

## Using classifiers to select objects for measuring brightness over time.

### Abstract

The capabilities of object classification can be used to select objects before measuring their brightness over time. This technique saves you the effort of manually outlining the objects you want to monitor.

### Instructions

#### Requirements

MetaMorph with the dropins IMA and traceobj installed.

#### Procedure

Note this procedure can be made into a journal for easier operation.

1. Set up measurement preferences for this procedure.  
Choose Preferences (Edit Menu) and select Measure Objects.  
Refer to the manual to decide whether you will want to enable or disable the following options:  
Exclude objects if centroid is not in the current region, exclude objects that touch the edge of the image.
2. Set the threshold on the image of interest. Use the image window threshold tools or Threshold Image (Process menu).
3. Set up the classifiers to include the objects of interest using Integrated Morphometry Analysis (Measure menu). Set the classifier parameters using the interactive Teach mode (set at the bottom of the dialog). Measure the image.
4. Create regions for each of the objects of interest. Use Create Regions Around Objects (Measure Menu).

**Document ID**  
T10125

**Product**  
MetaMorph  
Discovery-1

**Created**  
01-Jul-1998

**Last Reviewed**  
02-Oct-2003

**Keywords:** metamorph techniques  
**Issue Type:** Application