



MetaXpress[®] 6 Software Guide

Setting Up Batch Analysis

Date Revised 06/24/15 Version B



Chapter Purpose

The purpose of this chapter is to guide the user through setting up analysis of multiple plates (batch analysis) through the **Plate Data Utilities** dialog.

Batch analysis can be:

- Same analysis over multiple plates
- Different analyses over the same plate(s)
- Different analysis for different plate(s)



A Note About Batch Analysis

Batch analysis can be run on the acquisition computer, offline analysis workstation(s), or through MetaXpress PowerCore (requires licenses, contact your local sales account manager for more information).

NOTE Molecular Devices may not recommend to run analysis on the acquisition computer depending on the computer specifications. Some computers may not have the recommended specs for analysis.

During the set up process, you will be given the option to **Run on this computer** or **Add to auto run list** (described later). If you select the latter option and you have MetaXpress PowerCore software, PowerCore will automatically detect the plates in the auto run list and begin analysis. Otherwise, the current computer or another offline workstation(s) will need to be in Auto Run Mode (described later) to run the analysis.

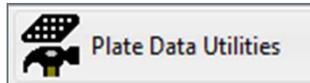
It is suggested to increase Auto run timeout in MDCStore Tools to at least 300 seconds (available on the **MX Settings** tab in MDCStore Tools).



Setting Up a Batch Analysis

1. Open **Plate Data Utilities**

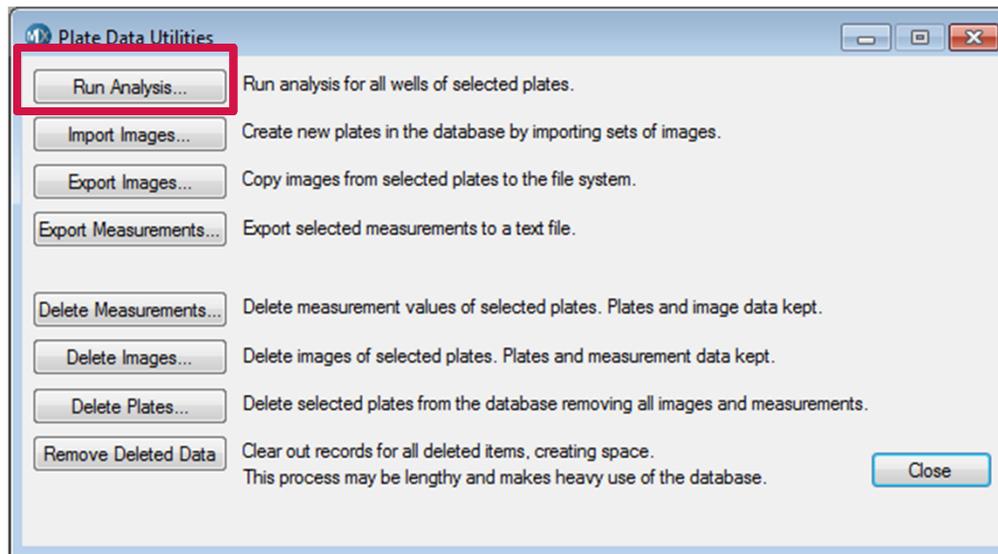
- In the main toolbar click on



OR

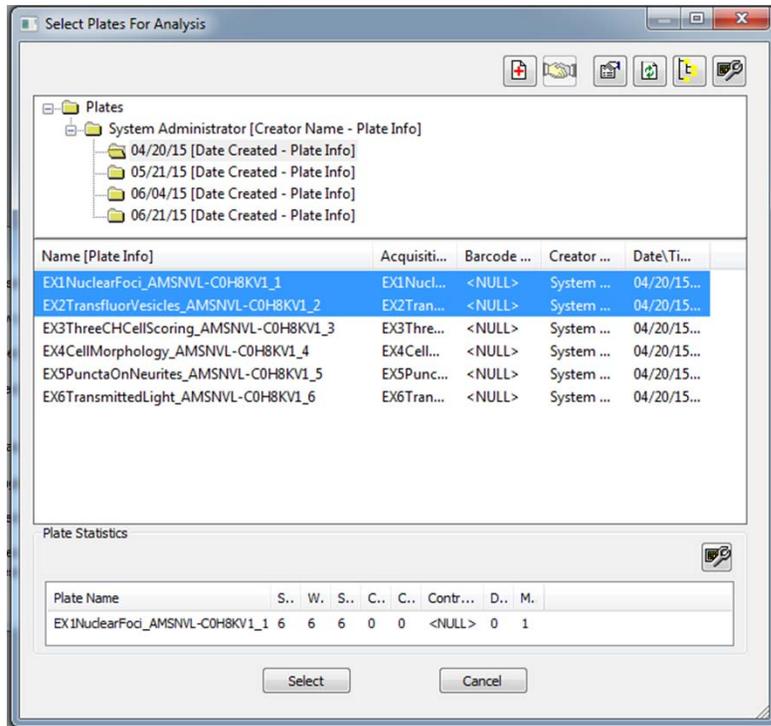
- Under the **Screening** menu, select **Plate Data Utilities**

2. Click on the **Run Analysis** button



Setting Up a Batch Analysis

3. Select the plate(s) of interest to be analyzed
 - Use the Shift or Ctrl keys on the keyboard for multiple selections
 - Click on the **Select** button

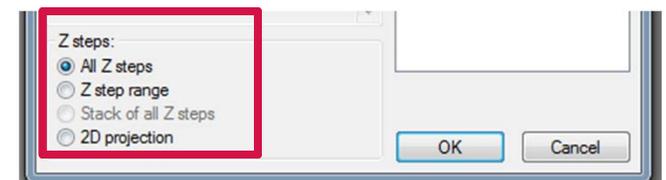
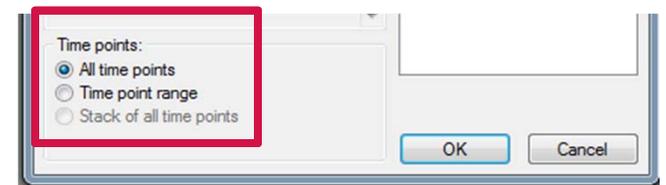
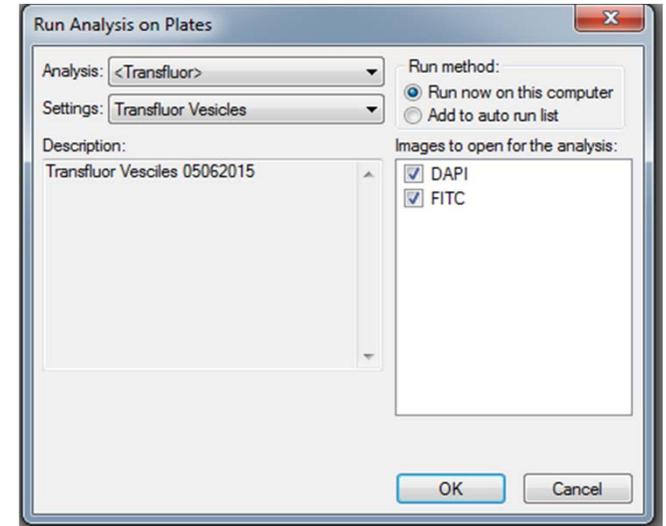


NOTE When selecting multiple plates, ensure that the wavelengths in the plate match those in the analysis. Otherwise, you will get an error at the end of the process.



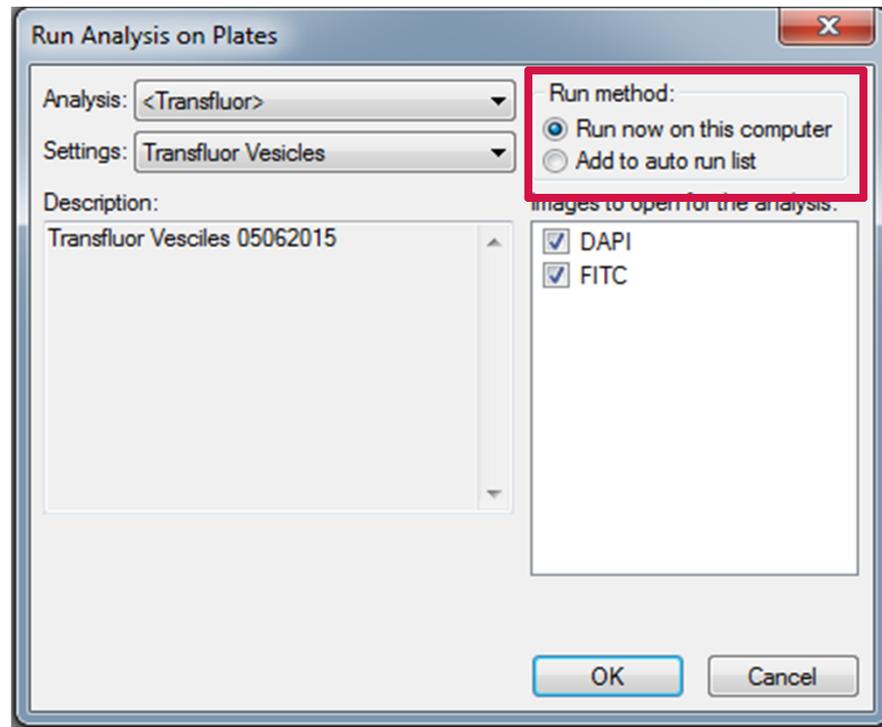
Setting Up a Batch Analysis

4. Select the desired **Analysis** and **Settings** from the drop-down menus
5. Select the **Run Method**:
 - **Run now on this computer**: analysis will begin immediately on the current computer
 - **Add to auto run list**: the analysis will be added to an auto run queue (Refer to section on Auto Run List). Use this option if adding additional analyses, using PowerCore or delaying analysis to a later time.
6. Enable the appropriate wavelengths for analysis
7. If the plate contains Time lapse or Z Series data, select which images to analyze
- *NOTE* Batch analysis will analyze all sites/wells in the plate.
8. Click on the **OK** button



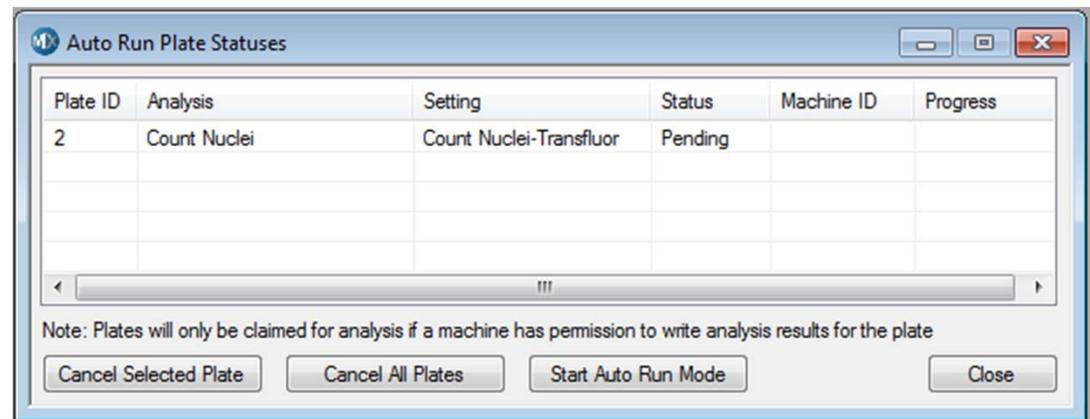
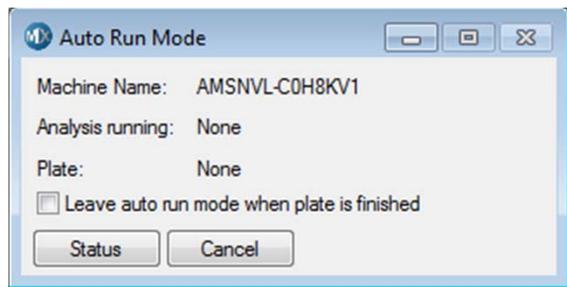
Setting Up a Batch Analysis for Additional Plates

9. To add additional analyses (for the same plate(s) or additional plate(s))
- Repeat steps 2-8
 - Always select **Add to auto run list** under **Run method**



Auto Run List

- If you have purchased MetaXpress PowerCore, PowerCore will automatically detect the plates in the Auto Run List and begin analysis (make sure that PowerCore is signed into the correct database)
- Otherwise, you will need to activate **Auto Run Mode**. This can be done on the acquisition computer or an offline workstation
 - In the main menu, select **Screening > Start Auto Run Mode**
 - Deselect **Leave auto run mode when plate is finished** to allow analysis of all plates in the queue
 - *NOTE* While the computer is in **Auto Run Mode**, other functions are not possible
 - Click on the **Status** button to display the status of the current plate, see a list of plates in the auto run list, and cancel analysis of plates
 - *NOTE* If you would like to see the list of plates and analyses before starting Auto Run Mode, go to **Screening > Auto Run Plate Statuses**. From this dialog you can also start Auto Run Mode.



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