



MetaXpress[®] 6 Software Guide

Creating Color Image Overlays in Review Plate Data

Date Revised 06/05/15 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 guide the user through creating up to a 3 color (RGB) overlay post-acquisition in both thumbnail montage and full resolution formats on the **Display** tab in **Review Plate Data**.

Note that this chapter covers this process within **Review Plate Data ONLY**. For additional color overlay options refer corresponding chapter for details.



Creating Image Overlays in Review Plate Data

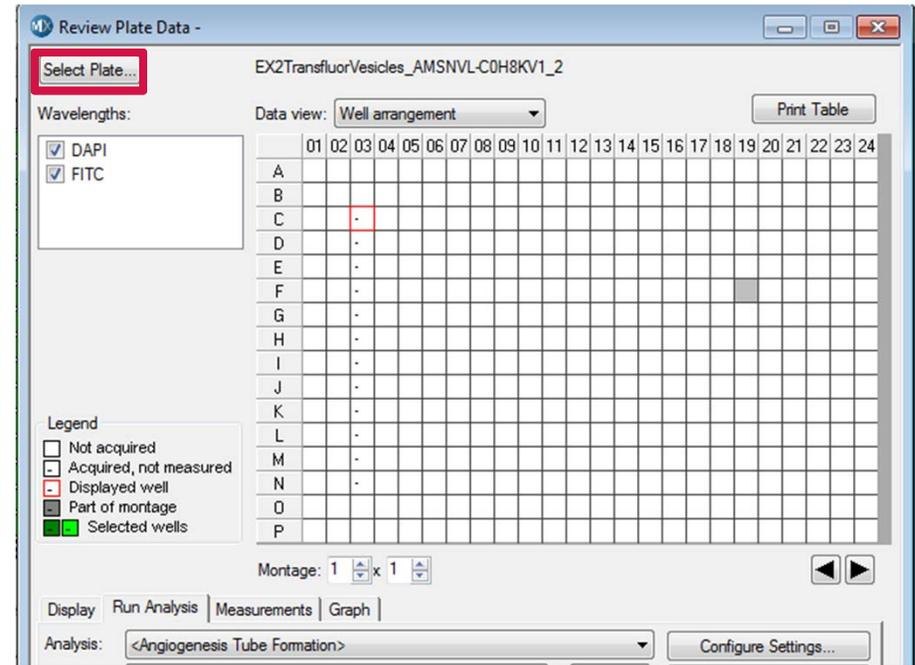
1. Open **Review Plate Data**

- In the main toolbar click on



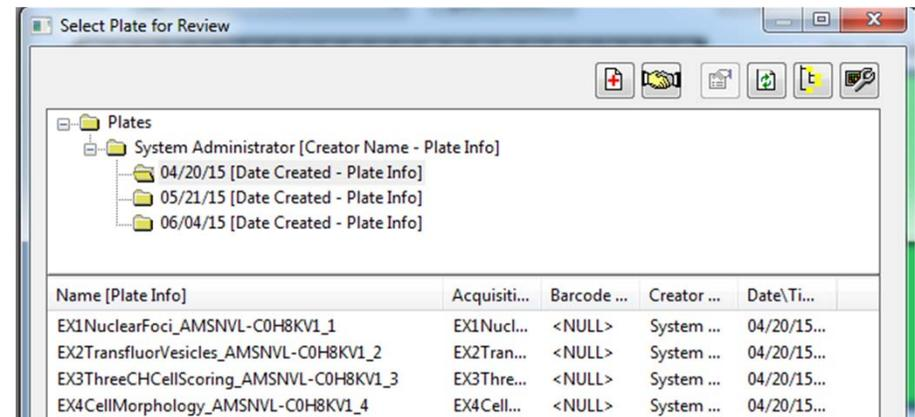
OR

- Under the **Screening** menu, select **Review Plate Data**



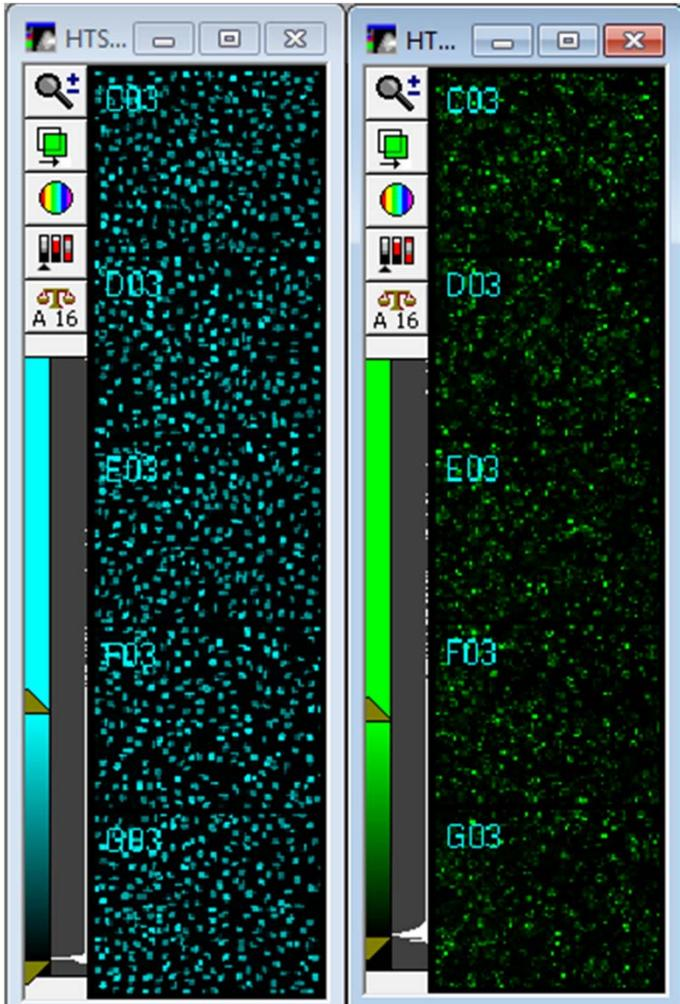
2. Click on the **Select Plate** button

3. Browse through the folders to open the plate of interest

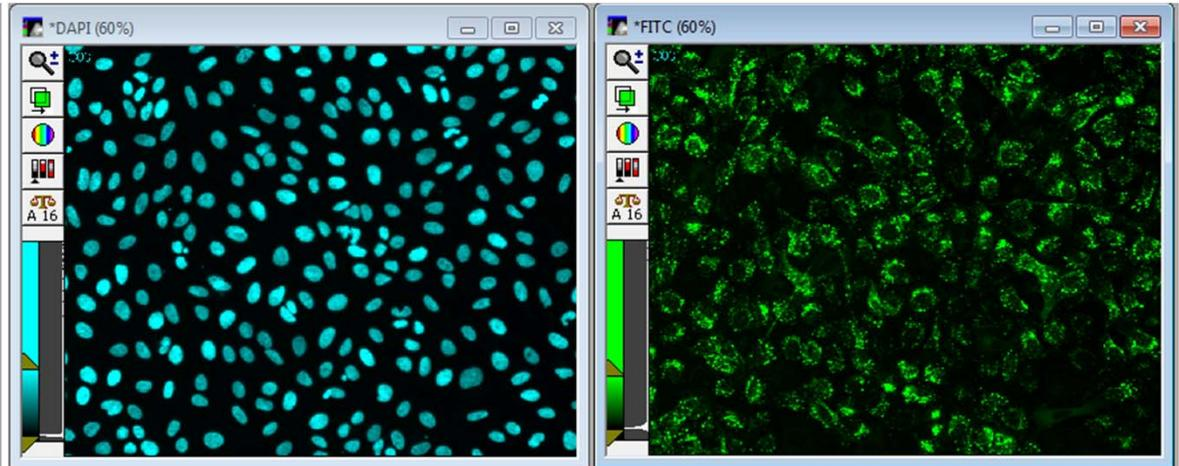


Creating Image Overlays in Review Plate Data

Thumbnail Montage (“HTS”) Images



High Resolution Images



4. In plate view, highlight a consecutive range of wells to display the **HTS-Thumbnail Montage**
 - Refer to the chapters on viewing images in **Review Plate Data** for more details



Creating Image Overlays in Review Plate Data

The screenshot shows the 'Review Plate Data' window for 'EX1TransfluorVesicles_AMSNVL-FDJRFV1_26'. The 'Wavelengths' section has 'DAPI' and 'FITC' checked. The 'Data view' is set to 'Well arrangement'. A grid shows well C03 highlighted in red. The 'Legend' includes 'Displayed well' (red square). The 'Display' tab is selected, and the 'Source RGB' dropdowns are set to 'None', 'FITC', and 'DAPI' respectively. The 'Image Overlay' is set to 'Show well information' and 'Col' is 'Cyan'.

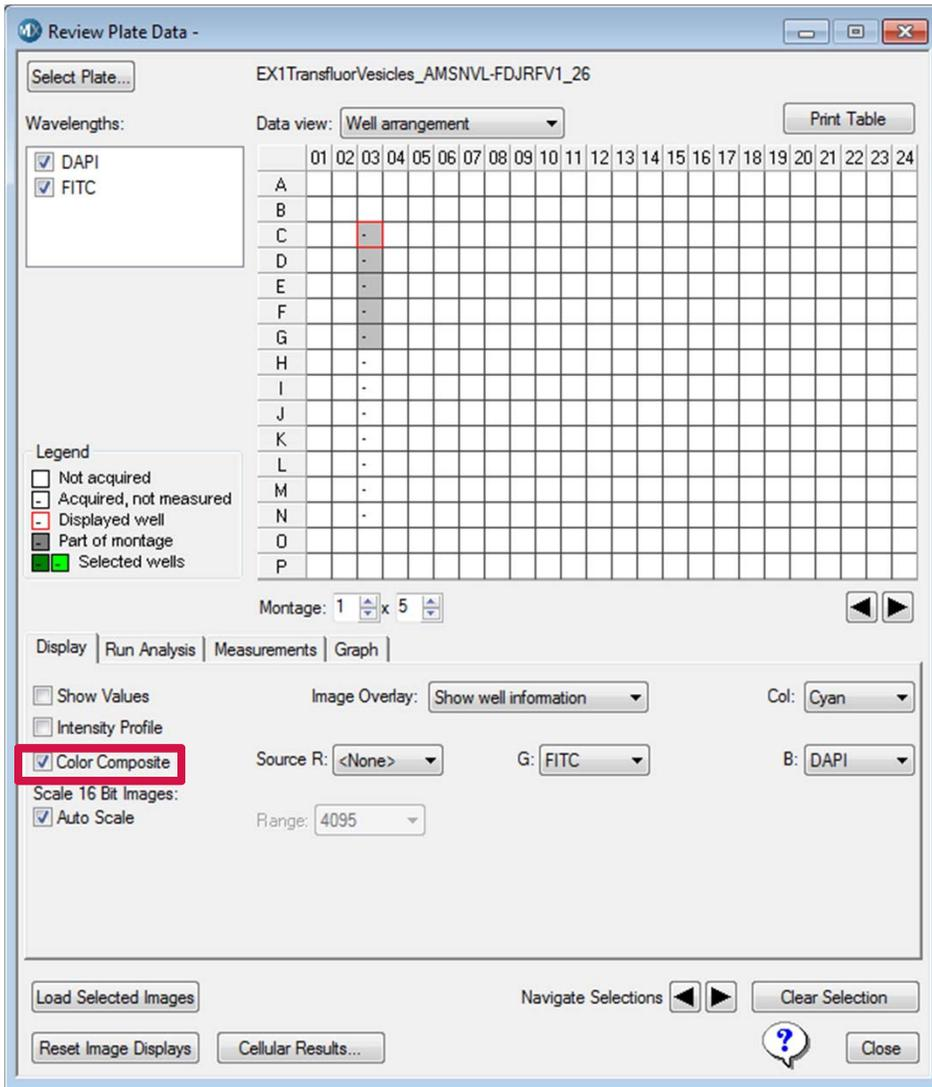
	01	02	03	04	05	06	07	08	09	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A																								
B																								
C			-																					
D																								
E																								
F																								
G																								
H																								
I																								
J																								
K																								
L																								
M																								
N																								
O																								
P																								

5. Select the **Display** tab
6. Select the appropriate wavelengths in the **Source RGB** drop-down menus

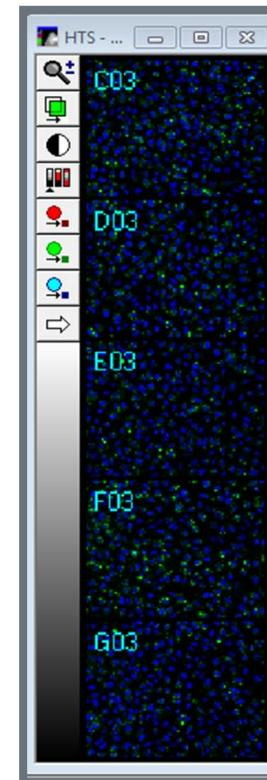
NOTE You are combining three monochrome images into a single full-color image. The “R,” “G,” and “B” annotations in this dialog stand for Red, Green, and Blue, respectively. These are the colors that will be assigned to your selected images in the resulting color overlay.



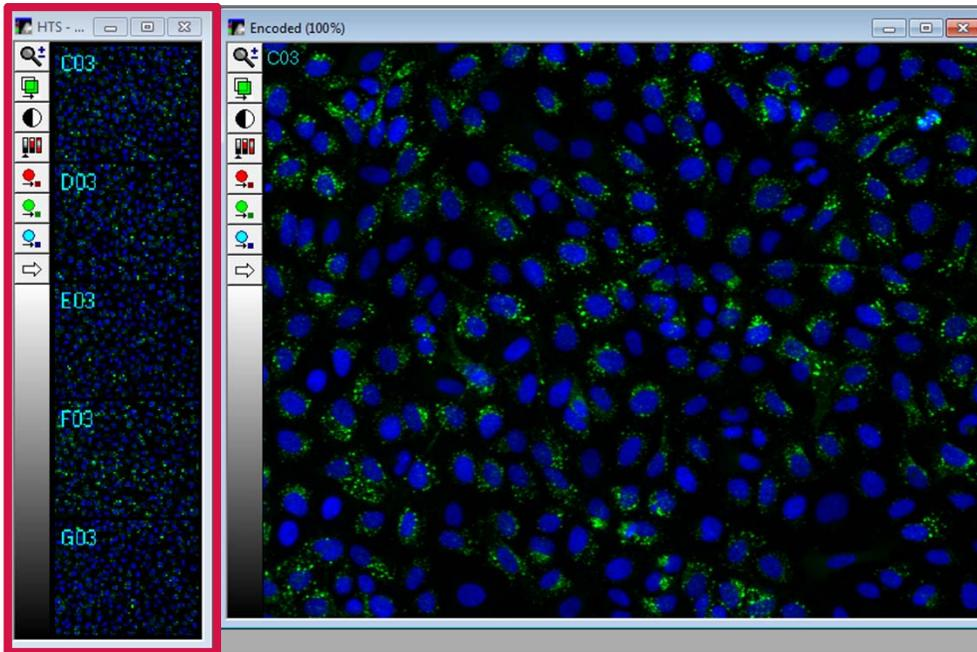
Creating Image Overlays in Review Plate Data



7. Enable **Color Composite** on the **Display** tab
 - A Thumbnail Montage will appear as a color overlay named **HTS-Encoded** in place of the individual thumbnail montages



Displaying High Resolution Color Overlay Images



8. Click in the thumbnail montage to display a high resolution color overlay image named **Encoded**
 - The high resolution image will update automatically if you click on an alternate thumbnail in the montage

NOTE To display a high resolution image, click on the thumbnail montage **ONLY**, do **NOT** click on the well arrangement within **Review Plata Data** as this will only the images displayed in the thumbnail montage.



Adjusting Display of Color Overlays

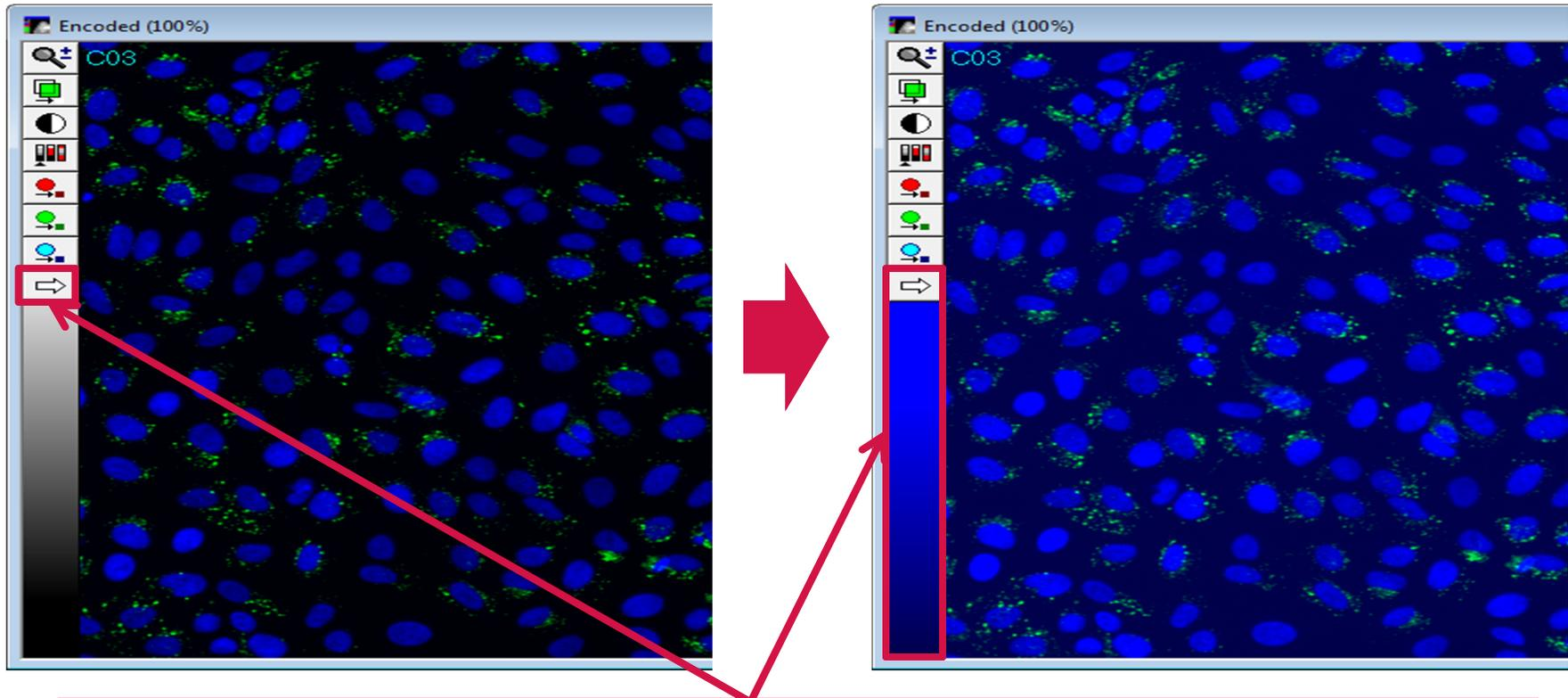
The screenshot shows the 'Review Plate Data' software interface. On the left, there is a 'Well arrangement' grid with columns 01-24 and rows A-P. Below the grid is a legend with options: Not acquired, Acquired, not measured, Displayed well, Part of montage, and Selected wells. The 'Color Composite' section is highlighted with a red box, showing 'Auto Scale' checked and a 'Range' dropdown menu open with values: 255, 1023, 4095, 16383, and 65535. On the right, there are two image windows: 'HTS - ...' and 'Encoded (100%)'. Both windows have a grayscale bar on the left side. Two orange arrows point from these grayscale bars to a text box at the bottom right.

Overlay images are auto scaled by default. To select an alternate intensity range to display in your images, disable **Auto Scale**, then select a desired intensity **Range** from the drop down menu.

To adjust brightness and contrast on an overlay image, click and drag within the grayscale bar on each image window. Dragging up increases brightness, dragging down increases contrast.



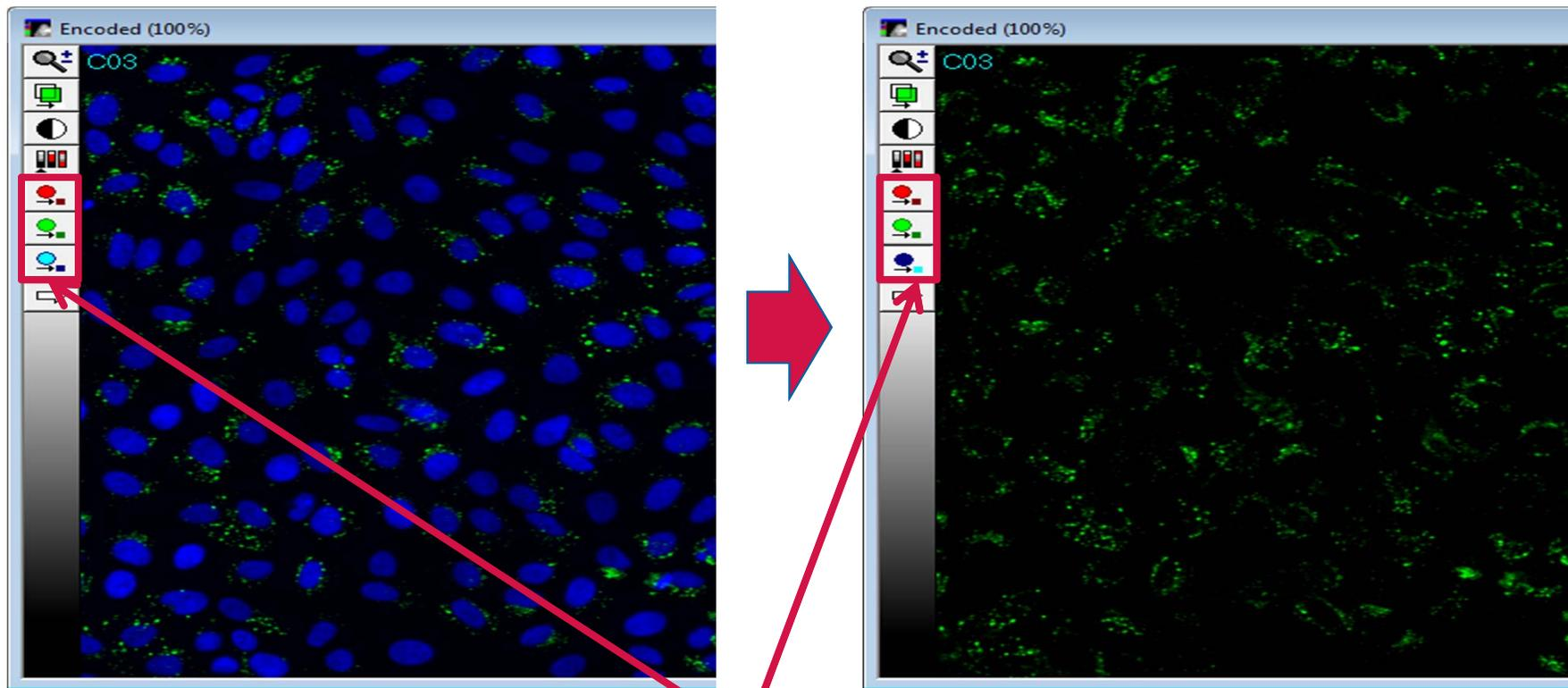
Adjusting Individual Color Channels



On the **Encoded** image window, click the **White Arrow** icon located on the left side to select the color channel to adjust independently. Click and drag the mouse within the section just below to adjust brightness and contrast. Click the **White Arrow** icon repeatedly to change between Red-Green-Blue.



Turning Individual Color Channels On or Off



Click on the individual **Red-Green-Blue** icons to toggle that particular color on or off.



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

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