



MetaXpress® 6 Software Guide

Logical Operations and Marked Objects

Date Revised 05/26/2020 Version B

Chapter Purpose

The purpose of this chapter is to illustrate the **Logical Operations**, **Remove Marked Objects**, and **Keep Marked Objects** options in the **Custom Module Editor (CME)** plugin.

These tools are found in the CME **Modify Objects** menu. They are used to combine two masks together as appropriate for the analysis.

Overview

Logical Operations

- Performs operations on two masks on a pixel-by-pixel basis
 - AND (intersection)
 - OR (union)
 - ANDNOT (subtract)
 - XOR (union – intersection)

Marked Objects

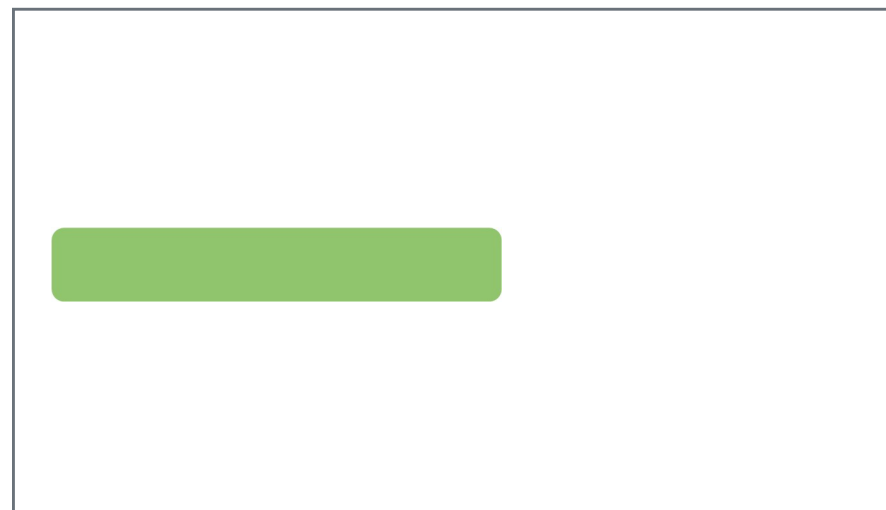
- Performs operations on two masks on an object-by-object basis
 - Remove Marked Objects
 - Keep Marked Objects



Example: Original Masks

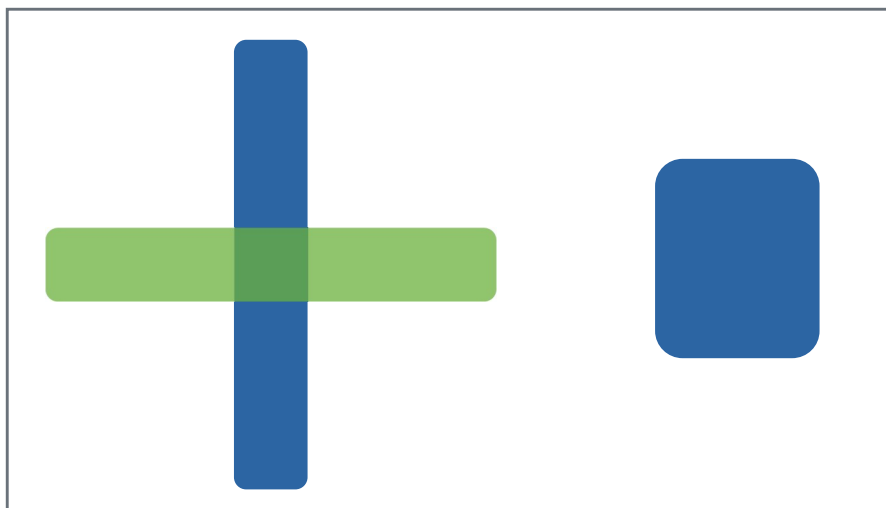


Mask 1 (blue)

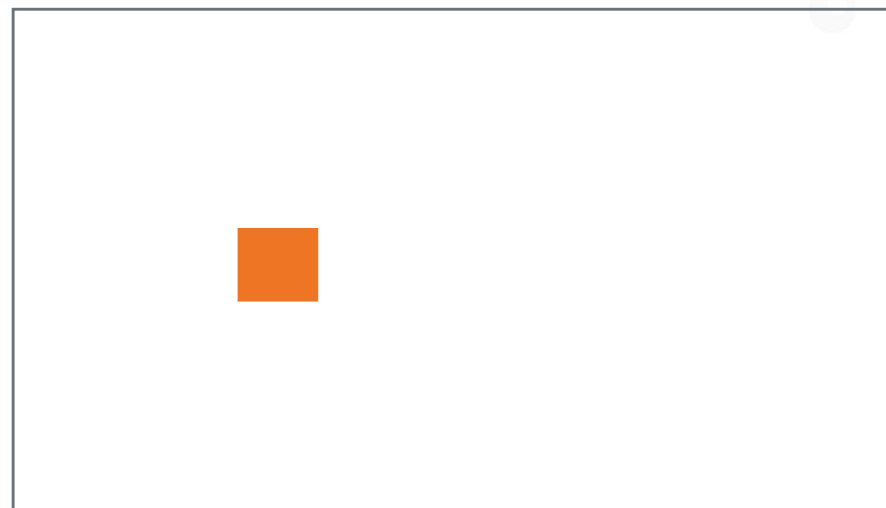


Mask 2 (green)

Logical Operations – AND

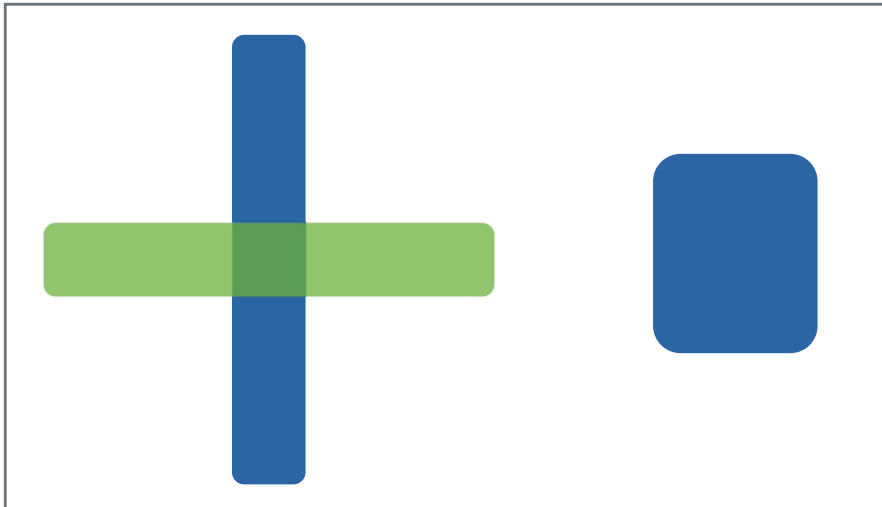


Original Masks:
Mask 1 (blue)
Mask 2 (green)

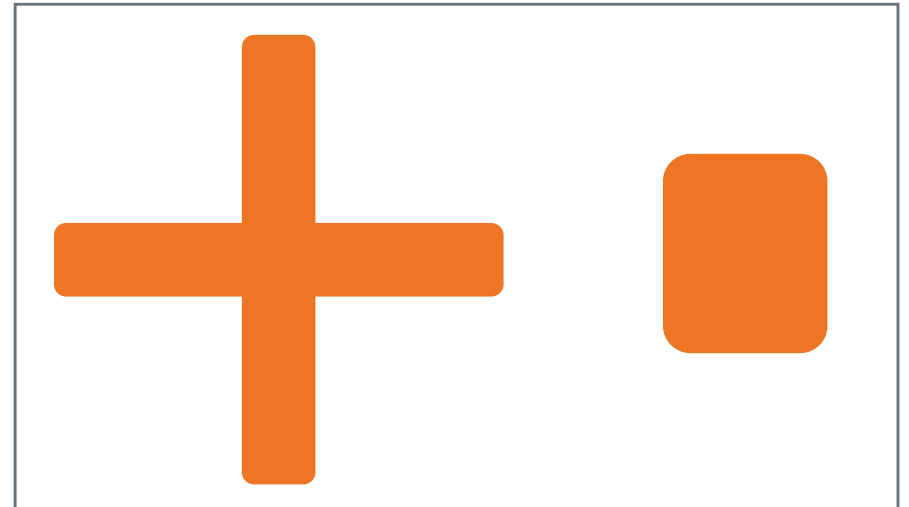


Logical Operations AND result:
Intersection of pixels from
Mask 1 (blue)
Mask 2 (green)

Logical Operations – OR

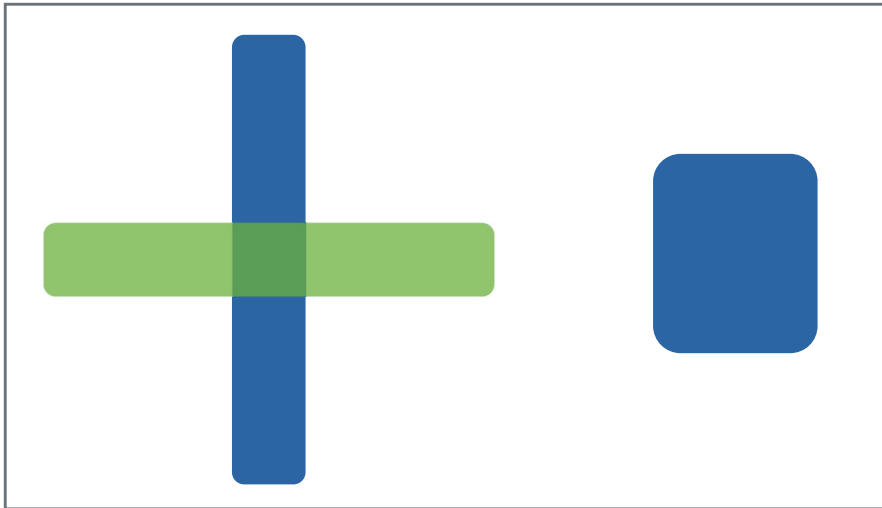


Original Masks:
Mask 1 (blue)
Mask 2 (green)



Logical Operations OR result:
Union of pixels from
Mask 1 (blue)
Mask 2 (green)

Logical Operations – ANDNOT

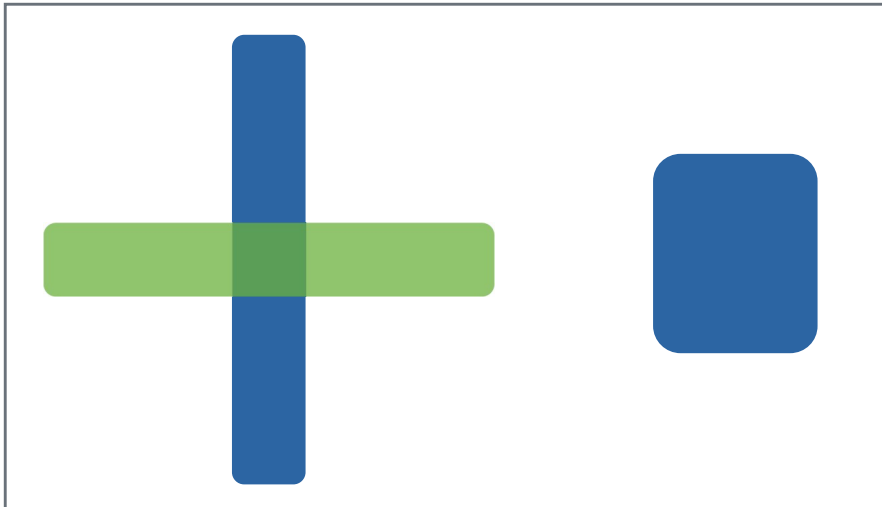


Original Masks:
Mask 1 (blue)
Mask 2 (green)

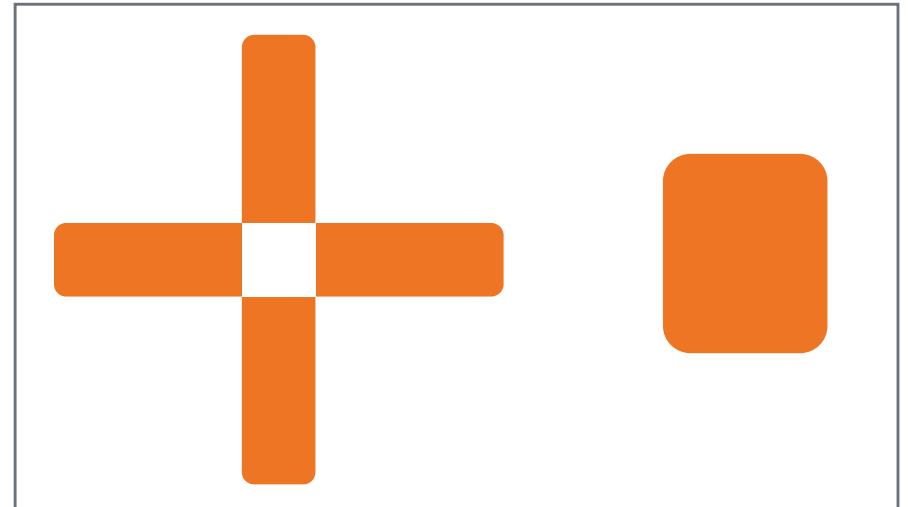


Logical Operations ANDNOT result:
Subtract
Mask 2 (green)
pixels from
Mask 1 (blue)

Logical Operations – XOR

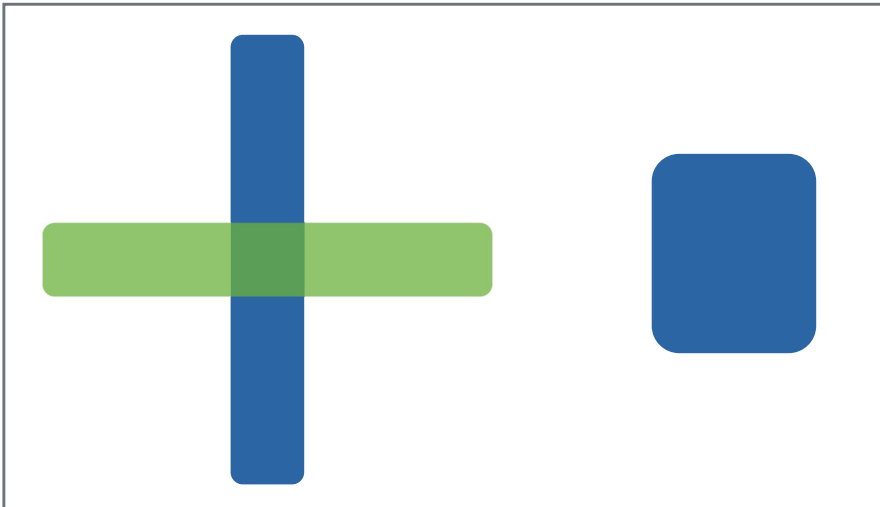


Original Masks:
Mask 1 (blue)
Mask 2 (green)

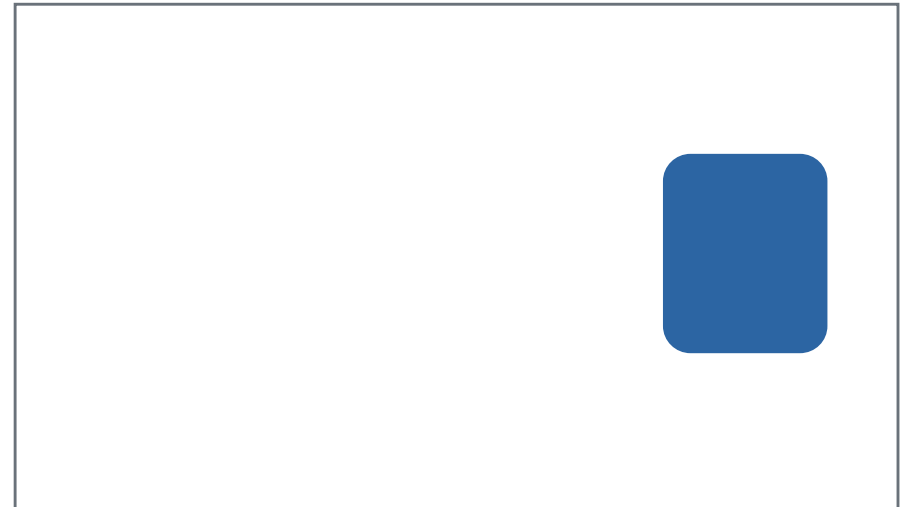


Logical Operations XOR result:
Union minus Intersection of
Mask 1 (blue)
Mask 2 (green)

Remove Marked Objects

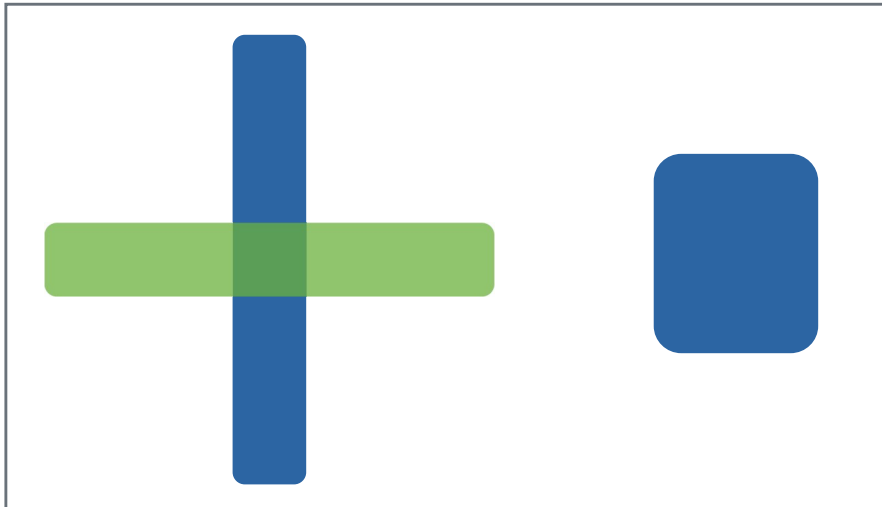


Original Masks:
Objects (blue)
Marker (green)

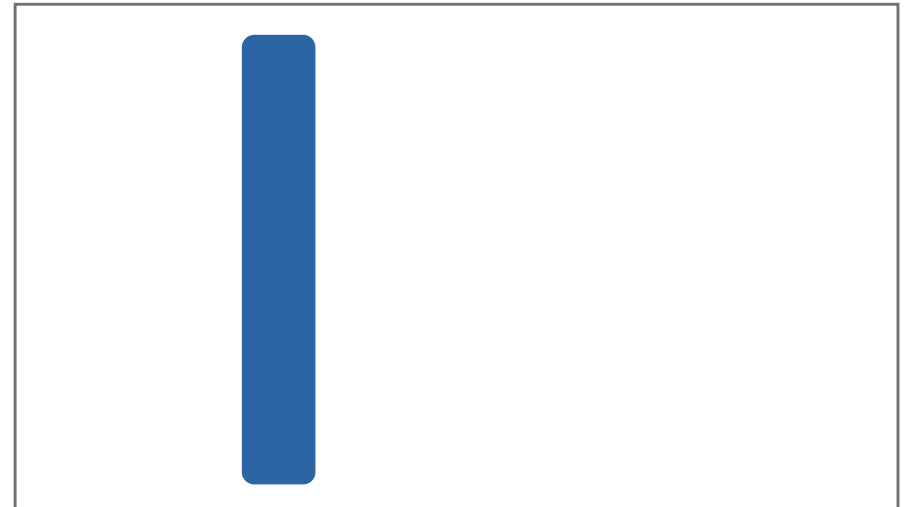


Remove Marked Objects result:
Objects (blue) not touching
Marker (green) remain

Keep Marked Objects



Original Masks:
Objects (blue)
Marker (green)



Keep Marked Objects result:
Objects (blue) touching
Marker (green) remain




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

