



# MetaXpress® 6.5 Software Guide

Custom Module Analysis in a journal

Date Revised 04/11/2019 Version B



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

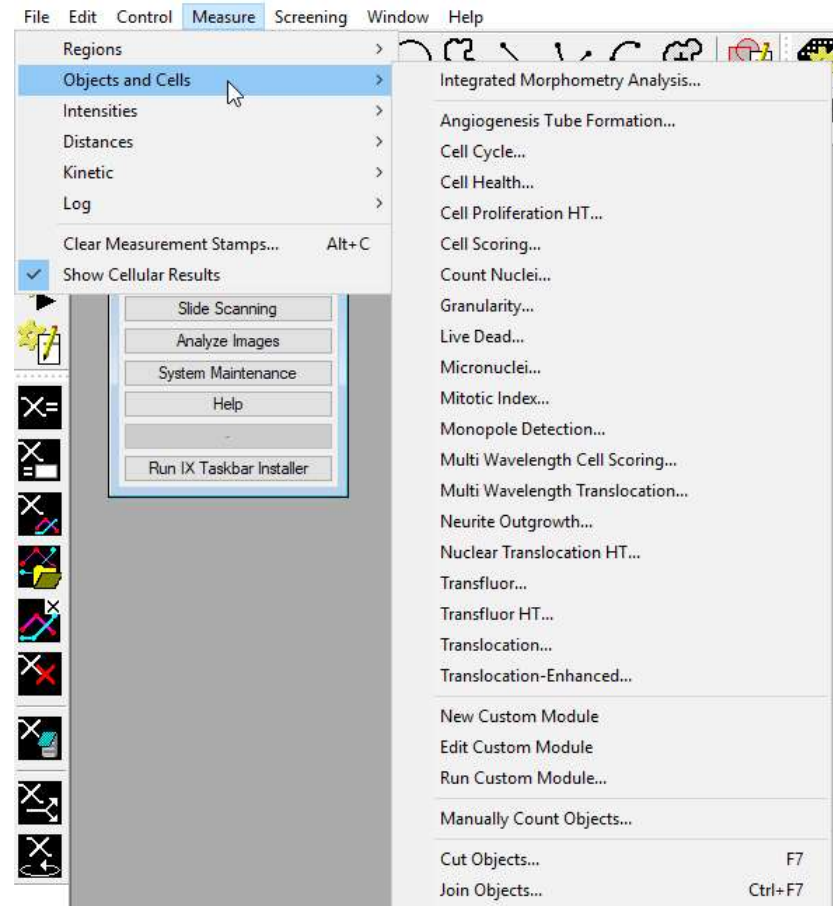
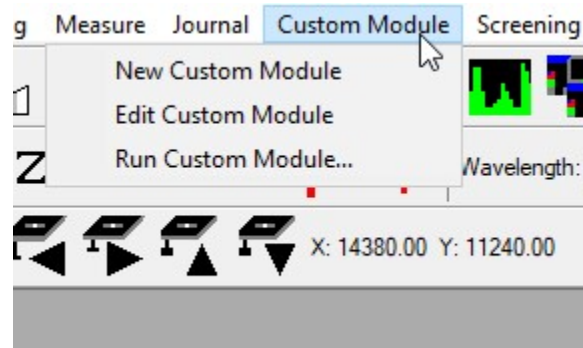
# Chapter Purpose

The purpose of this chapter is to guide the user through using the new Custom Module menu items added in MetaXpress 6.5.

With these new functions, custom module analysis can be designed and run on images outside of the database. Custom modules can also be incorporated within journals for further analysis customization.



# New Custom Module menu



# Enable Custom Module menu

- To access the new menu items, enable the “**custommodule**” drop-in via Meta Imaging Series Administrator
- Repeat this step for all Groups (profiles) that need to run this function.
- See [https://mdc.custhelp.com/app/answers/detail/a\\_id/19850/](https://mdc.custhelp.com/app/answers/detail/a_id/19850/) for more details on enabling drop-ins.

The screenshot shows two windows from the Meta Imaging Series Administrator. The left window, titled "Meta Imaging Series Administrator: Single User Configuration", displays a "List of Groups" table with columns for Group Name, Hardware Setting, and File Association. The "MX offline" group is selected, and the "Drop-ins/Toolbars" button is highlighted. The right window, titled "Configure Drop-ins/Toolbars for MX offline", shows the configuration for this group. In the "Drop-ins" list, the "custommodule" drop-in is checked and highlighted. In the "Menus" list, the "Custom Module" folder is expanded, showing three sub-items: "New Custom Module <custommodule drop-in>", "Edit Custom Module <custommodule drop-in>", and "Run Custom Module <custommodule drop-in>". The "CUSTOMMODULE Description" field contains "Custom Module for MetaMorph", and the "CUSTOMMODULE adds these Menu Items:" field lists "Custom Module: New Custom Module, Edit Custom Module, Run Custom Module".

Group Name	Hardware Setting	File Association
MetaFluor	Default	
MetaFluor Offline	Offline	
MetaMorph	Default	
MetaMorph Offline	Offline	
MetaVue	Default	
MetaXpress	Default	
MX offline	Offline	



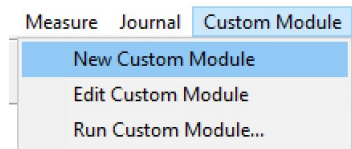
# Custom Module Editor Menu in MetaXpress

- Allows you to
  - set up a CME analysis on any individual images in MetaXpress software.
    - Open the images(s) via File > Open
    - Open image(s) via Review Plate data
  - Use CME analysis in journals
- Not yet compatible with stacks



# Create a new custom module

- Open the image(s).
- Make sure any images not being used for the analysis are closed.
- Go to **Custom Module > New Custom Module**.

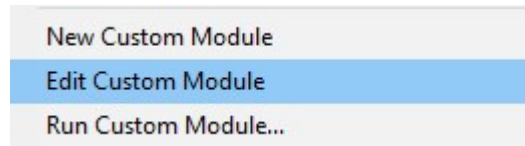


- The Custom Module Editor window opens.
- Develop an analysis routine and select measurement outputs.
- Click **Save** to save as an .xml file.



# Edit an existing custom module

- Open the image(s).
- Make sure any images not being used for the analysis are closed.
- Go to **Custom Module > Edit Custom Module**.

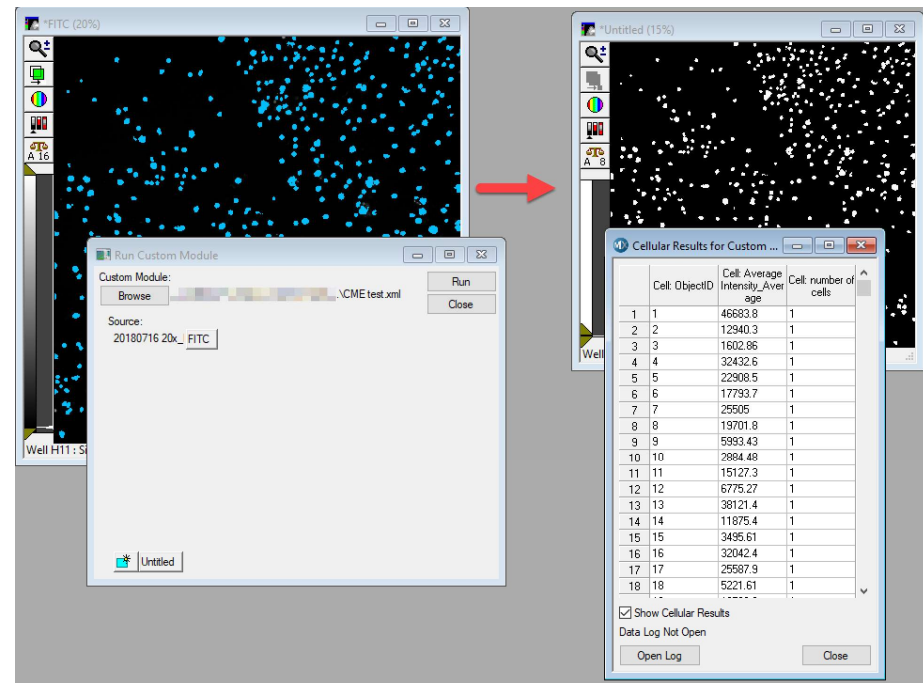


- When prompted, navigate and select the .xml file.
- The Custom Module Editor window opens.
- Modify the settings as needed.
- Click **Save** to overwrite or save as a new .xml file.



# Run Custom Module

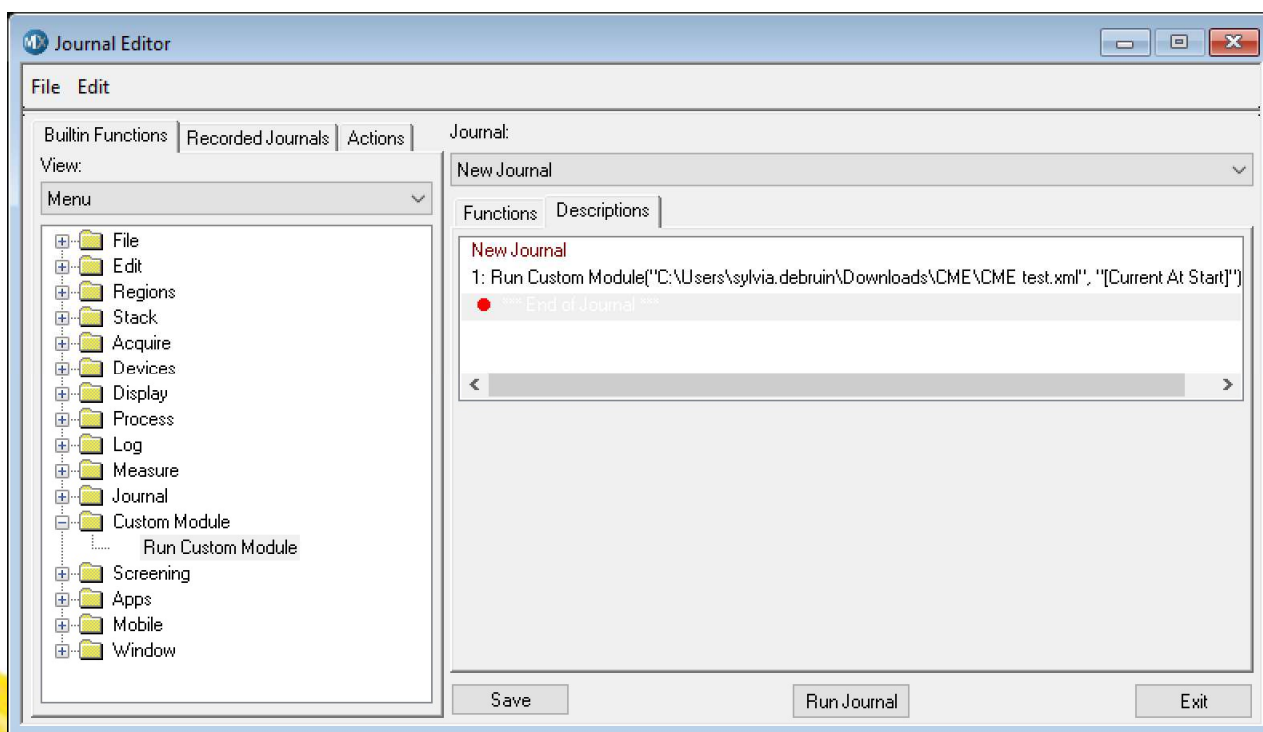
- Open the image(s).
- Go to **Custom Module > Run Custom Module**.
- Browse to the saved .xml file.
- Select the source image(s).
- Adjust the image result name.
- Click **Run**.
- To save the data:
  - Click **Open Log**.
  - Select DDE (Excel) or text file.
  - Click **Log Data**.





# Add CME to journal

- Create the CME analysis via the steps in the previous slides
- Open the journal editor
- Add the **Run Custom Module** step to the journal.
- To run over a plate via Review Plate Data, follow the steps in
- [https://mdc.custhelp.com/app/answers/detail/a\\_id/19370](https://mdc.custhelp.com/app/answers/detail/a_id/19370)



# 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





# MOLECULAR DEVICES

ADVANCING PROTEIN AND CELL BIOLOGY