



MetaXpress® 6 Software Guide

Selecting Masks for Measurement

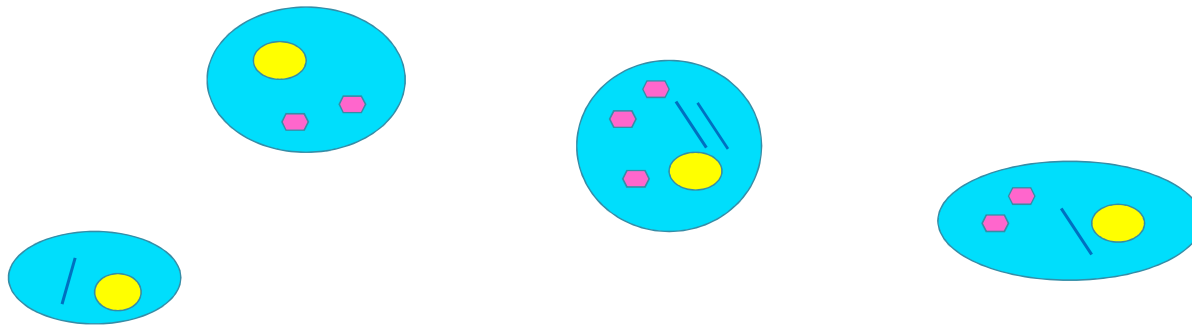
Date Revised 05/26/2020 Version B

Chapter Purpose

The purpose of this chapter is to describe the **Measure** tab in the **Custom Module Editor** (CME) plugin.

The **Measure** tab is used to select and measure the masks created in the **Segmentation** tab. Appropriate selection of masks is important for generating relevant measurements.

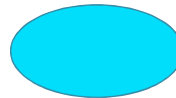
Masks for Measurement



Objects to Measure:

This is your primary mask.

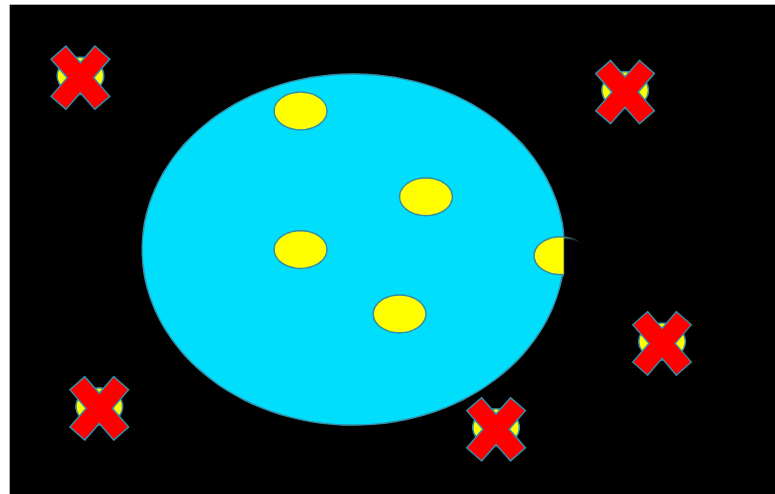
One row of data is generated for each of these objects.



Feature Group masks:

These masks are measured relative to the Objects to Measure mask.

Feature Groups measured inside main Objects



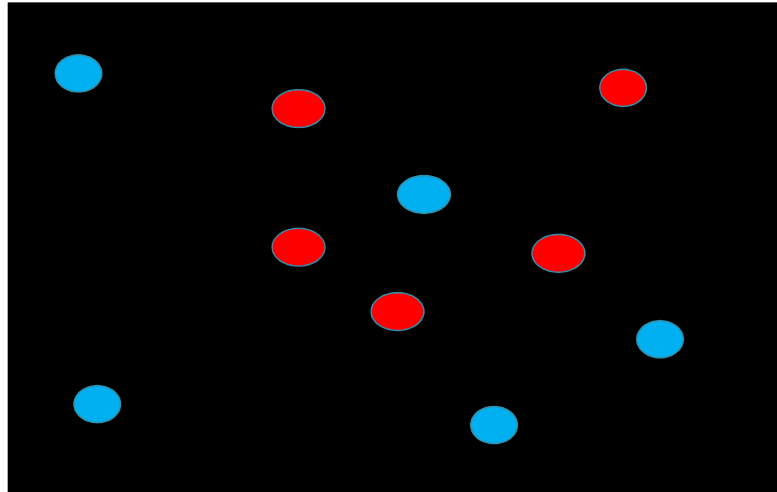
Objects to Measure

Feature Group

Feature Group objects outside of **main objects** are **not measured**.

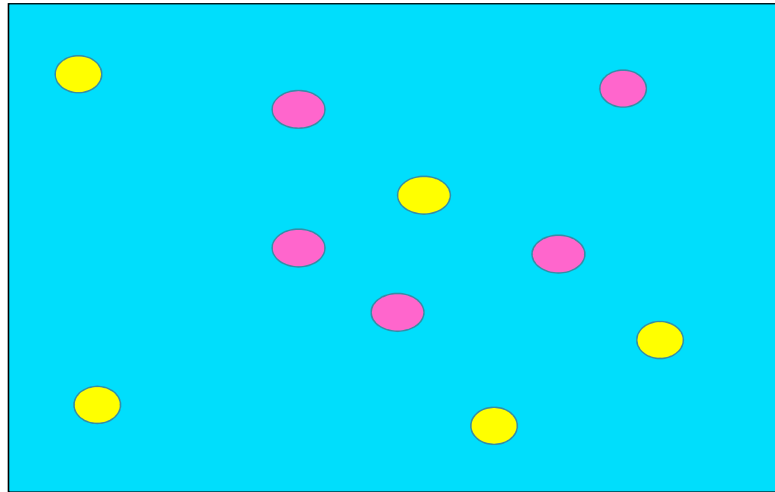
Feature Group objects partially inside **main objects** are **cropped**.

Example 1: Measure 2 distinct populations



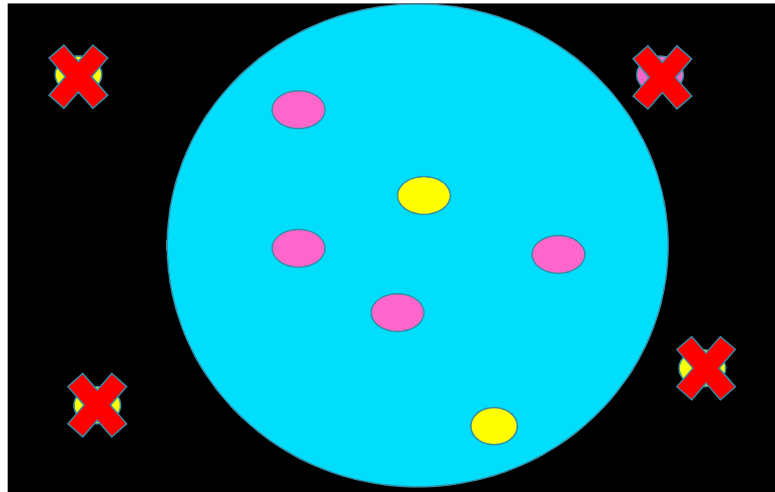
1. Create a mask for each population.
2. Use **Logical Operations > OR** to create a mask that is a union
3. Set **Objects to Measure** to be the union mask.
4. Add **Feature Groups** for each individual mask.

Example 2: Summary Measurements



1. Create a mask for each population.
2. Use **Simple Threshold 0-65535** to create a mask that represents the entire image.
3. Set **Objects to Measure** to be the simple threshold
4. Add **Feature Groups** for each individual mask.

Example 3: Cells in a well



1. Create a mask for the cells.
2. Create a mask for the well.
3. Set **Objects to Measure** to be the well mask.
4. Add **Feature Group** for the cell mask.




Support Resources


- F1 / HELP within MetaXpress® Software
- Support and Knowledge Base: <http://mdc.custhelp.com/>
- Request Support: <http://mdc.custhelp.com/app/ask>
- Technical Support can also be reached by telephone:
 - 1 (800) 635-5577 (North America)
 - <https://www.moleculardevices.com/contact> for other regions
 - Select options for Tech Support → Cellular Imaging Products → ImageXpress Instruments

ADVANCING DISCOVERY

Molecular Devices, LLC
3860 N First Street
San Jose, CA 95134 USA

 +1 408.747.1700

 support@moldev.com

 moleculardevices.com

