



MetaXpress[®] 6 Software Guide

Viewing Images in Review Plate Data

Date Revised 06/25/15 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 viewing images post-acquisition, both in thumbnail montage and full resolution formats, through **Review Plate Data**.

Viewing Time point or Z Series images will not be covered, refer to corresponding chapters for details.



Viewing Images in Review Plate Data

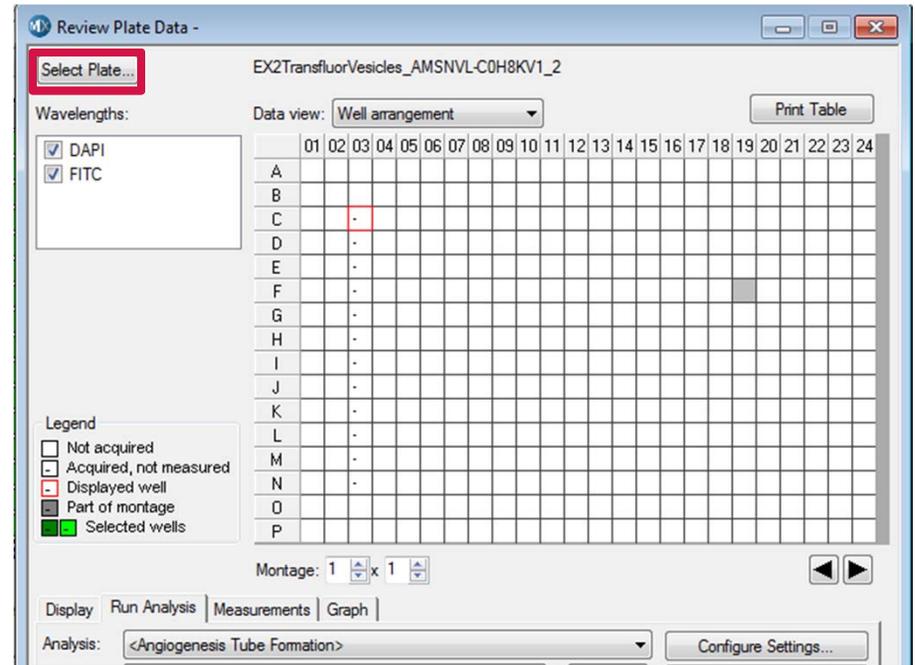
1. Open **Review Plate Data**

- In the main toolbar click on



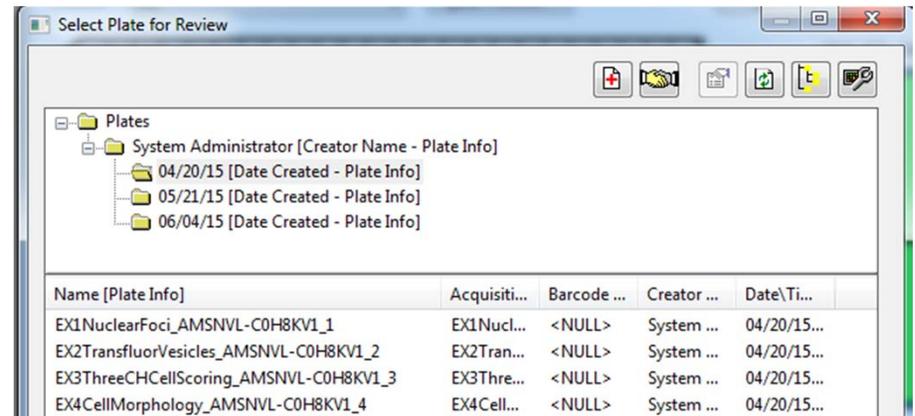
OR

- 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



Viewing Images in Review Plate Data

Review Plate Data - Transfluo-10X-Demo-19Nov14_AMSNVL-FDJRFV1_1

Select Plate...

Wavelengths: DAPI FITC

Data view: Well arrangement Print Table

	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																								

Montage: 9 x 11

Display | Run Analysis | Measurements | Graph

Show Values Intensity Profile Color Composite

Image Overlay: Show cell segmentation Col: Cyan

Source R: <None> G: <None> B: <None>

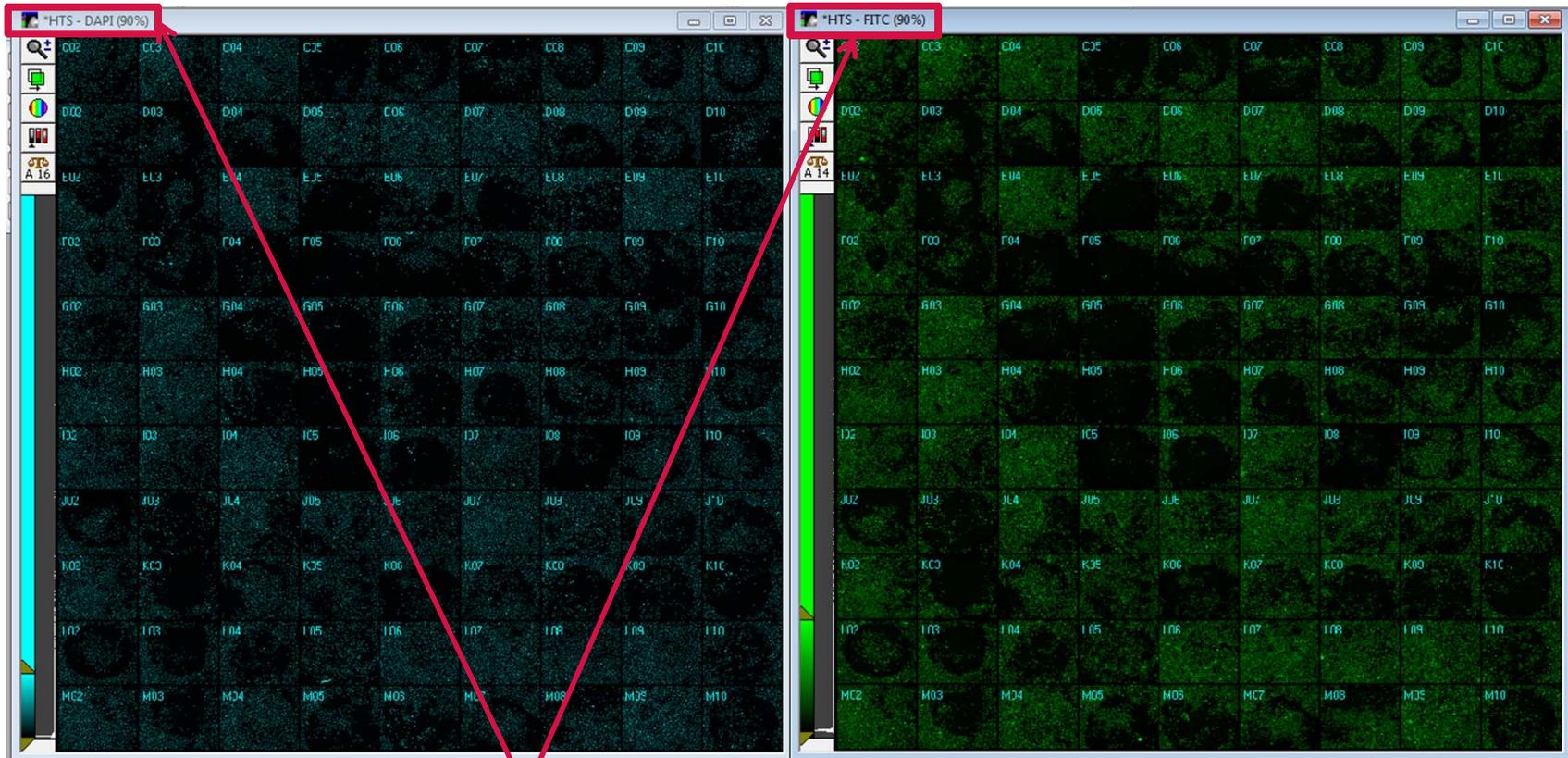
Load Selected Images Navigate Selections Clear Selection

Reset Image Displays Cellular Results... ? Close

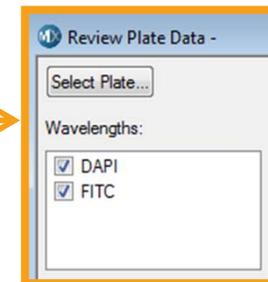
4. Left-click and drag over contiguous wells of interest (highlighted in gray) to display images in a thumbnail montage (see section on Thumbnail Montage)



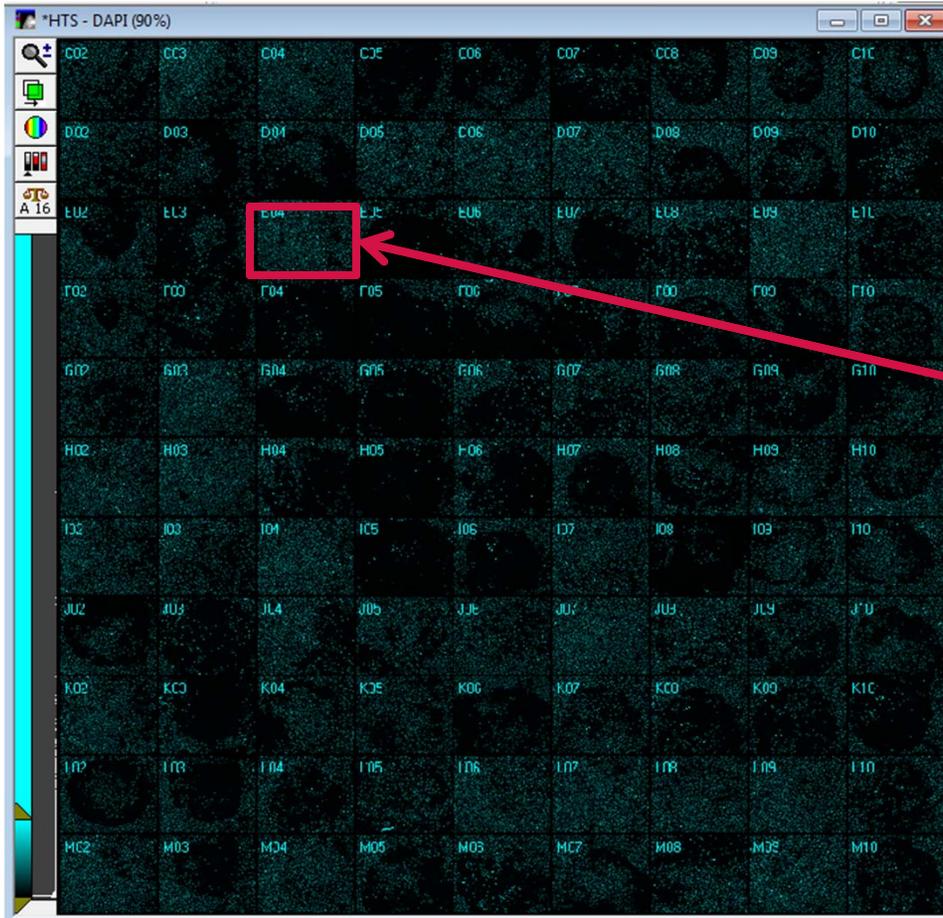
Viewing the Thumbnail Montage



Thumbnail montage(s) appear, one for each wavelength selected in the **Wavelengths** section of the **Review Plate Data** dialog. These montages will have “HTS” as the prefix in their image names.



Viewing High Resolution Images

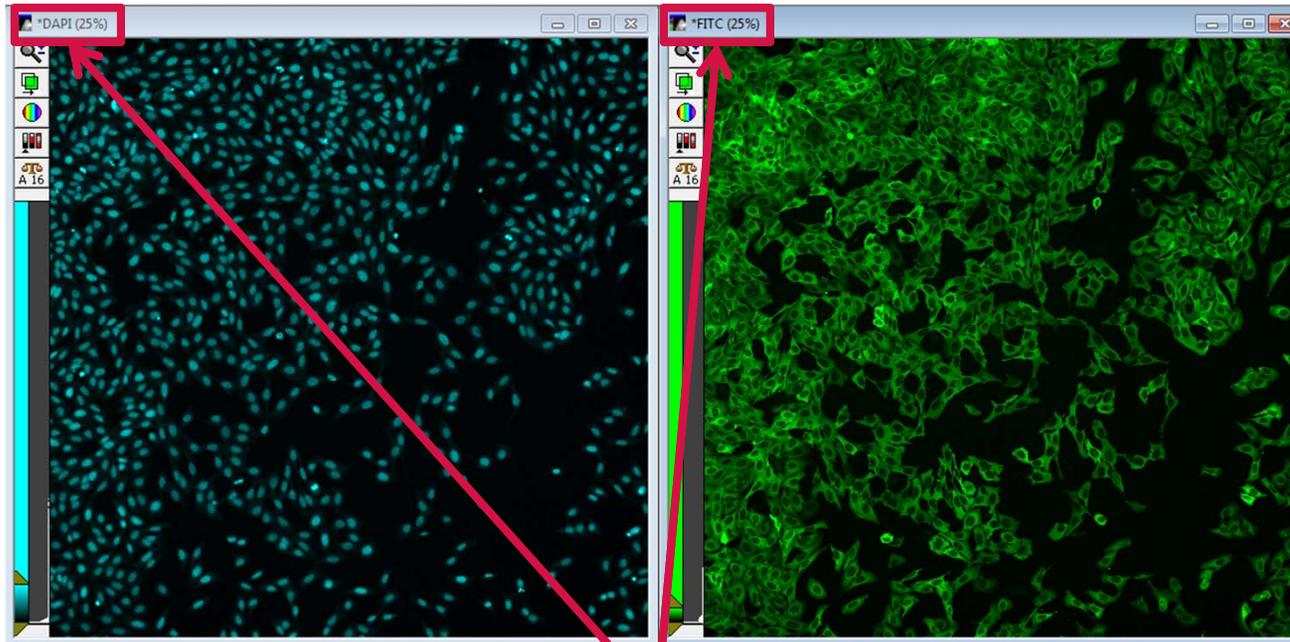


Click on a thumbnail image for the well of interest within the montage to open its corresponding high resolution images.

NOTE Do NOT click on the plate grid within **Review Plate Data** at this point, as this will only result in changing the wells that appear in the thumbnail montage.



Viewing High Resolution Images

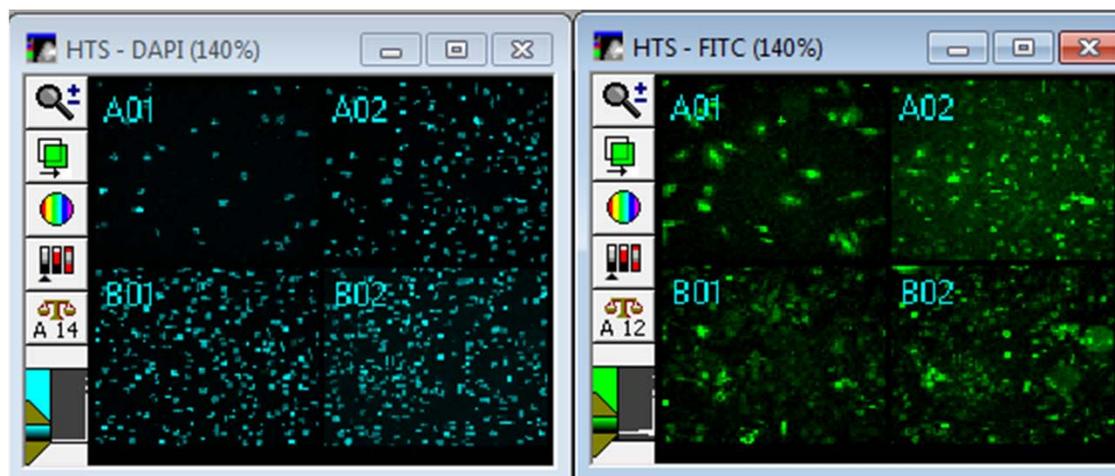
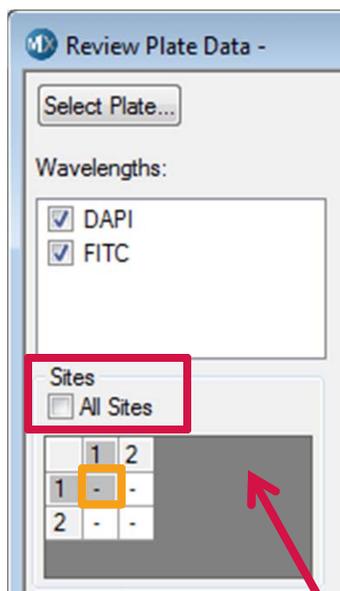


Full resolution copies for all wavelengths for the corresponding well/site will now appear. They will be titled with the wavelength name ONLY (no “HTS” prefix).

NOTE To view a different site/well in high resolution format, do not close the high resolution images. click on the thumbnail image for that well WITHIN THE MONTAGE and the images will automatically update.



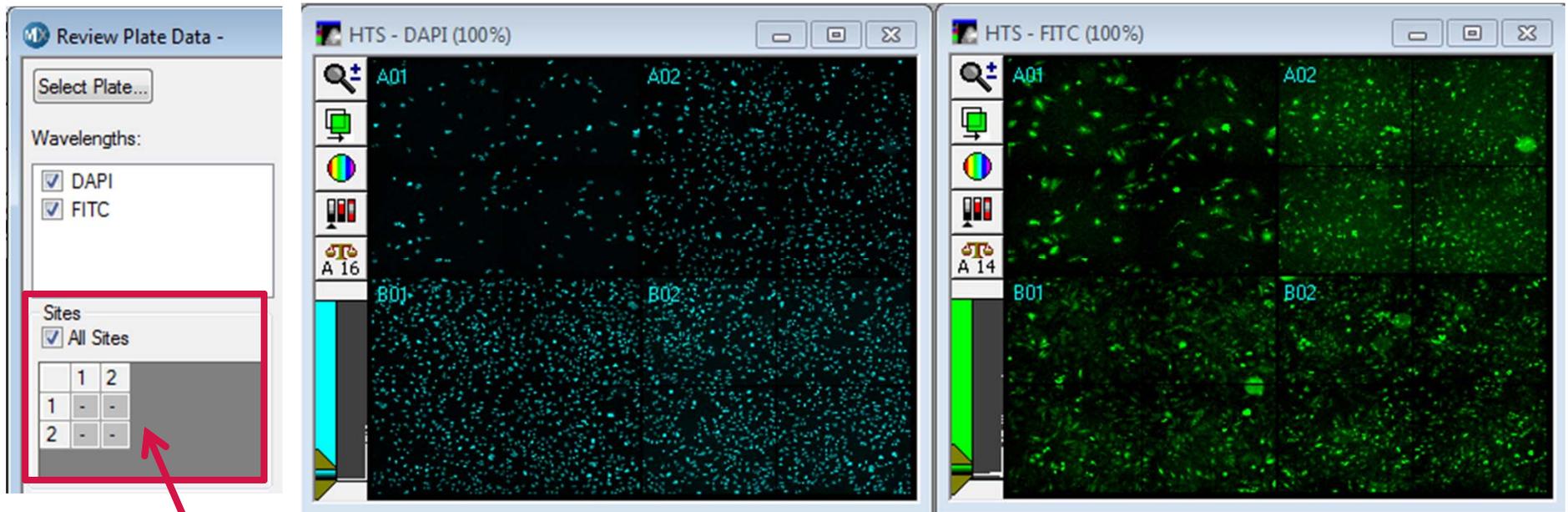
View Thumbnail Montage for a Single Site in Each Well



- To display a thumbnail montage showing a single site in each well:
- Disable **All Sites**
 - Select the site of interest within the site map (highlighted in gray)
 - Click and drag to select your wells of interest in plate view
 - To display a different site in the thumbnail montage, left-click on that site



Displaying Thumbnail Montage for All Sites-Multiple Wells



- To display a thumbnail montage showing all sites in each well:
- Enable **All Sites**
 - Click and drag to select your wells of interest in plate view

NOTE The exact distribution of sites within the well will be reflected in both the site map AND the resulting thumbnail montage. Empty sites will be represented by blank images in the montage.



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.