



MetaXpress[®] 6 Software Guide

Opening and Editing a Custom Module

Date Revised 07/13/15 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 opening and editing a custom module.

The user may want to:

- Test other sites/wells in the plate using the current custom module
- Edit the custom module to work on plates with different wavelengths

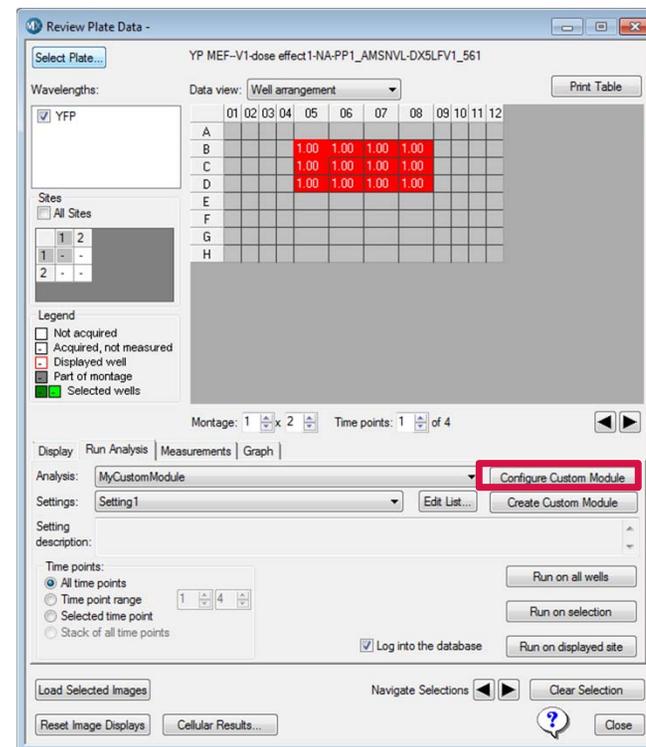
Refer to corresponding chapters on importing, exporting, and saving custom modules.



Testing Another Site / Well

Molecular Devices recommends testing any analysis with both positive and negative control wells. To do this in **Custom Module Editor**:

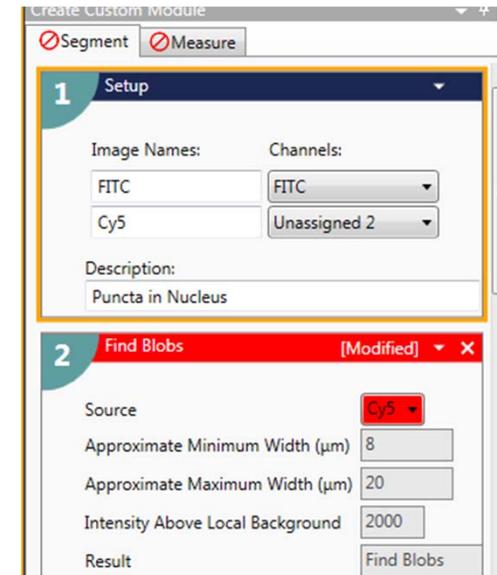
1. Save the custom module
2. Minimize the **Custom Module Editor** interface
3. Click on the desired site/well in the thumbnail montage to display the high resolution images
4. Click on the **Configure Custom Module** button on the **Run Analysis** tab
5. Maximize the **Custom Module Editor** interface. The images are now updated.



Editing a Custom Module for Different Wavelengths

The user may want to run a custom module that was originally compiled using wavelengths that are different than in the current plate. To change the wavelengths:

1. In the **Review Plate Data** dialog, open the new plate of interest and select a site/well to test settings
2. Select the **Run Analysis** tab and the desired custom module from the **Analysis** and **Settings** drop-down menus
3. Click on the **Configure Custom Module** button
4. In **Custom Module Editor**, the **Segment** and **Measure** tabs will have red warning icons indicating that the current wavelengths do not match what is in the custom module. In addition, any cards that use wavelengths not in the current plate will also be highlighted in red.
5. In the first card (**Setup**), select the new corresponding wavelength from the **Channel** drop-down menu. Type in the correct **Image Name** as selected in the drop-down menu. This will automatically change the **Source** images in each step card
6. Step through the custom module to ensure all the cards correspond to the correct source image(s)
7. Save the custom module as a new name



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