



# MetaXpress<sup>®</sup> 6 Software Guide

Using Modify Image Tools in CME

Date Revised 07/13/15 Version B

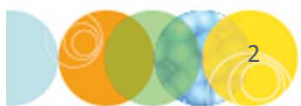


© 2012-2015. Trademarks property of Molecular Devices, LLC or their respective owners.  
For research use only. Not for use in diagnostic procedures.

# Chapter Purpose

The purpose of this chapter is to describe the **Modify Image** tools available in the **Custom Module Editor** (CME) plugin.

**Modify Image** tools can be used to modify a grayscale (8 or 16-Bit) image. The newly created grayscale image can be modified further or used to detect objects of interest using a **Find Objects** or **Application Modules Objects** step.



# Ribbon: Modify Image Tools

The screenshot displays the software interface with the 'Custom Module' ribbon selected. The 'Modify Image' group is highlighted, showing icons for 'Add', 'Subtract', 'Multiply', 'Divide', 'Maximum', 'Minimum', 'Morphology', 'Invert', 'Top Hat', 'Bottom Hat', 'HDome', 'HBasin', 'Regional Maximum', and 'Regional Minimum'. A red box highlights the 'Add' icon, with a red arrow pointing to it. A red callout box says 'Click on the appropriate icon'. A red callout box says 'Click here for a drop-down menu of Modify Image tools'. An orange box highlights the 'Add' dialog box, which has a title bar '10 Add [Modified] X'. The dialog contains fields for 'Source 1' (Invert), 'Source 2' (Invert), 'Constant' (0), and 'Result' (Add). The 'Description' field says 'Combines the intensity values of two images.' and there is an 'Apply' button. An orange arrow points from the 'Add' icon to the dialog box. The main workspace shows a binary image of cells. A 'Custom Module Step Images' panel at the bottom shows three steps: 1 (original image), 2 (thresholded image), and 3 (result of the 'Add' operation).

Click here for a drop-down menu of **Modify Image** tools

Click on the appropriate icon

Clicking on a **Modify Image** tool icon will add a card on the panel to the left



# Modify Image Tools Descriptions

There are 3 different classes of modify image tools: **Arithmetic**, **Morphological**, and **Special**

## ARITHMETIC

<b>ADD</b>	Combines the intensity values of each corresponding pixel in two images
<b>ADD CONSTANT</b>	Adds a user-specified intensity value to each pixel in the image
<b>SUBTRACT</b>	Subtracts the intensity values of each corresponding pixel in one image from the other
<b>MULTIPLY</b>	Multiplies intensity values of each corresponding pixel in two images
<b>MULTIPLY CONSTANT</b>	Multiplies a user-specified intensity value to each pixel in the image
<b>DIVIDE</b>	Divides the intensity values of each corresponding pixel in one image by the other
<b>MAXIMUM</b>	Compares the intensity values of corresponding pixels in two images and displays the higher of the two values in the resulting image
<b>MINIMUM</b>	Compares the intensity values of corresponding pixels in two images and displays the lower of the two values in the resulting image



# Modify Image Tools Descriptions

There are 3 different classes of modify image tools: **Arithmetic**, **Morphological**, and **Special**

## MORPHOLOGICAL

<b>ERODE</b>	Grows dark areas and shrinks bright areas
<b>DILATE</b>	Grows bright areas and shrinks dark areas
<b>OPEN</b>	Removes parts of bright areas where the filter shape and size does not fit
<b>CLOSE</b>	Removes parts of dark areas where the filter shape and size does not fit
<b>OPEN-CLOSE / CLOSE-OPEN</b>	Combines Open and Close filters in sequential order to correct for both bright and dark noise spots
<b>CENTER FILTER</b>	Combines Open-Close and Close-Open filters for more robust noise reduction
<b>GRADIENT</b>	Highlights the boundaries, or outlines, between the bright and dark regions in an image
<b>INVERT</b>	Inverses the image histogram so dark pixels become bright and bright pixels become dark
<b>TOP HAT</b>	Finds small bright spots based on a filter shape and size
<b>BOTTOM HAT</b>	Finds small dark spots based on a filter shape and size



# Modify Image Tools Descriptions

There are 3 different classes of modify image tools: **Arithmetic**, **Morphological**, and **Special**

## MORPHOLOGICAL CONT.

### HDOME

Finds bright spots where the relative intensity values are equal to or higher than the threshold value specified

### HBASIN

Finds dark spots where the relative intensity values are equal to or lower than the threshold value specified

### REGIONAL MAXIMUM

Displays a bright spot for all peaks in intensity (a pixel or group of pixels of a constant intensity value completely surrounded by lower intensity pixels)

### REGIONAL MINIMUM

Displays a bright spot for all valleys in intensity (a pixel or group of pixels of a constant intensity value completely surrounded by higher intensity pixels)

### HOLES

Finds and, optionally, fills light or dark holes in an image and displays them as bright areas

### BORDER OBJECTS

Intensifies or suppresses bright or dark objects that touch the edge of the image



# Modify Image Tools Descriptions

There are 3 different classes of modify image tools: **Arithmetic**, **Morphological**, and **Special**

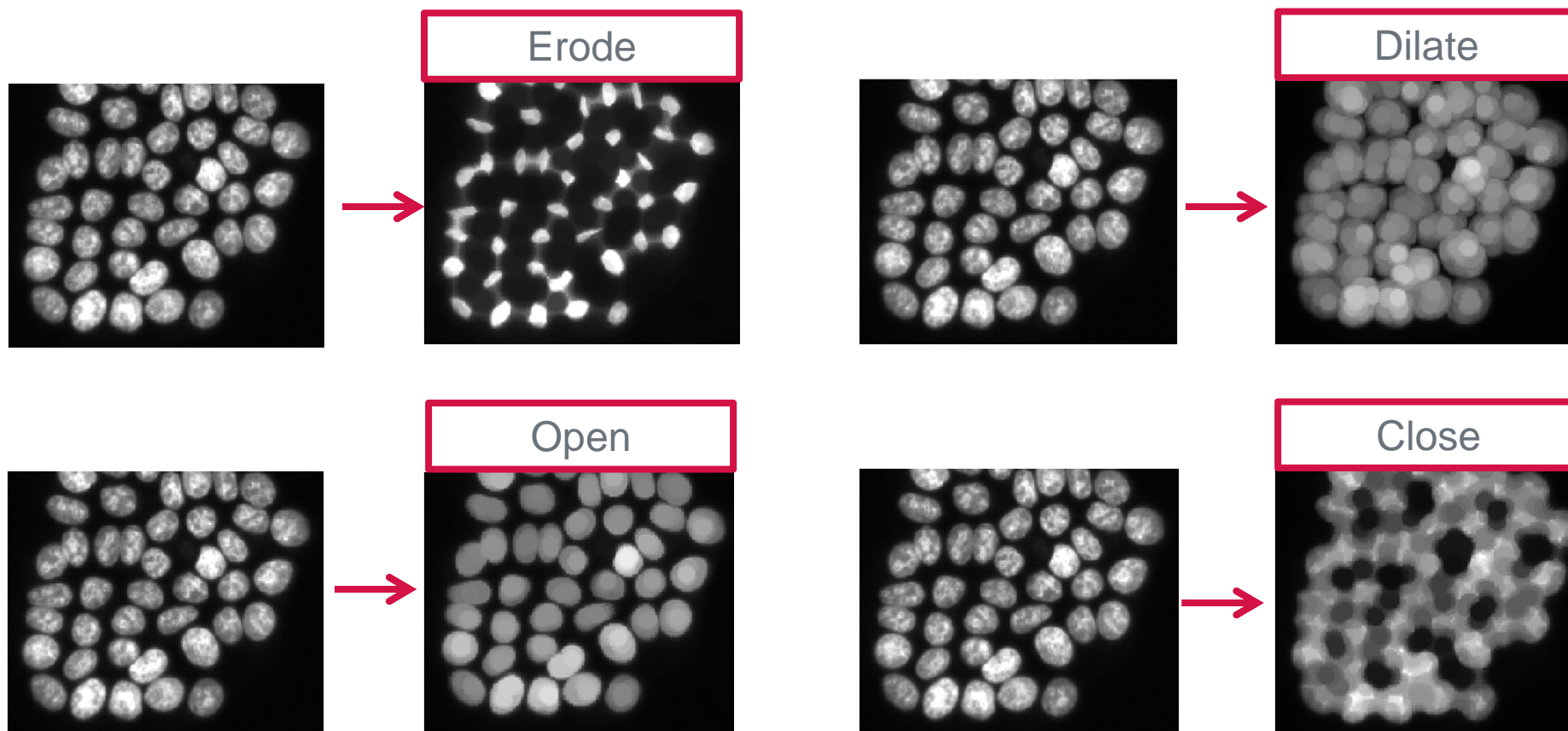
## SPECIAL

<b>AVERAGE FILTER</b>	Reduces the amount of variation between intensity values in adjacent pixels
<b>COLOR SEPARATE RGB</b>	Separates a color image into red, green, and blue components
<b>COLOR SEPARATE HSL</b>	Separates a color image into hue, saturations, and luminosity components
<b>COLOR SEPARATE CEI-LAB</b>	A non-linear color separation technique designed to better mimic human vision
<b>GAUSSIAN FILTER</b>	Smooths and blurs an image by a user-defined factor (sigma)
<b>LOG FILTER</b>	Laplacian of Gaussian, useful for finding areas or objects of a specific size
<b>DISTANCE</b>	Creates a 16-bit grayscale image from a binary mask image where the intensity value indicates the distance, measured in pixels, to the nearest white pixel in the original binary image
<b>MASK TO IMAGE</b>	Convert a segmentation mask to an image so that it can be used with the Modify Image tools



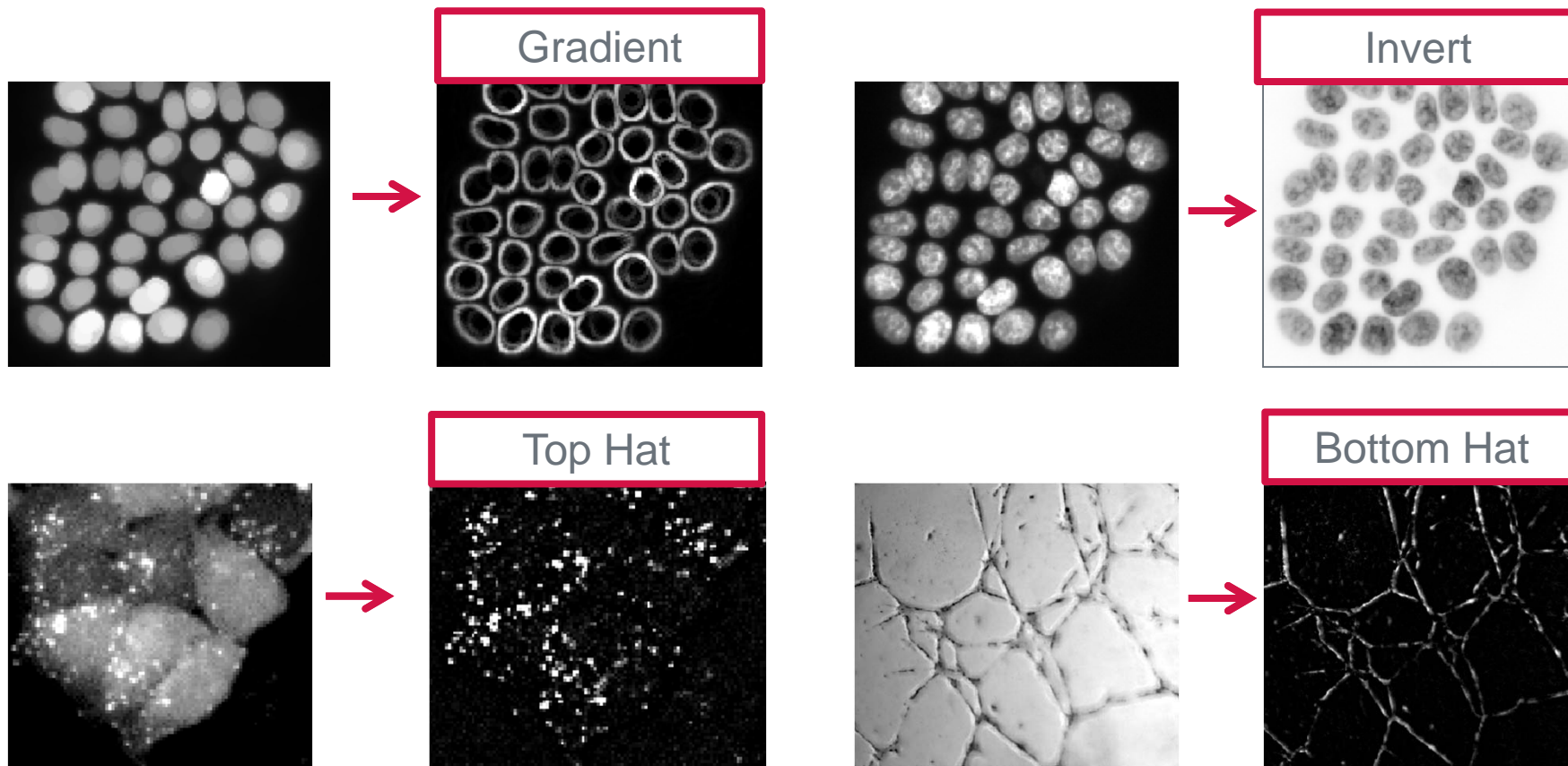


# Modify Image: Examples of Morphology Filters

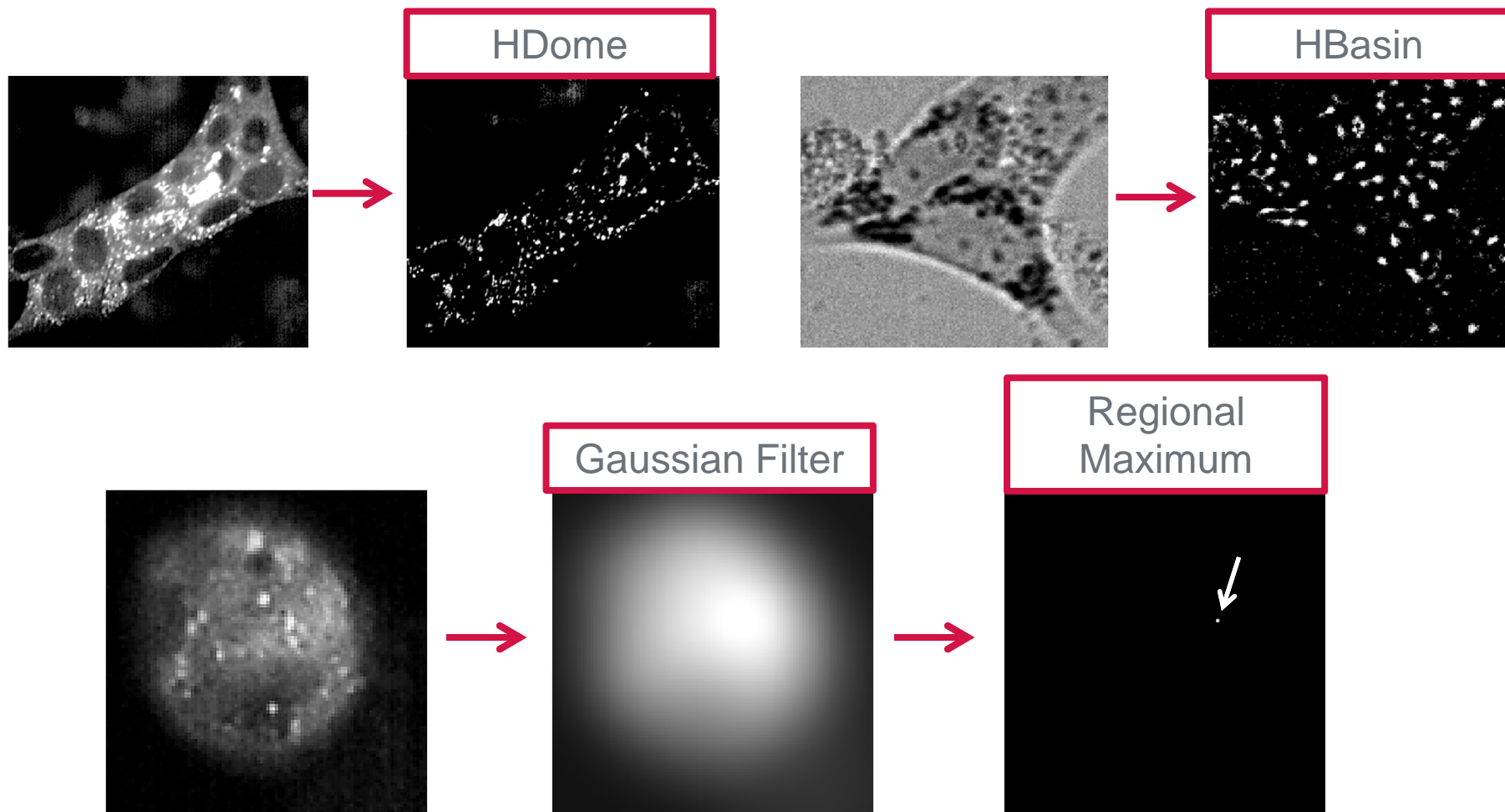




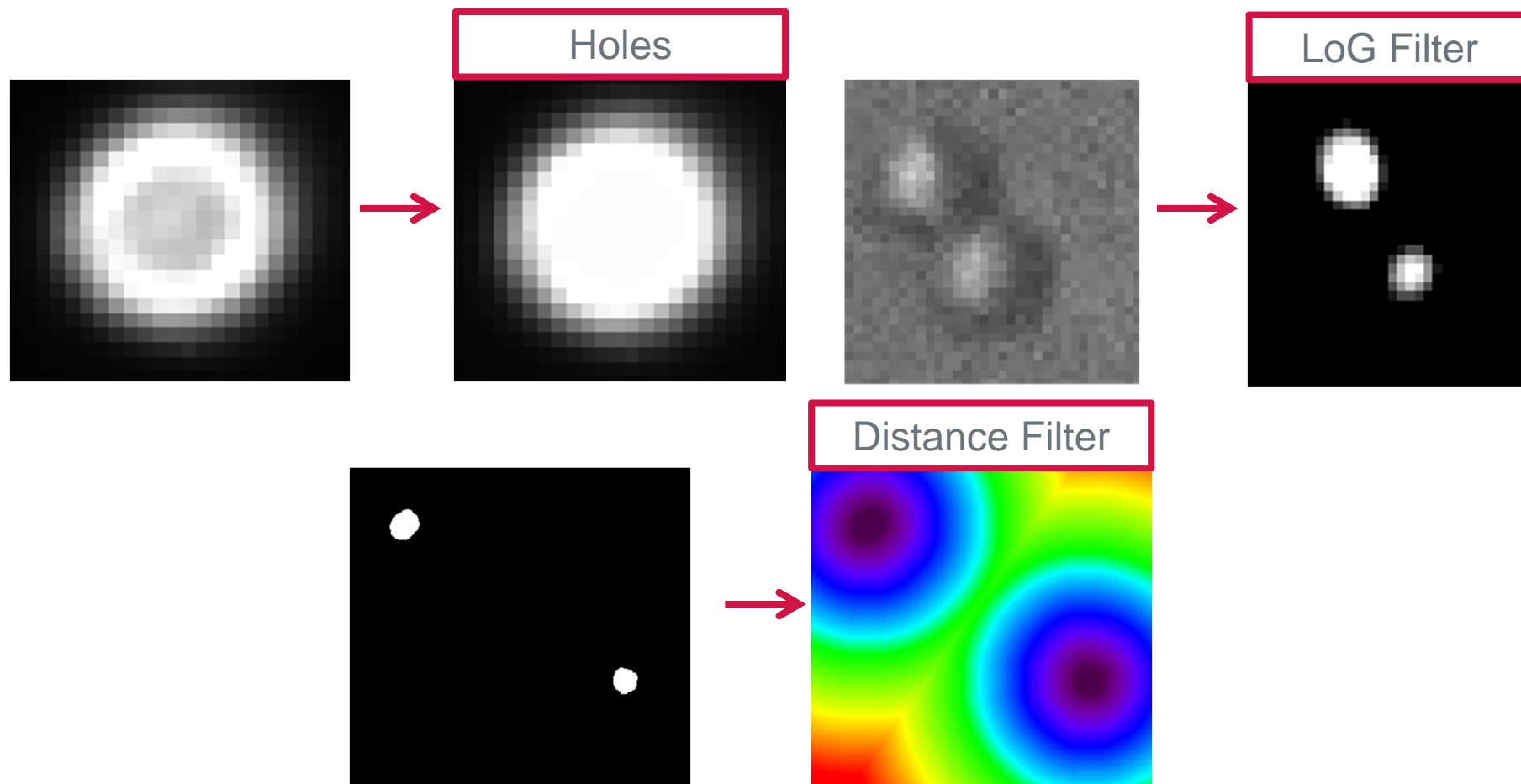
# Modify Image: Examples of Morphology Filters



# Modify Image: Examples of Morphology Filters



# Modify Image: Examples of Morphology Filters



# Support Resources

- F1 / HELP within MetaXpress® Software
- Support and Knowledge Base: <http://mdc.custhelp.com/>
- User Forum: <http://metamorph.moleculardevices.com/forum/>
- Request Support: <http://mdc.custhelp.com/app/ask>
- Technical Support can also be reached by telephone:
  - 1 (800) 635-5577
  - Select options for Tech Support → Cellular Imaging Products → ImageXpress Instruments





ADVANCING PROTEIN AND CELL BIOLOGY