

## MetaMorph NX Software Launches!!

Molecular Devices, Inc., is proud to announce the launch of the next generation of its MetaMorph® Software, streamlining the imaging workflow while preserving the acquisition capabilities and extensive hardware support that has made the software the best in its class.

The new MetaMorph® NX Software presents an entirely redesigned user-centered interface to simplify even the most complex imaging tasks. With one-click access to features, integrated hardware setup, and synchronized, unobstructed views of data, the new software modernizes the imaging process. Instead of spending time learning complicated software, scientists can spend their time on what is most important—their research.

The MetaMorph NX Software continues the strong market leadership position long held by the MetaMorph® Software Suite of products as the imaging software that supports hundreds of microscope devices and components. As such, it is the most flexible and powerful foundation available for a cohesive and customizable imaging acquisition and analysis system. Hardware configurations have never been as easy as they are now – with just a few clicks on the Hardware Setup tab, devices and components can be quickly configured for use with the MetaMorph NX Software.

To celebrate the launch of MetaMorph NX Software we are offering the following great promotions!

### Half price upgrades

Upgrade MetaMorph Software version 6 or 7 to MetaMorph NX Software for only \$1000.

### Half price crossgrade from Nikon NIS Elements

Move over to MetaMorph NX Software the best microscopy automation and image analysis package for half price. Any existing NIS Elements user who purchases MetaMorph NX Software will receive a 50% discount from list price. To qualify customers must surrender their NIS Elements license dongle within 10 days after receipt of ordered product. Any order not paired with returned NIS Elements license dongle will be invoiced at normal price.

Both these offers are valid February 1 to April 30, 2011.



## MetaMorph NX

Inside:

MetaMorph NX Software Launches	1
Focus: MetaMorph® NX Software	2-3
General information	4
Upcoming events	4

# Focus: MetaMorph® NX Software

## MetaMorph® NX Software Makes Imaging Easier

For nearly 25 years, MetaMorph® Software has given microscopists the means to gain extraordinary insight into cell morphology, function, and behavior. Researchers have relied on MetaMorph Software's myriad options for microscope component integration, task automation, image acquisition using a variety of illumination and collection methods, and thorough data analysis. MetaMorph Software pioneered the bio research imaging industry and, with unsurpassed power and flexibility, it continues to be the best in its class.

## Introducing MetaMorph® NX Software – The Best Gets Better

MetaMorph NX Software, the next generation of MetaMorph Software, reflects our commitment to ensuring that our customers' experience is the best it can be. Without diminishing any of the features that facilitate comprehensive image acquisition and analysis, MetaMorph NX Software streamlines the workflow for all tasks and provides an entirely new user-centered interface. With one-click access to features, integrated hardware setup, and synchronized, unobstructed views of your data, you can become an imaging expert in minutes.

## Organized Workspace Coupled With A Faster Workflow

In the new patent-pending Dataset View, all of the images and data that belong to a dataset stay together in one organized workspace. The Dataset View displays images in a grid using a layout that you specify – select a view mode to organize any number of images by time point, channel (wavelength), stage position, Z position, and so on. Or, view a composite overlay of channels on one image and turn on or off the display of each channel. When your image looks the way you want it, export it to PowerPoint with one click.

Display large amounts of data without sacrificing clarity. Multiple Dataset Views may be positioned side-by-side horizontally or vertically to compare various aspects of the same dataset or different datasets. A filmstrip at the bottom of the Dataset View displays thumbnail images, augmenting the image navigation options. Zoom and pan all displayed images at once, scale images of any bit depth, and assign colors to channels to further investigate even the most subtle objects in your images.

The tab-based “ribbon” interface provides convenient access to the most commonly used options. Select an acquisition mode and only the options that are relevant to the mode appear, guiding you quickly through the

task by providing just the right tools at the right time. The ribbon's left-to-right workflow sets the path for hardware configuration, experiment setup, image acquisition, and data analysis, while extensive tooltips and online help offer assistance along the way. To make things even easier, you can create your own shortcuts to the options you use most.

## Built-in Hardware Setup and Device Control

Installing and configuring microscope devices and cameras takes only a few steps using the options on the Hardware Setup tab. MetaMorph NX Software does not impose limits on how hardware is configured – any supported component can be configured as a magnification or illumination component. And, because MetaMorph NX Software leverages the continuous development of drivers throughout its history, it has hundreds of high quality drivers for devices and components on the market today.

## Exceptional Display Performance and Multi-Tasking Capability

Optimum image and data display performance is ensured through configurable cache and display memory settings which let you determine how memory is allocated to the MetaMorph NX Software. And with support for 64-bit Microsoft Windows 7, more memory is available, making data analysis and manipulation even faster.

With its new multi-threading capability, MetaMorph NX Software remains responsive during even the longest acquisitions. and run multiple measurements at the same time. Multi-threading also allows for asynchronous control of hardware peripherals, improving the speed of acquisitions.

## Comprehensive Uncomplicated Image Acquisition

The numerous options for image acquisition continue in MetaMorph NX Software but the new interface and workflow simplify the configuration process and eliminate the need for journals or macros. In most cases, you can select an acquisition mode and configure a single or multidimensional image acquisition using

## Benefits

- Reduce training and support with new interface.
- Faster results with clear view and comparison of images.
- More productive with simultaneous image acquisition and analysis.
- Enables unique applications through customization.

# Focus: MetaMorph® NX Software (cont.)

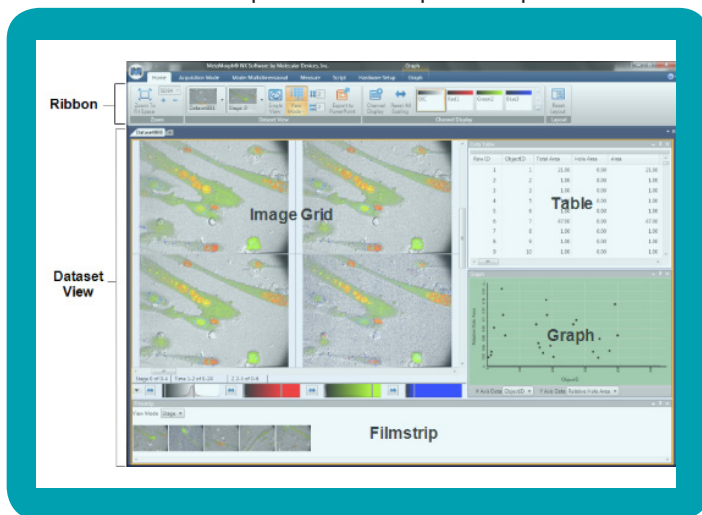
just the settings on the ribbon. Options are available for acquiring images at various sites, at certain times, at multiple focal planes, with a variety of illumination conditions using built-in image autofocus or microscope focus stabilization. You can hide the options for the parts of the acquisition configuration that you do not need. Advanced settings (for example, the order in which images are acquired with multiple channels and a Z series) are a click away in the configuration dialog box.

In multidimensional acquisition mode, MetaMorph NX Software saves images as they are acquired and automatically stores acquisition settings. You will never lose your work, and you can quickly repeat an acquisition at a later date using the stored settings.

## Coherent Image Review, Analysis, and Data Management

MetaMorph NX Software provides several ways to measure and evaluate objects in images. Measurements can include the area, dimensions, shape, orientation, and intensity of objects and regions, as well as a statistical summary of the measurements made. You can measure all the images from one sample at once without needing to write a journal or script. Once you measure images, you can sort and graph the results. Each set of measurement results and the settings used to make the measurement are automatically saved with the dataset and can be accessed, reconfigured, and filtered using MetaMorph NX Software at any time. In the Dataset View, you can arrange images next to tables and graphs to explore how measurements made on some or all of

**MetaMorph NX Desktop Workspace**



Two main areas comprise the MetaMorph NX Software workspace:

- The ribbon across the top, which contains tabs and options for performing tasks.
- The Dataset View, which contains images and can optionally include a filmstrip, the data table, and one or more graphs that belong to a dataset.

the images relate to the dataset as a whole, to other measurements, and to the visual information in the images. When you have sorted your analysis data according to your criteria, export your data to Excel with one click.

## Features

- Easy-to-Find Tools
- Workflow Oriented
- Clean Desktop
- Experiment History
- Hundreds of Hardware Drivers
- Multi-Threaded Application

Every time you create a new experiment, MetaMorph NX Software creates a folder for the experiment on the hard drive of your computer. An experiment folder contains a subfolder for each dataset in the experiment. Each dataset folder contains images (single image .tif files), image metadata, measurement sets, and, if applicable, the acquisition settings used to acquire the images. You can add other files to the folder that support your experiment, such as emails, presentations, documents, and notes.

## Full Support For Extendibility and Task Automation Through Scripting

The Script tab provides an environment for writing scripts using IronPython, an implementation of the Python scripting language, which provides native support for all standard .NET tools. With IronPython, you can extend the MetaMorph NX Software to solve virtually any experimental problem and you can automate tasks within MetaMorph NX Software. The scripting environment uses auto-completion, which presents a list of available successive functions as you write a script.

## Modular Architecture Means Faster Delivery of Features and Customization Opportunities

The MetaMorph NX Software is built upon a new modular architecture that assures rapid addition of subsequent features, applications, and support for new hardware. Enhancements will be frequent, readily available, and easy to install.

## Contact Us

Molecular Devices Inc.  
402 Boot Road  
Downingtown, PA 19335  
USA

Phone Toll Free: (800) 635-5577  
Phone Intl.: (610) 873-5610  
Fax: (610) 873-5492  
Support: (800) 635-5577 x1820

On the web: [www.moleculardevices.com](http://www.moleculardevices.com)

Sales: [meta.admin@moldev.com](mailto:meta.admin@moldev.com)  
Orders: [om-meta@moldev.com](mailto:om-meta@moldev.com)  
Support: [support.dtn@moldev.com](mailto:support.dtn@moldev.com)  
Training: [training.dtn@moldev.com](mailto:training.dtn@moldev.com)

**MetaMorph® Software...**  
**making imaging easy!**

<b>Product Information:</b>	
Overview of available MetaMorph Software products	<a href="http://www.moleculardevices.com/Products/Software/Meta-Imaging-Series.html">http://www.moleculardevices.com/Products/Software/Meta-Imaging-Series.html</a>
Comparison of available MetaMorph Software products	<a href="http://www.moleculardevices.com/Products/Software/Meta-Imaging-Series/Comparison-Table.html">http://www.moleculardevices.com/Products/Software/Meta-Imaging-Series/Comparison-Table.html</a>
Application modules available for MetaMorph Software	<a href="http://www.moleculardevices.com/Products/Software/Meta-Imaging-Series/Application-Modules.html">http://www.moleculardevices.com/Products/Software/Meta-Imaging-Series/Application-Modules.html</a>
List of distribution partners	<a href="http://www.moleculardevices.com/Company/Who-We-Are/Contact-Us/Distributors/Meta-Morph.html">http://www.moleculardevices.com/Company/Who-We-Are/Contact-Us/Distributors/Meta-Morph.html</a>
<b>Support:</b>	
MetaMorph Software updates and upgrades	<a href="http://www.meta.moleculardevices.com/updates/">http://www.meta.moleculardevices.com/updates/</a>
MetaMorph Software installation guides	<a href="http://support.meta.moleculardevices.com/install/index.php">http://support.meta.moleculardevices.com/install/index.php</a>
MetaMorph Software supported hardware database	<a href="http://support.meta.moleculardevices.com/hardware/hardware.php">http://support.meta.moleculardevices.com/hardware/hardware.php</a>
MetaMorph Software supported operating systems	<a href="http://support.meta.moleculardevices.com/info/os-compatibility.php">http://support.meta.moleculardevices.com/info/os-compatibility.php</a>
Technical and application notes for MetaMorph Software	<a href="http://support.meta.moleculardevices.com/docs/">http://support.meta.moleculardevices.com/docs/</a>
Back issues of MetaMatters newsletter	<a href="http://support.moleculardevices.com/pages/software/meta_matters.html">http://support.moleculardevices.com/pages/software/meta_matters.html</a>
Information and registration for MetaMorph Software training courses and webinars	<a href="http://www.moleculardevices.com/Support/Product-Training/Research-Imaging.html">http://www.moleculardevices.com/Support/Product-Training/Research-Imaging.html</a>

# Upcoming Training, Courses, and Exhibitions

## Training

March 25, 2011  
Webinar: Introduction to MetaMorph NX Software

March 28, 2011  
MetaMorph NX Software Introductory Training Course  
Downingtown, PA

March 29-30, 2011  
MetaMorph® Software Basic Training Course  
Downingtown, PA

March 31-April 1, 2011  
Metamorph Software Advanced Training Course  
Downingtown, PA

May 17, 2011  
Webinar: MetaMorph Software in Working Lab Environment

June 6, 2011  
MetaMorph NX Software Introductory Training Course  
Sunnyvale, CA

June 7-8, 2011  
Metamorph® Software Basic Training Course  
Sunnyvale, CA

June 9-10, 2011  
Metamorph Software Advanced Training Course  
Sunnyvale, CA

July 15, 2011  
Webinar: TBD

September 19, 2011  
MetaMorph NX Software Introductory Training Course  
Downingtown, PA

September 20-21, 2011  
Metamorph® Software Basic Training Course  
Downingtown, PA

September 22-23, 2011  
Metamorph Software Advanced Training Course  
Downingtown, PA

September 30, 2011  
Webinar: TBD

November 18, 2011  
Webinar: TBD

## Supported Courses

May 4-13, 2011  
AQLM  
MBL, Woods Hole, MA

May 20-27, 2011  
QFM  
MDIBL, Mount Desert Island, ME

October 11-21, 2011  
OMIBS  
MBL, Woods Hole, MA

October 19-November 1, 2011  
In-situ Hybridization & Live Cell Imaging  
CSHL, Cold Spring Harbor, NY

## Exhibitions

November 12-16, 2011  
Society for Neuroscience Annual Meeting  
Washington, DC

December 3-7, 2011  
ASCB Annual Meeting  
Denver, CO