



MetaXpress® 6 Software Guide

Viewing Z Series Image Sets in Review Plate Data

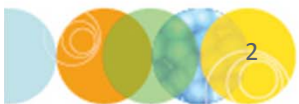
Date Revised 06/09/15 Version B



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

Chapter Purpose

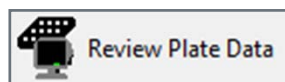
The purpose of this chapter is to guide the user through viewing Z Series images in **Review Plate Data** post-acquisition in both thumbnail montage and full resolution formats. Z Series images can be viewed as a thumbnail montage (filmstrip-like view), a series of individual high resolution images, or in a single stack (.stk) file.



Viewing Z Series Images

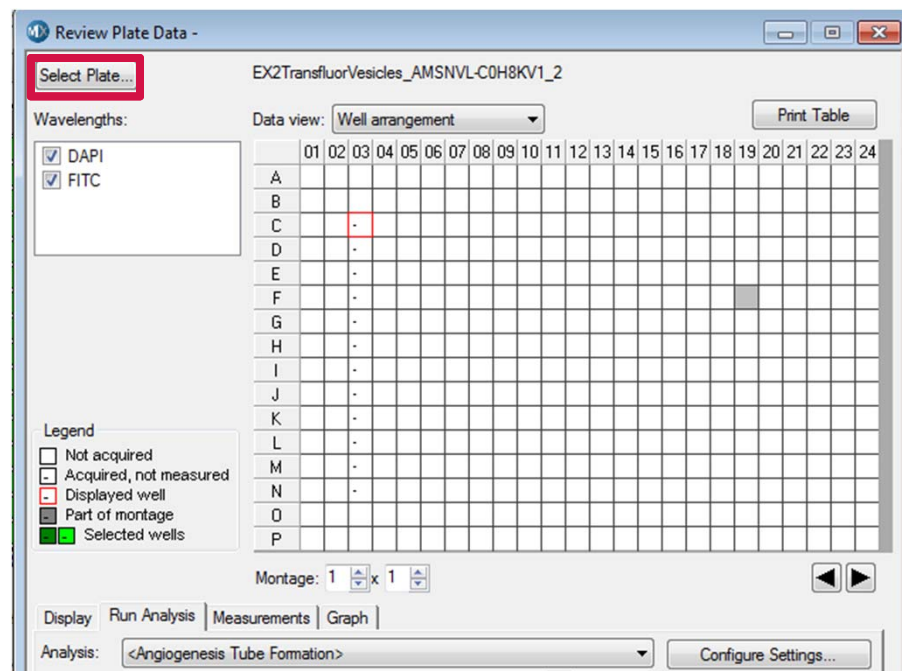
1. Open **Review Plate Data**

- In the main toolbar click on



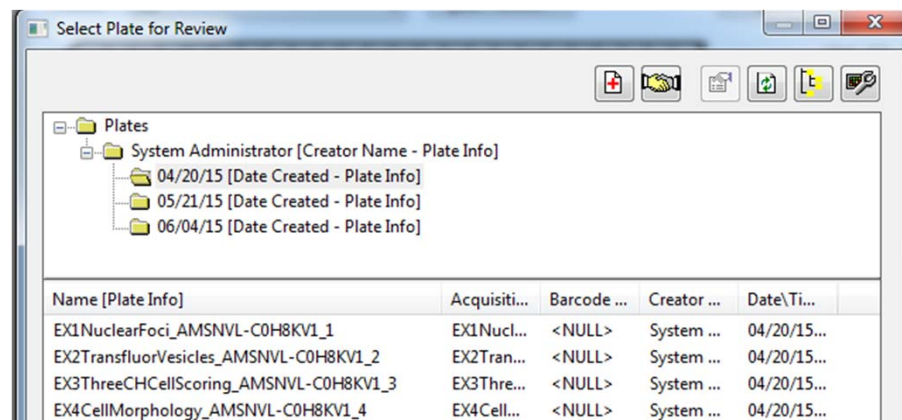
OR

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



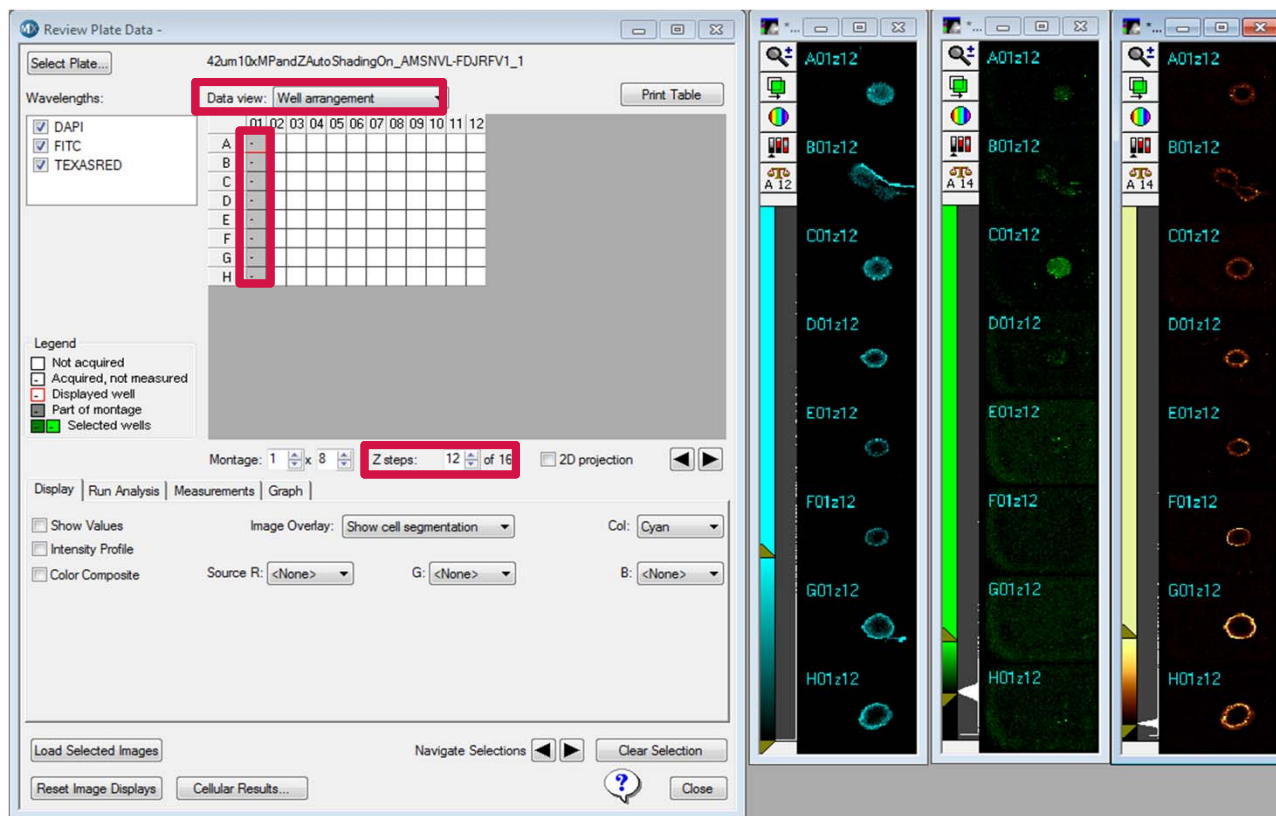
2. Click on the **Select Plate** button

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

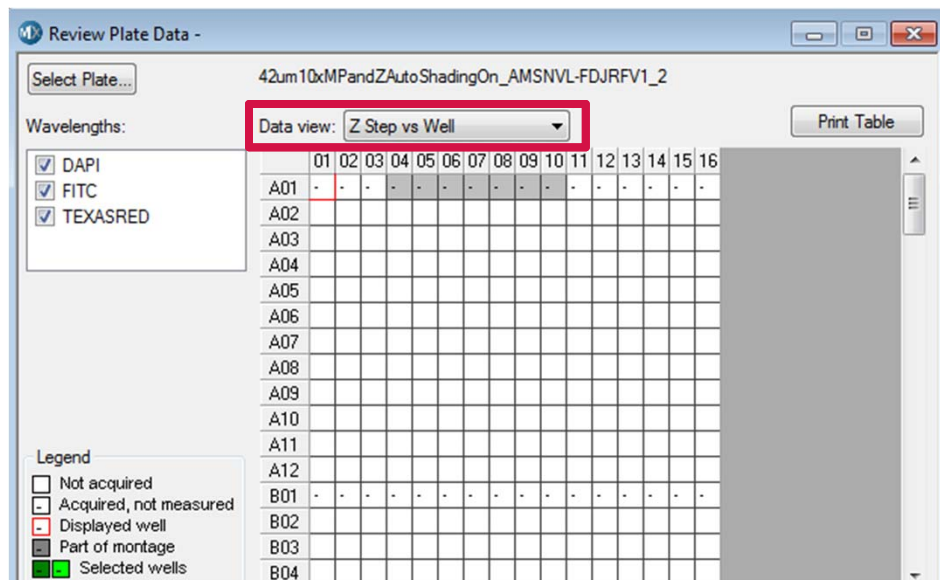


Viewing Single Z Step from Multiple Wells

4. In order to view a thumbnail montage of multiple wells at a single Z step, change the **Data View** to **Well Arrangement**
- Select desired **Z Step** in the Z Step Section
 - Left-click to select a range of wells to display in thumbnail montage



Viewing Multiple Z Steps from Multiple Wells



5. In order to view a thumbnail montage (filmstrip) of multiple Z Series, change the **Data View** to **Z Steps vs Well**

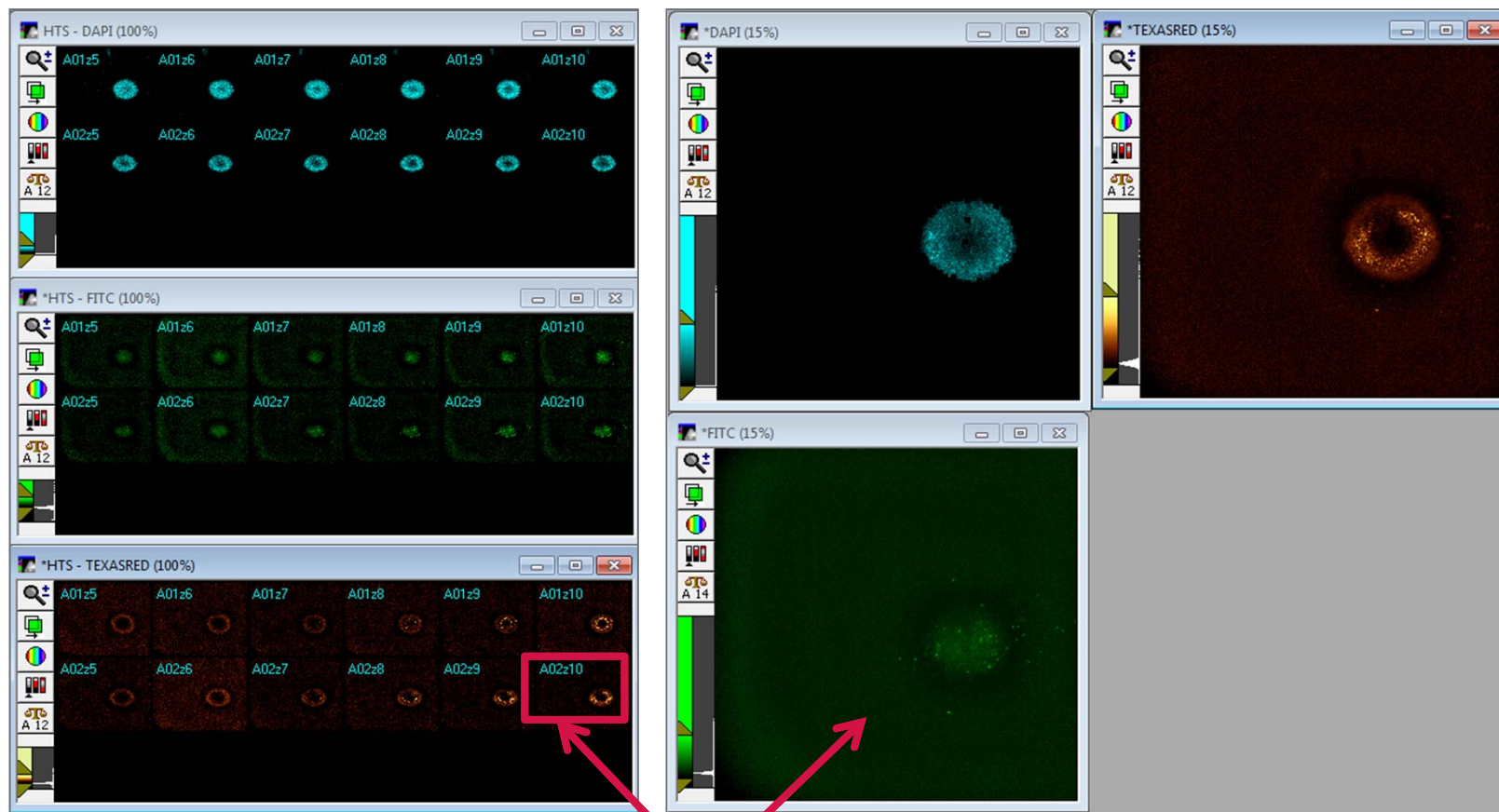
- Plate View now changes such that Z steps are arrayed in the X direction and the Well ID arrayed in the Y direction
- Click and drag to select a range of Z steps or select a range of adjacent wells to create a Z Series thumbnail montage



Z step montage for your selected well(s). The annotation indicates both well ID and Z step (z1, z2, z3, etc.) in the upper left corner of each thumbnail image.



Viewing High Resolution Images



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

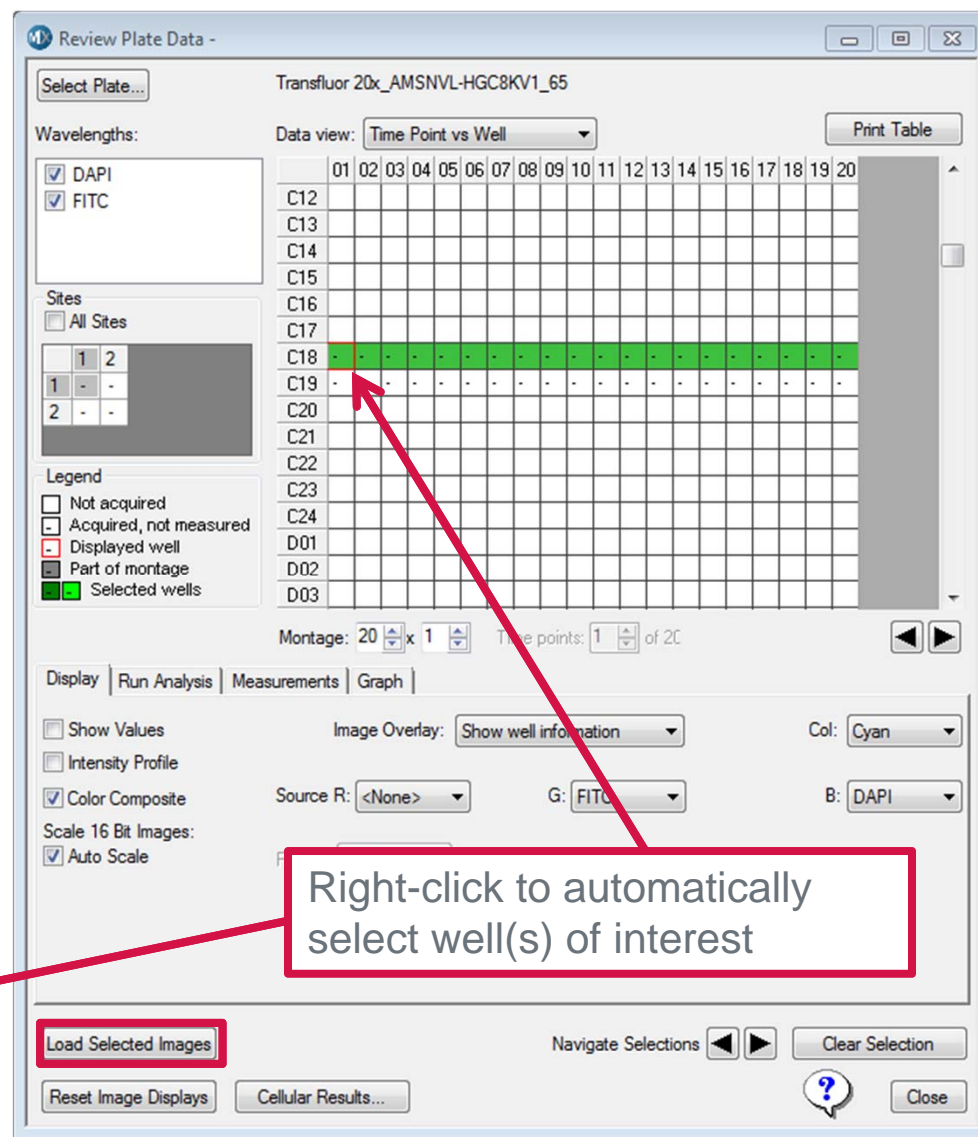
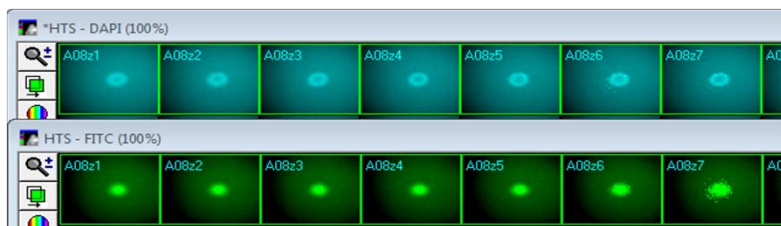


Viewing Z Series as a Stack

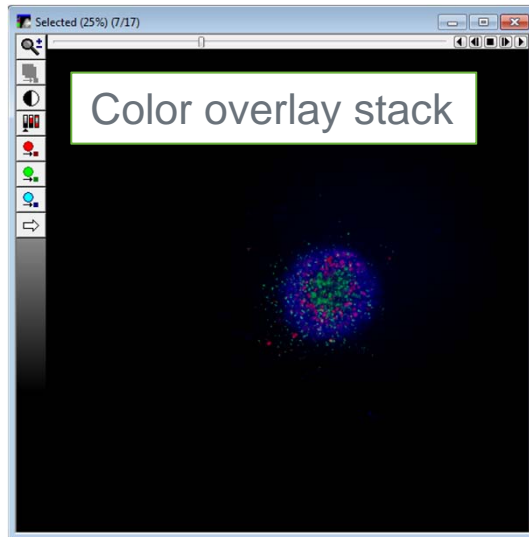
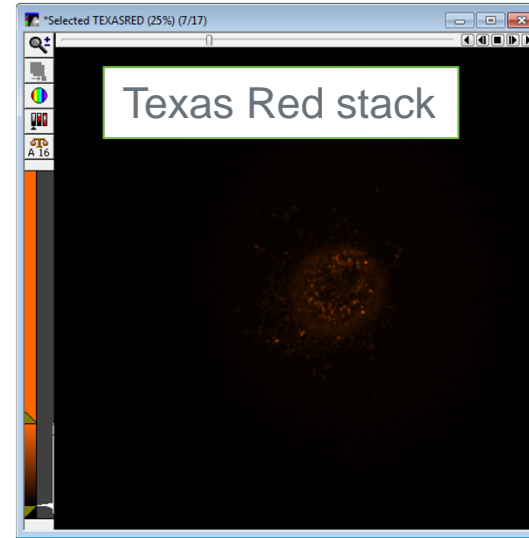
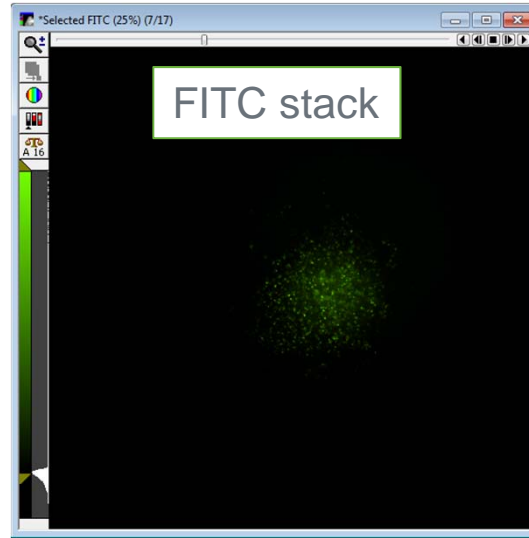
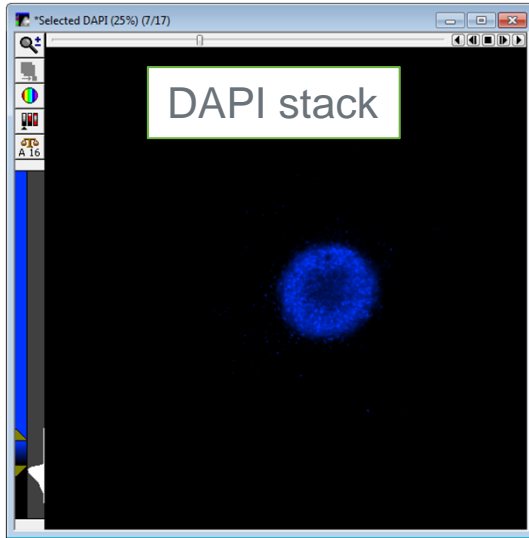
6. To create a stack (.stk) file of a Z Series of images from one site/well

- 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

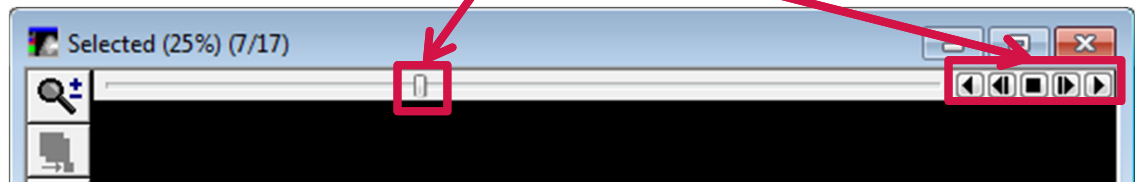
NOTE If you have more than one site per well, then deselect **All Sites** and highlight only one site to view.



Scrolling Through Z Series Images in a Stack

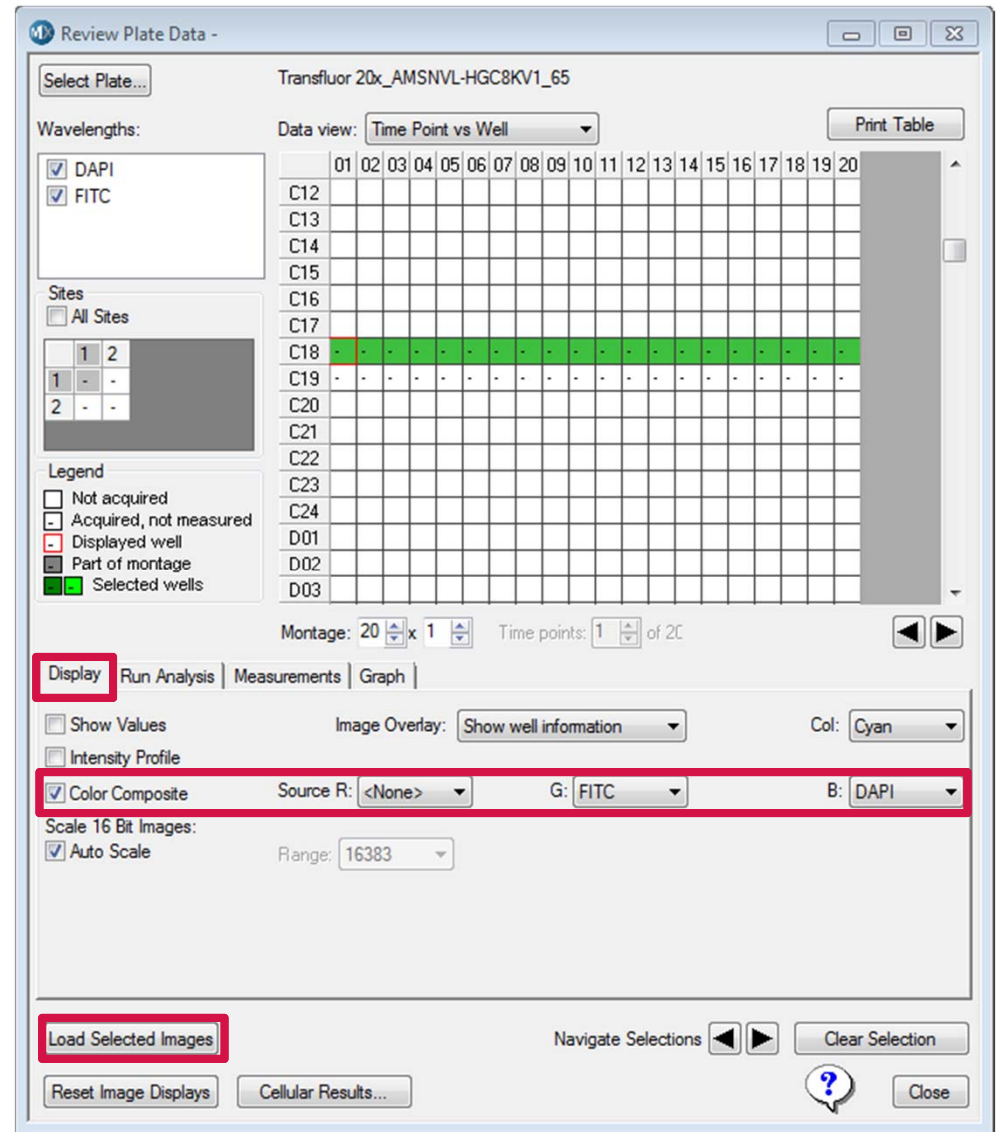


To scroll through the Z stack, click and drag the image slider or use the playback controls.



Viewing Individual Wavelengths or Color Overlays

7. Select 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 corresponding chapters on creating overlay images
6. Click on **Load Select Images** button to load the Z Series into a stack (.stk) file



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