

ImageXpress[®] Micro XLS & MetaXpress[®] 6 for systems with Sutter XL Light Source



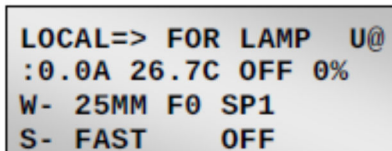
The purpose of this guide is to briefly describe:

- I. Turn on system and acquire plate with saved settings
- II. Test acquisition settings
- III. Define new acquisition settings
- IV. View images and run an analysis

I. Turn on System and Acquire Plate with Saved Settings

Turn on the system:

- Light source (if not already on)
 - a. * Always power on the light source before launching MetaXpress software. Please consult the Lambda XL Light Source quickstart guide and manual for more information on operation.
 - Turn on the power switch in the back of the light source close to the power cord.
 - The Lambda XL light source powers on and 1 minute later, the LED display should look similar to this:



- Press **Local** on the XL light source keypad.
- Press **1** to start the lamp.
- Please make sure “U@” is showing at the upper right hand corner. If “U@” is not showing at the upper right hand corner, press Local key on the keypad until @ shows up in the upper right hand corner, then press Online key to switch to U mode.
- The XL light source is powered up and ready to use. Please see page 14 on proper procedure for powering down the light source. (WARNING: DO NOT STOP the lamp within 1 minute of starting the lamp to avoid damaging the lamp and reduce life time of the lamp)

- IXM power supply controller box
- IXM options controller box (for Transmitted light, Environmental control or Fluidics modules)
- Computer and Monitor

2. Go to the MetaXpress folder and double-click on the appropriate hardware profile shortcut

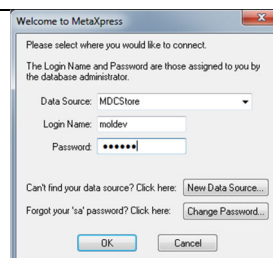


Login to MDCStore database with username and password

3.

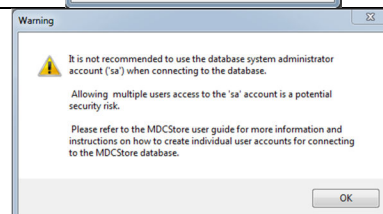
Username	moldev
Password	moldev

NOTE Your database, username, and password maybe different.
Refer to your administrator for this information



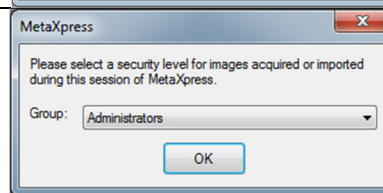
4.

If you log in as system administrator (sa), the next window is a warning regarding security risks; click **OK**

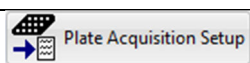


5.

Select Group (security level) and click **OK**



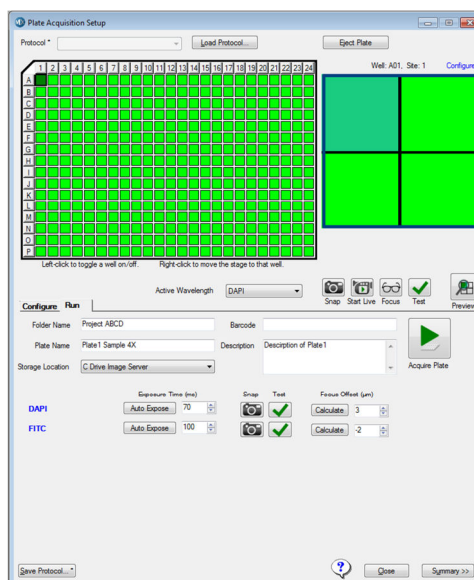
In the main toolbar, click



or in the main menu select **Screening > Plate Acquisition Setup**

Acquisition Setup

6.

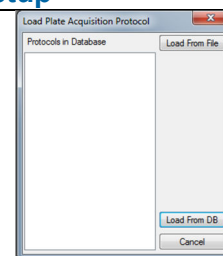




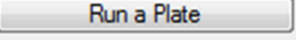
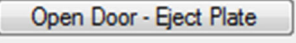
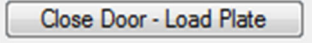
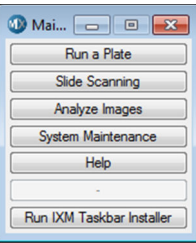
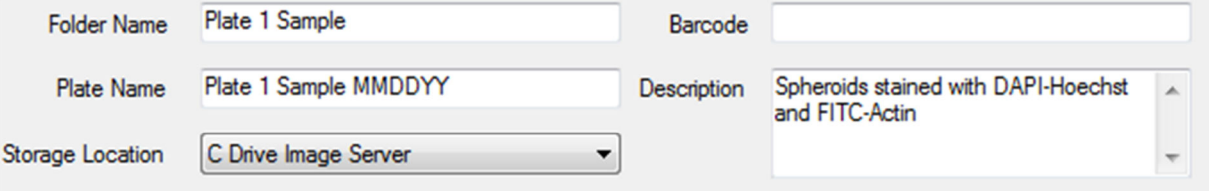


7.

To load a previous saved protocol, click on **Load Protocol...** in **Plate Acquisition Setup**

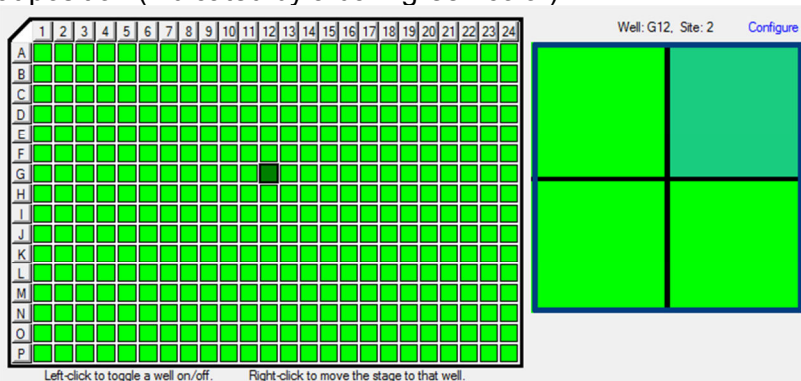
8.




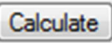
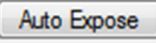
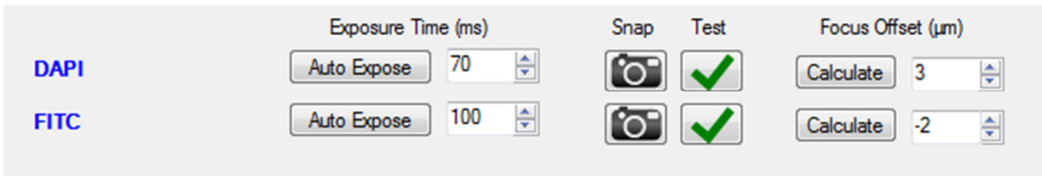
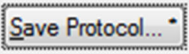

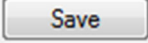

- Click **Load From File** to search windows for the appropriate .hts file.
- If the settings file is saved to the database, highlight the protocol and click **Load From DB**



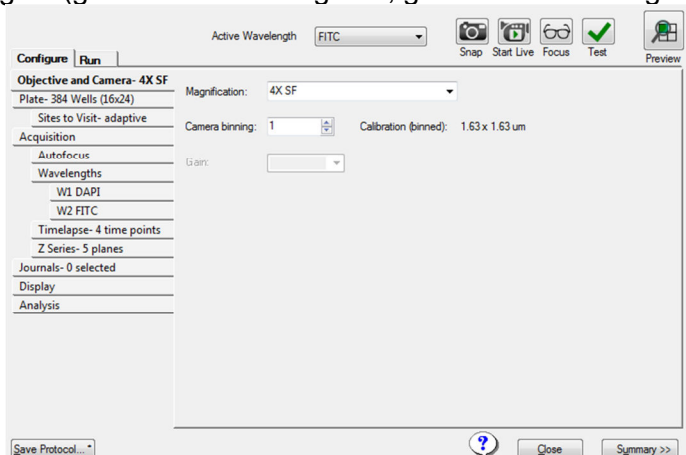
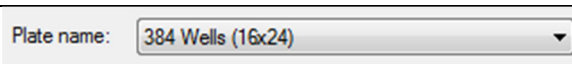
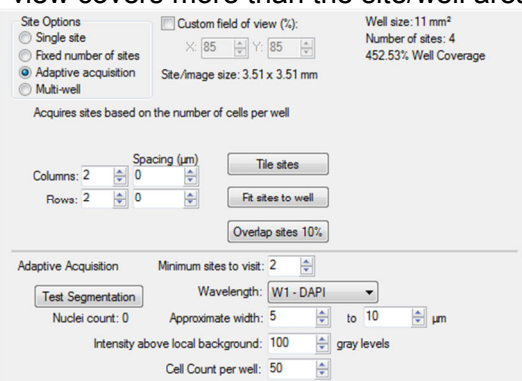
9. • Click  to open the door and place the plate in the in the system
- Click  to close the door
- Alternatively, you can use the **Main Taskbar** to open and close the door.
10. • Click 
- Click  or 
- 
- On the **Run** tab, update the folder name, plate name, and description as desired
11. 
12.  Click  to begin acquiring the plate

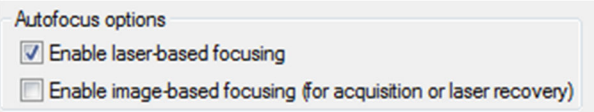
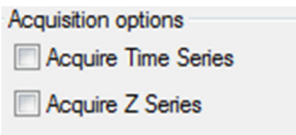
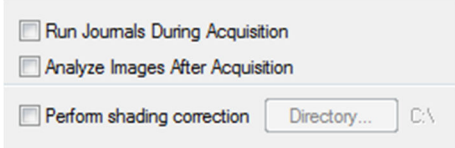
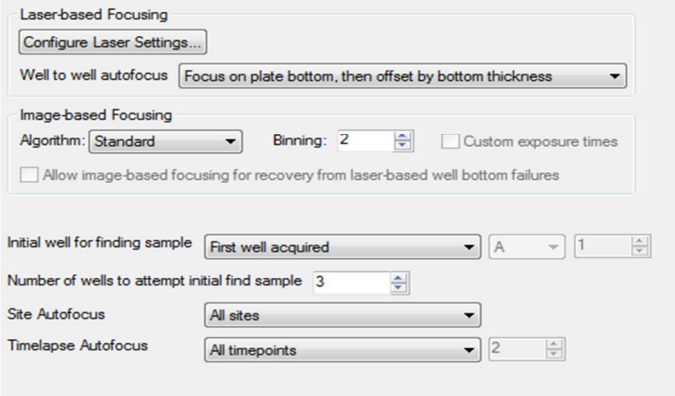
II. Test Acquisition Settings

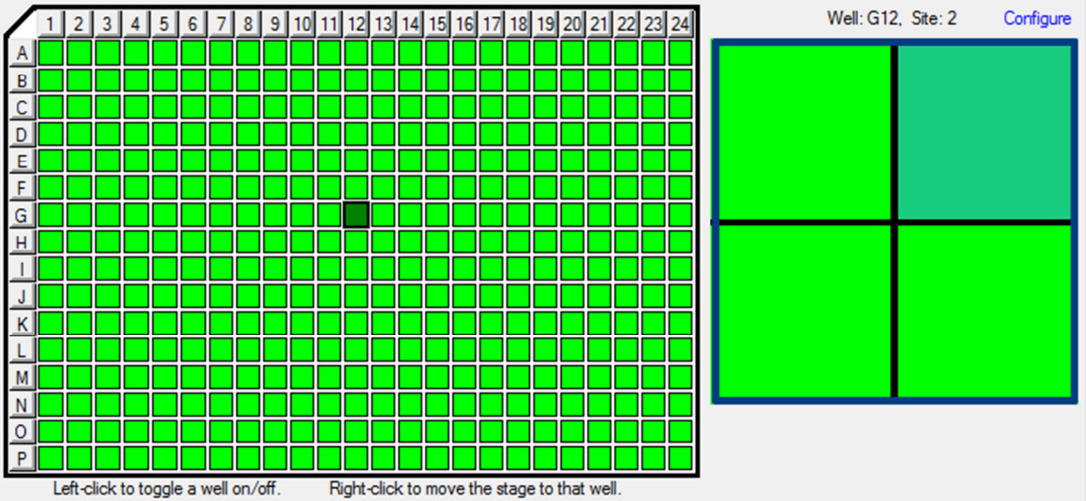
1. Open **Plate Acquisition Setup**
2. In the plate and site section of **Plate Acquisition Setup**, right-click on the desired well and/or site to move the plate to that position (indicated by a dark green color)
- 

3.	<p>Test the acquisition settings by clicking</p> <ul style="list-style-type: none">  Focus to perform a large range autofocus and snap image routine  Test to perform a focus and snap image routine (if Z series has been activated, all planes will be acquired)  Preview to perform an autofocus and snap image routine all for all wavelengths (if Z series has been activated, all planes will be acquired)
4.	<p>Adjust the acquisition settings, if necessary:</p> <ul style="list-style-type: none"> Adjust the focus offset by clicking  or adjusting the number manually Adjust the exposure time by clicking  or changing the number manually  <p><i>*NOTE*</i> Click on the wavelength name to open the corresponding wavelength tab for advanced options</p>
5.	<p>When you have optimized settings, click </p> <ul style="list-style-type: none"> Molecular Devices recommends enabling  Click  to search for a location on the hard drive.
6.	<p>Click  to begin acquiring the plate</p>

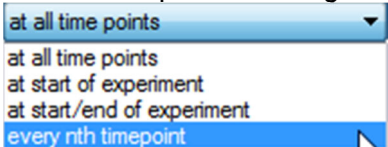
III. Define New Acquisition Settings

1.	Open Plate Acquisition Setup
2.	Select the Configure tab
3.	<p>Select the Objective and Camera tab</p> <ol style="list-style-type: none"> Select the appropriate magnification from the drop-down menu Set binning (2 for cell counting and cell scoring; 1 for fine sub-cellular detail) If necessary set gain (gain of 2 for binning of 1, gain of 1 for binning of 2 or higher) 
4.	Adjust the objective correction collar if necessary (setting on objective should match physical plate bottom thickness). On the Run a Plate Taskbar , click on Adjust Correction Collar to step through the process.
5.	<p>Select the Plate tab and select the appropriate plate type from the drop-down list</p> 
6.	<p>Select the Sites to Visit tab and select the appropriate number of sites</p> <ul style="list-style-type: none"> Single Site: image one site per well in the center Fixed number of sites: image the number of selected sites for every well. Adjust number and spacing of sites. Left-click on sites to select (green) and deselect (grey). Right-click on any site to move the plate to that site position (dark green) Adaptive acquisition: collect the minimum number of sites to image at least the cell count indicated by the user. The Adaptive Acquisition section will appear allowing the user to choose wavelength, size and threshold settings, and desired minimum count for cells Multi-well: collect multiple wells within one image which is then cropped to define single wells automatically Custom field of view (%): reduce the size of each image by the percentage entered. This is useful when the field of view covers more than the site/well area desired 
7.	Select the Acquisition tab to select Autofocus and Acquisition options

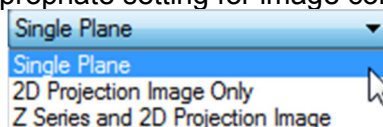
8.	<p>Autofocus options:</p> <ul style="list-style-type: none"> Always select Enable laser-based focusing Enable image-based focusing for thick samples or those with different focal planes from site-to-site or well-to-well 	
9.	<p>Acquisition options:</p> <ul style="list-style-type: none"> Enable Acquire Time series for timelapse experiments Enable Acquire Z series for Z step acquisition 	
10.	<p>Other options:</p> <ul style="list-style-type: none"> If running a journal during acquisition, enable this option to activate the Journals tab If an analysis has already been setup, enable Analyze Images After Acquisition <i>*NOTE* this requires an offline computer to be in Auto-run mode or running PowerCore software</i> To correct for uneven background, enable Perform shading correction and select the appropriate directory where shading correction images are saved 	
11.	<p>Select the Autofocus tab:</p> <ol style="list-style-type: none"> Set Well to well autofocus to Focus on well bottom. This is the default acquisition setup, however when imaging thin-bottom plates with low magnification objectives (4x and below) or microscope slides, select Focus on plate bottom, then offset by bottom thickness For Image-based Focusing refer to corresponding MetaXpress 6 Software Guide modules for suggested settings Set Initial well for finding sample to First well acquired Set Number of wells to attempt initial find sample to 3 If more than one site is acquired, set Site Autofocus to All sites If timelapse is enabled, set Timelapse Autofocus to All timepoints for long term timelapse, and First timepoint only for fast kinetic experiments 	

12. Select the **Wavelengths** tab and select the number of wavelengths (colors) including transmitted light that you would like to acquire
- Number of wavelengths:
13. In the plate and site section of **Plate Acquisition Setup**, right-click on the desired well (typically a control well) and/or site to move the plate to that position (indicated by a dark green color)
- 
14. Select the **W1** (wavelength) tab
- Select the desired filter set from the drop-down menu under **Illumination setting**
 - Click **Focus**
 - Examine the image
 - If the image appears to be dim or saturated, first adjust the image scaling, then adjust exposure time if necessary
 - If a blank or snowy image appears, this can indicate that a plate is not in the system or laser autofocus settings are incorrect
 - Click the **Calculate Offset** button to perform an automatic focus determination
 - For more control, enable ☐ **Use Z stack** and follow the prompts
 - If necessary, enable ☐ **Custom Range**

Range (um) Step (um)

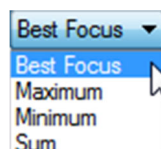
312.5 12.5
 - Click **Focus** again to test the new post-laser offset. Image should now be in focus.
 - Examine the image for brightness
 - If necessary, click **Auto Expose** with **Target max intensity: 45000** set to **33000 – 45000**
 - You can also increase or decrease exposure manually
15. If acquiring a Timelapse, select how often to acquire this image from the drop down menu
- 

If acquiring a Z Stack, select the appropriate setting for image collection



NOTE Z Series and 2D Projection Image is not available when acquiring a Timelapse

16. If saving the 2D Projection Image, select the appropriate projection method (press F1 for more information)

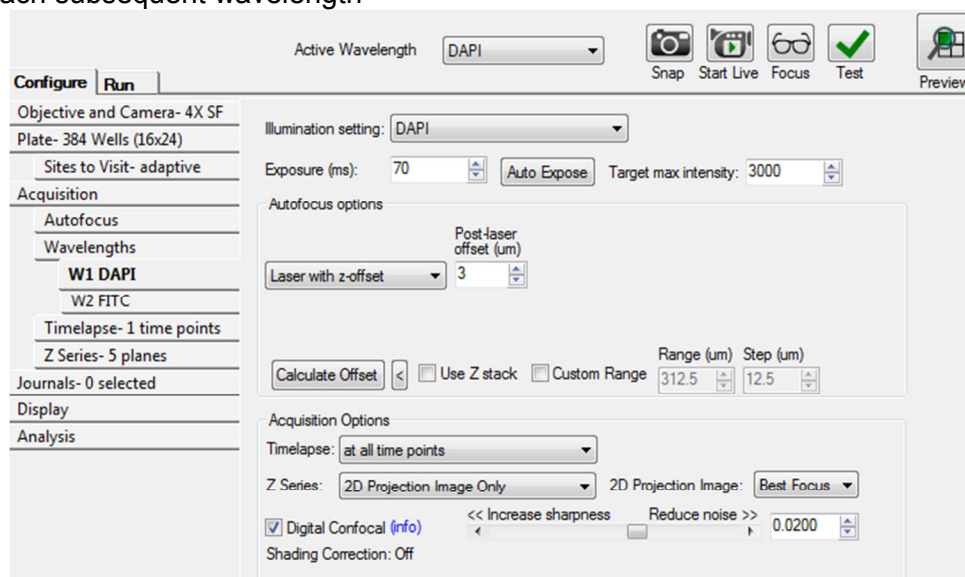


NOTE Best Focus is not recommended for comparison of intensity measurements

17. If the option is available, you can enable **Digital confocal** and select the appropriate K value using the slider bar (press F1 for more information)

Repeat for each subsequent wavelength

18.

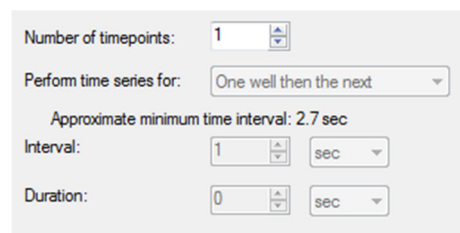


If acquiring with Timelapse, select the **Timelapse** tab



- i. Enter the number of **Time points** desired
- ii. Set **Interval** as the time between each time point
- iii. Set **Duration** as the total time of the experiment
- iv. Set **Perform time series** for:

19.



- **One well then the next:** entire timelapse is run for one well before acquiring next well
- **One column then the next:** entire timelapse is run for one column before acquiring next column
- **One row then the next:** entire timelapse is run for one row before acquiring next row
- **All selected wells:** all wells are imaged before continuing with next time point




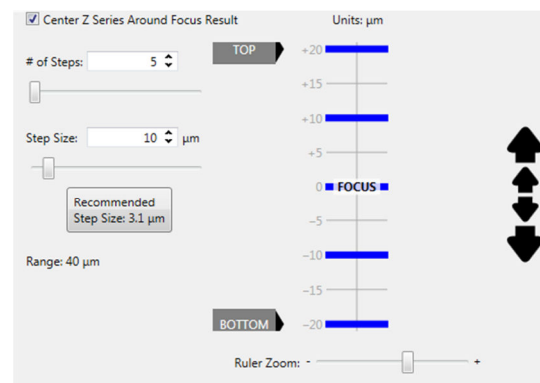
If acquiring a Z Series, select the **Z Series** tab

- i. Deselect ☐ Center Z Series Around Focus Result
- ii. Adjust **Step size** for spacing between each Z plane
- iii. Click  to determine the Z start position
- iv. Click  to start **Live Mode**
- v. Use the large and small arrows to move to the top of the focus range for the sample and click **Set T**
- vi. Use the large and small arrows to move to the bottom of the focus range for the sample and click **Set B**

20.

- vii. Click  to stop **Live Mode**
- Click  to perform focus and snap image routine to acquire all Z steps for the active wavelength. The last image in the stack will be the selected 2D projection image.

- Click  to perform focus and snap image routine to acquire all Z steps for all wavelengths. The last image in the stack will be the selected 2D projection image.



Current 2912.4 μm

Set T
Set B

Select the **Journal** tab (enabled on acquisition tab) and activate journals where appropriate. It might also be necessary to enable

21.


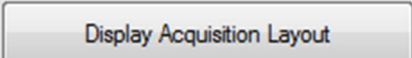
- ☐ Prevent asynchronous hardware moves (recommended if any journals are dependent on hardware positioning).

for certain journals (refer to documentation accompanying journals for details)

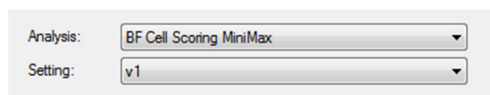
Acquisition Step	Journal
<input type="checkbox"/> Before each image	[None]
<input type="checkbox"/> After each image	[None]
<input type="checkbox"/> Before focusing	[None]
<input type="checkbox"/> Start of z	[None]
<input type="checkbox"/> End of z	[None]
<input type="checkbox"/> Start of site	[None]
<input type="checkbox"/> End of site	[None]
<input type="checkbox"/> Start of well	[None]
<input type="checkbox"/> End of well	[None]
<input type="checkbox"/> Start of time point	[None]
<input type="checkbox"/> End of time point	[None]
<input type="checkbox"/> Start of plate	[None]
<input type="checkbox"/> End of plate	[None]
<input type="checkbox"/> Prevent asynchronous hardware moves (recommended if any journals are dependent on hardware positioning).	

Select the **Display** tab to:

22.

-  will use default settings to arrange displayed images
-  manually adjust image display prior to acquisition
- Enable ☒ Display images during autofocus
- Enable ☒ Display images during acquisition
- (Optional) Enable ☒ Display a color overlay of wavelength images during acquisition

23. Select the **Analysis** tab (enabled on the acquisition tab) to specify the appropriate optimized **Analysis** routine and **Settings** from the drop down-menus



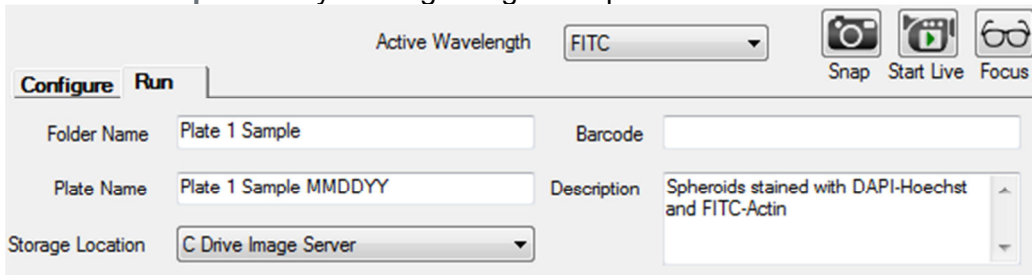
Analysis: BF Cell Scoring MiniMax
Setting: v1

**NOTE* This requires an offline computer set in Auto-run mode or running PowerCore software*

Under the **Run** tab, enter:

- **Folder Name:** Project name, your name, PI, etc. All your plates will go under this name.
- **Plate Name:** Name of this particular experiment
- **Storage location:** Select appropriate server for image storage.
**NOTE* There may only be one choice.*
- **Barcode:** Enter a barcode if desired
- **Description:** Any text regarding the experiment

24.



Active Wavelength: FITC

Snap Start Live Focus

Configure Run

Folder Name: Plate 1 Sample Barcode:


Plate Name: Plate 1 Sample MMDDYY Description: Spheroids stained with DAPI-Hoechst and FITC-Actin


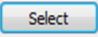
Storage Location: C Drive Image Server

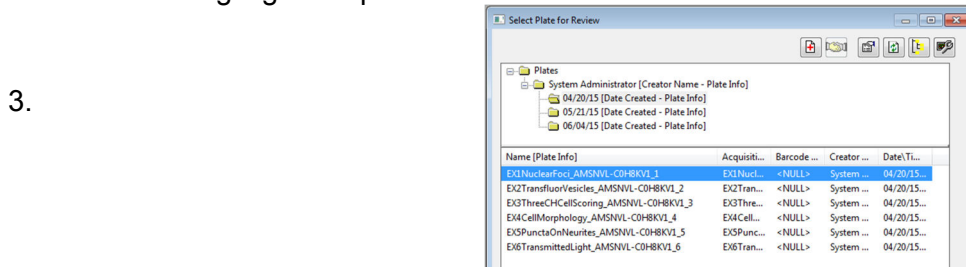
25. When you have optimized settings, click
- Molecular Devices recommends enabling ☐ Save to file rather than database
 - Click to search for a location on the hard drive.

26. Click on  to begin acquiring the plate

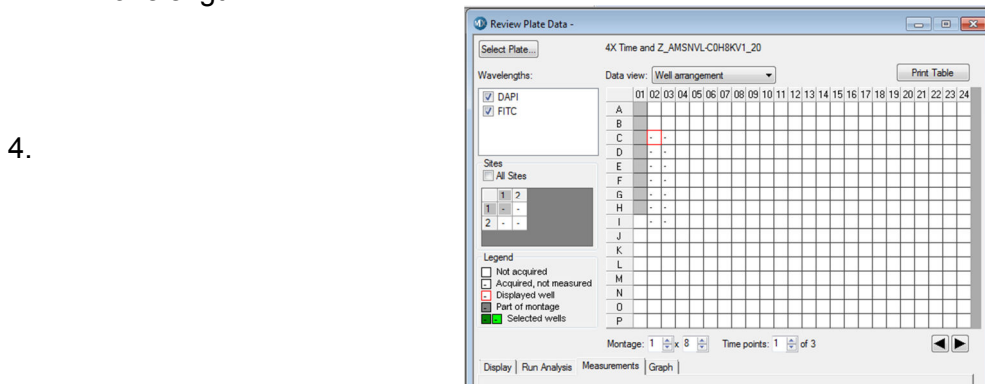
IV. Review Images and Run an Analysis


1. In the main toolbar, click  **Plate Acquisition Setup** or in the main menu select **Screening > Plate Acquisition Setup**

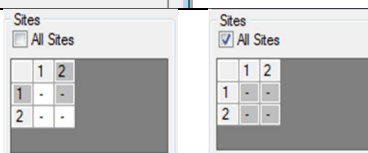
2. On the **Review Plate Data** dialog, click 
 - Navigate through the folders to find the plate of interest.
 - Highlight the plate and click .

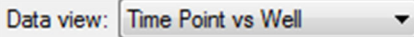
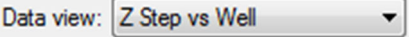


In the Plate View section, you will see a “-” in each well that was imaged. Left-click and drag across the wells you want to view. A thumbnail montage of these wells will open for each wavelength

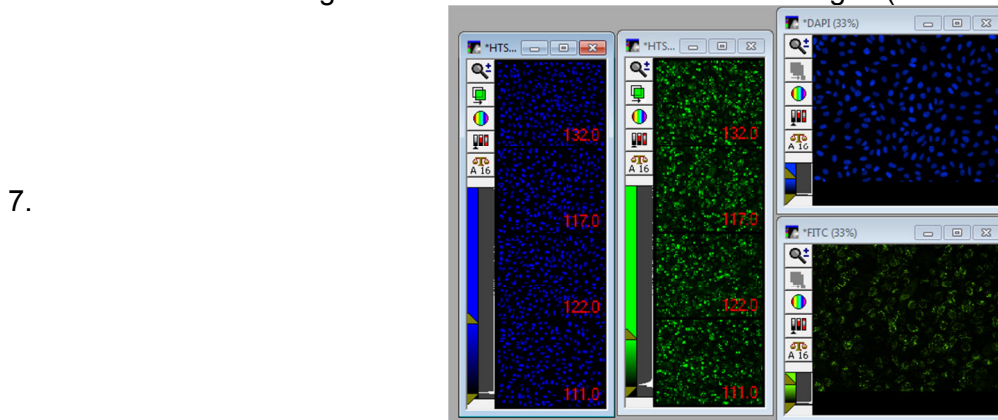


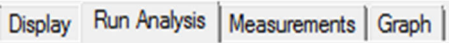
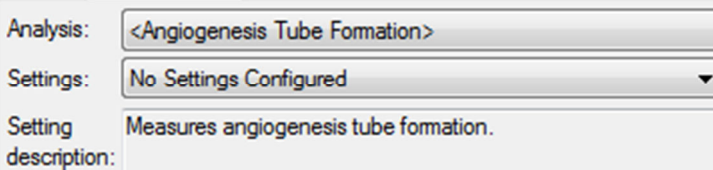


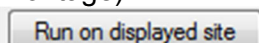
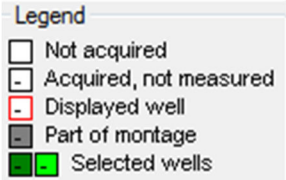
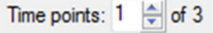
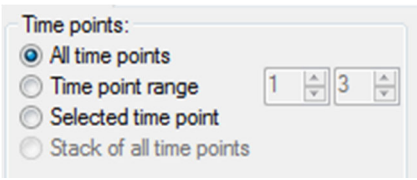

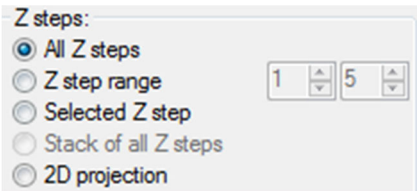
5. If there are multiple sites per well, select an appropriate site to view, or enable  **All Sites**. The image montages will automatically adjust.



6. To view all Timelapse or Z Series images at once, change the **Data view** to  or , respectively.

Left-click on a single thumbnail to view full resolution images (all wavelengths)

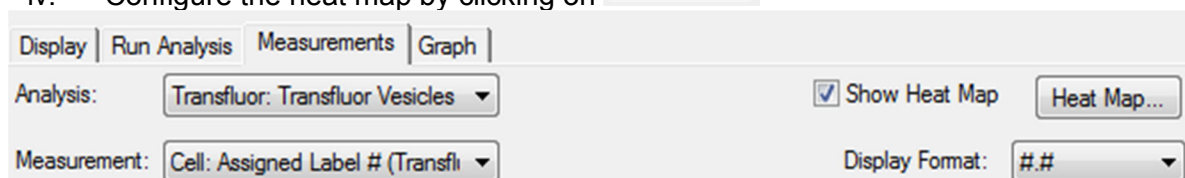


8.	To run or set up an analysis, select the the Run Analysis tab	
9.	If analysis settings have already been optimized, select the analysis routine (application module, custom module, or journal) and settings from the drop-down menus	
10.	<p>Under the Run Analysis tab, select the appropriate button to run the analysis:</p> <ul style="list-style-type: none">  analysis we be run on all acquired images  analysis will be run on selected wells (selected wells are indicated in green; to select wells, right click well(s) in the plate section or image montage)  analysis will be run only on the currently displayed site 	
11.	<p>For a Timelapse data set, select the appropriate option for analysis under the Time points section</p> <ul style="list-style-type: none"> All time points: run analysis on all time points in the data set Time point range: run analysis on a consecutive range of time points Selected time point: run analysis on only one time point that is select in the Time point section below the plate layout  Stack of all time points: use if, in the Analysis field, you select a legacy timelapse journal which analyzes the planes in a stack 	
12.	<p>For a Z Series data set where all Z planes were saved, select the appropriate option in the Z steps section</p> <ul style="list-style-type: none"> All Z Steps: run analysis on all Z planes Z Step range: run analysis on consecutive range of Z planes Selected Z step: run analysis on only one Z plane that is selected in the Z step section below the plate layout  Stack of all Z steps: run an analysis with a journal that requires a stack of images 2D projection: only run analysis on the saved 2D projection image 	
13.	If the selected analysis has already been run on the plate, a warning will appear asking to overwrite the data. If you are not sure, save the analysis settings with a new name before analyzing your plate.	

To view analysis results, select the **Measurements** tab

- i. Select the **Analysis** (module and settings name) from the drop down menu
- ii. Select a measurement from the drop-down menu. The values will be shown in the plate layout.
 - a. Measurements starting with a "Cell" are cell-by-cell data and will give the average of all cells in the well

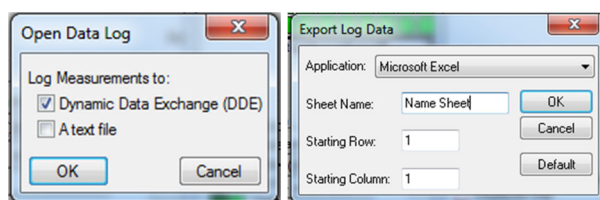
14. iii. Activate the heat map by enabling ☒ Show Heat Map
- iv. Configure the heat map by clicking on **Heat Map...**



15. To view the cell-by-cell data, click **Cellular Results...** at the bottom of the **Review Plate Data** dialog. Data will be automatically updated based on the well and site selected in the montage view

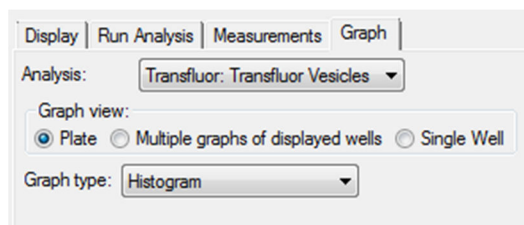
To export data to Excel:

16. i. On the **Measurements** tab, click on **Open Log**
- ii. Select only **Dynamic Data Exchange**
- iii. Select **Microsoft Excel** and name worksheet as desired. This opens an empty worksheet.
- iv. Click **Log Data**. Currently viewed data will be logged into the Excel sheet.



To create simple graphs in MetaXpress:

17. i. Go to **Graph** tab
- ii. From Graph Type, select:
- iii. Select measurements to plot from the drop-down menu
- iv. Click **Show Graph**
- v. Right-click on the graph for more options



Shutting down the system after use

1. Turning off the system:

- b. Save settings as needed and exit MetaXpress software ("X" in upper right corner or File>Exit). Be sure to exit the software before turning off the hardware components.
 - Turn off the IXM Power Supply
 - Turn off the Options Controller
 - Power down the lamp:
 - Press Local key on the Lambda XL light source keypad. This should give you options then to stop the lamp. If you do not see, then select Local again.
 - Press 2 to stop the lamp (WARNING: DO NOT STOP the lamp within 1 minute of starting the lamp to avoid damaging the lamp and reduce life time of the lamp)
 - Press 1 to confirm stop the lamp.
 - The lamp will turn off, wait until the temperature of the lamp drop below **40° C** before turning the lamp power off.
 - Turn off Computer/monitor (optional)