



MetaXpress[®] 6 Software Guide

Viewing Timelapse Image Sets in Review Plate Data

Date Revised 06/09/15 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

This purpose of this chapter is to guide the user through viewing Timelapse image sets in **Review Plate Data** post-acquisition. Timelapse images can be viewed as a thumbnail montage (filmstrip-like view), a series of individual high resolution images, or as a single stack (.stk) file.



Viewing Timelapse Images

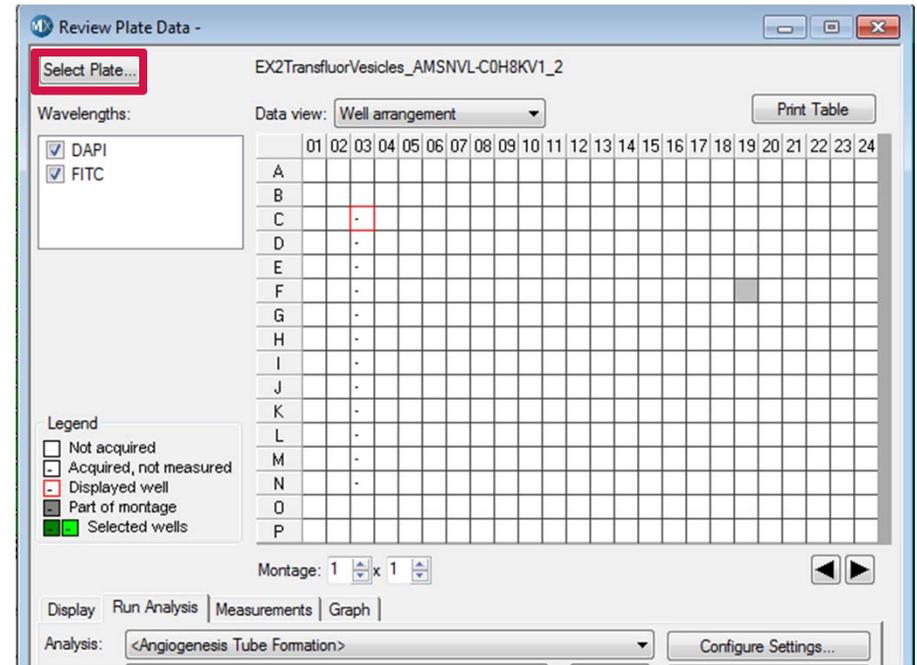
1. Open **Review Plate Data**

- In the main toolbar click on



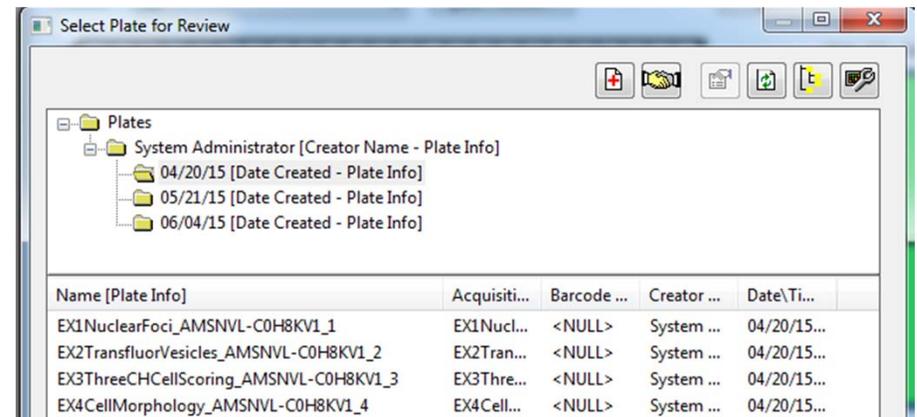
OR

- 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



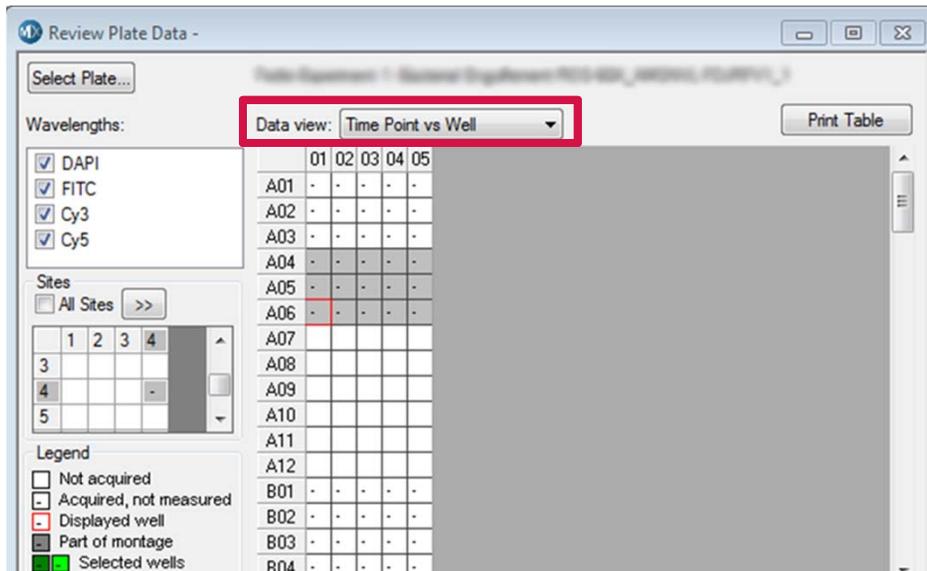
Viewing Multiple Wells – Single Time point

- To view a thumbnail montage of multiple wells at a single time point, change the **Data View** to **Well Arrangement**
 - Select desired **time point** in the Time point section
 - Left-click and drag to select a range of wells to display as a thumbnail montage

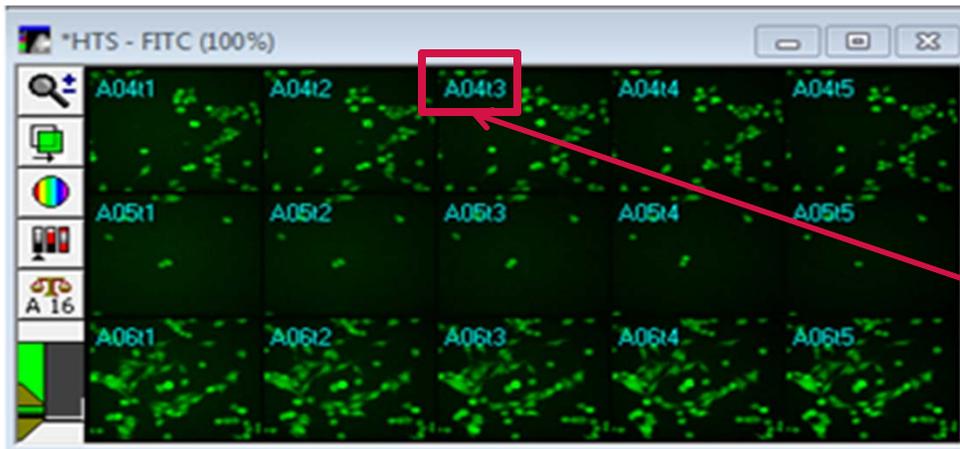
The screenshot displays the 'Review Plate Data' software interface. The main window is titled 'Fadie-Experiment 1- Bacterial Engulfment-ROS-60X_AMSNVL-FDJRFV1_1'. The 'Data view' dropdown is set to 'Well arrangement', and a red box highlights a 4x4 grid of wells (A1-D4) in the well arrangement grid. The 'Time points' dropdown is set to '3 of 5'. The interface shows four fluorescence images: 'HTS - DAPI (80%)', 'HTS - Cy3 (80%)', 'HTS - FITC (80%)', and 'HTS - Cy5 (80%)'. The DAPI image shows blue fluorescence, Cy3 shows red, FITC shows green, and Cy5 shows yellow. The well arrangement grid is also visible in the bottom right corner of the main window.



Viewing Multiple Wells – Multiple Time Points



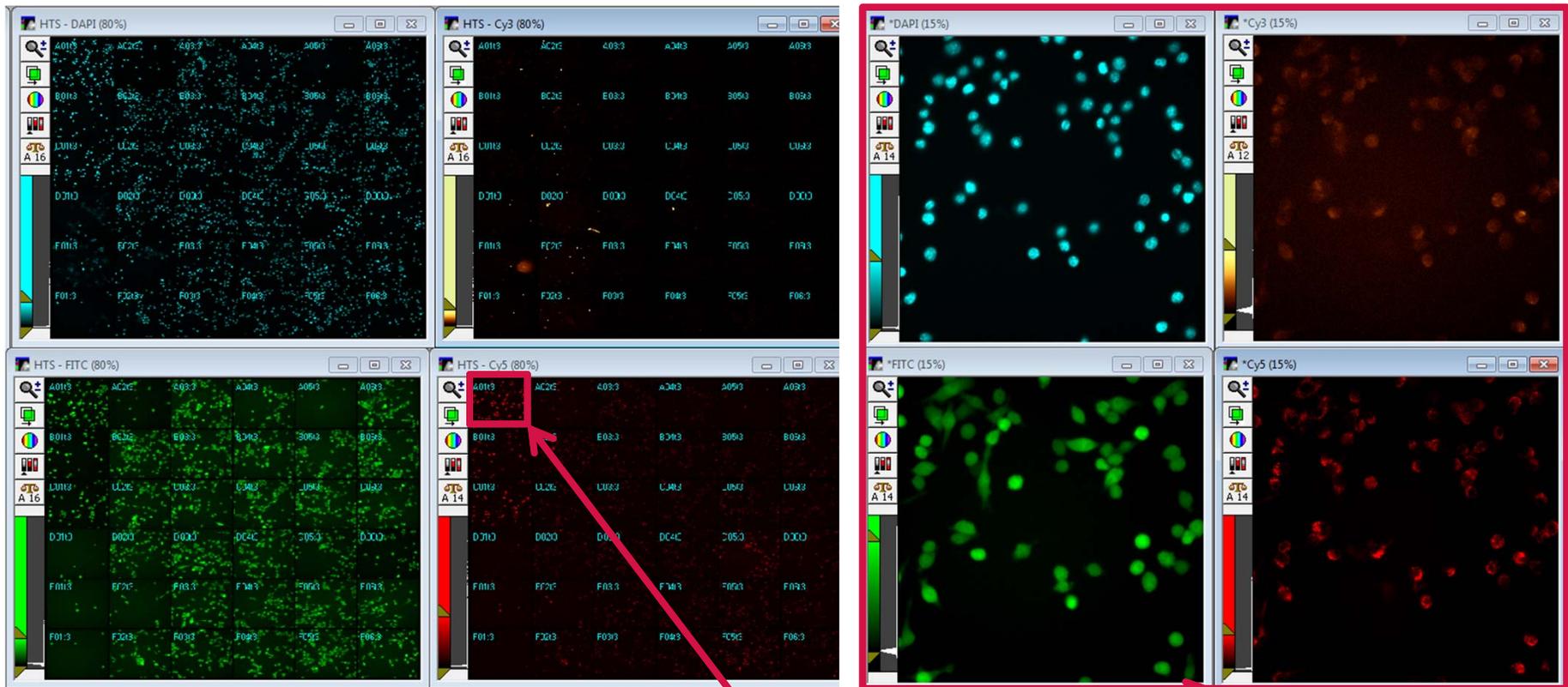
- To view a thumbnail montage (filmstrip) of multiple time points, change the **Data View** to **Time Point vs Well**
 - Plate View now changes such that time points are arrayed in the X direction and the Well ID arrayed in the Y direction
 - Click and drag to select a range of time points or select a range of adjacent wells to create a time point thumbnail montage



Time lapse montage(s) for the selected well(s). Image annotation indicates both well ID and time point (t1, t2, t3, etc.) in the upper left corner of each thumbnail image.



Viewing High Resolution Images



To review high resolution images from selected wells, simply click on any thumbnail image in any of the montages to open the corresponding high resolution image set.



Viewing Time Points as a Stack

- To create a stack (.stk) file of a series of time points from one site/well
 - Right-click on a well of interest either in the Plate View or in the thumbnail montage
 - Corresponding images and well will be highlighted in green
 - Click on the **Load Selected Images** button

NOTE If you have more than one site per well, then deselect **All Sites** and highlight only one site to view.

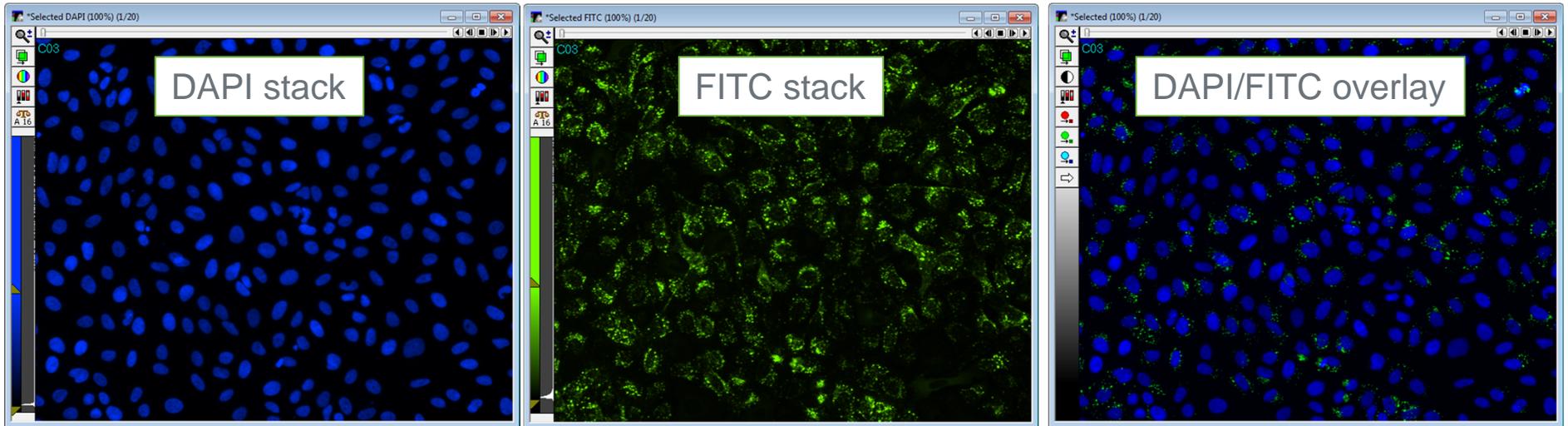


The screenshot shows the 'Review Plate Data' window for 'Transfluor 20x_AMSNVL-HGC8KV1_65'. The 'Data view' is set to 'Time Point vs Well'. The grid shows wells C01 through C24 and D01 through D03. Row C18 is highlighted in green. A red arrow points from a text box 'Right-click to select well(s) of interest' to the C18 row. Another red arrow points from the 'Load Selected Images' button to the thumbnail montage.

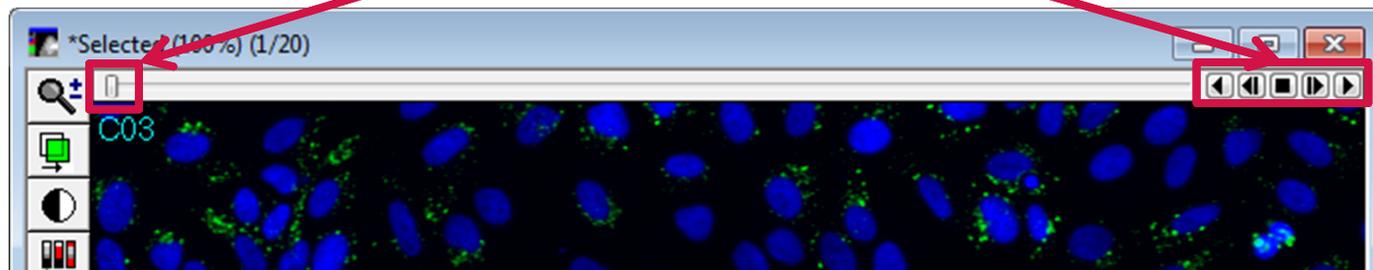
	01	02	03	04	05	06	07	08	09	10	11	12	13	14	15	16	17	18	19	20
C12																				
C13																				
C14																				
C15																				
C16																				
C17																				
C18																				
C19																				
C20																				
C21																				
C22																				
C23																				
C24																				
D01																				
D02																				
D03																				



Scrolling Through Time Lapse Images



To scroll through the timelapse, click and drag the image slider or use the playback controls.



Viewing Individual Wavelengths or Color Overlays

- Click on the **Display** tab to turn color composite on or off
 - To view channels separately uncheck **Color Composite**
 - To view a color overlay of channels, enable **Color Composite** and select the corresponding R, G, and B images from the source drop-down menus
 - Refer to chapters on overlaying images for more details
- Click on **Load Select Images** to load the Z Series into a stack (.stk) file

Review Plate Data - Transflur 20x_AMSNVL-HGC8KV1_65

Select Plate...

Wavelengths: DAPI FITC

Sites All Sites

	1	2
1	-	-
2	-	-

Legend Not acquired Acquired, not measured Displayed well Part of montage Selected wells

	01	02	03	04	05	06	07	08	09	10	11	12	13	14	15	16	17	18	19	20
C12																				
C13																				
C14																				
C15																				
C16																				
C17																				
C18																				
C19																				
C20																				
C21																				
C22																				
C23																				
C24																				
D01																				
D02																				
D03																				

Montage: 20 x 1 Time points: 1 of 20

Display Run Analysis Measurements Graph

Show Values Image Overlay: Show well information Col: Cyan

Intensity Profile

Color Composite Source R: <None> G: FITC B: DAPI

Scale 16 Bit Images: Auto Scale Range: 16383

Load Selected Images Navigate Selections Clear Selection

Reset Image Displays Cellular Results... ? Close



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