



MetaXpress[®] 6 Software Guide

Importing and Exporting 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.

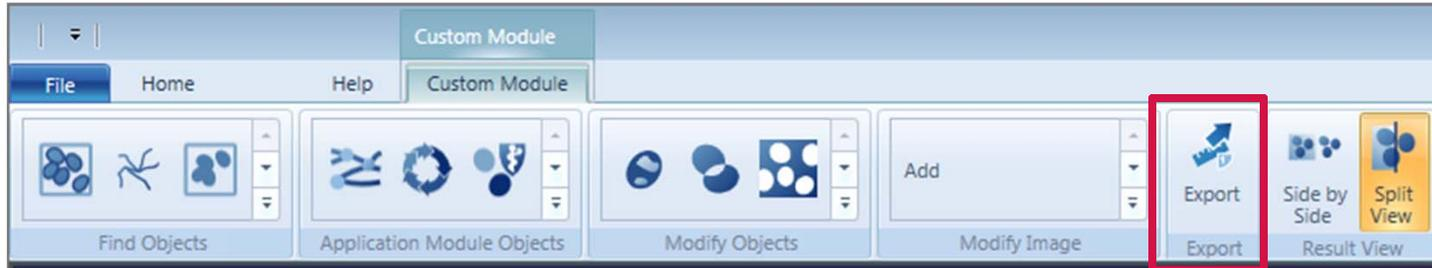
Chapter Purpose

The purpose of this chapter is to guide the user in importing and exporting a custom module.

Importing and exporting custom modules allow for easy collaboration or for troubleshooting with technical support.



Exporting a Custom Module in CME

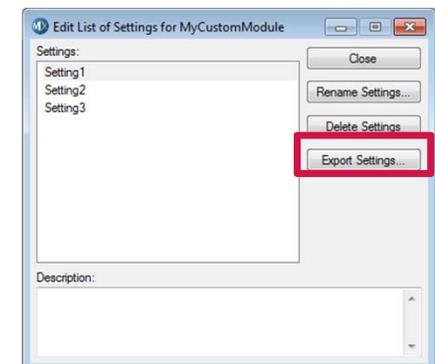
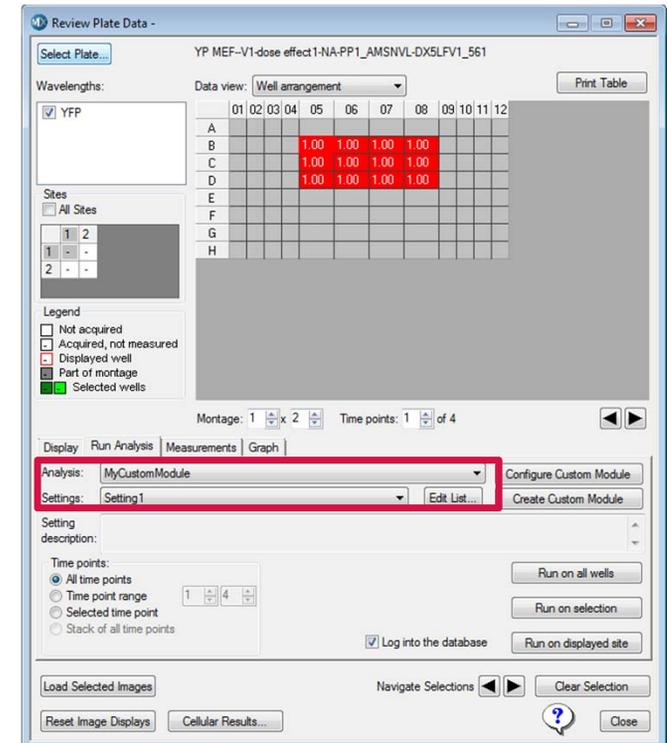


1. Click on the **Custom Module** tab in **Custom Module Editor**
2. Click on the **Export** button
3. Navigate through Windows for a save location
4. Name the file. This will appear as the **Measurement Name** in Review Plate Data when the custom module is imported
5. Click on the **Save** button
6. Custom Modules have a file extension **.xml**



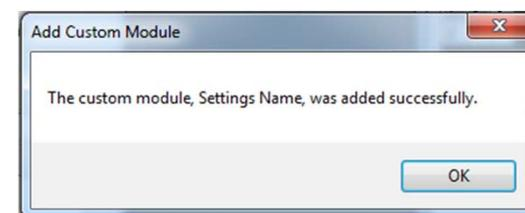
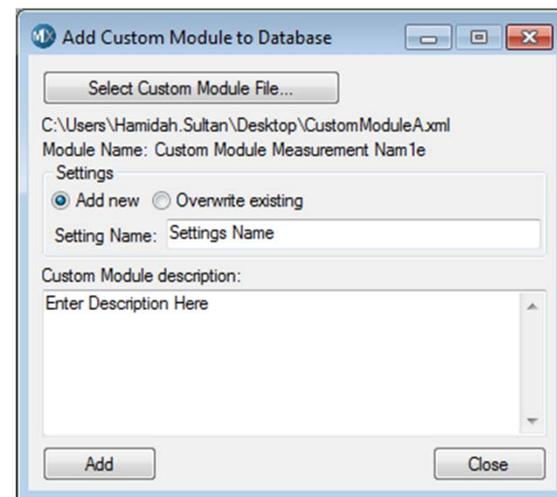
Exporting a Custom Module in Review Plate Data

1. In the **Review Plate Data** dialog, select the desired custom module from the **Analysis** and **Settings** drop-down menus
2. Click on the **Edit List** button
3. Highlight the appropriate settings in the list and click on the **Export Settings** button
4. Navigate Windows to the save location
5. Name the file. This will appear as the **Measurement Name** in Review Plate Data when the custom module is imported
6. Click on the **Save** button
7. Custom Modules have a file extension .xml



Importing a Custom Module

1. In the main menu, select **Screening > Add Custom Module to Database**
2. Click on the **Select Custom Module File** button
3. Navigates Windows to locate the .xml file
4. Select **Add New** and enter the **Settings Name** to add a new custom module to the database
5. Select **Overwrite existing** and a drop-down menu of exiting settings in the database will appear. Select the appropriate settings.
6. (optional) Enter and/or modify the custom module description
7. Click on the **Add** button to add the custom module to the database
8. A message will appear confirming that the custom module was successfully added to the database. Click **OK**



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