



# MetaXpress® 6 Software Guide

Creating a Movie from Timelapse or Z Series Images

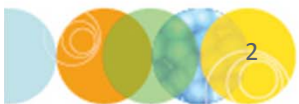
Date Revised 06/09/15 Version B



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

# Chapter Purpose

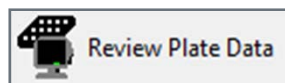
The purpose of this chapter is to guide the user in creating movie files (AVI) post-acquisition from a Timelapse or Z Series set of images through **Review Plate Data**.



# Creating Movies

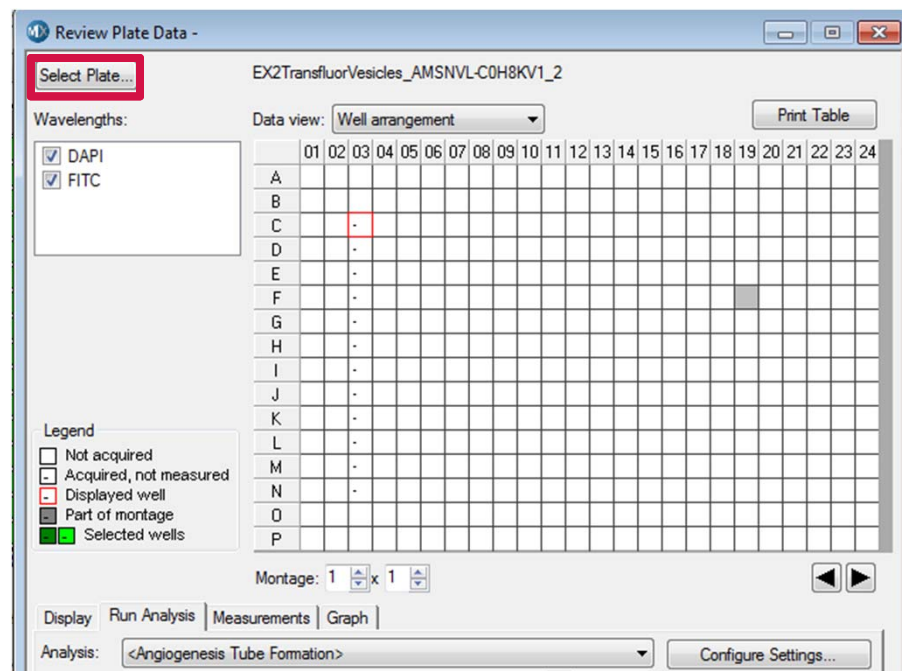
## 1. Open **Review Plate Data**

- In the main toolbar click on



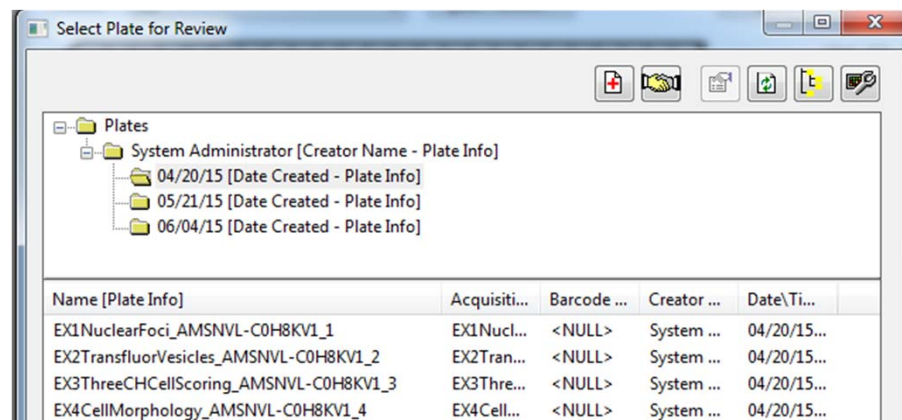
OR

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



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

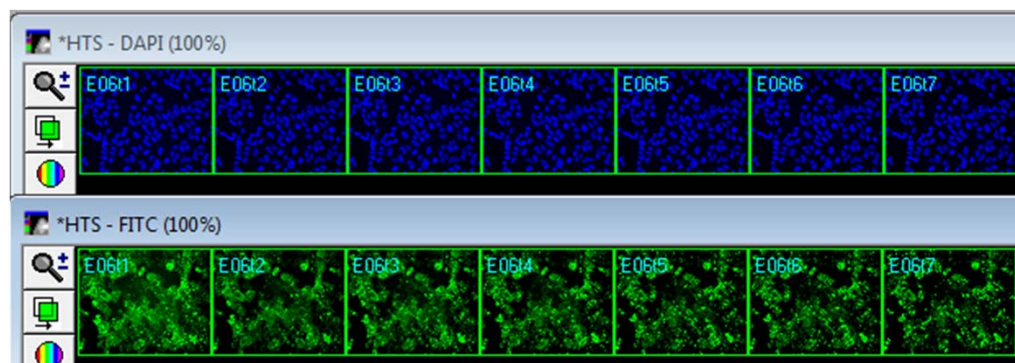
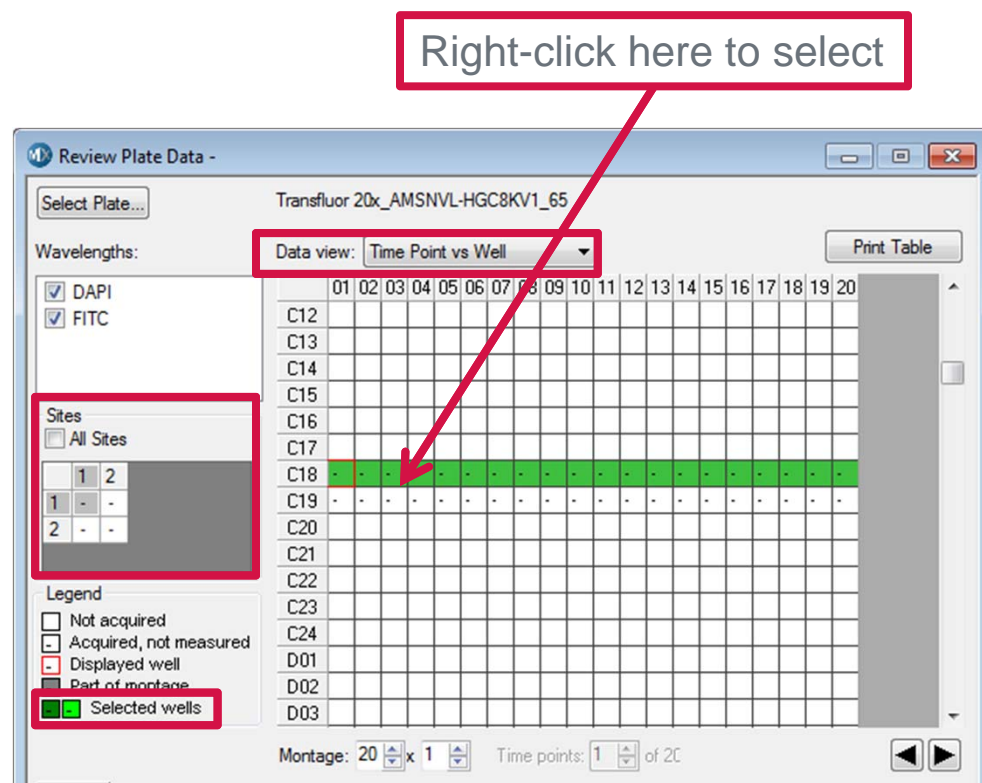
## 3. Browse through the folders to open the plate of interest



# Creating Movies

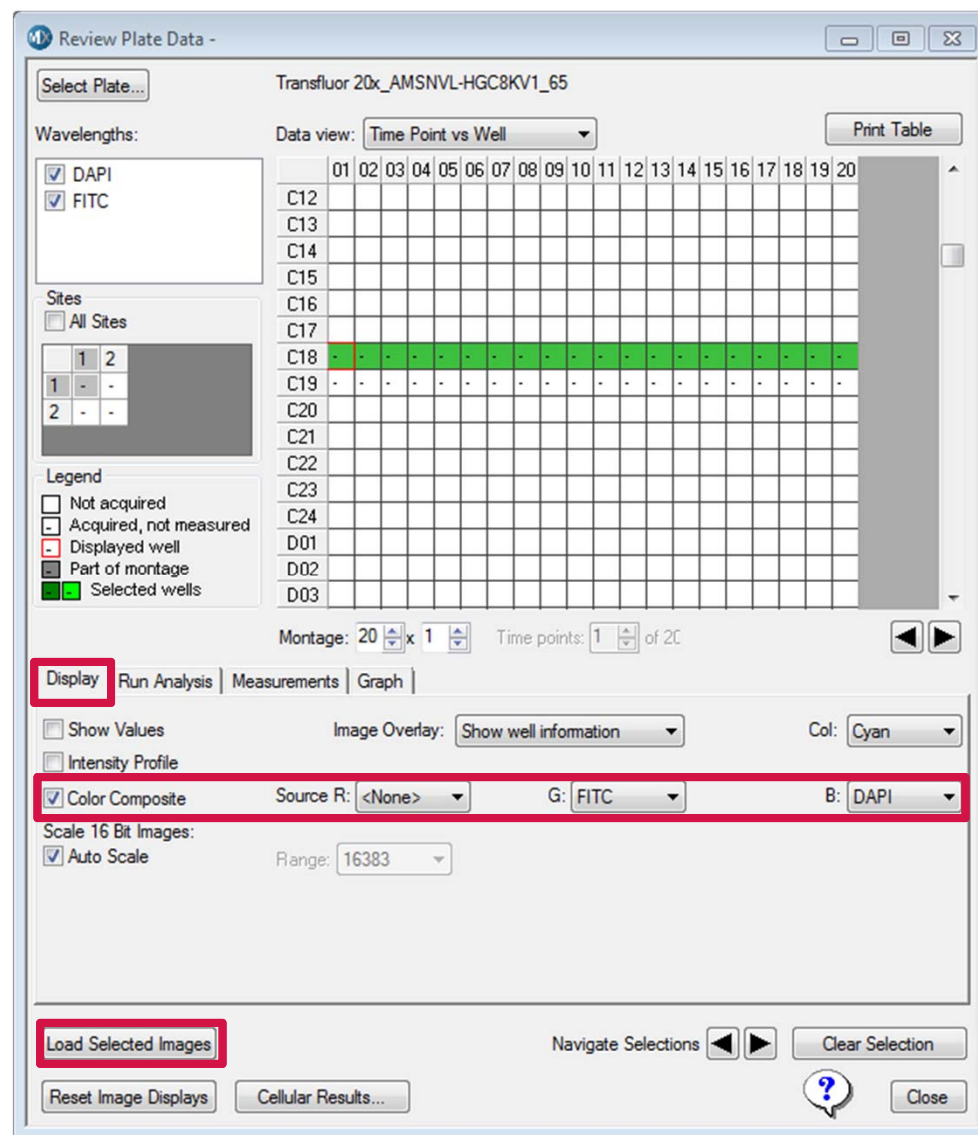
3. Change the **Data View** to:
  - Z Step vs Well
  - Time Point vs Well
4. Right-click on any Time point or Z Series in the Plate View or thumbnail montage to select all Time points or Z Series for the well of interest
  - Selected well will be highlighted in green
  - Corresponding images in thumbnail montage will also be highlighted in green in the thumbnail montage

**\*NOTE\*** If you have more than site per well, then deselect **All Sites** and click on the site of interest.



# Creating Movies

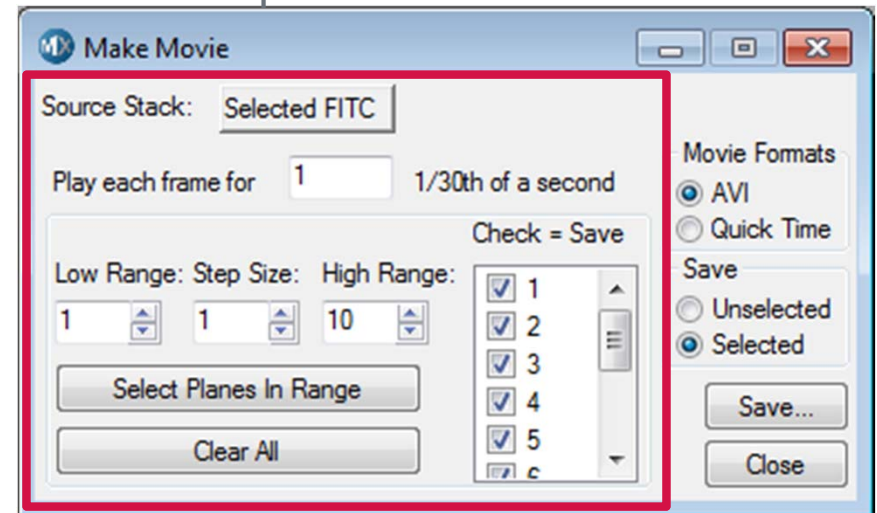
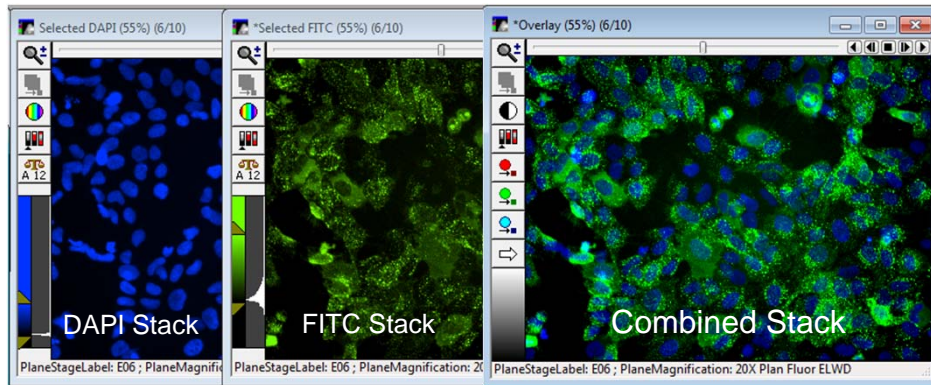
5. Select the **Display** tab to turn **Color Composite** on or off
  - To view channels separately deselect **Color Composite**
  - To view a color overlay of channels, enable **Color Composite** and select the corresponding R, G, and B images from the source drop-down menus
  - See corresponding chapter for creating overlay images
6. Click on **Load Select Images** to load the Time points or Z Series into a stack (.stk) file





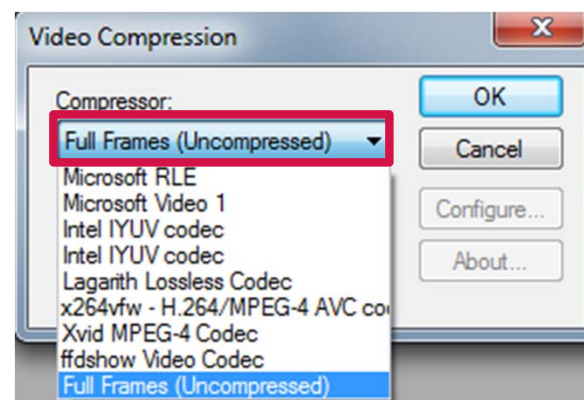
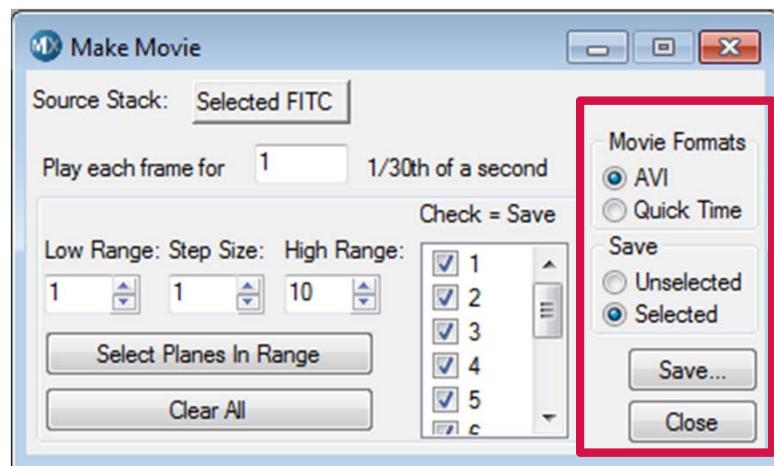
# Creating Movies

7. Adjust image scaling (brightness and contrast) as desired
8. In the main menu select **Edit > Stack > Make Movie** to open the **Make Movie** dialog
  - Click on the button next to **Source Stack** and select the .stk file you want to make a movie from
  - Enter a value for **Play each frame for**: enter 30 to play each frame for 1 second in the movie (typical film is 30 frames per second)
  - Select Planes in Range: select/deselect planes to include in the movie or click on the **Select Planes in Range** button to select all planes



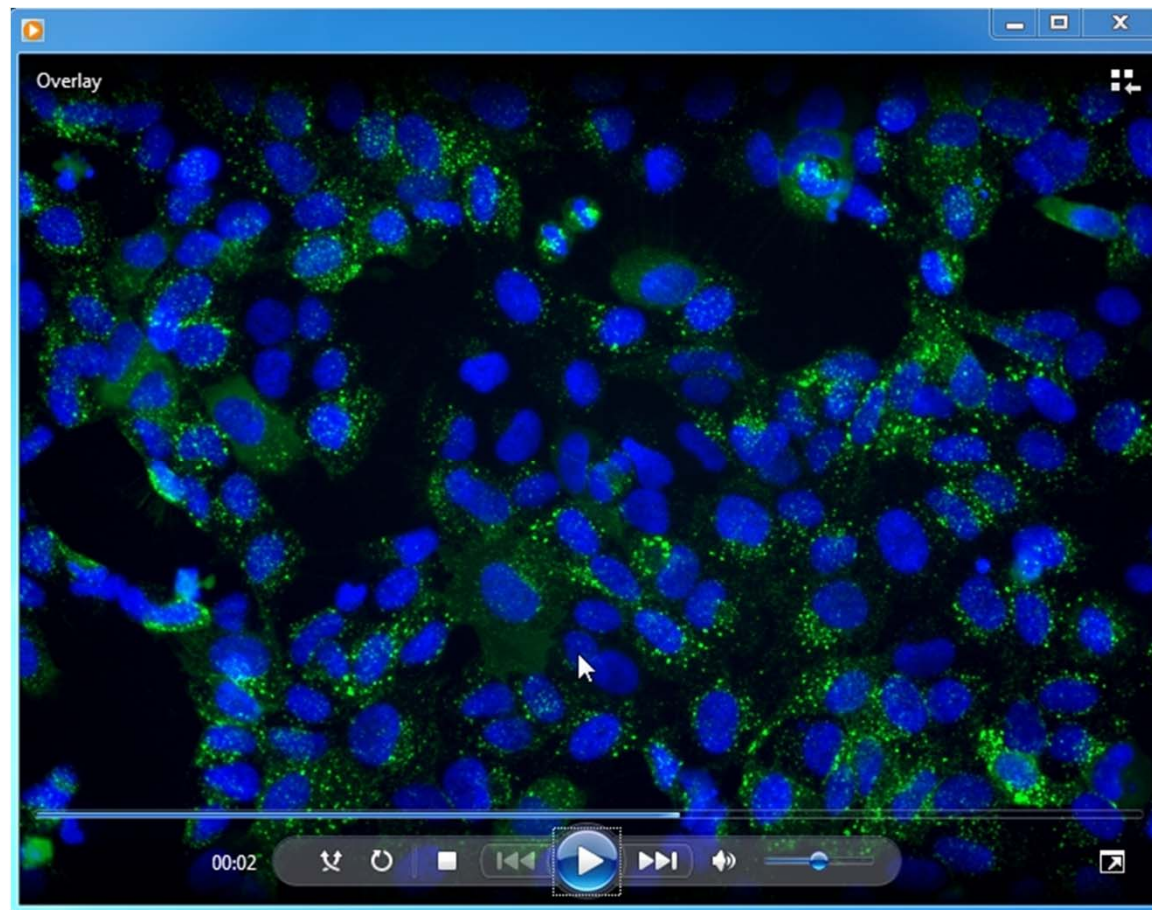
# Configuring Movie Format

9. Under **Movie Formats**, select **AVI**
  10. Under the **Save** section, select **Selected** and click on the **Save** button
    - Navigate to a directory and name the movie file
  11. In the **Video Compression** dialog, select **Full Frames (Uncompressed)**
    - This will create an uncompressed movie file
    - To ensure wide compatibility, Molecular Devices does not recommend using the compression choices in the drop down menu.
- \*NOTE\* Third party software can be used to compress the movie if desired*



# Playing the Movie

11. AVI movie files can be opened with Windows Media Player or most other software media players.



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



# 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