



# MetaXpress<sup>®</sup> 6 Software Guide

## Saving 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 saving a custom module so that it can be ran on plate(s) through **Review Plate Data** or **Plate Data Utilities**.

Refer to corresponding chapters on running a custom module.

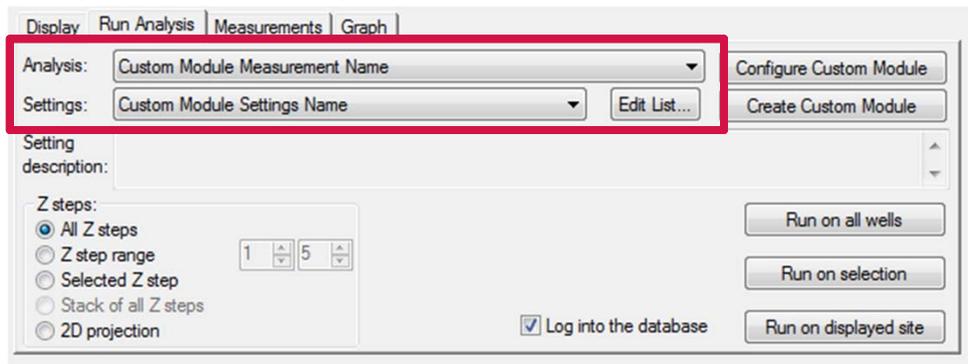


# Custom Module Naming Convention

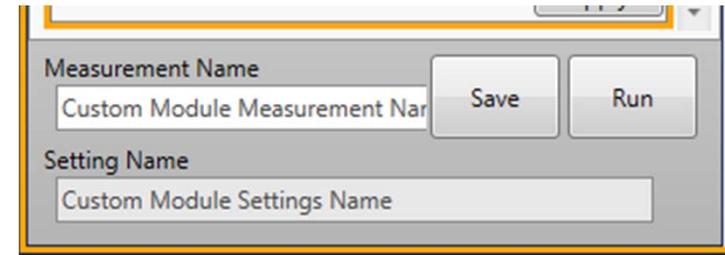
Custom modules have a two-part name

- **Measurement Name:** appears under the **Analysis** drop-down menu
- **Settings Name:** appears under the **Settings** drop-down menu

Similar custom modules can be saved with different settings names so they are grouped together in **Review Plate Data**.



The screenshot shows a software interface with a menu bar at the top containing 'Display', 'Run Analysis', 'Measurements', and 'Graph'. Below the menu bar, there are two dropdown menus: 'Analysis:' and 'Settings:'. The 'Analysis:' dropdown is set to 'Custom Module Measurement Name' and the 'Settings:' dropdown is set to 'Custom Module Settings Name'. A red box highlights these two dropdown menus. To the right of the dropdowns are buttons for 'Configure Custom Module' and 'Create Custom Module'. Below the dropdowns is a section for 'Setting description:' with a scrollable area. Underneath, there are radio buttons for 'Z steps' options: 'All Z steps' (selected), 'Z step range' (with input fields for 1 and 5), 'Selected Z step', 'Stack of all Z steps', and '2D projection'. There is also a checkbox for 'Log into the database' and buttons for 'Run on all wells', 'Run on selection', and 'Run on displayed site'.



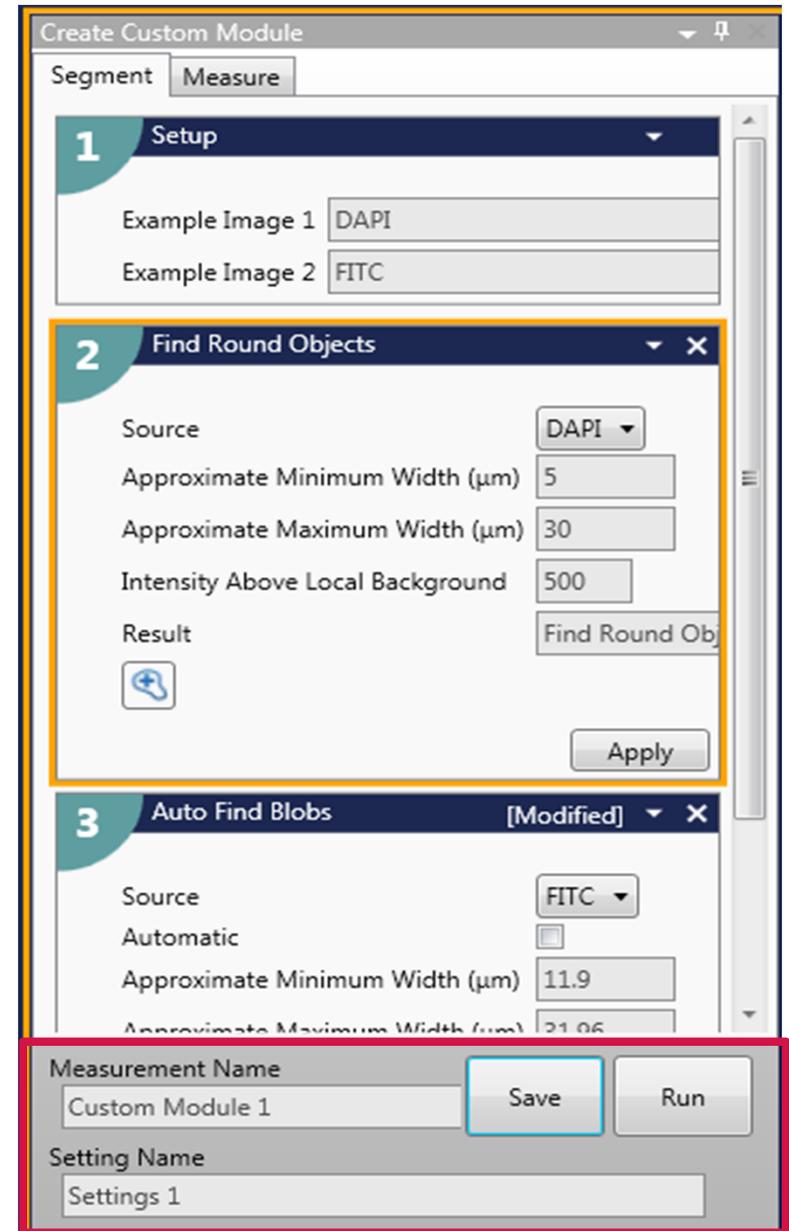
The screenshot shows a software interface with two input fields. The first is labeled 'Measurement Name' and contains the text 'Custom Module Measurement Name'. To its right are 'Save' and 'Run' buttons. The second is labeled 'Setting Name' and contains the text 'Custom Module Settings Name'.



# Saving a Custom Module

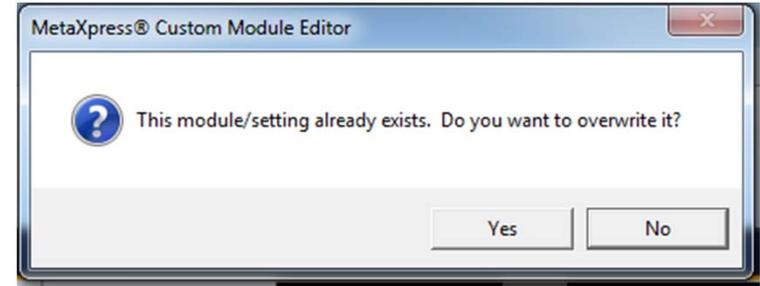
Molecular Devices recommends saving custom modules periodically in order to prevent loss of data in case of a computer crash.

- Save a custom module:
  - Enter a name in **Measurement Name** and **Settings Name**
  - Click the **Save** button
- Overwrite current custom module:
  - Click the **Save** button without changing the **Measurement Name** and **Setting Name**
- Save current custom module as a new custom module:
  - Change the Measurement Name and/or Setting Name
  - Click the **Save** button



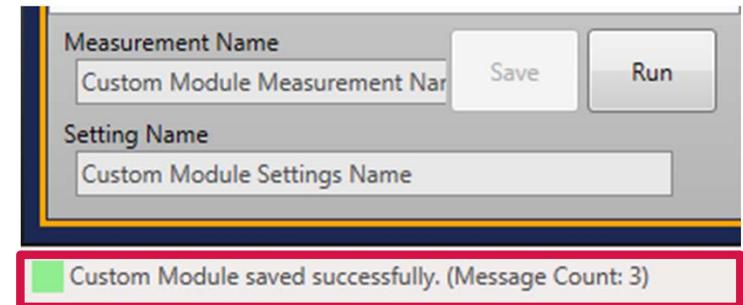
# Confirmation of Saved Custom Module

If you are overwriting settings, a dialog box will appear asking if you want to replace settings, click **Yes** to save, click **No** to cancel



Before closing the **Custom Module Editor**

- Verify that the custom module has been saved by examining the message bar at the bottom of the **Custom Module Editor** dialog box.
- If the message reads "**Custom Module saved successful**", you may now exit the **Custom Module Editor** interface.



\*NOTE\* Molecular Devices always recommends exporting the module in case the saving procedure was not completed successfully. Refer to the corresponding chapters on exporting custom modules.



# 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

For research use only. Not for use in diagnostic procedures.