		
Parameters	hImage Specifies the handle of the image to be closed.		
Example	' Close the current desktop image Dim <i>src</i> As Long <i>MM</i> .GetCurrentImage <i>src</i> <i>MM</i> .CloseImage <i>src</i>		
See also:	ForceCloseImage, GetCurrentImage (Section 4.3), GetImage (Section 4.3)		

CopyImage	
Description	Copies a source image, hSourceImage, over a destination image, hDestImage.
Syntax	CopyImage(hSourceImage As Long, hDestImage As Long) As Long
Remarks	This function differs from CloneImage in that CopyImage requires the existence of a destination image, <i>hDestImage</i> , that will be overwritten. Only the image data, annotation, Z-position, wavelength, and image times are copied.
Parameters	<i>hSourceImage</i> Specifies the handle of the source image to be copied.<i>hDestImage</i> Specifies the handle of the destination image to be overwritten.
Example	' Copy the current desktop image to an existing image named ' "Temp" Dim src As Long, dest As Long MM.GetCurrentImage src MM.GetNamedImage "Temp", dest MM.CopyImage src, dest
See also:	CloneImage, GetCurrentImage (Section 4.3), GetImage (Section 4.3)

CopyImagePlane

Description	Copies a specified source image plane to the current plane in a specified destination image.	
Syntax	CopyImagePlane (<i>hSourceImage</i> As Long, <i>hDestImage</i> As Long, <i>nSrcPlane</i> As Integer) As Long	
Remarks	Plane numbering starts at 0. The information that will be copied includes the image data, the annotation, zoom factor, calibration setting, thresholding, position, wavelength, scaling, and image times.	
Parameters	hSourceImage Specifies the handle of the source image.	
	<i>hDestImage</i> Specifies the handle of the destination image into which the plane will be copied.	
	<i>nSrcPlane</i> Specifies the number of the plane that will be copied.	
Example	' Copy the first plane of the current image to the current ' plane of an existing image named "Stack". Dim <i>src</i> As Long, <i>dest</i> As Long <i>MM</i> .GetCurrentImage <i>src</i> <i>MM</i> .GetNamedImage "Stack", <i>dest</i> <i>MM</i> .CopyImagePlane <i>src</i> , <i>dest</i> , 0	
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3)	

CreateImage

Description	Creates a new image, using a specified name, size, and bit-depth.	
Syntax	CreateImage (<i>width</i> As Integer, <i>height</i> As Integer, <i>depth</i> As Integer, <i>name</i> As String, <i>hRetNewImage</i> As Long) As Long	
Remarks	The SetDisplayedImagesWhenCreated function sets whether or not images will be automatically drawn on the screen when they are created. If you set its <i>bState</i> parameter to FALSE, the image will not appear on the screen when CreateImage is called, although the image will still be created. This may be useful for performing intermediate work on a temporary work image. In this case, you will need to use ShowImage to force the undisplayed image to be displayed.	

4.2 Loading, Creating, Copying, and Closing Images, continued

CreateImage (continued)		
Parameters	<i>width</i> Specifies a width (X-axis size), in pixels, for the new image.	
	<i>depth</i> Specifies a bit-depth for the new image. This must be 1, 8, 16, or 24.	
	<i>name</i> Specifies a name for the new image. If an image with that name already exists, "-1" will be appended to the name of the new image.	
Return values	hRetNewImage Returns the handle of the new image.	
Example	' Create a 512 x 480 pixel 8-bit image named "myimage" and ' place it in 'im'. Dim <i>im</i> As Long <i>MM</i> . CreateImage 512, 480, 8, "myimage", <i>im</i>	
See also:	SetDisplayImagesWhenCreated, ShowImage	

ForceCloseImage

Description	Closes a specified image and removes it from the desktop.	
Syntax	ForceCloseImage(hImage As Long) As Long	
Remarks	If the image has not been saved, it will be deleted anyway, and you will lose any unsaved information. In this regard, ForceCloseImage differs from CloseImage , which will prompt you to save a modified image before closing it. This function does not affect any disk-based copies of the image.	
Parameters	<i>hImage</i> Specifies the handle of the image to be closed.	
Example	<pre>' Close the current desktop image even if it has not yet been ' saved to disk. Dim im As Long MM.GetCurrentImage im MM.ForceCloseImage im</pre>	
See also:	CloseImage, GetCurrentImage (Section 4.3), GetImage (Section 4.3)	

LoadImage

Description	Loads a specified image file.	
Syntax	LoadImage(sFileName As String, hRetImage As Long) As Long	
Parameters	<i>sFileName</i> Specifies the name of the image to be loaded. This should contain the complete path of the image.	
Return values	hRetImage Returns the handle of the loaded image.	
Example	<pre>' Load the image stack Nerve.stk which is stored in the ' directory C:\MM\Images, and place it in nerveIm. MM.LoadImage "c:\mm\images\nerve.stk", nerveIm</pre>	
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3)	

SaveImage

Description	Saves the specif	fied image to a file on the hard disk.
Syntax	SaveImage(<i>hIm</i> As Integer) As I	nage As Long, sFileName As String, bPartial As Boolean, nFormat Long
Remarks	<i>nFormat</i> is a nusaved. This value	Immeric value which denotes the format in which the image is to be be may be any of the following: Image-1 File Type (*.img) Windows File Type (*.bmp) TIFF File Type (*.tif) Stack File Type (*.stk)
	5	MRC-500 File Type (*.ptc) Photometrics CC200 File Type (*.cc2)
	7	Hamamatsu File Type (*.ham)
	8	RGB TIFF File Type (*.tif)
	9	WinView File Type (*.spe)
	10	Argus File Type (*.ham)

4.2 Loading, Creating, Copying, and Closing Images, continued

SaveImage (continued)	
Parameters	hImage Specifies the handle of the image to be saved.
	<i>sFileName</i> Gives the name to be used for the image being saved. This should include the full path.
	<i>bPartial</i> For an image with an active region of interest, this Boolean logic variable specifies whether only the image inside the region is to be saved (TRUE) or if the whole image is to be saved regardless of the presence of an active region (FALSE). If there is no active region, the whole image will be saved.
	<i>nFormat</i> Supplies a numeric value that indicates what format the image is to be saved. (See Remarks.)
Example	<pre>' Save the current desktop image as a TIFF file named ' Newimage.tif in the directory C:\MM\Images. Dim currentIm As Long MM.GetCurrentImage currentIm MM.SaveImage currentIm, "c:\mm\images\newimage.tif", FALSE, 3</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3)

SetDisplayImagesWhenCreated

Description	Configures whether or not images are to be drawn automatically on the screen when they are created.
Syntax	SetDisplayImagesWhenCreated(bState As Boolean) As Long
Remarks	SetDisplayedImagesWhenCreated sets Use ShowImage to force an undisplayed image to be displayed.
Parameters	<i>bState</i> This Boolean logic variable specifies whether created images are to be displayed (TRUE) or kept hidden (FALSE). If <i>bState</i> is set to FALSE when CreateImage or other functions are called which create an image, the image will be created, but will not appear on the screen. This is useful for performing intermediate work on a temporary work image which you don't necessarily want to see. If <i>bState</i> is set to TRUE, new images will be displayed upon creation.

4.2 Loading, Creating, Copying, and Closing Images, continued

```
SetDisplayImagesWhenCreated
```

(continued)

Example	<pre>' Create an image and do some operations on it. Don't let the ' user see the image until all the operations are done. MM.SetDisplayImagesWhenCreated FALSE Dim im As Long MM.CreateImage 512, 480, 8, "temp", im</pre>
	' Do some operations on 'im' ' Cause 'im' to appear on the screen <i>MM.ShowImage im,</i> 236
	' Reset state so images are once again displayed as soon as ' they are created <i>MM</i> . SetDisplayImagesWhenCreated TRUE

See also: CreateImage

ShowImage

Description	Causes the specified image to be displayed, if it is not already being displayed.
Syntax	ShowImage(hImage As Long, nPalEntries As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image to be displayed. <i>nPalEntries</i> Gives the number of palette entries to use in displaying the image: 2, 4, 8, 16, 32, 64, 128, or 236.
Example	' Create an image and do some operations on it. Don't let the ' user see the image until all the operations are done. MM.SetDisplayImagesWhenCreated FALSE Dim <i>im</i> As Long MM.CreateImage 512, 480, 8, "temp", <i>im</i>
	' Do some operations on 'im' ' Cause 'im' to appear on the screen MM.ShowImage im, 236
	' Reset state so images are once again displayed as soon as ' they are created MM.SetDisplayImagesWhenCreated TRUE
See also:	GetCurrentImage, GetImage

4.3 Finding Loaded Images

Introduction	The next set of functions are crucial for control of the procedures being carried out by
	your user program. These functions are used for finding an image's handle, the
	"indexing tag" that both Visual Basic and MetaMorph use to keep track of the image,
	as well as for determining how many images are loaded and whether a handle is
	currently in use. The GetCurrentImage and GetImage functions, which obtain an
	image's handle, are particularly important, and you will see these two functions
	referenced frequently in the "See Also" lists throughout this manual.

GetCurrentImage

Description	Returns the handle of the currently active image.
Syntax	GetCurrentImage(hRetImage As Long) As Long
Remarks	A handle is a unique number used by MetaMorph to keep track of the appropriate image while working with it. This number is used extensively as a variable by most of the MetaMorph Visual Basic functions in this manual. The current (active) image is the one at the front of all the loaded images.
Return values	<i>hRetImage</i> Returns the handle of the current image.
Example	' Get the current desktop image and put it in 'im' Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i>
See also:	GetImage

GetImage

Description	All images loaded in MetaMorph have a unique index number, from 0 to $(n - 1)$, where <i>n</i> is the number of loaded images. GetImage returns the handle of the image denoted by <i>nIndex</i> .
Syntax	GetImage(nIndex As Integer, hRetImage As Long) As Long
Remarks	The image handle is returned in <i>nRetImage</i> . If an error occurs (for example, if <i>nIndex</i> is not a valid number), <i>nRetImage</i> will contain 0.
Parameters	<i>nIndex</i> Gives the index number of the desired image. (See Description.) If an error occurs (for example, if <i>nIndex</i> is not a valid number), MetaMorph will return a handle of "0".

4.3 Finding Loaded Images, continued

See also:	GetCurrentImage
	<pre>For i = 0 To MM.GetNumberOfImages MM.GetImage i, im MM.GetImageName im, name MM.PrintMsg name Next i</pre>
Example	' Print out the names of all loaded images in the message ' window Dim <i>im</i> As Long Dim <i>i</i> As Integer Dim <i>nam</i> e As String
Return values	<i>hRetImage</i> Returns the handle of the specified image. If an error occurs (for example, if <i>hRetImage</i> is not a valid number), MetaMorph will return a handle of "0".
GetImage (continued)	

GetNumberOfImages

Description	Returns the number of images loaded in MetaMorph.
Syntax	GetNumberOfImages() As Long
Remarks	A value of 0 will be returned if there are no images or if an error occurred.
Return values	This function returns a value corresponding to the number of images that are currently loaded. If an error has occurred, or if there are no images currently loaded, a value of 0 will be returned.
Example	' Print out the names of all loaded images in the message ' window Dim <i>im</i> As Long Dim <i>i</i> As Integer Dim <i>name</i> As String
	<pre>For i = 0 To MM.GetNumberOfImages MM.GetImage i, im MM.GetImageName im, name MM.PrintMsg name Next i</pre>

4.3 Finding Loaded Images, continued

IsValidImage

Description	Tests whether or not the image handle you pass is valid.
Syntax	IsValidImage(hImage As Long) As Long
Parameters	<i>hImage</i> Gives the handle of the image being tested. If the handle value that you pass is not valid, IsValidImage will return a value of "0".
Return values	hImage Returns a value of 0 if the handle you pass is not valid.
Example	<pre>' Determine if the image handle 'im' denotes a valid image. ' 'im' is an image handle that was initialized previously. If MM.IsValidImage(im) Then MM.PrintMsg "image is valid" Else MM.PrintMsg "Image is not valid" End If</pre>
See also:	GetCurrentImage, GetImage

4.4 Manipulating Image Windows

Introduction This section describes the functions that are used for controlling the appearance of your image windows. The following functions are used for changing the size and position of image windows, as well as minimizing (shrinking to a desktop icon) and maximizing (restoring to full size).

GetImageWindowPosition

Description	Obtains the on-screen position of the specified image.
Syntax	GetImageWindowPosition(hImage As Long, nRetXPos As Integer, nRetYPos As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose position you want to know.
Return values	<i>nRetXPos</i> Gives the X-coordinate of the upper left corner of the image.<i>nRetYPos</i> Gives the Y-coordinate of the upper left corner of the image.
Example	' Print the size and position of the window holding the current image Dim <i>x</i> As Integer, <i>y</i> As Integer, <i>dx</i> As Integer, <i>dy</i> As Integer Dim <i>im</i> As Long
	<pre>MM.GetCurrentImage im MM.GetImageWindowSize im, dx, dy MM.GetImageWindowPosition im, x, y MM.PrintMsg "The current image is at " + Str(x) + ", " + Str(y) MM.PrintMsg "The size of the current image is " + Str(dx) + " X " + Str(dy)</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetImageWindowSize, SetImageWindowPosition

GetImageWindowSize

Description	Obtains the width and height of the specified image.
Syntax	GetImageWindowSize(hImage As Long, nRetXSize As Integer, nRetYSize As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose size you want to know.
	Continued on next page

MetaMorph

4.4 Manipulating Image Windows, continued

GetImageWindowSi (continued)	ze
Return values	<i>nRetXSize</i> Gives the width, in pixels, of the image.
	<i>nRetYSize</i> Gives the height, in pixels, of the image.
Example	' Print the size and position of the window holding the current ' image
	Dim x As Integer, y As Integer, dx As Integer, dy As Integer Dim im As Long
	MM.GetCurrentImage im MM.GetImageWindowSize im, dx, dy
	MM.GetImageWindowPosition im, x, y
	<pre>MM.PrintMsg "The current image is at " + Str(x) + ", " + Str(y) MM.PrintMsg "The size of the current image is " + Str(dx) + " X " + Str(dy)</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetImageWindowPosition, SetImageWindowSize

Description	Expands the window of the specified image to full-size. If the image has been minimized (shrunk to a desktop icon), it will expand the window to its normal size.
Syntax	MaximizeImageWindow(hImage As Long) As Long
Parameters	<i>hImage</i> Specifies the handle of the image you want to maximize.
Example	' Maximize the current image Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i> MM.MaximizeImageWindow <i>im</i>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), MinimizeImageWindow

Continued on next page

MaximizeImageWindow

MinimizeImageWindow

Description	Shrinks the window of the specified image to a desktop icon.
Syntax	MinimizeImageWindow(hImage As Long) As Long
Parameters	hImage Specifies the handle of the image you want to minimize.
Example	' Minimize the current image Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i> MM.MinimizeImageWindow <i>im</i>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), MaximizeImageWindow

SetImageWindowPosition

Description	Sets the position of the window for the specified image.
Syntax	SetImageWindowPosition(<i>hImage</i> As Long, <i>nXPos</i> As Integer, <i>nYPos</i> As Integer)
Remarks	This function dictates the position of the upper left corner of the image window.
Parameters	<i>hImage</i> Specifies the handle of the image whose position you want to set. <i>nXPos</i> Specifies the X-coordinate of the upper left corner of the image window. <i>nYPos</i> Specifies the Y-coordinate of the upper left corner of the image window.
Example	<pre>' Make the window holding the current image 100 X 100 and put ' it in the upper left corner of the screen Dim im As Long MM.GetCurrentImage im MM.SetImageWindowPosition im, 0, 0 MM.SetImageWindowSize im, 100, 100</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetImageWindowSize

SetImageWindowSize

Description	Sets the size of the window for the specified image.
Syntax	SetImageWindowSize(hImage As Long, nXSize As Integer, nYSize As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose size you want to set.<i>nXSize</i> Specifies the width, in pixels, of the image.<i>nYSize</i> Specifies the height, in pixels, of the image.
Example	' Make the window holding the current image 100 X 100 and put ' it in the upper left corner of the screen Dim <i>im</i> As Long <i>MM</i> .GetCurrentImage <i>im</i> <i>MM</i> .SetImageWindowPosition <i>im</i> , 0, 0 <i>MM</i> .SetImageWindowSize <i>im</i> , 100, 100
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetImageWindowPosition

4.5 Reading and Using Image Properties

Introduction It is frequently necessary or useful to change certain characteristics about an image, such as its dimensions, bit-depth, or name. The functions that follow can be used to read or configure such image properties, as well as to read or create an annotation, change zoom factor, or switch to a different plane in a stack. One programming example is given for all of the image "Get" functions (see Figure 4.1 on page 56).

GetActivePlane

Description	Obtains the number of the currently active plane in the specified image.
Syntax	GetActivePlane(hImage Long, nRetPlaneNumber As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image in which the plane resides.
Return values	<i>nRetPlaneNumber</i> Returns the number of the active plane.
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetNumberOfPlanes SetActivePlane

GetDepth

Description	Obtains the bit-depth of a specified image.
Syntax	GetDepth(hImage As Long, nRetDepth As Integer) As Long
Parameters	hImage Specifies the handle of the image.
Return values	<i>nRetDepth</i> Returns the bit-depth—1, 8, 16, or 24.
See also:	GetCurrentImage (Section 4.3), GetHeight, GetImage (Section 4.3), GetWidth

GetHeight

Description	Obtains the height (Y-axis size) of a specified image. (Note: This may differ from the height of the image <i>window</i> .)
Syntax	GetHeight(hImage As Long, nRetHeight As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image.

4.5 Reading and Using Image Properties, continued

GetHeight (continued)	
Return values	<i>nRetHeight</i> Returns the Y-axis size, in pixels, of the image.
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetImageWindowSize , GetWidth

GetImageAnnotation

Description	Obtains the annotation of the specified image.
Syntax	GetImageAnnotation(hImage As Long, planeNumber As Integer, sRetAnnotation As String) As Long
Remarks	Returns the annotation of the given plane of the given image. The annotation is returned in <i>sRetAnnotation</i> .
Parameters	<i>hImage</i> Specifies the handle of the image. <i>planeNumber</i> Specifies the number of the plane in a stack whose annotation you want to obtain. If the image is a single plane, this variable must be 0.
Return values	sRetAnnotation Returns the annotation text.
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetImageAnnotation

GetImageName

Description	Obtains the name of a specified image.
Syntax	GetImageName(hImage As Long, sRetName As String) As Long
Parameters	hImage Specifies the handle of the image.
Return values	sRetName Returns the name of the image.
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetImageName

GetNumberOfPlanes

Description	Obtains the number of planes in a specified image or stack of images.
Syntax	GetNumberOfPlanes(hImage As Long, nRetNumberOfPlanes As Integer) As Long
Parameters	hImage Specifies the handle of the image.
Return values	<i>nRetNumberOfPlanes</i> Returns the number of planes
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetActivePlane

GetWidth

Description	Obtains the width (X-axis size) of the specified image. (Note: This may differ from the width of the image <i>window</i> .)
Syntax	GetWidth(hImage As Long, nRetWidth As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image.
Return values	nRetWidth Returns the X-axis size, in pixels, of the image.
See also:	GetCurrentImage (Section 4.3), GetHeight, GetImage (Section 4.3), GetImageWindowSize

GetZoom

Description	Obtains the zoom factor level of the specified image.
Syntax	GetZoom(hImage As Long, nRetZoom As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image.
Return values	nRetZoom Returns the zoom level, expressed as a percentage of the full size (100 = normal size).
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetZoom

4.5 Reading and Using Image Properties, continued

Figure 4.1 Image "Get" Function Programming Example

```
' Print information about the current image
Dim im As Long
MM.GetCurrentImage im
Dim name As String
MM.GetImageName im, name
Dim w As Integer, h As Integer, d As Integer
MM.GetWidth im, w
MM.GetHeight im, h
MM.GetDepth im, d
MM.PrintMsg "The image " + name + " is " + Str(w)
 + " X " + Str(h)
MM.PrintMsg "and has a depth of " + Str(d)
Dim nPlanes As Integer
Dim curPlane As Integer
MM.GetNumberOfPlanes im, nPlanes
MM.GetActivePlane im, curPlane
MM.PrintMsg "It has " + Str(nPlanes) +
 " planes, and the current plane is " + Str(curPlane)
Dim ann As String
MM.GetImageAnnotation im, curPlane, ann
MM. PrintMsg "The annotation of the current plane is "
  + ann
Dim z As Integer
MM.GetZoom im, z
MM.PrintMsg "and the zoom is " + Str(z)
```

SetActivePlane

GetActivePlane, GetCurrentImage (Section 4.3), GetImage (Section 4.3)
End If
PrintMsg "Error! Image only has " + Str(<i>nPlanes</i>) + " planes"
MM.SetACtivePiane 10, 3
II nPlanes > 3 Then
MM.GetNumberOfPlanes im, nPlanes
Dim <i>nPlanes</i> As Integer
MM.GetCurrentImage im
Dim <i>im</i> As Long
' that many planes
' Set the active plane of the current image to 3 if it has
r
<i>planeNumber</i> Specifies the number of the plane to be made active.
hImage Specifies the handle of the image.
Planes are numbered starting at 0.
Serverver lane (nimage 113 Long, planetvanber 113 integer) 113 Long
Set Active Plane (hImage As Long plane Number As Integer) As Long
Selects a plane in a specified image stack and makes it the active plane.

SetImageAnnotation

Description	Assigns an annotation to an image or stack plane.
Syntax	SetImageAnnotation(hImage As Long, planeNumber As Integer, annotation As String) As Long
Parameters	hImage Specifies the handle of the image.
	<i>planeNumber</i> Specifies the number of the plane to be made active.
	annotation Provides the annotation text to be assigned to the selected image or plane.
Example	' Set the annotation of plane 0 of the current image Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i>
	MM.SetImageAnnotation im, 0, "This is the first plane"
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetImageAnnotation

SetImageName

Description	Assigns a new name to a selected image.
Syntax	SetImageName(hImage As Long, name As Long) As Long
Parameters	<i>hImage</i> Specifies the handle of the image. <i>name</i> Specifies the name for the image.
Example	' Set the name of the current image to "Fred" Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i> MM.SetImageName <i>im</i> , "Fred"
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetImageName

SetImageTimestamp

Description	Sets the timestamp (either creation time or last saved time) for a given image to the specified date and time.
Syntax	SetImageTimestamp (<i>hImage</i> As Long, <i>nStampType</i> As Integer, <i>nMonth</i> As Integer, <i>nDay</i> As Integer, <i>nYear</i> As Integer, <i>nHours</i> As Integer, <i>nMinutes</i> As Integer, <i>nSeconds</i> As Integer, <i>nMilliseconds</i> As Integer)
Parameters	hImage Specifies the handle of the image.
	<i>nStampType</i> Specifies which timestamp will be set, 0 for creation time and non-0 for last saved time.
	<i>nMonth</i> Specifies the month, with 1 signifying January, 6 signifying June, etc.
	<i>nDay</i> Specifies the date, from 1 to 31.
	<i>nYear</i> Specifies the year, using four digits.
	<i>nHours</i> Specifies the hour, from 0 to 23, counting forward from midnight.
	<i>nMinutes</i> Specifies the minute, from 0 to 59.
	<i>nSeconds</i> Specifies the seconds, from 0 to 59.
	<i>nMilliseconds</i> Specifies the milliseconds, from 0 to 999.
	Continued on next page

Reading and Using Image Properties, continued 4.5

SetImageTimestamp (continued)

Example	' Set the creation timestamp of the current image to 2:33:20.68 ' PM on November 3, 2001.
	Dim <i>im</i> As Long MM. GetCurrentImage <i>im</i> MM. SetImageTimestamp <i>im</i> , 0, 11, 3, 2001, 14, 33, 20, 680
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3)

SetZoom

Description	Sets the zoom factor of a specified image.
Syntax	SetZoom(<i>hImage</i> As Long, <i>zoom</i> As Integer, <i>nXCenter</i> As Integer, <i>nYCenter</i> As Integer) As Long
Remarks	Full size is assigned a numeric value of 100. Numbers smaller than 100 will make the image smaller, and those greater than 100 will make the image larger. If the image becomes larger than its display window, <i>nXCenter</i> and <i>nYCenter</i> give the pixel coordinates at which to center the image.
Parameters	hImage Specifies the handle of the image.
	zoom Specifies the zoom level, as a percent of the full-sized image.
	<i>nXCenter</i> Specifies the X-coordinate for the pixel at which the zoomed image is to be centered.
	<i>nYCenter</i> Specifies the Y-coordinate for the pixel at which the zoomed image is to be centered.
Example	' Set the zoom of the current image to 200 and center the image ' at 100, 100 Dim <i>im</i> As Long <i>MM</i> .GetCurrentImage <i>im</i> <i>MM</i> .SetZoom <i>im</i> , 200, 100, 100
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetZoom

Chapter 5 – Adjusting Image Display

5.1 Overview

Introduction	Image display can be modified in a number of ways—you contrast, you can brighten or darken the image, you can adj resolution by altering the number of entries in the image's p the values in the image's look-up table. This chapter covers need to use to accomplish these tasks.	can increase image ust the grayscale or color palette, or you can modify the functions you will
In this chapter	This chapter contains the following topics:	
	Торіс	See Page
	Updating the Image After Changing the Display	61
	Adjusting Brightness and Contrast	62
	Autoscaling 16-Bit Images	66
	Working with Look-up Tables and Palettes	71

5.2 Updating the Image After Changing the Display

Introduction	The first function in this chapter is of such vital importance that we feel it warrants a separate section of its own. Any time you execute a function that makes a change to
	an image, such as when you create a region of interest, alter the brightness or contrast, switch look-up tables, or assign a set of intensity values to a row or column of pixels, you will need to update the image display. The UpdateDisplay function carries out this operation.

UpdateDisplay

Description	Redraws any parts of the image that need updating. This function should be called after pixel modification operations to cause those changes to be drawn.
Syntax	UpdateDisplay(hImage As Long) As Long
Parameters	<i>hImage</i> Specifies the handle of the image to be updated.
Example	' Draw a white box on the current image and then update the ' display so the changed pixels are shown
	Dim <i>i</i> As Integer Dim <i>im</i> As Long Dim <i>dat</i> a(20) As Byte
	For <i>i</i> = 1 To 20 <i>data</i> (<i>i</i>) = 255 Next <i>i</i>
	MM.GetCurrentImage im
	For <i>i</i> = 1 To 20 <i>MM.</i> WriteRow <i>im</i> , 0, <i>i</i> , 20, 8, <i>data</i> Next <i>i</i>
	MM.UpdateDisplay im
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3)

5.3 Adjusting Brightness and Contrast

Introduction	The image display adjustments you are most often likely to perform are those which modify an image's brightness and contrast. Brightness settings can range from 0 to
	100, with 50 representing the brightness value of the original, unaltered image. Contrast settings range from 50 to 100 (MetaMorph can only <i>increase</i> image contrast
	from the original.) The following functions are used for reading the current settings
	and changing them to suit your needs. If you need to adjust the brightness or contrast of a 16-bit image, you will need to use the functions described in Section 5.4.

AutoEnhance

Description	Automatically enhances the brightness and contrast of a specified image.
Syntax	AutoEnhance(hImage As Long) As Long
Parameters	<i>hImage</i> Specifies the handle of the image to be autoenhanced.
Example	' Autoenhance the image contrast then fix it Dim <i>im</i> As Long <i>MM</i> .GetCurrentImage <i>im</i> <i>MM</i> .AutoEnhance <i>im</i> <i>MM</i> .FixImage <i>im</i>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetBrightness, SetContrast

FixImage

Description	Makes changes to the brightness and/or contrast of an image permanent.
Syntax	FixImage(hImage As Long) As Long
Parameters	<i>hImage</i> Specifies the handle of the image in which brightness and/or contrast changes are to be made permanent.
Example	' Autoenhance the image contrast then fix it Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i> MM.AutoEnhance <i>im</i> MM.FixImage <i>im</i>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetBrightness, SetContrast

5.3 Adjusting Brightness and Contrast, continued

GetBrightness

Description	Obtains the brightness setting of a specified image.
Syntax	GetBrightness(hImage As Long, nRetBrightness As Double) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose brightness setting you want to obtain.
Example	<pre>' Double the brightness and contrast of the current image Dim im As Long Dim bright As Double Dim contrast As Long MM.GetCurrentImage im MM.GetBrightness im, bright MM.GetContrast im, contrast MM.SetBrightness im, bright * 2 MM.SetContrast im, contrast * 2</pre>
Return values	<i>nRetBrightness</i> Returns the brightness setting value.
See also:	GetContrast, GetCurrentImage (Section 4.3), GetImage (Section 4.3)

GetCo	ntrast
-------	--------

Description	Obtains the contrast setting of a specified image.
Syntax	GetContrast(hImage As Long, nRetContrast As Double) As Long
Parameters	hImage Specifies the handle of the image whose contrast setting you want to obtain.
Return values	nRetContrast Returns the contrast setting value.
Example	<pre>' Double the brightness and contrast of the current image Dim im As Long Dim bright As Double Dim contrast As Long MM.GetCurrentImage im MM.GetBrightness im, bright MM.GetContrast im, contrast MM.SetBrightness im, bright * 2 MM.SetContrast im, contrast * 2</pre>
See also:	GetBrightness, GetCurrentImage (Section 4.3), GetImage (Section 4.3)

ResetContrast

Description	Resets the contrast and brightness of a specified image to the settings it had when last saved or "fixed."
Syntax	ResetContrast(hImage As Long) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose brightness and/or contrast settings you want to reset.
Example	' Reset the contrast of the current image Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i> MM.ResetContrast <i>im</i>
See also:	FixImage, GetCurrentImage (Section 4.3), GetImage (Section 4.3)

SetBrightness

Description	Sets the brightness of an image to a specified intensity value.
Syntax	SetBrightness(hImage As Long, brightness As Double) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose brightness you want to set. <i>brightness</i> Specifies a new intensity setting.
Example	<pre>' Double the brightness and contrast of the current image Dim im As Long Dim bright As Double Dim contrast As Long MM.GetCurrentImage im MM.GetBrightness im, bright MM.GetContrast im, contrast MM.SetBrightness im, bright * 2 MM.SetContrast im, contrast * 2</pre>
See also:	AutoEnhance, FixImage, GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetContrast

SetContrast

Description	Sets the contrast of an image to a specified setting.
Syntax	SetContrast(hImage As Long, contrast As Double) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose contrast you want to set. <i>contrast</i> Specifies a new contrast setting.
Example	<pre>' Double the brightness and contrast of the current image Dim im As Long Dim bright As Double Dim contrast As Long MM.GetCurrentImage im MM.GetBrightness im, bright MM.GetContrast im, contrast MM.SetBrightness im, bright * 2 MM.SetContrast im, contrast * 2</pre>
See also:	AutoEnhance, FixImage, GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetBrightness

5.4 Autoscaling 16-Bit Images

Introduction Sixteen-bit images require a different approach when it comes to adjusting brightness and contrast. Because of the greater range of intensity values that are involved, autoscaling a 16-bit image relies on a different set of functions from those described in the preceding section.

GetAutoScale

Description	Obtains the current autoscaling state of a specified 16-bit image.
Syntax	GetAutoScale(hImage As Long, bRetOnOff As Boolean) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose autoscaling state you want to obtain.
Return values	<i>bRetOnOff</i> Returns the current autoscaling state. A value of TRUE means it is currently enabled, and FALSE means it is currently disabled.
Example	' Get the autoscale properties of the current image
	Dim <i>im</i> As Long Dim <i>onoff</i> As Boolean Dim <i>min</i> As Integer Dim <i>max</i> As Integer
	MM.GetCurrentImage im MM.GetAutoScale im, onoff MM.GetMinScale im, min MM.GetMaxScale im, max
	<pre>MM.PrintMsg "The min and max scale of the current image is " + Str(min) + " and " + Str(max) If onoff = TRUE Then MM PrintMag "The outcompling of the surrent image is ON"</pre>
	MMM.FrintMsg "The autoscaling of the current image is ON" Else MM.PrintMsg "The autoscaling of the current image is OFF" End If
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetAutoScale

5.4 Autoscaling 16-Bit Images, continued

GetMaxScale

Description	Obtains the maximum scaling value of a specified 16-bit image.
Syntax	GetMaxScale(hImage As Long, nRetMaxScale As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose maximum scaling value you want to obtain.
Return values	<i>nRetMaxScale</i> Returns the maximum scaling grayscale value.
Example	' Get the autoscale properties of the current image
	Dim <i>im</i> As Long Dim <i>onoff</i> As Boolean Dim <i>min</i> As Integer Dim <i>max</i> As Integer
	MM.GetCurrentImage im MM.GetAutoScale im, onoff MM.GetMinScale im, min MM.GetMaxScale im, max
	<pre>MM.PrintMsg "The min and max scale of the current image is " + Str(min) + " and " + Str(max) If onoff = TRUE Then MM.PrintMsg "The autoscaling of the current image is ON" Else MM.PrintMsg "The autoscaling of the current image is OFF" End If</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetMinScale
GetMinScale

Description	Obtains the minimum scaling value of a specified 16-bit image.
Syntax	GetMinScale(hImage As Long, nRetMinScale As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose minimum scaling value you want to obtain.
Return values	<i>nRetMinScale</i> Returns the minimum scaling grayscale value.
Example	' Get the autoscale properties of the current image
	Dim <i>im</i> As Long Dim <i>onoff</i> As Boolean Dim <i>min</i> As Integer Dim <i>max</i> As Integer <i>MM</i> .GetCurrentImage <i>im</i> <i>MM</i> .GetAutoScale <i>im</i> , <i>onoff</i>
	MM.GetMinScale im, min MM.GetMaxScale im, max
	<pre>MM.PrintMsg "The min and max scale of the current image is " + Str(min) + " and " + Str(max) If onoff = TRUE Then MM.PrintMsg "The autoscaling of the current image is ON" Else MM.PrintMsg "The autoscaling of the current image is OFF"</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetMaxScale

SetAutoScale

Description	Enables or disables autoscaling for the specified 16-bit image.
Syntax	SetAutoScale(hImage as Long, bOnOff As Boolean) As Long
Parameters	<i>hImage</i> Specifies the handle of the image for which you want to enable or disable autoscaling.
	<i>bOnOff</i> Sets the scaling state. If you set <i>bOnOff</i> to TRUE, autoscaling will be enabled. To disable autoscaling, set <i>bOnOff</i> to FALSE.

5.4 Autoscaling 16-Bit Images, continued

SetAutoScale

(continued)

Example	<pre>' Enable autoscaling for the current image and set the ' range from 0 to 4095 Dim im As Long MM.GetCurrentImage im MM.SetAutoScale im, TRUE MM.SetMinScale im, 0 MM.SetMaxScale im, 4095</pre>
See also:	GetAutoScale, GetCurrentImage (Section 4.3), GetImage (Section 4.3)

SetMaxScale

Description	Sets the maximum value for the grayscale range being used to autoscale a specified 16-bit image.
Syntax	SetMaxScale(hImage As Long, nMaxScale As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image for which you want to set the maximum autoscaling value.
	<i>nMaxScale</i> Defines the maximum grayscale value for the autoscaling range.
Example	<pre>' Enable autoscaling for the current image and set the ' range from 0 to 4095 Dim im As Long MM.GetCurrentImage im MM.SetAutoScale im, TRUE MM.SetMinScale im, 0 MM.SetMaxScale im, 4095</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetAutoScale, SetMinScale

SetMinScale

Description	Sets the minimum value for the grayscale range being used to autoscale a specified 16-bit image.
Syntax	SetMinScale(hImage As Long, nMinScale As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image for which you want to set the minimum autoscaling value.
	<i>nMinScale</i> Defines the minimum grayscale value for the autoscaling range.
Example	<pre>' Enable autoscaling for the current image and set the ' range from 0 to 4095 Dim im As Long MM.GetCurrentImage im MM.SetAutoScale im, TRUE MM.SetMinScale im, 0 MM.SetMaxScale im, 4095</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetAutoScale, SetMaxScale

5.5 Working with Look-up Tables and Palettes

Introduction A *look-up table*, or LUT, is a "table" of values that translates each original grayscale value in an image to another intensity or to a color. This allows you to change the display of an image without affecting its data. The functions in this section can be used for making changes to the current LUT or for switching to another LUT. In addition, you can change the number of grayscale levels or colors that are displayed by changing the number of *palette* entries with the **SetNumPaletteEntries** function.

GetLut – for Visual Basic 6 & earlier

GetLutEx2 – for Visual Basic .NET (2002 – 2008)

Description	Reads the elements from the given look-up table (LUT) into the red, green, and blue arrays that are passed.
Syntax	GetLut(<i>hImage</i> As Long, <i>lutNumber</i> As Integer, <i>startElement</i> As Integer, <i>nElements</i> As Integer, <i>red()</i> As Byte, <i>green()</i> As Byte, <i>blue()</i> As Byte) As Long GetLutEx2(<i>hImage</i> As Long, <i>lutNumber</i> As Integer, <i>startElement</i> As Integer, <i>nElements</i> As Integer, <i>red()</i> As Byte, <i>green()</i> As Byte, <i>blue()</i> As Byte, <i>blue()</i> As Byte) As Long
Remarks	The LUT that you want to read is specified by <i>lutNumber</i> . Use one of the following constants:
	0 = Monochrome 1 = Pseudocolor 2 = Red 3 = Green 4 = Blue
	or use a user LUT number from 5 to 15.
Parameters	<i>hImage</i> Specifies the handle of the image for which you want to read the LUT values.
	<i>lutNumber</i> Specifies the LUT to be read. (See Remarks.)
	<i>startElement</i> Indicates the number of the element in the LUT where reading is to start.
	<i>nElements</i> Specifies the number of LUT elements to read. This will determine the size of the <i>red()</i> , <i>green()</i> , and <i>blue()</i> arrays (see Return values).
Return values	<i>red()</i> Passes the values of the elements from the red channel of the LUT. The size of this array will be determined by the number of elements in the look-up table.
	<i>green()</i> Passes the values of the elements from the green channel of the LUT. The size of this array will be determined by the number of elements in the look-up table.
	<i>blue()</i> Passes the values of the elements from the blue channel of the LUT. The size of this array will be determined by the number of elements in the look-up table.

GetLut	
GeiLui	

(continued)

Example	<pre>' Read the first user lut of the current image into memory. ' Since LUTs 0 through 4 are the fixed LUTs, the first user ' LUT is LUT number 5 Dim im As Long Dim r(256) As Byte Dim g(256) As Byte Dim b(256) As Byte MM.GetCurrentImage im MM.GetLut im, 5, 0, 256, r, g, b</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetLutModel, SetLut

GetLutModel

Description	Obtains a value that indicates the look-up table (LUT) currently being used to display the given image.
Syntax	GetLutModel(hImage As Long, nRetLutModel As Integer) As Long
Remarks	The LUT that is in use is indicated by a numeric value. The values returned in <i>nRetLutModel</i> are
	0 = Monochrome 1 = Pseudocolor 2 = Red 3 = Green 4 = Blue or a user-defined LUT, with a number between 5 and 15
Demonstrations	or a user-defined LOT, with a number between 5 and 15.
Parameters	<i>hImage</i> Specifies the handle of the image for which you want to determine the LUT.
Return values	<i>nRetLutModel</i> Returns the LUT value. (See Remarks.)

GetLutModel

(continued)

Example	' Determine which LUT is currently being used on the current ' image
	Dim <i>im</i> As Long Dim <i>lutnum</i> As Integer
	MM.GetCurrentImage im MM.GetLutModel im, lutnum
	<pre>If lutnum = 0 Then MM.PrintMsg "LUT is Monochrome"</pre>
	Elself <i>lutnum</i> = 1 Then <i>MM</i> . PrintMsg "LUT is Pseudocolor" Elself <i>lutnum</i> = 2 Then
	MM.PrintMsg "LUT is Red" ElseIf lutnum = 3 Then
	MM.PrintMsg "LUT is Green" ElseIf <i>lutnum</i> = 4 Then MM.PrintMsg "LUT is Blue"
	<pre>Else MM.PrintMsg "LUT is user LUT number " + Str(lutnum = 5) End If</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetLut, SetLutModel

SetLut	
Description	Writes the contents of the red, green, and blue arrays passed into the elements of the given look-up table (LUT).
Syntax	SetLut (<i>hImage</i> As Long, <i>lutNumber</i> As Integer, <i>startElement</i> As Integer, <i>nElements</i> As Integer, <i>red()</i> As Byte, <i>green()</i> As Byte, <i>blue()</i> As Byte) As Long
Parameters	<i>hImage</i> Specifies the handle of the image for which you want to set the LUT.
	<i>lutNumber</i> Denotes the user-defined LUT to which you want to write (5 to 15). You cannot write to the predefined LUTs (Monochrome, etc).
	<i>startElement</i> Indicates the number of the element in the LUT where writing is to start.
	<i>nElements</i> Specifies the number of LUT elements to write.
	<i>red()</i> Indicates the array for the red channel of the LUT. LUTs have a maximum possible total of 256 elements. The red array must contain at least <i>nElements</i> elements.
	<i>green()</i> Indicates the array for the green channel of the LUT. LUTs have a maximum possible total of 256 elements. The green array must contain at least <i>nElements</i> elements.
	<i>blue()</i> Indicates the array for the blue channel of the LUT. LUTs have a maximum possible total of 256 elements. The blue array must contain at least <i>nElements</i> elements.
Example	' Configure user LUT 0 on the current image as an inverted ' contrast monochrome LUT and then use it
	Dim <i>im</i> As Long Dim <i>i</i> As Integer Dim <i>lut</i> (256) As Byte
	For i = 0 To 255 lut(i) = 255 - i Next i
	MM.GetCurrentImage im MM.SetLut im, 5, 0, 256, lut, lut, lut MM.SetLutModel im, 5
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetLutModel, SetNumPaletteEntries

SetLutModel

Description	Specifies a look-up table (LUT) to use to display the given image.
Syntax	SetLutModel(hImage As Long, nLutModel As Integer) As Long
Remarks	The model of LUT to use is specified with <i>nLutModel</i> . Values you can use are
	0 = Monochrome 1 = Pseudocolor 2 = Red 3 = Green 4 = Blue or a user-defined LUT, with a number between 5 and 15.
Parameters	<i>hImage</i> Specifies the handle of the image for which you want to set the LUT model.
	<i>nLutModel</i> Specifies the value of the LUT you want to use. (See Remarks.)
Example	' Configure user LUT 0 on the current image as an inverted ' contrast monochrome LUT and then use it
	Dim <i>im</i> As Long Dim <i>i</i> As Integer Dim <i>lut</i> (256) As Byte
	For <i>i</i> = 0 To 255 <i>lut</i> (<i>i</i>) = 255 - <i>i</i> Next <i>i</i>
	MM.GetCurrentImage im MM.SetLut im, 5, 0, 256, lut, lut, lut MM.SetLutModel im, 5
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetLut, SetNumPaletteEntries

SetNumPaletteEntries

Description	Specifies the number of palette entries to use for displaying an image.
Syntax	SetNumPaletteEntries(hImage As Long, newNumEntries As Integer) As Long
Remarks	The maximum possible number of palette entries is 236. This is because the Windows operating system reserves 20 palette colors for use by such interface elements as dialog box buttons, title bars, and the like.
Parameters	<i>hImage</i> Specifies the handle of the image for which you want to set the number of palette entries.
	<i>newNumEntries</i> Specifies the number of palette entries: 2, 4, 8, 16, 32, 64, 128, or 236.
Example	' Set the number of palette entries on the current image to 8. ' This will make the image appear as if it is quantized. Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i> MM.SetNumPaletteEntries <i>im</i> , 8
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetLut, SetLutModel

Chapter 6 – Reading and Using Image Pixel Data

6.1 Overview

Introduction	Digital images are constructed of individual points, or <i>pixels</i> , each of specific, quantifiable intensity value or color. Because each pixel call in memory by a set of numeric values that encode its location (X and brightness or color (grayscale or color value), a variety of mani applied that analyze and process the information contained in an impleals with the functions that read and manipulate image pixel data.	of which has a an be represented ad Y coordinates) pulations can be lage. This chapter
In this chapter	This chapter contains the following topics:	
	Торіс	See Page
	Applying Thresholding	78
	Reading and Manipulating Image Data	82

6.2 Applying Thresholding

Introduction

Most procedures in image analysis require that a distinction be made between the objects to be measured and the rest of the image. MetaMorph makes this distinction on the basis of the intensity values of pixels in the object vs. those in the background regions. This procedure, which is called *segmentation* or *thresholding*, involves defining a range of intensity values that will be considered as belonging to the objects being measured. A threshold range can be inclusive or exclusive. An *inclusive* threshold is one which considers all grayscale values between the upper and lower limits of the threshold considers all grayscale values that are equal to, or outside of, the upper and lower limits as belonging to the objects being measured. The four functions described in this section are used for determining the current threshold settings and for defining a new set of threshold settings.

GetThresholdRange

Description	Obtains the current threshold range (maximum and minimum grayscale settings) of the specified image.
Syntax	GetThresholdRange(hImage As Long, nRetLow As Integer, nRetHigh As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose threshold range you want to obtain.
Return values	<i>nRetLow</i> Returns the grayscale value of the lower threshold limit.<i>nRetHigh</i> Returns the grayscale value of the upper threshold limit.
Example	<pre>' Get the threshold range and state of the current image Dim im As Long Dim low As Integer, high As Integer, state As Integer MM.GetCurrentImage im MM.GetThresholdRange im, low, high MM.GetThresholdState im, state MM.PrintMsg "The low and high threshold values of the current image are " + Str(low) + " and " + Str(high)</pre>
	<pre>Image are " + Str(Iow) + " and " + Str(IIgh) If state = 0 Then MM.PrintMsg "Thresholding is currently off" ElseIf state = 1 Then MM.PrintMsg "Thresholding is set to 'On-Inclusive' mode" ElseIf state = 2 Then MM.PrintMsg "Thresholding is set to 'On-Exclusive' mode" End If</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetThresholdState, SetThresholdRange

GetThresholdState

Description	Obtains the threshold state of the specified image
Syntax	GetThresholdState(hImage As Long, nRetState As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose thresholding state you want to obtain.
Return values	<i>nRetState</i> Returns the current thresholding state of the given image. If thresholding is disabled, <i>nRetState</i> will return a value of 0. If the threshold state is set to On-Inclusive (pixels with grayscale values between the upper and lower threshold limits are included), <i>nRetState</i> will be set to 1. If the threshold state is set to On-Exclusive (pixels with grayscale values equal or lower than the lower threshold limit, and those with values equal or higher than the upper limit, are included), <i>nRetState</i> will return a value of 2.
Example	<pre>' Get the threshold range and state of the current image Dim im As Long Dim low As Integer, high As Integer, state As Integer MM.GetCurrentImage im MM.GetThresholdRange im, low, high MM.GetThresholdState im, state</pre>
	<pre>MM.PrintMsg "The low and high threshold values of the current image are " + Str(low) + " and " + Str(high) If state = 0 Then MM.PrintMsg "Thresholding is currently off" ElseIf state = 1 Then MM.PrintMsg "Thresholding is set to 'On-Inclusive' mode" ElseIf state = 2 Then MM.PrintMsg "Thresholding is set to 'On-Exclusive' mode" End If</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetThresholdRange, SetThresholdState

SetThresholdRange

Description	Sets the upper and lower limit of a threshold range for a specified image.
Syntax	SetThresholdRange(hImage As Long, low As Integer, high As Integer) As Long
Parameters	<i>hImage</i> Specifies the handle of the image whose thresholding range you want to set.<i>low</i> Specifies the grayscale value of the lower threshold limit.<i>high</i> Specifies the grayscale value of the upper threshold limit.
Example	' Enable thresholding in the current image, and highlight pixels ' between grayscale values 50 and 187 Dim <i>im</i> As Long MM.GetCurrentImage <i>im</i> MM.SetThresholdRange <i>im</i> , 50, 187 MM.SetThresholdState <i>im</i> , 1
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetThresholdState

SetThresholdState

Description	Sets the thresh	hold state of the specified image.
Syntax	SetThreshold	IState(hImage As Long, state As Integer) As Long
Remarks	The threshold state is set with the <i>state</i> variable. Possible values are	
	0:	Disables thresholding (Off).
	1:	Enables inclusive thresholding, in which all pixels with grayscale values between the <i>high</i> and <i>low</i> values (see SetThresholdRange) are highlighted.
	2:	Enables exclusive thresholding, in which all pixels with grayscale values equal to or outside the <i>high</i> and <i>low</i> values are highlighted.
Parameters	<i>hImage</i> Specifies the handle of the image whose thresholding state you want to set. <i>state</i> Indicates the thresholding state for the image. (See Remarks.)	

6.2 Applying Thresholding, continued

SetThresholdState (continued)	
Example	<pre>' Enable thresholding in the current image, and highlight pixels ' between grayscale values 50 and 187 Dim im As Long MM.GetCurrentImage im MM.SetThresholdRange im, 50, 187 MM.SetThresholdState im, 1</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetThresholdRange

6.3 Reading and Manipulating Image Data

Introduction	Pixel intensities are represented by numeric values in computer memory, and can be
	easily read and manipulated. This section describes the functions that perform such
	"read" and "write" procedures. The BinarizeImage and WriteText functions are
	included here because the process of converting all pixels to black or white and the
	process of adding a text label to an image are little more than a reassignment of the
	intensity values of the underlying pixels.

BinarizeImage

Description	Creates a binarized (1-bit) version of a selected image.
Syntax	BinarizeImage(hImage As Long, hDest As Long) As Long
Remarks	This function places a binarized version of <i>hImage</i> in the image given by <i>hDest</i> . <i>hImage</i> should be thresholded before this function is called, or you will get a blank result image.
Parameters	hImage Specifies the handle of the image you want to binarize.
	hDest Specifies the handle of a destination image that will be overwritten with the binarized version of $hImage$. The destination image must exist before you can apply this function. Typically, you will create an image with the CreateImage command.
Example	' Binarize the current image and place it in "Binarized"
	Dim <i>curIm</i> As Long Dim <i>binIm</i> As Long
	MM.GetCurrentImage curIm MM.CreateImage 512, 512, 1, "Binarized", binIm
	' <i>curIm</i> must be thresholded before it can be binarized MM.SetThresholdState <i>curIm</i> , 1 MM.SetThresholdRange <i>im</i> , 50, 150
	MM.BinarizeImage curIm, binIm
See also:	GetCurrentImage (Section 4.3), CreateImage (Section 4.2), GetImage (Section 4.3)

ReadColumn

Description	Reads the intensity values of a vertical line of pixels from a selected image and writes them into an array.
Syntax	ReadColumn (<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nPixels</i> As Integer, <i>depth</i> As Integer, <i>pixelBuffer()</i> As Variant) As Long
Parameters	<i>hImage</i> Specifies the handle of the image to be read.
	<i>xPos</i> Specifies the X-coordinate of the pixel where reading should start.
	<i>yPos</i> Specifies the Y-coordinate of the pixel where reading should start.
	<i>nPixels</i> Specifies the number of pixels to be read, starting at the position given by $xPos$ and $yPos$. The size of the array will depend on the image bit-depth and the data type used for the array buffer. (See Section 1.6, <i>Data Types and Arrays.</i>)
	<i>depth</i> Specifies what the bit-depth of the pixel data values should be. (See Section 1.6, <i>Data Types and Arrays.</i>)
	<i>pixelBuffer()</i> Defines a buffer into which the pixel values will be read. If <i>depth</i> is 8, <i>pixelBuffer</i> should be defined as <i>Byte</i> . If <i>depth</i> is 16, <i>pixelBuffer</i> should be <i>Integer</i> . If <i>depth</i> is 24, <i>pixelBuffer</i> should be <i>Byte</i> .
Return values	<i>pixelBuffer()</i> The pixel values will be read into this predefined buffer. If the source image is 24-bit, the array will be read out in the format B, G, R, B, G, R, etc. Thus, array(0) will be the first blue pixel, array(1) will be the first green pixel, and so forth. (See Parameters.)
Example	<pre>' Read a column of 50 pixels starting at location (10, 0) from ' the current image Dim im As Long Dim buf(50) As Byte MM.GetCurrentImage im MM.ReadColumn im, 10, 0, 50, 8, buf</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), ReadColumnEx, ReadPixel, ReadRow

ReadColumnEx – for Visual Basic 6 & earlier ReadColumnEx2 – for Visual Basic .NET (2002 – 2008)

Description	Reads the intensity values of a vertical line of pixels from a selected image and writes the values into an array, starting at a specified starting location in the array.
Syntax	ReadColumnEx (<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nPixels</i> As Integer, <i>depth</i> As Integer, <i>xStart</i> As Integer, <i>yStart</i> As Integer, <i>pixelBuffer()</i> As Variant) As Long ReadColumnEx2 (<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nPixels</i> As Integer, <i>depth</i> As Integer, <i>xStart</i> As Integer, <i>yStart</i> As Integer, <i>pixelBuffer()</i> As Variant) As Long
Parameters	<i>hImage</i> Specifies the handle of the image to be read.
	<i>xPos</i> Specifies the X-coordinate of the pixel where reading should start.
	<i>yPos</i> Specifies the Y-coordinate of the pixel where reading should start.
	<i>nPixels</i> Specifies the number of pixels to be read, starting at the position given by $xPos$ and $yPos$. The size of the array will depend on the image bit-depth and the data type used for the array buffer. (See Section 1.6, <i>Data Types and Arrays.</i>)
	<i>depth</i> Specifies what the bit-depth of the pixel data values should be. (See Section 1.6, <i>Data Types and Arrays.</i>)
	<i>xStart</i> Specifies the X-coordinate of the position in the array where writing should start.
	<i>yStart</i> Specifies the Y-coordinate of the start position in the array for writing.
	<i>pixelBuffer()</i> Defines a buffer into which the pixel values will be read. This is a two- dimensional array. Data are always placed in the array in "column-major" fashion. If <i>depth</i> is 8, <i>pixelBuffer</i> should be defined as <i>Byte</i> . If <i>depth</i> is 16, <i>pixelBuffer</i> should be <i>Integer</i> . If <i>depth</i> is 24, <i>pixelBuffer</i> should be <i>Byte</i> .
Return values	<i>pixelBuffer()</i> The pixel values will be read into this predefined buffer. If the source image is 24-bit, the array will be read out in the format B, G, R, B, G, R, etc. Thus, array(0) will be the first blue pixel, array(1) will be the first green pixel, and so forth. (See Parameters.)
Example	' Read a 50 x 50 block of pixels from the current image ' and write it back rotated 90 degrees counterclockwise Dim <i>im</i> As Long Dim <i>x</i> As Integer Dim <i>buf</i> (50, 50) As Byte <i>MM</i> . GetCurrentImage <i>im</i>
	<pre>For x = 0 To 49 MM.ReadColumnEx im, x, 0, 50, 8, 0, x, buf MM.WriteRowEx im, 0, x + 50, 50, 8, 0, x, buf Next x</pre>

ReadPixel

Description	Reads the intensity value of a specified pixel in a selected image.
Syntax	ReadPixel (<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nRetPixelValue</i> As Integer) As Long
Remarks	This function does not operate on 24-bit color images.
Parameters	<i>hImage</i> Specifies the handle of the image to be read.<i>xPos</i> Specifies the X-coordinate of the pixel to be read.<i>yPos</i> Specifies the Y-coordinate of the pixel to be read.
Return values	<i>nRetPixelValue</i> Returns the grayscale value of the pixel.
Example	' Read a pixel at location 43, 98 on the current image Dim <i>im</i> As Long Dim <i>pix</i> As Integer MM.GetCurrentImage <i>im</i> MM.ReadPixel <i>im</i> , 43, 98, <i>pix</i>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), ReadColumn, ReadRow

ReadRow – for Visual Basic 6 & earlier

Description	Reads the intensity values of a horizontal line of pixels from a selected image and writes them into an array.
Syntax	ReadRow (<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nPixels</i> As Integer, <i>depth</i> As Integer, <i>pixelBuffer()</i> As Variant) As Long
Parameters	<i>hImage</i> Specifies the handle of the image to be read.
	<i>xPos</i> Specifies the X-coordinate of the pixel where reading should start.
	<i>yPos</i> Specifies the Y-coordinate of the pixel where reading should start.
	<i>nPixels</i> Specifies the number of pixels to be read, starting at the position given by $xPos$ and $yPos$. The size of the array will depend on the image bit-depth and the data type used for the array buffer. (See Section 1.6, <i>Data Types and Arrays.</i>)

ReadRow (continued)	
	<i>depth</i> Specifies what the bit-depth of the pixel data values should be. (See Section 1.6, <i>Data Types and Arrays.</i>)
	<i>pixelBuffer()</i> This is a defined buffer into which the pixel values will be read. If <i>depth</i> is 8, <i>pixelBuffer</i> should be defined as <i>Byte</i> . If <i>depth</i> is 16, <i>pixelBuffer</i> should be <i>Integer</i> . If <i>depth</i> is 24, <i>pixelBuffer</i> should be <i>Byte</i> .
Return values	<i>pixelBuffer()</i> The pixel values will be read into this predefined buffer. If the source image is 24-bit, the array will be read out in the format B, G, R, B, G, R, etc. Thus, array(0) will be the first blue pixel, array(1) will be the first green pixel, and so forth. (See Parameters.)
Example	' Read a row of 50 pixels from the current image Dim <i>im</i> As Long Dim <i>buf</i> (50) As Byte <i>MM</i> .GetCurrentImage <i>im</i> <i>MM</i> .ReadRow <i>im</i> , 0, 0, 50, 8, <i>buf</i>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), ReadColumn, ReadPixel, ReadRowEx

ReadRowEx – for Visual Basic 6 & earlier ReadRowEx2 – for Visual Basic .NET (2002 – 2008)

Reads the intensity values of a horizontal line of pixels from a selected image and writes the values into an array, starting at a specified starting location in the array.
ReadRowEx (<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nPixels</i> As Integer, <i>depth</i> As Integer, <i>xStart</i> As Integer, <i>yStart</i> As Integer, <i>pixelBuffer()</i> As Variant) As Long
<i>hImage</i> Specifies the handle of the image to be read.
<i>xPos</i> Specifies the X-coordinate of the pixel where reading should start.
<i>yPos</i> Specifies the Y-coordinate of the pixel where reading should start.
<i>nPixels</i> Specifies the number of pixels to be read, starting at the position given by $xPos$ and $yPos$. The size of the array will depend on the image bit-depth and the data type used for the array buffer. (See Section 1.6, <i>Data Types and Arrays.</i>)
<i>depth</i> Specifies what the bit-depth of the pixel data values should be. (See Section 1.6, <i>Data Types and Arrays.</i>)
<i>xStart</i> Specifies the X-coordinate (row) of the position in the array where writing should start.

ReadRowEx 2 (continued)	
	<i>yStart</i> Specifies the Y-coordinate (column) of the position in the array where writing should start.
	<i>pixelBuffer()</i> This is a defined buffer into which the pixel values will be read. This is a two-dimensional array. Data are always placed in the array in "column-major" fashion. If <i>depth</i> is 8, <i>pixelBuffer</i> should be defined as <i>Byte</i> . If <i>depth</i> is 16, <i>pixelBuffer</i> should be <i>Integer</i> . If <i>depth</i> is 24, <i>pixelBuffer</i> should be <i>Byte</i> .
Return values	<i>pixelBuffer()</i> The pixel values will be read into this predefined buffer. If the source image is 24-bit, the array will be read out in the format B, G, R, B, G, R, etc. Thus, array(0) will be the first blue pixel, array(1) will be the first green pixel, and so forth. (See Parameters.)
Example	' Read a 50 x 50 block of pixels from the current image ' and write it back rotated 90 degrees counterclockwise Dim <i>im</i> As Long Dim <i>x</i> As Integer Dim <i>buf</i> (50, 50) As Byte <i>MM</i> . GetCurrentImage <i>im</i>
	<pre>For x = 0 To 49 MM.ReadColumnEx im, x, 0, 50, 8, 0, x, buf MM.WriteRowEx im, 0, x + 50, 50, 8, 0, x, buf Next x</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), ReadColumnEx, ReadPixel, ReadRow

WriteColumn – for Visual Basic 6 & earlier

Description	Reads a set of intensity values from an array and writes them to a vertical line of pixels in a selected image.
Syntax	WriteColumn(<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nPixels</i> As Integer, <i>depth</i> As Integer, <i>pixelBuffer()</i> As Variant) As Long
Parameters	hImage Specifies the handle of the image to be written to.
	<i>xPos</i> Specifies the X-coordinate of the pixel where writing should start.
	yPos Specifies the Y-coordinate of the pixel where writing should start.
	<i>nPixels</i> Specifies the number of pixels to be written, starting at the position given by $xPos$ and $yPos$. The size of the array will depend on the image bit-depth and the data type used for the array buffer. (See Section 1.6, <i>Data Types and Arrays.</i>)

WriteColumn (continued)	
	<i>depth</i> Specifies what the bit-depth of the pixel data values should be. (See Section 1.6, <i>Data Types and Arrays.</i>)
	<i>pixelBuffer()</i> This is a buffer from which the pixel values will be read. If <i>depth</i> is 8, <i>pixelBuffer</i> should be defined as <i>Byte</i> . If <i>depth</i> is 16, <i>pixelBuffer</i> should be <i>Integer</i> . If <i>depth</i> is 24, <i>pixelBuffer</i> should be <i>Byte</i> .
Example	' Write a white vertical line on the current image
	Dim <i>im</i> As Long Dim <i>buf</i> (50) As Byte Dim <i>x</i> As Integer
	For $x = 0$ To 49 buf(x) = 255 Next x
	MM. GetCurrentImage im MM. WriteColumn im, 50, 0, 50, 8, buf
WriteColum WriteColum Description	WritePixel, WriteRow nEx – for Visual Basic 6 & earlier nEx2 – for Visual Basic .NET (2002 – 2008) Reads a set of intensity values from a specified starting location in an array and writes
	them to a vertical line of pixels in a selected image
Syntax	 WriteColumnEx(<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nPixels</i> As Integer, <i>depth</i> As Integer, <i>xStart</i> As Integer, <i>yStart</i> As Integer, <i>pixelBuffer()</i> As Variant) As Long WriteColumnEx2(<i>hImage</i> As Long, <i>xPos</i> As Integer, <i>yPos</i> As Integer, <i>nPixels</i> As Integer, <i>depth</i> As Integer, <i>xStart</i> As Integer, <i>vStart</i> As Integer, <i>pixelBuffer()</i> As
	Variant) As Long
Parameters	<i>hImage</i> Specifies the handle of the image to be written to.
	<i>xPos</i> Specifies the X-coordinate of the pixel where writing should start.
	<i>yPos</i> Specifies the Y-coordinate of the pixel where writing should start.
	<i>nPixels</i> Specifies the number of pixels to be written, starting at the position given by $xPos$ and $yPos$. The size of the array will depend on the image bit-depth and the data type used for the array buffer. (See Section 1.6, <i>Data Types and Arrays.</i>)
	<i>depth</i> Specifies what the bit-depth of the pixel data values should be. (See Section 1.6, <i>Data Types and Arrays.</i>)

<i>xStart</i> Specifies the X-coordinate (row) of the position in the array where reading should start.
<i>yStart</i> Specifies the Y-coordinate (column) of the position in the array where reading should start.
<i>pixelBuffer()</i> This is a buffer from which the pixel values will be read. If <i>depth</i> is 8, <i>pixelBuffer</i> should be defined as <i>Byte</i> . If <i>depth</i> is 16, <i>pixelBuffer</i> should be <i>Integer</i> . If <i>depth</i> is 24, <i>pixelBuffer</i> should be <i>Byte</i> .
' Read a 50 x 50 block of pixels from the current image ' and write it back rotated 90 degrees clockwise Dim <i>im</i> As Long Dim <i>x</i> As Integer Dim <i>buf</i> (50, 50) As Byte
MM.GetCurrentImage im
<pre>For x = 0 To 49 MM.ReadRowEx im, 0, x, 50, 8, 0, x, buf MM.WriteColumnEx im, x + 50, 0, 50, 8, 0, x, buf Next x</pre>
GetCurrentImage (Section 4.3), GetImage (Section 4.3), WriteColumn, WritePixel, WriteRowEx
Writes a specified intensity value to a selected image pixel.
WritePixel(hImage As Long, xPos As Integer, yPos As Integer, pixelValue As Integer) As Long
No pixel conversion occurs in this function. This function does not operate on 24-bit color images.
<i>hImage</i> Specifies the handle of the image to be written to.
<i>xPos</i> Specifies the X-coordinate of the pixel where writing should start.
<i>yPos</i> Specifies the Y-coordinate of the pixel where writing should start.
<i>pixelValue</i> Provides the intensity value to be written to the selected pixel.

WritePixel (continued)	
Example	' Write a white pixel at location 50, 50 on the current image Dim <i>im</i> As Long <i>MM</i> .GetCurrentImage <i>im</i> <i>MM</i> .WritePixel <i>im</i> , 50, 50, 255
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), ReadPixel, WriteColumn, WriteRow

WriteRow - for Visual Basic 6 & earlier

Description Reads a set of intensity values from an array and writes them to a horizontal line of pixels in a selected image. Syntax WriteRow(hImage As Long, xPos As Integer, yPos As Integer, nPixels As Integer, depth As Integer, pixelBuffer() As Variant) As Long **Parameters** *hImage* Specifies the handle of the image to be written to. *xPos* Specifies the X-coordinate of the pixel where writing should start. *vPos* Specifies the Y-coordinate of the pixel where writing should start. *nPixels* Specifies the number of pixels to be written, starting at the position given by *xPos* and *yPos*. The size of the array will depend on the image bit-depth and the data type used for the array buffer. (See Section 1.6, Data Types and Arrays.) *depth* Specifies what the bit-depth of the pixel data values should be. (See Section 1.6, Data Types and Arrays.) *pixelBuffer()* This is a buffer from which the pixel values will be read. If *depth* is 8, pixelBuffer should be defined as Byte. If depth is 16, pixelBuffer should be Integer. If depth is 24, pixelBuffer should be Byte.

WriteRow (continued)	
Example	' Write a white horizontal line on the current image Dim <i>im</i> As Long Dim <i>buf</i> (50) As Byte Dim <i>x</i> As Integer
	For $x = 0$ To 49 buf(x) = 255 Next x
	MM. GetCurrentImage im MM. WriteRow im, 50, 0, 50, 8, buf
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), WriteColumn, WritePixel, WriteRowEx

WriteRowEx – for Visual Basic 6 & earlier WriteRowEx2 – for Visual Basic .NET (2002 – 2008)

Description	Reads a set of intensity values from a specified starting location in an array and writes them to a horizontal line of pixels in a selected image
Syntax	 WriteRowEx(hImage As Long, xPos As Integer, yPos As Integer, nPixels As Integer, depth As Integer, xStart As Integer, yStart As Integer, pixelBuffer() As Variant) As Long WriteRowEx2(hImage As Long, xPos As Integer, yPos As Integer, nPixels As Integer, depth As Integer, xStart As Integer, yStart As Integer, pixelBuffer() As Variant) As Variant) As Long

Parameters *hImage* Specifies the handle of the image to be written to.

xPos Specifies the X-coordinate of the pixel where writing should start.

yPos Specifies the Y-coordinate of the pixel where writing should start.

nPixels Specifies the number of pixels to be written, starting at the position given by xPos and yPos. The size of the array will depend on the image bit-depth and the data type used for the array buffer. (See Section 1.6, *Data Types and Arrays.*)

depth Specifies what the bit-depth of the pixel data values should be. (See Section 1.6, *Data Types and Arrays.*)

xStart Specifies the X-coordinate (row) of the position in the array where reading should start.

yStart Specifies the Y-coordinate (column) of the position in the array where reading should start.

pixelBuffer() This is a buffer from which the pixel values will be read. If *depth* is 8, *pixelBuffer* should be defined as *Byte*. If *depth* is 16, *pixelBuffer* should be *Integer*. If *depth* is 24, *pixelBuffer* should be *Byte*.

WriteRowEx2

(continued)

	' and write it back rotated 90 degrees clockwise Dim <i>im</i> As Long Dim x As Integer
	Dim <i>buf</i> (50, 50) As Byte
	MM.GetCurrentImage im
	<pre>For x = 0 To 49 MM.ReadRowEx im, 0, x, 50, 8, 0, x, buf MM.WriteColumnEx im, x + 50, 0, 50, 8, 0, x, buf Next x</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), WriteColumnEx, WritePixel, WriteRow

WriteText

Description	Writes a specified string of text onto an image.
Syntax	WriteText(hImage As Long, xPos As Integer, yPos As Integer, text As String, fillBackground As Boolean, foreColor As Integer, backColor As Integer) As Long
Parameters	hImage Specifies the handle of the destination image.
	<i>xPos</i> Specifies the X-coordinate where writing of the text is to start.
	<i>yPos</i> Specifies the Y-coordinate where writing of the text is to start.
	<i>text</i> Provides the string of text that will be written onto the image.
	<i>fillBackground</i> Selects whether the area behind the text is to be filled in with an opaque background of a grayscale value selected by <i>backColor</i> . If <i>fillBackground</i> is set to TRUE, a background "fill" will be used. If <i>fillBackground</i> is set to FALSE, no "fill" will be used, and <i>backColor</i> will be ignored.
	foreColor Specifies the grayscale value of the text.
	<i>backColor</i> Specifies the grayscale value of the "fill" behind the text on the image.
Example	' Write "Nerve Cell" at the top of the current image Dim <i>im</i> As Long <i>MM</i> .GetCurrentImage <i>im</i> <i>MM</i> .WriteText <i>im</i> , 50, 20, "Nerve Cell", TRUE, 255, 0
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3)

Chapter 7 – Working with Regions of Interest

7.1 Overview

Introduction	A <i>region of interest</i> , or <i>ROI</i> , is an area in an image window that you define and specify for subsequent processing and analysis. For the purposes of user programs, MetaMorph only uses rectangular regions. If you need to use elliptical, linear, or irregularly shaped regions, you will need to use a journal to manipulate them. This chapter presents the Visual Basic functions that you will need for creating and manipulating regions and for measuring and obtaining data from regions.	
	Торіс	See Page
	Creating and Removing Regions	97
	Finding Regions	99
	Reading and Manipulating Region Properties	103
	Reading Image Data from Regions	109

Introduction The following functions are used for creating and removing regions of interest. At the time you create an ROI, MetaMorph assigns it a handle, just as it assigns handles to functions and entire image windows. All of the remaining functions that involve regions in this chapter will need to be supplied with the region handle. If you need to obtain the handle after you have already created the region, you will need to use the **GetRegion** or **GetActiveRegion** functions which are covered in Section 7.3.

CreateRectRegion

Description	Creates a rectangular region of interest in a specified image. You can specify the size and placement of a region you create simply by dictating the locations of its four corners.
Syntax	CreateRectRegion (<i>hImage</i> As Long, <i>x1</i> As Integer, <i>y1</i> As Integer, <i>x2</i> As Integer, <i>y2</i> As Integer, <i>hRetRegion</i> As Long) As Long
Parameters	<i>hImage</i> Specifies the handle of the image in which the region is to be created.
	<i>x1</i> Specifies the X-coordinate of the upper left corner of the region.
	<i>y1</i> Specifies the Y-coordinate of the upper left corner of the region.
	x^2 Specifies the X-coordinate of the lower right corner of the region.
	y2 Specifies the Y-coordinate of the lower right corner of the region.
Return values	hRetRegion Returns the handle for the created region.
Example	' Create a recangular region on the current image Dim <i>im</i> As Long Dim r As Long <i>MM</i> . GetCurrentImage <i>im</i> <i>MM</i> . CreateRectRegion <i>im</i> , 10, 10, 50, 50, r
See also:	DestroyRegion , GetCurrentImage (Section 4.3), GetImage (Section 4.3), SetActiveRegion (Section 6.4)

DestroyRegion

Description	Removes the specified region of interest.
Syntax	DestroyRegion(hRegion As Long) As Long
Parameters	<i>hRegion</i> Specifies the handle of the region to be removed. The handle can be obtained with the GetRegion function.
Example	' Delete the active region on the current image, if there ' is one Dim <i>im</i> As Long Dim <i>r</i> As Long
	MM.GetCurrentImage im MM.GetActiveRegion im, r
	<pre>If r <> 0 Then MM.DestroyRegion r Else MM.PrintMsg "No active region on the current image" End If</pre>
See also:	GetRegion (Section 7.3), GetActive Region (Section 7.3)

Introduction	The functions in this section are essential for the control of procedures involving
	regions. These functions are used for finding a region's handle, the "indexing tag"
	that both Visual Basic and MetaMorph use to keep track of the region, as well as for
	determining how many regions have been defined and whether a handle is currently in
	use. The GetActiveRegion and GetRegion functions, which obtain a region's handle,
	are particularly important, and you will see these two functions referenced in the "See
	Also" lists throughout the rest of this chapter.

GetActiveRegion

Description	Obtains the handle number of the active region, if any, in a specified image.
Syntax	GetActiveRegion(hImage As Long, lRetActiveRegion As Long) As Long
Parameters	hImage Specifies the handle of the image.
Return values	<i>lRetActiveRegion</i> Returns the handle of the active region. If there is no active region, the value will be returned as 0. Note: An image can contain regions without any of them being active.
Example	' Delete the active region on the current image, if there ' is one Dim <i>im</i> As Long Dim <i>r</i> As Long
	MM.GetCurrentImage im MM.GetActiveRegion im, r
	<pre>If r <> 0 Then MM.DestroyRegion r Else MM.PrintMsg "No active region on the current image" End If</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetRegion, SetActiveRegion

GetNumberOfRegions

Description	Returns the number of regions that exist on a specified image.
Syntax	GetNumberOfRegions(hImage As Long, nRetNumberOfRegions As Integer) As Long
Parameters	hImage Specifies the handle of the image.
Return values	<i>nRetNumberOfRegions</i> Returns the number of regions that currently exist on the image.
Example	' Get the number of regions on the current image
	Dim <i>im</i> As Long Dim <i>rCount</i> As Integer
	MM.GetCurrentImage im
	<pre>MM.GetNumberOfRegions im, rCount MM.PrintMsg "There are " + Str(rCount) + " regions on the current image"</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3)

GetRegion

Description	Obtains the handle number of a specified region.
Syntax	GetRegion(hImage As Long, index As Integer, lRetRegion As Long) As Long
Remarks	All of the regions loaded in MetaMorph have an index number, from 0 to $(n - 1)$, where <i>n</i> is the number of loaded regions. Regions are numbered according to their order of creation; thus, the first region is 0, the next is 1, and so on.
Parameters	<i>hImage</i> Specifies the handle of the image. <i>index</i> Specifies the number of the region. (See Remarks.)
Return values	<i>lRetRegion</i> Returns the handle of the selected region. (Note: The handle is a number used by MetaMorph to carry out programmatic functions. This differs from the <i>index</i> number associated with the region.) If an error occurs, for example, if <i>index</i> is not a valid number, <i>lRetRegion</i> will contain 0.

7.3 Finding Regions, continued

GetRegion (continued)	
Example	' Get the first region on the current image Dim <i>im</i> As Long Dim r As Long <i>MM</i> . GetRegion <i>im</i> , 0, r
See also:	GetActiveRegion, GetCurrentImage (Section 4.3), GetImage (Section 4.3)

IsValidRegion

Description	Tests whether or not the region handle you pass is valid.
Syntax	IsValidRegion(hRegion As Long) As Long
Parameters	<i>hRegion</i> Specifies the handle of the region in question. If the handle value that you pass is not valid, IsValidRegion returns a value of "0". The function will return a nonzero value if the region handle is valid.
Example	' Check if some region is valid, and if it is, make it the ' active region. 'r' and 'im' are region and image handles ' that were declared and used previous to the code here.
	<pre>If MM.IsValidRegion(r) <> 0 Then MM.PrintMsg "Region is valid" MM.SetActiveRegion im, r Else MM.PrintMsg "Region is no longer valid" End If</pre>
See also:	GetRegion

SetActiveRegion

Description	Makes a selected region the active region.
Syntax	SetActiveRegion(hImage As Long, hRegion As Long) As Long
Parameters	<i>hImage</i> Specifies the handle of the image.<i>hRegion</i> Specifies the handle of the region to be made active.
Example	' Check if some region is valid, and if it is, make it the ' active region. 'r' and 'im' are region and image handles ' that were declared and used previous to the code here.
	<pre>If MM.IsValidRegion(r) <> 0 Then MM.PrintMsg "Region is valid" MM.SetActiveRegion im, r Else MM.PrintMsg "Region is no longer valid" End If</pre>
See also:	GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetRegion

7.4 Reading and Manipulating Region Properties

Introduction It is often necessary or useful to obtain information about a region, such as its width, height, area, perimeter, or location. The functions that follow can be used to read or configure such region properties, as well as to obtain the X and Y coordinates of the region's outline. One programming example is given for all of the region property "Get" functions (see Figure 7.1 on page 106).

GetRegionArea

Description	Obtains the area, expressed as the number of pixels, of a specified region of interest.
Syntax	GetRegionArea(hRegion As Long, lRetArea As Long) As Long
Parameters	<i>hRegion</i> Specifies the handle of the region in question.
Return values	<i>lRetArea</i> Returns the number of pixels in the region.
See also:	GetActiveRegion (Section 7.3), GetRegion (Section 7.3), GetRegionSize

GetRegionDistance

Description	Obtains the perimeter of a specified region of interest. The distance measurement will be reported in calibrated units.
Syntax	GetRegionDistance(hRegion As Long, dRetDistance As Double) As Long
Remarks	This function differs from RegionGetNumEdgePixels in its expression of the perimeter in calibrated units. RegionGetNumEdgePixels expresses the perimeter in terms of the number of pixels.
Parameters	<i>hRegion</i> Specifies the handle of the region in question.
Return values	<i>dRetDistance</i> Returns the perimeter of the region, in calibrated units.
See also:	GetActiveRegion (Section 7.3), GetRegion (Section 7.3), RegionGetNumEdgePixels
GetRegionPosition

Description	Obtains the on-screen position of a specified region of interest.
Syntax	GetRegionPosition(hRegion As Long, nRetXPos As Integer, nRetYPos As Integer) As Long
Parameters	<i>hRegion</i> Specifies the handle of the region in question.
Return values	<i>nRetXPos</i> Returns the starting (upper left) X-coordinate of the specified region. <i>nRetYPos</i> Returns the starting Y-coordinate of the specified region.
See also:	GetActiveRegion (Section 7.3), GetRegion (Section 7.3), GetRegionSize, SetRegionPosition

GetRegionSize

Description	Obtains the width and height of a specified region of interest.
Syntax	GetRegionSize(hRegion As Long, nRetXSize As Integer, nRetYSize As Integer) As Long
Parameters	<i>hRegion</i> Specifies the handle of the region in question.
Return values	<i>nRetXSize</i> Returns the width, in pixels, of the specified region.<i>nRetYSize</i> Returns the height, in pixels, of the specified region.
See also:	GetActiveRegion (Section 7.3), GetRegion (Section 7.3), GetRegionArea, GetRegionPosition, SetRegionSize

RegionGetNumEdgePixels

Description	Obtains the perimeter of a specified region of interest. The distance measurement will be reported in terms of the number of pixels.
Syntax	RegionGetNumEdgePixels(hRegion As Long, lNumPixels As Long) As Long
Remarks	RegionGetNumEdgePixels returns in <i>lNumPixels</i> the number of pixels comprising the edge of the given region. This function will usually be used in conjunction with RegionGetEdgePixelCoordinates . You would call RegionGetNumEdgePixels first to determine how big to make the arrays you pass to RegionGetEdgePixelCoordinates .
Parameters	<i>hRegion</i> Specifies the handle of the region in question.
Return values	<i>lNumPixels</i> Returns the perimeter of the region, in pixels.
See also:	GetActiveRegion (Section 7.3), GetRegion (Section 7.3), GetRegionDistance

RegionGetEdgePixelCoordinates – for Visual Basic 6 & earlier RegionGetEdgePixelCoordinatesEx2 – for Visual Basic .NET (2002 – 2008)

Description	Obtains the X and Y coordinates of all the pixels lying under the edge of a region.
Syntax	RegionGetEdgePixelCoordinates(hRegion As Long, lBufferSize As Long, aRetXBuffer() As Integer, aRetYBuffer() As Integer) As Long RegionGetEdgePixelCoordinatesEx2(hRegion As Long, lBufferSize As Long, aRetXBuffer() As Integer, aRetYBuffer() As Integer) As Long
Parameters	hRegion Specifies the handle of the region in question.
	<i>lBufferSize</i> Sets the size of the array into which the X and Y coordinates will be read. You can use RegionGetNumEdgePixels to determine how many pixels that should be.
Return values	<i>aRetXBuffer()</i> Returns the X-coordinates of the pixels lying under the edge of the region. The size of this array will be determined by the number of pixels in the edgelist.
	<i>aRetYBuffer()</i> Returns the Y-coordinates of the pixels lying under the edge of the region. The size of this array will be determined by the number of pixels in the edgelist.
See also:	GetActiveRegion (Section 7.3), GetRegion (Section 7.3), RegionGetNumEdgePixels
	Continued on next page

7.4 Reading and Manipulating Region Properties, continued

Figure 7.1 Region Property "Get" Function Programming Example

```
' Get property information about the active region
Dim im As Long
Dim r As region
MM.GetCurrentImage im
MM.GetActiveRegion im, r
Dim area As Long
MM.GetRegionArea r, area
MM.PrintMsg "Pixel area of the active region is " + Str(area)
Dim distance As Double
MM.GetRegionDistance r, distance
MM.PrintMsg "Calibrated perimeter of the active region is "
  + Str(distance)
Dim x As Integer, y As Integer
MM.GetRegionPosition r, x, y
Dim dx As Integer, dy As Integer
MM.GetRegionSize r, dx, dy
MM.PrintMsg "Position of the region is " + Str(x) + ", "
  + Str(y)
MM.PrintMsg "Size of the region is " + Str(dx) + ", "
  + Str(dy)
Dim numpix As Long
' First find out how many pixels are in the edge of
' the region
MM.RegionGetNumEdgePixels r, numpix
' Then get the coordinates of all the edge pixels
Dim x(numpix) As Integer, y(numpix) As Integer
MM.RegionGetEdgePixelCoordinates r, numpix, x, y
```

7.4 Reading and Manipulating Region Properties, continued

SetRegionPosition

Description	Moves a selected region of interest to a specified position.
Syntax	SetRegionPosition(hRegion As Long, xPos As Integer, yPos As Integer) As Long
Parameters	hRegion Specifies the handle of the region in question.
	xPos Specifies the X-coordinate of the new starting point (upper left corner) of the selected region.
	<i>yPos</i> Specifies the Y-coordinate of the new starting point of the selected region.
Example	' Move the active region to the upper left corner of the image ' and make it 100 x 100 pixels in size
	Dim <i>im</i> As Long Dim <i>r</i> As Long
	MM.GetCurrentImage im MM.GetActiveRegion im, r
	MM.SetRegionPosition r, 0, 0 MM.SetRegionSize r, 100, 100
See also:	GetRegion (Section 7.3), GetRegionPosition

SetRegionSize

Description	Resizes a selected region of interest to a specified width and height.
Syntax	SetRegionSize(hRegion As Long, xSize As Integer, ySize As Integer) As Long
Parameters	<i>hRegion</i> Specifies the handle of the region in question.<i>xSize</i> Specifies the new width, in pixels, of the selected region.<i>ySize</i> Specifies the new height, in pixels, of the selected region.

Reading and Manipulating Region Properties, continued 7.4

SetRegionSize (continued)

See also:	GetActiveRegion (Section 7.3), GetRegion (Section 7.3), GetRegionSize
	MM.SetRegionPosition r, 0, 0 MM.SetRegionSize r, 100, 100
	MM.GetCurrentImage im MM.GetActiveRegion im, r
	Dim <i>im</i> As Long Dim <i>r</i> As Long
Example	' Move the active region to the upper left corner of the image ' and make it 100 x 100 pixels in size

SwapRegionNumbers

Description	Exchanges the numbers associated with two selected regions.
Syntax	SwapRegionNumbers(rhRegion1 As Long, rhRegion2 As Long)
Parameters	<i>rhRegion1</i> Specifies the handle of one of the two selected regions.<i>rhRegion2</i> Specifies the handle of the other selected region.
See also:	GetActiveRegion (Section 7.3), GetRegion (Section 7.3)

7.5 Reading Image Data from Regions

Introduction This section describes a number of functions which are invaluable for obtaining grayscale information from a region of interest. You can derive the minimum, maximum, and average intensity value in a region, as well as the standard deviation around the mean. One programming example is given for all of the region data "Get" functions (see Figure 7.2 on page 112). You will need to apply the **MeasureRegion** function as a preliminary step to obtaining these data. If your images have been calibrated in MetaMorph with the Calibrate Distances or Calibrate Gray Levels commands, the "read" functions in this section will return data expressed in calibrated units, rather than as numbers of pixels or as grayscale levels.

GetRegionAverageValue

See also:	GetActive Region (Section 7.3), GetRegion (Section 7.3), GetRegionMaximumValue, GetRegionMinimumValue, GetRegionStdDeviation, MeasureRegion
Return values	<i>dRetAverageValue</i> Returns the average intensity value from the region.
Parameters	hRegion Specifies the handle of the region.
Remarks	Before you apply this function, the region must first be measured with the MeasureRegion function.
Syntax	GetRegionAverageValue(hRegion As Long, dRetAverageValue As Double) As Long
Description	Obtains the average of all of the grayscale values in a specified region of interest.

GetRegionMinimumValue

Description	Obtains the lowest grayscale value in a specified region of interest.
Syntax	GetRegionMinimumValue(hRegion As Long, dRetMinValue As Double) As Long
Remarks	Before you apply this function, the region must first be measured with the MeasureRegion function.
Parameters	hRegion Specifies the handle of the region.
Return values	<i>dRetMinValue</i> Returns the minimum intensity value in the region.
See also:	GetActive Region (Section 7.3), GetRegion (Section 7.3), GetRegionAverageValue, GetRegionMaximumValue, GetRegionStdDeviation, MeasureRegion
	- · · ·

GetRegionMaximumValue

See also:	GetActive Region (Section 7.3), GetRegion (Section 7.3), GetRegionAverageValue, GetRegionMinimumValue, GetRegionStdDeviation, MeasureRegion
Return values	<i>dRetMaxValue</i> Returns the maximum intensity value in the region.
Parameters	hRegion Specifies the handle of the region.
Remarks	Before you apply this function, the region must first be measured with the MeasureRegion function.
Syntax	GetRegionMaximumValue(hRegion As Long, dRetMaxValue As Double) As Long
Description	Obtains the highest grayscale value in a specified region of interest.

GetRegionStdDeviation

Description	Obtains the standard deviation of the grayscale values in a specified region of interest.
Syntax	GetRegionStdDeviation(hRegion As Long, dRetStdDeviation As Double) As Long
Remarks	Before you apply this function, the region must first be measured with the MeasureRegion function.
Parameters	<i>hRegion</i> Specifies the handle of the region.
Return values	<i>dRetStdDeviation</i> Returns the standard deviation of the intensity values in the region.
See also:	GetActive Region (Section 7.3), GetRegion (Section 7.3), GetRegionAverageValue, GetRegionMaximumValue, GetRegionMinimumValue, MeasureRegion

GetRegionThresholdArea

Description	Obtains the number of pixels in a specified region that are inside the current threshold.
Syntax	GetRegionThresholdArea(hRegion As Long, lRetArea As Long) As Long
Remarks	Before you apply this function, the region must first be measured with the MeasureRegion function.
Parameters	<i>hRegion</i> Specifies the handle of the region.
Return values	<i>lRetArea</i> Returns the number of thresholded pixels in the region.
See also:	GetActive Region (Section 7.3), GetRegion (Section 7.3), MeasureRegion

MeasureRegion

Description	Measures the grayscale data in a specified region of interest of an image.
Syntax	MeasureRegion (<i>hRegion</i> As Long, <i>hImage</i> As Long, <i>bUseThreshold</i> As Boolean) As Long
Parameters	hRegion Specifies the handle of the region.
	hImage Specifies the handle of the image.
	<i>bUseThreshold</i> Determines whether the measurement will be made of just those pixels that are thresholded, or if all pixels are to be measured. If only thresholded pixels are to be measured, <i>bUseThreshold</i> should be set to TRUE. If all pixels are to be measured, set <i>bUseThreshold</i> to FALSE.
See also:	GetActive Region (Section 7.3), GetCurrentImage (Section 4.3), GetImage (Section 4.3), GetRegion (Section 7.3)

7.5 Reading Image Data from Regions, continued

Figure 7.2Region Data "Get" Function Programming Example

```
' Measure the active region on the current image and print
' out the information that was measured
Dim im As Long
Dim r As region
MM.GetCurrentImage im
MM.GetActiveRegion r
MM.MeasureRegion r, im, FALSE ' Do not use threshold
Dim average As Double
Dim min As Double
Dim max As Double
Dim stddev As Double
Dim area As Long
MM.GetRegionAverageValue r, average
MM.GetRegionMinimumValue r, min
MM.GetRegionMaximumValue r, max
MM.GetRegionStdDeviation r, stddev
MM.GetRegionThresholdArea r, area
MM.PrintMsg "Average is " + Str(average)
MM.PrintMsg "Minimum is " + Str(min)
MM.PrintMsg "Maximum is " + Str(max)
MM.PrintMsg "Standard deviation is " + Str(stddev)
MM. PrintMsg "Threshold area should be 0, because the
  region was"
MM. PrintMsg "Measured without thresholding. The area is: "
  + Str(area)
```

Chapter 8 – Performing Morphometry

8.1 Overview

Introduction	As the premier morphometric analysis system, MetaMorph offers an extensive array of functions for measurement and analysis of image objects. This chapter discusses the Visual Basic functions that you will need for configuring your morphometric measurements, creating classifier filters, and measuring objects. The operative function that measures your image, MorphMeasureObjects , is covered in Section 8.5. Before you can apply this function, however, please remember to define a threshold range, as detailed in Section 6.2, <i>Applying Thresholding</i> .		
In this chapter	This chapter contains the following topics:		
	Торіс	See Page	
	Configuring Measurement Preferences	114	
	Configuring Object Measurements	115	
	Configuring Classifier Filters	119	
	Measuring All Objects in an Image	121	
	Measuring Single Objects	126	

8.2 Configuring Measurement Preferences

Introduction Before you measure your images, you may want to configure MetaMorph's behavior during the measurement procedure. The settings you use for the two functions that follow will determine two types of behavior: whether objects that "fail" a classifier filter (see Section 8.4) are redrawn in the result image, and whether "holes" (unthresholded areas in the middle of an object) are "filled" prior to measurement.

MorphSetDrawFailedObjectsFlag

DescriptionSets whether objects that fail to pass any classifiers should be drawn in the resultant
measurement image.SyntaxMorphSetDrawFailedObjectsFlag(bRedraw As Boolean) As LongParametersbRedraw
Determines whether failed objects will be drawn. If bRedraw is set to
TRUE, failed objects will be drawn. If bRedraw is set to FALSE, failed objects will
not be drawn.Example' The next time measurements are performed, don't draw objects
' that don't pass any classifiers and fill holes in objects
MM.MorphSetDrawFailedObjectsFlag FALSE
MM.MorphSetFillHoleFlag TRUE

MorphSetFillHoleFlag

Description	Sets whether "holes" (unthresholded areas enclosed within thresholded areas) will be filled in (treated as thresholded) when measurement occurs.
Syntax	MorphSetFillHoleFlag(bRedraw As Boolean) As Long
Remarks	If <i>bRedraw</i> is TRUE, they will be filled in. If FALSE, they will not be.
Parameters	<i>bRedraw</i> Determines whether or not holes are to be filled. If <i>bRedraw</i> is set to TRUE, holes will be filled and subsequent measurements will include their areas. If <i>bRedraw</i> is set to FALSE, holes will not be filled.
Example	' The next time measurements are performed, don't draw objects ' that don't pass any classifiers and fill holes in objects MM.MorphSetDrawFailedObjectsFlag FALSE MM.MorphSetFillHoleFlag TRUE

Introduction Just as image windows, functions, and regions are manipulated by their handles, so too are object parameters manipulated by using an index number, the *parameter number*. Several configuration and measurement functions depend on the correct parameter number being passed to them. In this section, we discuss several functions that allow you to obtain a parameter's number, or to obtain other information about a parameter based on the number you pass. In addition, another measurement configuration function, **MorphSetupMeasurements**, is used for specifying the image to be measured and a second image, the mask image, which will be used to define the distribution of the thresholding overlay.

MorphFindParmIndex

Description	Obtains the number of a specified parameter.	
Syntax	MorphFindParmIndex(sParmName As String, nRetParmNumber As Integer) As Long	
Parameters	<i>sParmName</i> Specifies the name of the parameter for which you want to obtain a parameter number. The parameter names that you can use consist of those used for the MetaMorph morphometry functions. You can see these names in the Configure Object Classifiers, Configure Object Measurements, or Integrated Morphometry Analysis dialog boxes.	
Return values	<i>nRetParmNumber</i> Returns the number of the parameter specified by <i>sParmName</i> . If no parameter by that name exists, <i>nRetParmNumber</i> will be returned as -1.	
Example	' Find the width of object 5. This code assumes that ' measurements have already been made. Dim I As Integer Dim <i>nWidth</i> As Single <i>MM</i> .MorphFindParmIndex "Width", <i>n</i> <i>MM</i> .MorphGetParmMeasurement 5, <i>n</i> , <i>nWidth</i> <i>MM</i> .PrintMsg "The width of object 5 is " + Str(<i>nWidth</i>)	
See also:	MorphGetParmDescription, MorphGetParmMeasurement (Section 8.6), MorphGetParmName	

MorphGetNumberOfParms

Description	Obtains the total number of measurable parameters.
Syntax	MorphGetNumberOfParms(nRetParms As Integer) As Long
Return values	<i>nRetParms</i> Returns the total number of measurable parameters.
Example	' Print out the names and descriptions of all the measurement ' parameters
	Dim <i>nParms</i> As Integer Dim <i>i</i> As Integer Dim <i>parmName</i> As String Dim <i>desc</i> As String
	<pre>MM.MorphGetNumberOfParms nParms For i = 0 To nParms - 1 MM.MorphGetParmName i, parmName MM.MorphGetParmDescription i, desc MM.PrintMsg parmName + ": " + desc Next i</pre>

MorphGetParmDescription

Description	Obtains a description of a parameter for which you have a parameter number.
Syntax	MorphGetParmDescription(<i>nParmNumber</i> As Integer, <i>sRetParmDescription</i> As String) As Long
Parameters	<i>nParmNumber</i> Specifies the number of the parameter for which you want to obtain a description. This number can be obtained with the MorphFindParmIndex function.
Return values	sRetParmDescription Returns a textual description of the parameter in question.
	Continued on next page

8.3 Configuring Object Measurements, continued

MorphGetParmDescription

(continued)

See also:	MorphFindParmIndex, MorphGetParmName
	<pre>MM.MorphGetNumberOfParms nParms For i = 0 To nParms - 1 MM.MorphGetParmName i, parmName MM.MorphGetParmDescription i, desc MM.PrintMsg parmName + ": " + desc Next i</pre>
	Dim <i>nParms</i> As Integer Dim <i>i</i> As Integer Dim <i>parmName</i> As String Dim <i>desc</i> As String
Example	' Print out the names and descriptions of all the measurement ' parameters

MorphGetParmName

Description	Obtains the name of a parameter for which you have a parameter number.
Syntax	MorphGetParmName(nParmNumber As Integer, sRetParmName As String) As Long
Parameters	<i>nParmNumber</i> Specifies the number of the parameter for which you want to obtain a description. This number can be obtained with the MorphFindParmIndex function.
Return values	<i>sRetParmName</i> Returns the name of the parameter in question.
Example	' Print out the names and descriptions of all the measurement ' parameters
	Dim <i>nParms</i> As Integer Dim <i>i</i> As Integer Dim <i>parmName</i> As String Dim <i>desc</i> As String
	<pre>MM.MorphGetNumberOfParms nParms For i = 0 To nParms - 1 MM.MorphGetParmName i, parmName MM.MorphGetParmDescription i, desc MM.PrintMsg parmName + ": " + desc Next i</pre>
See also:	MorphFindParmIndex, MorphGetParmDescription

MorphSetupMeasurements

Description	Selects images to be used as measurement and mask images for subsequent measurement with MorphMeasureObjects .
Syntax	MorphSetupMeasurements(hGrayImage As Long, hMaskImage As Long, fStdArea As Single) As Long
Remarks	This function selects the measurement image and mask image for use in subsequent measurements with MorphMeasureObjects . Also, if you want to count objects, this function allows you to define and make measurements with a <i>standard area</i> . Counting objects can sometimes be difficult because they may overlap or are poorly defined in the image. A standard area is a value that you believe represents the area of a standard object, based on the assumption that the objects being measured are of a fairly uniform size. The clumps of objects will be counted by dividing the area of the clump by the standard area.
	If you are not interested in using a standard area, you can set <i>fStdArea</i> to 1.
Parameters	<i>hGrayImage</i> Specifies the handle of the image to be used for measurement. Image handles can be obtained with GetImage .
	<i>hMaskImage</i> Specifies the handle of the image to be used as a mask image. Typically, <i>hMaskImage</i> is an image created by setting a threshold on <i>hGrayImage</i> and then using BinarizeImage to create a binary image from that.
	<i>fStdArea</i> Defines the size of the standard area. The units used depends on whether you have calibrated the image for distance.
Example	' Create a binary mask of the current image and then measure ' the image Dim <i>im</i> As Long
	' Put a threshold on the current image MM.GetCurrentImage im MM.SetThresholdState im, 1 MM.SetThresholdRange im, 50, 150
	' Create the binarized image Dim <i>bin</i> As Long <i>MM.</i> CreateImage 512, 512, 1, "binary image", <i>bin</i> <i>MM.</i> BinarizeImage <i>im, bin</i>
	' Measure MM.MorphSetupMeasurements im, bin, 1# MM.MorphMeasureObjects TRUE
See also:	BinarizeImage (Section 6.3), GetCurrentImage (Section 4.3), GetImage (Section 4.3), MorphMeasureObjects (Section 8.5)

Introduction Classifier filters are morphometric measurement ranges through which an object's measurements must "pass" to be included in the final set. Using a classifier filter, you can restrict your measurements to just those objects that meet your set criteria, while excluding other classes of objects. The two functions which follow are used for reading the current settings of a classifier filter for a specified parameter and for configuring those settings.

MorphGetFilter

Description	Obtains the settings of the classifier filter for a specified parameter.
Syntax	MorphGetFilter (<i>nClassifierNumber</i> As Integer, <i>nParmNumber</i> As Integer, <i>fRetMinVal</i> As Single, <i>fRetMaxVal</i> As Single, <i>bRetInclusive</i> As Boolean, <i>bRetEnableFilter</i> As Boolean) As Long
Parameters	<i>nClassifierNumber</i> Specifies the number of the classifier filter in question $(0 - 7)$. This classifier is one you will have defined, and the number is one you will have assigned with the MorphSetFilter function.
	<i>nParmNumber</i> Specifies the parameter number of the classifier for which the filter is being configured. This number can be obtained with the MorphFindParmIndex function. These predefined classifiers are those that correspond to the parameters used for the MetaMorph morphometry functions. You can see these parameters in the Configure Object Classifiers, Configure Object Measurements, or Integrated Morphometry Analysis dialog boxes.
Return values	<i>fRetMinVal</i> Returns the minimum value of the filter.
	<i>fRetMaxVal</i> Returns the maximum value of the filter.
	<i>bRetInclusive</i> Indicates whether the range minimum and maximum are inclusive (that is, the filter "passes" objects if they have values between the minimum and maximum) or exclusive (that is, the filter "passes" objects if they have values equal to or outside the minimum and maximum). If the filter has been set to the inclusive state, a <i>bRetInclusive</i> value of TRUE will be returned. If the filter is exclusive, a value of FALSE will be returned.
	<i>bRetEnableFilter</i> Indicates whether the filter is active or not. If the filter is active, <i>bRetEnableFilter</i> will return with a value of TRUE. If the filter is inactive, a value of FALSE will be returned.

MorphGetFilter (continued)	
Example	<pre>' Tell the first filter to stop filtering on total area, while ' leaving the filtering parameters unchanged. Dim min As Single, max As Single Dim inclusive As Boolean, enable As Boolean MM.MorphGetFilter 0, 0, min, max, inclusive, enable MM.MorphSetFilter 0, 0, min, max, inclusive, FALSE</pre>
See also:	MorphFindParmIndex (Section 8.3), MorphSetFilter

MorphSetFilt	er
--------------	----

Description	Configures the settings of a classifier filter for a specified parameter.
Syntax	MorphSetFilter (<i>nClassifierNumber</i> As Integer, <i>nParmNumber</i> As Integer, <i>fMinVal</i> As Single, <i>fMaxVal</i> As Single, <i>bInclusive</i> As Boolean, <i>bEnableFilter</i> As Boolean) As Long
Parameters	<i>nClassifierNumber</i> Specifies a number for the classifier $(0 - 7)$. Typically, this number will be based on the order of configuration.
	<i>nParmNumber</i> Specifies the parameter number of the classifier for which the filter is being configured. This number can be obtained with the MorphFindParmIndex function.
	fMinVal Specifies the minimum value of the filter.
	fMaxVal Specifies the maximum value of the filter.
	<i>bInclusive</i> Specifies whether the range minimum and maximum are to be inclusive (that is, the filter will "pass" objects that have values between the minimum and maximum) or exclusive (that is, the filter will "pass" objects that have values equal to or outside the minimum and maximum). To set the filter to the inclusive state, assign a value of TRUE to <i>bInclusive</i> . To make the filter exclusive, assign a value of FALSE.
	<i>bEnableFilter</i> Specifies whether the filter will be active or inactive. To activate the filter, set <i>bEnableFilter</i> to TRUE. To deactivate the filter, assign a value of FALSE.
Example	' Set the first filter to only pass objects whose total area ' is between 10 and 50. MM.MorphSetFilter 0, 0, 10, 50, TRUE, TRUE
See also:	MorphFindParmIndex (Section 8.3), MorphGetFilter

8.5 Measuring All Objects in an Image

Introduction MetaMorph automatically measures all objects in an image, but the resulting data are managed either as a group or on an object-by-object basis. Data for the entire group are presented as a statistical summary, and can be saved in a summary log. The functions in this section deal with obtaining and managing group measurement data. This includes the important **MorphMeasureObjects** function. (For a discussion of single object measurement functions, see Section 8.6.)

MorphDeleteObject

See also:	MorphRecalc
Example	' Delete object 4 from the morph summary data. This code ' assumes measurements have already been performed. MM.MorphDeleteObject 4 MM.MorphRecalc
Parameters	<i>nObjectID</i> Specifies the number of the object to be deleted. Objects are numbered from 0 to $(n - 1)$, where <i>n</i> is the total number of objects.
Remarks	You can prevent a specific object from having a contribution to the summary data by using this function in conjunction with MorphRecalc .
Syntax	MorphDeleteObject(nObjectID As Integer) As Long
Description	Removes a specified object from the list of measured objects.

MorphGetNumberOfObjects

Description	Obtains the number of objects that were last measured by MorphMeasureObjects.
Syntax	MorphGetNumberOfObjects(nRetObjects As Integer) As Integer
Return values	<i>nRetObjects</i> Returns the number of objects that were measured.
See also:	MorphMeasureObjects

MorphGetParmAccumSummary

Description	Obtains the cumulative statistics for all objects that passed a specified classifier filter for a selected parameter. This function returns the same information as MorphGetParmSummary , but combines the information with all previous measurements that have been made using the selected classifier filter.
Syntax	MorphGetParmAccumSummary (<i>nClassifierNumber</i> As Integer, <i>nParmNumber</i> As Integer, <i>lRetCount</i> As Long, <i>fRetAverage</i> As Single, <i>fRetMinVal</i> As Single, <i>fRetMaxVal</i> As Single, <i>fRetStdDeviation</i> As Single, <i>fRetTotal</i> As Single) As Long
Parameters	<i>nClassifierNumber</i> Specifies the classifier number $(0 - 7)$. This number is one that you will have assigned with the MorphSetFilter function.
	<i>nParmNumber</i> Specifies the parameter number of the classifier filter. This number can be obtained with the MorphFindParmIndex function.
Return values	<i>lRetCount</i> Returns the total number of objects which passed the classifier.
	<i>fRetAverage</i> Returns the average value of the given parameter for all passed objects.
	<i>fRetMinVal</i> Returns the lowest value of the parameter for all objects.
	<i>fRetMaxVal</i> Returns the highest value of the parameter for all objects.
	<i>fRetStdDeviation</i> Returns the standard deviation of the parameter for all objects.
	<i>fRetTotal</i> Returns the sum of the values for all objects.
Example	<pre>' Get the parameter summary and the accumulated summary data of ' the first filter for total area MM.MorphGetParmSummary 0, 0, count, average, min, max, stddev, total</pre>
	Dim <i>count</i> As Long Dim <i>average</i> As Single, <i>min</i> As Single, <i>max</i> As Single Dim <i>stddev</i> As Single, <i>total</i> As Single
	<pre>MM.PrintMsg "The count is " + Str(count) MM.PrintMsg "The average is " + Str(average) MM.PrintMsg "The min is " + Str(min) MM.PrintMsg "The max is " + Str(max) MM.PrintMsg "The standard deviation is " + Str(stddev) MM.PrintMsg "The total is " + Str(total)</pre>
	MM.MorphGetParmAccumSummary 0, 0, count, average, min, max, stddev, total

8.5 Measuring All Objects in an Image, continued

MorphGetParmAccumSummary (continued) MM.PrintMsg "The accumulated count is " + Str(count) MM.PrintMsg MM.PrintMsg "The accumulated average is " + Str(average) MM.PrintMsg MM.PrintMsg "The accumulated min is " + Str(min) MM.PrintMsg MM.PrintMsg "The accumulated max is " + Str(max) MM.PrintMsg MM.PrintMsg "The accumulated standard deviation is " + Str(stddev) MM.PrintMsg See also: MorphFindParmIndex (Section 8.3), MorphGetParmSummary, MorphSetFilter (Section 8.4)

MorphGetParmSummary

Description	Obtains the statistics for all objects that passed a specified classifier filter for a selected parameter.
Syntax	MorphGetParmSummary (<i>nClassifierNumber</i> As Integer, <i>nParmNumber</i> As Integer, <i>lRetCount</i> As Long, <i>fRetAverage</i> As Single, <i>fRetMinVal</i> As Single, <i>fRetMaxVal</i> As Single, <i>fRetStdDeviation</i> As Single, <i>fRetTotal</i> As Single) As Long
Parameters	<i>nClassifierNumber</i> Specifies the classifier number $(0 - 7)$. This number is one that you will have assigned with the MorphSetFilter function.
	<i>nParmNumber</i> Specifies the parameter number of the classifier filter. This number can be obtained with the MorphFindParmIndex function.
Return values	 <i>lRetCount</i> Returns the total number of objects which passed the classifier. <i>fRetAverage</i> Returns the average value of the given parameter for all passed objects. <i>fRetMinVal</i> Returns the lowest value of the parameter for all objects. <i>fRetMaxVal</i> Returns the highest value of the parameter for all objects. <i>fRetStdDeviation</i> Returns the standard deviation of the parameter for all objects. <i>fRetTotal</i> Returns the sum of the values for all objects.
	<i>fRetTotal</i> Returns the sum of the values for all objects.

8.5 Measuring All Objects in an Image, continued

MorphGetParmSummary (continued) Example ' Get the parameter summary and the accumulated summary data of ' the first filter for total area MM.MorphGetParmSummary 0, 0, count, average, min, max, stddev, total Dim count As Long Dim average As Single, min As Single, max As Single Dim stddev As Single, total As Single MM. PrintMsg "The count is " + Str(count) MM.PrintMsg "The average is " + Str(average) MM.PrintMsg "The min is " + Str(min) MM.PrintMsg "The max is " + Str(max) MM. PrintMsg "The standard deviation is " + Str(stddev) MM. PrintMsg "The total is " + Str(total) MM.MorphGetParmAccumSummary 0, 0, count, average, min, max, stddev, total MM. **PrintMsg** "The accumulated count is " + Str(count) MM.PrintMsg "The accumulated average is " + Str(average) MM. **PrintMsg** "The accumulated min is " + Str(min) MM. PrintMsg "The accumulated max is " + Str(max) MM.PrintMsg "The accumulated standard deviation is " + Str(stddev) MM. PrintMsg "The accumulated total is " + Str(total) See also: MorphFindParmIndex (Section 8.3), MorphGetParmAccumSummary, **MorphSetFilter** (Section 8.4)

MorphMeasureObjects

DescriptionMeasures the hGrayImage image that was last specified by
MorphSetupMeasurements.SyntaxMorphMeasureObjects(bMeasureAll As Boolean) As LongParametersbMeasureAll
basis only or as a whole. If you set bMeasureAll to TRUE, all of the objects in the
image will be measured and their data made available to the various summary data
functions, as well as to the object data functions. If you set this variable to FALSE,
data will be passed only to the object data functions.

8.5 Measuring All Objects in an Image, continued

MorphMeasureObjects (continued) Example ' Create a binary mask of the current image and then measure ' the image Dim im As Long ' Put a threshold on the current image MM.GetCurrentImage im MM.SetThresholdState im, 1 MM.SetThresholdRange im, 50, 150 ' Create the binarized image Dim *bin* As Long MM.CreateImage 512, 512, 1, "binary image", bin MM.BinarizeImage im, bin ' Measure MM.MorphSetupMeasurements im, bin, 1# MM.MorphMeasureObjects TRUE See also: MorphSetupMeasurements (Section 8.3)

MorphRecalc	
Description	Recomputes the statistical summary data for the current list of objects to be measured.
Syntax	MorphRecalc() As Long
Remarks	If any objects have been deleted (using MorphDeleteObject), those objects will be eliminated from the measurement of summary data.
Example	' Delete object 4 from the morph summary data. This code ' assumes measurements have already been performed. MM.MorphDeleteObject 4 MM.MorphRecalc
See also:	MorphDeleteObject

8.6 Measuring Single Objects

Introduction	Data for individual objects are measured for all objects in an image simultaneously,
	but can be obtained and used on an object-by-object basis. This section describes the
	functions that you can use to obtain measurement data for an individual object. (For a
	discussion of group object measurement functions, see Section 8.5.) One
	programming example is given for all of the single object data "Get" functions (see
	Figure 8.1 on page 132).

MorphGetCentroid

Description	Obtains the coordinates of the centroid (point that represents the center of mass) of a specified object. The centroid will be given in fractional pixel coordinates.
Syntax	MorphGetCentroid (<i>nObjectID</i> As Integer, <i>fRetX</i> As Single, <i>fRetY</i> As Single) As Long
Remarks	This function differs from MorphGetCentroidPixel in its use of fractional coordinates. MorphGetCentroidPixel expresses coordinates as integer values by locating the pixel nearest to the "fractional" centroid.
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
Return values	<i>fRetX</i> Returns the X-coordinate of the object centroid.<i>fRetY</i> Returns the Y-coordinate of the object centroid.
See also:	MorphGetCentroidPixel

MorphGetCentroidPixel

Description	Obtains the coordinates of the pixel nearest to the centroid (the point that represents the center of mass) of a specified object.
Syntax	MorphGetCentroidPixel (<i>nObjectID</i> As Integer, <i>nRetX</i> As Integer, <i>nRetY</i> As Integer) As Long
Remarks	This function differs from MorphGetCentroid in that it expresses coordinates as integer values by locating the pixel nearest to the "fractional" centroid. MorphGetCentroid expresses the exact location of the centroid using fractional coordinates.
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.

8.6 Measuring Single Objects, continued

MorphGetCentroidPixel (continued)	
Return values	<i>fRetX</i> Returns the X-coordinate of the pixel nearest to the object centroid.<i>fRetY</i> Returns the Y-coordinate of the pixel nearest to the object centroid.
See also:	MorphGetCentroid

MorphGetClassifiersPassed

Description	Obtains the number of classifier filters that a specified object has passed.
Syntax	MorphGetClassifiersPassed(nObjectID As Integer, nRetClassifiersPassed As Integer) As Long
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
Return values	<i>nRetClassifiersPassed</i> Returns the number of classifier filters that the object has passed.
Example	' Find out how many classifiers object 9 passed Dim <i>n</i> As Integer <i>MM</i> . MorphGetClassifiersPassed 9, <i>n</i>

MorphGetClosestObject

Description	Obtains the number of the object closest to a specified set of X and Y coordinates.
Syntax	MorphGetClosestObject (<i>nX</i> As Integer, <i>nY</i> As Integer, <i>nRetObjectID</i> As Integer) As Long
Parameters	<i>nX</i> Specifies the X-coordinate.<i>nY</i> Specifies the Y-coordinate.
Return values	<i>nRetObjectID</i> Returns the object number of the object closest to the specified coordinates.
Example	' Get the object number of the object nearest to coordinate ' 100, 100 Dim <i>obj</i> As Integer MM.MorphGetClosestObject 100, 100, <i>obj</i>

MorphGetInternalPoint

Description	Obtains the set of coordinates of some pixel within a specified object.
Syntax	MorphGetInternalPoint (<i>nObjectID</i> As Integer, <i>nRetX</i> As Integer, <i>nRetY</i> As Integer) As Long
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
Return values	<i>nRetX</i> Returns the X-coordinate of a randomly selected pixel in the object.<i>nRetY</i> Returns the Y-coordinate of the same randomly selected pixel.

MorphGetMeasurements – for Visual Basic 6 & earlier MorphGetMeasurementsEx2 – for Visual Basic .NET (2002 – 2008)

Description	Fills a defined array with all measurement data for a specified object.
Syntax	MorphGetMeasurements(<i>nObjectID</i> As Integer, <i>aMeasurements()</i> As Single) As Long MorphGetMeasurementsEx2(<i>nObjectID</i> As Integer, <i>aMeasurements()</i> As Single) As Long
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
Return values	<i>aMeasurements()</i> Defines an array into which the object's measurement data will be read. The array should have at least the number of elements returned by MorphGetNumberOfParms .
Example	' Get all the measurements for object number 12. This code ' assumes measurements have already been made. Dim <i>n</i> As Integer <i>MM</i> . MorphGetNumberOfParms <i>nParms</i> Dim <i>measurements(nParms)</i> As Single <i>MM</i> . MorphGetMeasurements 12, <i>measurements</i>
See also:	MorphGetNumberOfParms (Section 8.3), MorphMeasureObjects (Section 8.5)

MorphGetNumberOfRuns

Description	Obtains the number of horizontal rows that compose a specified object.
Syntax	MorphGetNumberOfRuns (<i>nObjectID</i> As Integer, <i>nRetNumberOfRuns</i> As Integer) As Long
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
Return values	<i>nRetNumberOfRuns</i> Returns the number of horizontal rows in the object.

MorphGetNumberOfVertices

Description	Obtains the number of vertices that make up the edgelist of a specified object.
Syntax	MorphGetNumberOfVertices (<i>nObjectID</i> As Integer, <i>nRetNumVertices</i> As Integer) As Long
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
Return values	<i>nRetNumVertices</i> Returns the number of vertices in the object's edgelist.
See also:	MorphGetVertexList

MorphGetObjectBoundingRect

Description	Obtains the coordinates of the upper left and lower right corners of a specified object's bounding rectangle.
Syntax	MorphGetObjectBoundingRect (<i>nObjectID</i> As Integer, <i>nRetX1</i> As Integer, <i>nRetY1</i> As Integer, <i>nRetX2</i> As Integer, <i>nRetY2</i> As Integer) As Long
Remarks	A <i>bounding rectangle</i> is a device used by MetaMorph to work with irregularly shaped objects. This contrivance is created by placing an imaginary rectangle over the object's outline. The sides of the rectangle will be perfectly horizontal and vertical, and the smallest rectangle possible will be used. The MorphGetObjectBounding-Rect function provides the coordinates of the starting point (upper left corner) and ending point (lower right corner).

Measuring Single Objects, continued 8.6

MorphGetObjectBoundingRect (continued)

Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
Return values	<i>nRetX1</i> Returns the X-coordinate of the upper left corner of the bounding rectangle.
	<i>nRetY1</i> Returns the Y-coordinate of the upper left corner of the bounding rectangle.
	<i>nRetX2</i> Returns the X-coordinate of the lower right corner of the bounding rectangle.
	<i>nRetY2</i> Returns the Y-coordinate of the lower right corner of the bounding rectangle.
See also:	MorphGetVertexList

MorphGetParmMeasurement

Description	Obtains the measured value of a selected parameter for a specified object.
Syntax	MorphGetParmMeasurement (<i>nObjectID</i> As Integer, <i>nParmNumber</i> As Integer, <i>fRetMeasurement</i> As Single) As Long
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
	<i>nParmNumber</i> Specifies the parameter number. This number can be obtained with the MorphFindParmIndex function.
Example	' Find the width of object 5. This code assumes that ' measurements have already been made. Dim <i>n</i> As Integer Dim <i>nWidth</i> As Single <i>MM</i> .MorphFindParmIndex "Width", <i>n</i> <i>MM</i> .MorphGetParmMeasurement 5, <i>n</i> , <i>nWidth</i> <i>MM</i> .PrintMsg "The width of object 5 is " + Str(<i>nWidth</i>)
Return values	<i>fRetMeasurement</i> Returns the measured value of the selected parameter.
See also:	MorphFindParmIndex (Section 8.3)

MorphGetPixelArea

Description	Obtains the number of pixels in a specified object.
Syntax	MorphGetPixelArea(nObjectID As Integer, lRetNumberOfPixels As Long) As Long
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
Return values	<i>lRetNumberOfPixels</i> Returns the number of pixels in the object.

MorphGetVertexList – for Visual Basic 6 & earlier MorphGetVertexListEx2 – for Visual Basic .NET (2002 – 2008)

Description	Obtains the X and Y coordinates of all vertices in the edgelist of a specified object.
Syntax	MorphGetVertexList(<i>nObjectID</i> As Integer, <i>aX</i> () As Integer, <i>aY</i> () As Integer) As Long MorphGetVertexListEx2(<i>nObjectID</i> As Integer, <i>aX</i> () As Integer, <i>aY</i> () As Integer) As Long
Remarks	returns the coordinates of the vertices making up the edge list of the given object. The number of elements in the arrays aX and aY should at least the number of vertices in the object. use MorphGetNumberOfVertices to obtain this number.
Parameters	<i>nObjectID</i> Specifies the number of the object. Objects are numbered from 0 to $(n-1)$, where <i>n</i> is the total number of objects.
	aX() Defines a buffer into which the X-coordinates of the vertices will be read.
	aY() Defines a buffer into which the Y-coordinates of the vertices will be read.
Return values	aX() The X-coordinates of the vertices will be read into this predefined buffer. (See Parameters.)
	aY() The Y-coordinates of the vertices will be read into this predefined buffer. (See Parameters.)
See also:	MorphGetNumberOfVertices
	Continued on next name

8.6 Measuring Single Objects, continued

Figure 8.1 Single Object Data "Get" Function Programming Example

```
' For all the measured objects, print out some information
' obtained during measurement about each one. This code
' assumes object measurements were previously performed.
Dim nObjects As Integer
MM.MorphGetNumberOfObjects nObjects
Dim i As Integer
For i = 0 To nObjects - 1
 MM.PrintMsg "Object " + Str(i) + " information:"
 Dim area As Long
 MM.MorphGetPixelArea i, area
 MM. PrintMsg " Pixel area is " + Str(area)
 Dim x As Single, y As Single
 MM.MorphGetCentroid i, x, y
 MM.PrintMsg " Centroid is at " + Str(x) + ", " + Str(y)
 Dim nx As Integer, ny As Integer
 MM.MorphGetCentroidPixel i, nx, ny
 MM.PrintMsg " Centroid pixel is at " + Str(nx) + ", "
    + Str(y)
 MM.MorphGetInternalPoint i, nx, ny
 MM. PrintMsg " The coordinates of a point inside the object
   are " + Str(nx) + ", " + Str(ny)
 Dim n As Integer
 MM.MorphGetNumberOfVertices i, n
 MM.PrintMsg " The object has " + Str(n) + " vertices"
 Dim ax(n) As Integer, ay(n) As Integer
 MM.MorphGetVertexList i, ax, ay
 Dim j As Integer
  For j = 0 To n
 MM.PrintMsg " Vertex " + Str(j) + " is " + Str(ax(j))
    + ", " + Str(ay(j))
  Next j
```

8.6 Measuring Single Objects, continued

Single Object Data "Get" Function Programming Example (continued)

```
MM.MorphGetNumberOfRuns i, n
MM.PrintMsg " There are " + Str(n) + " runs in the object"
Dim x1 As Integer, y1 As Integer, x2 As Integer,
y2 As Integer
MM.MorphGetObjectBoundingRect i, x1, y1, x2, y2
MM.PrintMsg " The bounding rectangle coordinates are "
+ Str(x1) + ", " + Str(y1) + " and " + Str(x2) + ", "
+ Str(y2)
Next i
```

Index

—A—

Annotations Creating and assigning 54 Reading 51 Area, measuring thresholded 106 Arrays 14 ASCII control codes 21 AutoEnhance 59 Autoscaling Current setting, reading 63 Enabling and disabling 65 Range maximum Configuring 66 Reading 64 Range minimum Configuring 67 Reading 65

—B—

BinarizeImage 81 Bit-depth 51 Bounding rectangle 126 Brightness Autoenhancing 60 Configuring 62 Making changes permanent 60 Reading the current setting 61 Reverting to the previous setting 62

Centroids, obtaining coordinates 123 Class Modules 9 Classifiers Failed objects, drawing 111 Measurement data, reading 119, 120 Number of filters passed, determining for an object 124 Settings Configuring 117 Reading 116 CloneImage 37 CloseImage 38 Command Line text box 5 Contrast Autoenhancing 60 Configuring 63 Making changes permanent 60 Reading the current setting 61 Reverting to the previous setting 62 Coordinates

Bounding rectangle, determining for 126 Internal point of an object 125 Nearest object, finding 124 Vertices in an object, reading 128 CopyImage 38 CopyImagePlane 39 Copying images 37 CreateImage 39 CreateRectRegion 93

—D—

DestroyRegion 94 DIGetFirst 23 DIGetIOStatus 23 DIGetLineCount 24 DIGetLineState 25 DIGetName 25 DIGetNext 26 Digital I/O High vs. Low state, reading 25 Input vs. output status of a line, reading 23 Number of lines, reading 24 Receiving signals from a device 27 Sending signals to a device 26 DoCommand 6, 8

—Е—

Edgelists Coordinates for a region, reading 101 Vertices Coordinates, reading for an object 128 Number of, reading 126 Exclusive thresholding 78, 79

—F—

FixImage 60 ForceCloseImage 40 Functions required by MetaMorph 7

—G—

GetActivePlane 51 GetActiveRegion 95 GetAutoScale 64 GetBrightness 61 GetContrast 61 GetCurrentImage 44 GetDepth 51 GetFunctionHandle 30 GetHeight 51 GetImage 44 GetImageAnnotation 52 GetImageName 52 GetImageWindowPosition 47 GetImageWindowSize 47 GetLut 69 GetLutModel 70 GetMaxScale 65 GetMinScale 66 GetNumberOfImages 45 GetNumberOfPlanes 53 GetNumberOfRegions 96 GetRegion 96 GetRegionArea 99 GetRegionAverageValue 105 GetRegionDistance 99 GetRegionMaximumValue 106 GetRegionMinimumValue 105 GetRegionPosition 100 GetRegionSize 100 GetRegionStdDeviation 106 GetRegionThresholdArea 107 GetThresholdRange 77 GetThresholdState 78 GetWidth 53 GetZoom 53 gParentWnd variable 7 gUserID variable 7

—H—

Handles Obtaining Digital I/O MetaDevice 23, 26 Function 30 Image 44 Region of interest 95, 96 Testing validity Image 46 Region 97 Height, image 51

I

Image windows Maximizing 48 Minimizing 49 Position Configuring 49 Reading 47 Size Configuring 50 Reading 47 Images Binarizing 81 Closing 38, 40 Copying 37 Copying a plane 39

Handles, finding 44 Height, reading 51 Loading 41 Name, reading 52 New images, creating 39 Number of loaded images, finding 45 Overwriting 38 Renaming 56 Saving 41 Selecting for measurement 115 Setting the timestamp 56 Updating display 59 Width, reading 53 Inclusive thresholding 78, 79 Intensity Average 105 Maximum 106 Measuring 107 Minimum 105 Reading From a column of pixels 82, 83 From a row of pixels 84, 85 From a single pixel 84 Standard deviation 106 Writing values To a column of pixels 86, 87 To a row of pixels 89, 90 To a single pixel 88 IsValidImage 46

___J___

Journals, running 33

IsValidRegion 97

—K—

Keep Program in Memory After Execution check box 6

—L—

Labeling an image 91 LoadImage 41 Look-up tables Assigning a LUT model 73 Current LUT model in use, determining 70 Reading a table's elements 69 Writing to a table's elements 72

—M—

Mask images 115 MaximizeImageWindow 48 Measurement Centroids 123 Data

Filling an array 125 Reading 119, 120 Measuring objects in an image 121 Number of pixels in an object, measuring 128 Parameter data, reading for an object 127 Recalculating 122 Vertices in an object, reading 128 Measurement images 115 MeasureRegion 107 Message windows Configuring size and position 34 Displaying 30 **MetaDevices** Obtaining a handle 23 Obtaining a name from a handle 25 MinimizeImageWindow 49 MM variable 7 MorphDeleteObject 118 MorphFindParmIndex 112 MorphGetCentroid 123 MorphGetCentroidPixel 123 MorphGetClassifiersPassed 124 MorphGetClosestObject 124 MorphGetFilter 116 MorphGetInternalPoint 125 MorphGetMeasurements 125 MorphGetNumberOfObjects 118 MorphGetNumberOfParms 113 MorphGetNumberOfRuns 126 MorphGetNumberOfVertices 126 MorphGetObjectBoundingRect 126 MorphGetParmAccumSummary 119 MorphGetParmDescription 113 MorphGetParmMeasurement 127 MorphGetParmName 114 MorphGetParmSummary 120 MorphGetPixelArea 128 MorphGetVertexList 128 MorphMeasureObjects 121 MorphRecalc 122 MorphSetDrawFailedObjectsFlag 111 MorphSetFillHoleFlag 111 MorphSetFilter 117 MorphSetupMeasurements 115

__N__

New images, creating 39

-0--

Object linking and embedding 7 Objects Measuring 121 Nearest object, finding 124 Number last measured, finding 118 Parameter data, reading 127 Pixels in an object, measuring 128 Remeasuring 122 Removing from a measurement list 118 Rows in an object, measuring 126 Vertices, determining number of 126 OLE 7

—P—

Palettes, configuring number of entries 74 Parameters Description, obtaining 113 Name, obtaining 114 Number of measurable parameters, determining 113 Parameter number, determining from a parameter's name 112 Perimeter, determining for a region 99 Planes Active plane, setting 55 Copying 39 Number of planes in stack, reading 53 Plane number, reading 51 Preferences Failed objects, drawing 111 Holes, filling 111 PrintMsg 30 Program Name drop-down list 5 Public variants 9

—R—

ReadColumn 82 ReadColumnEx 83 ReadPixel 84 ReadRow 84 ReadRowEx 85 RegionGetEdgePixelCoordinates 101 RegionGetNumEdgePixels 101 Regions Active region, setting 98 Area, determining 99 Creating 93 Handles, finding 95, 96 Intensity, measuring 107 Number of regions, determining 96 Numbers, switching 104 Perimeter, determining Calibrated units 99 Pixel units 101 Position Configuring 103 Reading 100 Removing 94 Size Configuring 103 Reading 100 Thresholded area, measuring 107 Remove command button 6

ResetContrast 62 Run User Program Dialog box 5 Options 5 RunFunction 31 RunFunctionEx 32 RunJournal 33 Running a function 31 Running a journal 33 RUNUSER drop-in 5

SaveImage 41 SendSerialData 18 Serial communication ASCII control codes 21 Receiving a data stream from a device 19 Sending a data stream to a device 18 Syntax rules 20 Set 9 SetActivePlane 55 SetActiveRegion 98 SetAutoScale 66 SetBrightness 62 SetContrast 63 SetDigitalIO 26 SetDisplayImagesWhenCreated 42 SetFunctionVariable 33 SetImageAnnotation 55 SetImageName 56 SetImageTimestamp 56 SetImageWindowPosition 49 SetImageWindowSize 50 SetLut 72 SetLutModel 73 SetMaxScale 67 SetMinScale 68 SetNumPaletteEntries 74 SetPrintMsgWindowPositionAndSize 34 SetRegionPosition 103 SetRegionSize 103 SetThresholdRange 79 SetThresholdState 79 SetZoom 57 ShowImage 43 Shutdown 6,8 Standard area 115 Startup 6, 8 Summary data, recalculating 122 SwapRegionNumbers 104

—T—

Text, writing on an image 91 Thresholding Area, measuring thresholded 107 Holes, filling 111 Range Configuring 79 Reading 77 State Configuring 79 Reading 78

—U—

UpdateDisplay 59 User programs Compiling 12 Creating With Visual Basic 4.0 10 With Visual Basic 5.0 11 Overview 7 Registering and unregistering 6 UserMethods 9

__V__

Variables required by MetaMorph 7 Variables, setting values for 33 Vertices Determining number in an object 126 Reading coordinates from an object 128

WaitForDigitalIO 27 WaitForSerialData 19 Width, image 53 WriteColumn 86 WriteColumnEx 87 WritePixel 88 WriteRow 89 WriteRowEx 90 WriteText 91

—Z—

Zoom factor Configuring 57 Reading 53