

# MetaXpress Custom Module Editor

## Background Subtraction for Complex Images

Rev A 2018-08-21



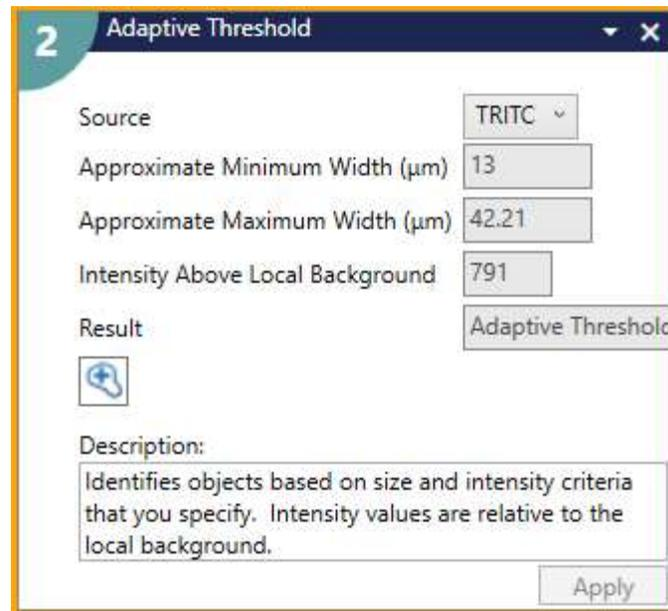
# Chapter Purpose

- This guide explains a method for doing background subtraction within the Custom Module Editor. This method is suitable for more complex images/objects, such as cytoplasmic or mitochondrial stains where cells are touching one another.
- The resulting object intensities will have the background subtracted. This may help in comparing results from plate to plate, where the fluorescent staining may vary.



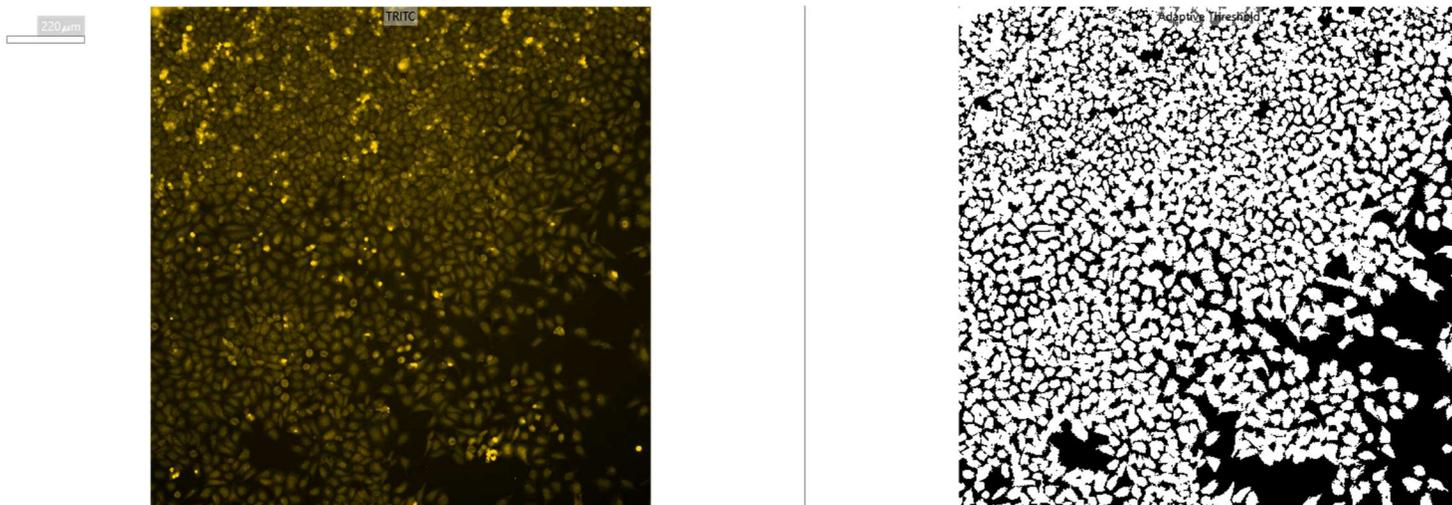
# Step 1: Find Objects > Adaptive Threshold

- Use the Click to Find tool to define settings.
- This should find all objects, even irregularly-shaped ones.



# Step 1: Find Objects > Adaptive Threshold

- Use the Click to Find tool to define settings.
- This should find all objects, even irregularly-shaped ones.



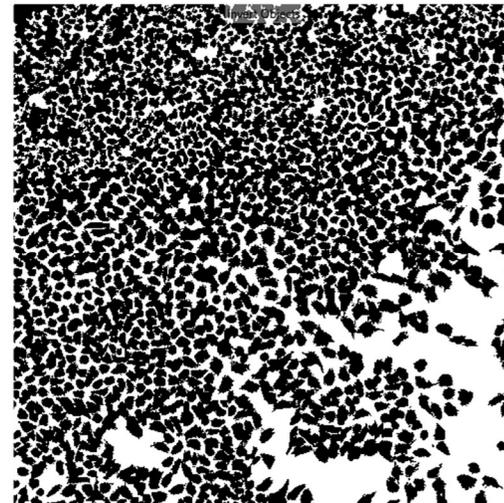
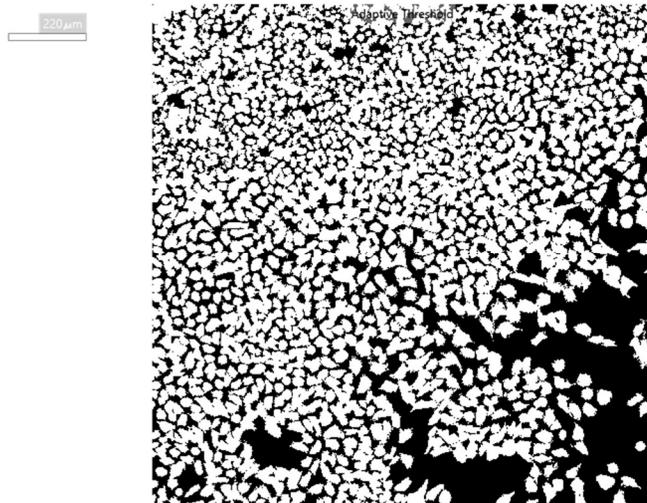
## Step 2: Modify Objects > Invert Objects

- Invert the objects mask, so that all of the background area is selected.



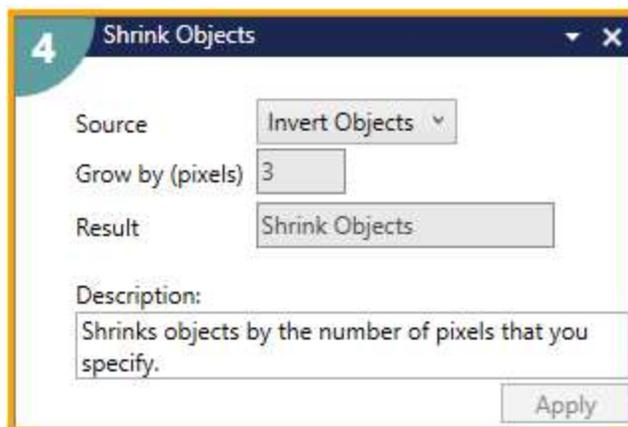
## Step 2: Modify Objects > Invert Objects

- Invert the objects mask, so that all of the background area is selected.



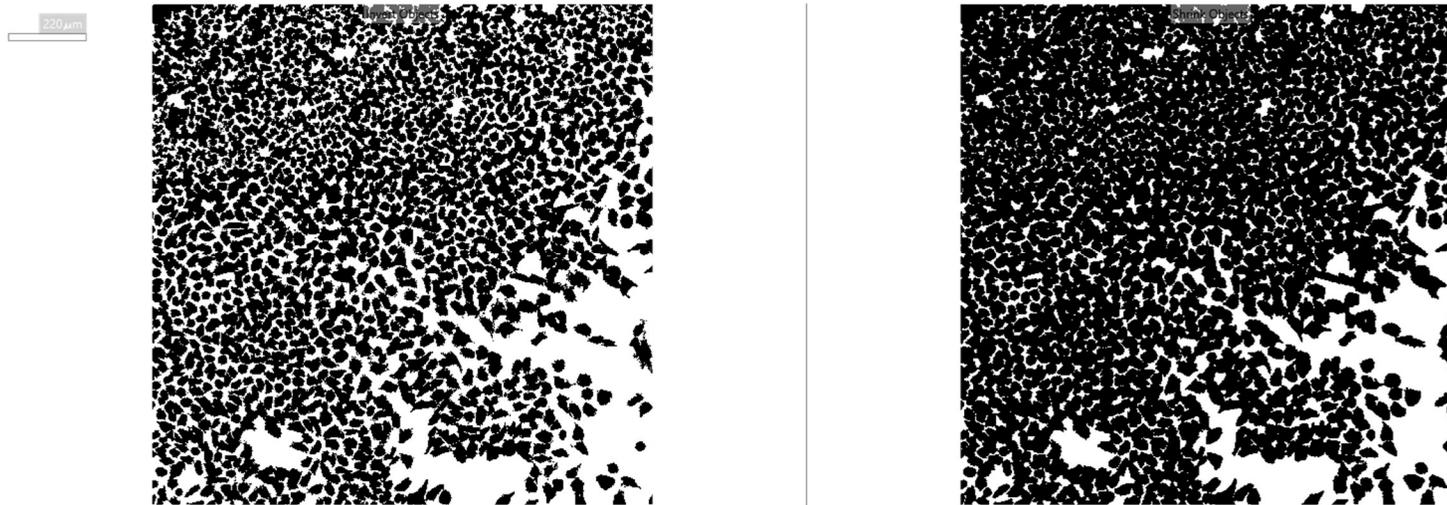
## Step 3: Modify Objects > Shrink Objects

- Shrink the background area by a few pixels, to avoid edge effects.



# Step 3: Modify Objects > Shrink Objects

- Shrink the background area by a few pixels, to avoid edge effects.



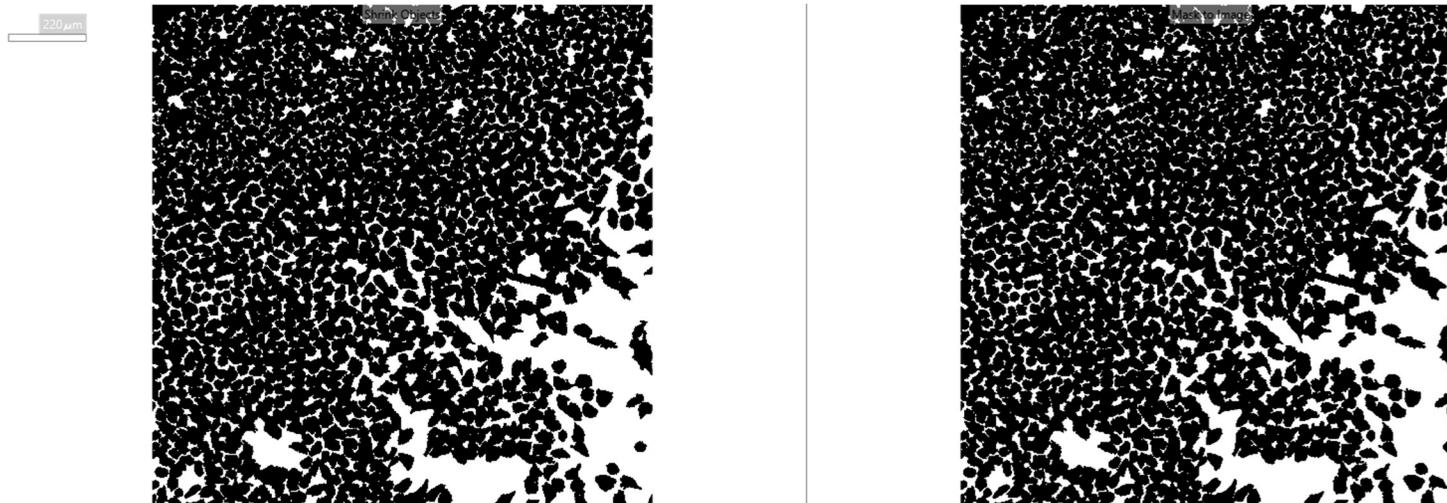
## Step 4: Modify Image > Mask to Image

- Convert the Shrink Objects from a binary mask to a 16-bit image so that it can be used for further image calculations



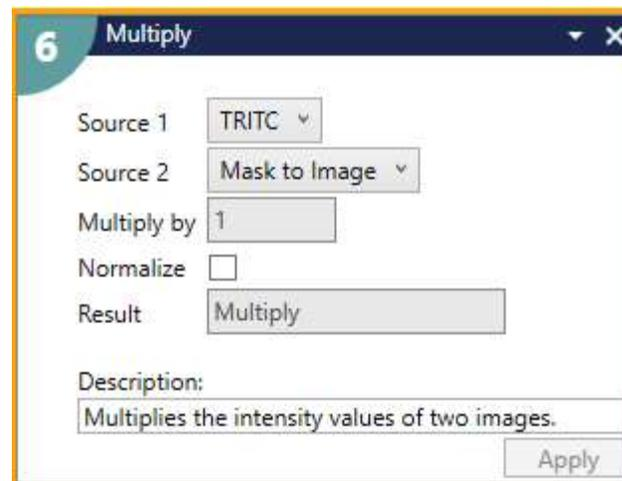
## Step 4: Modify Image > Mask to Image

- Convert the Shrink Objects from a binary mask to a 16-bit image so that it can be used for further image calculations



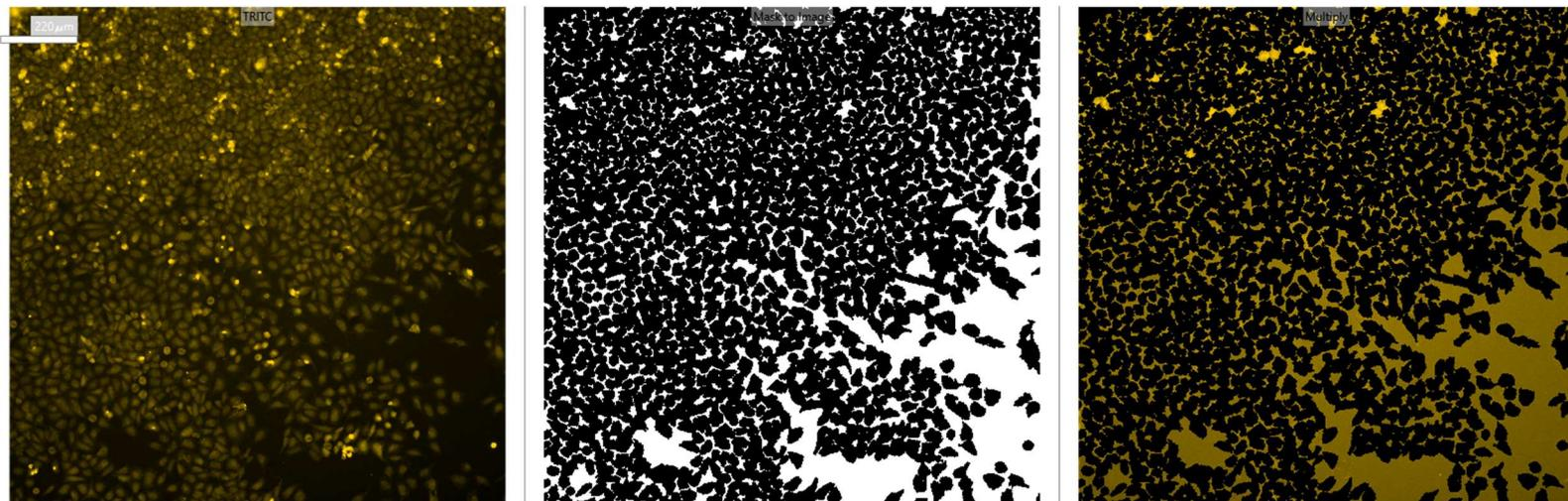
## Step 5: Modify Image > Multiply

- Multiply the original image by the Mask to Image. The background pixels are kept, and the object pixels are changed to zero.



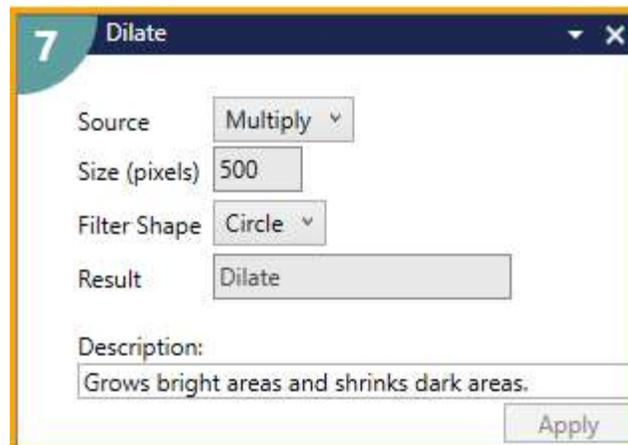
# Step 5: Modify Image > Multiply

- Multiply the original image by the Mask to Image. The background pixels are kept, and the object pixels are changed to zero.



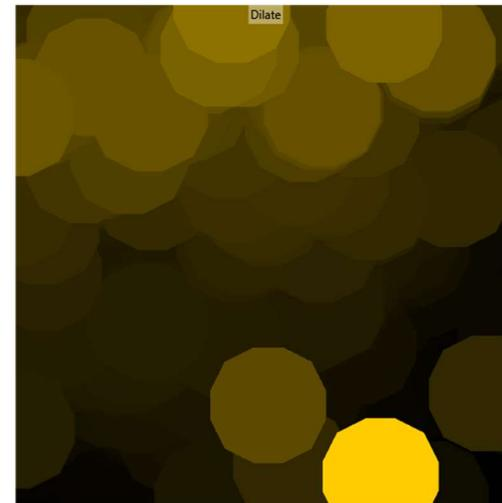
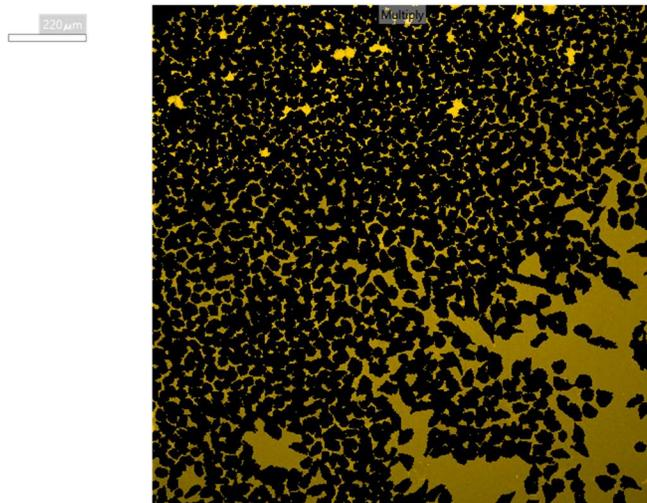
# Step 6: Modify Image > Dilate

- Dilate the Multiply result by a large number of pixels, so that the background area is expanded to cover the object area.



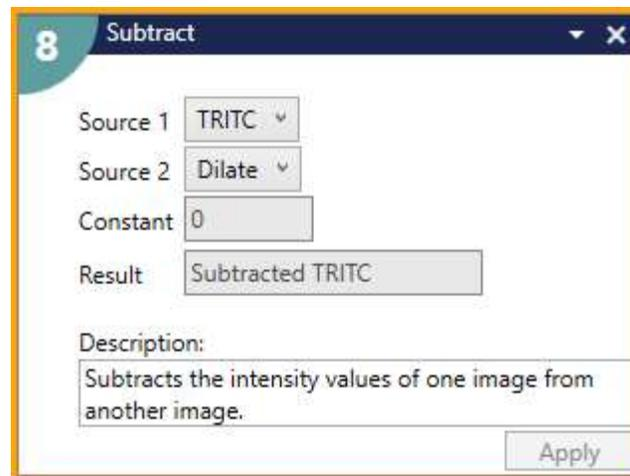
## Step 6: Modify Image > Dilate

- Dilate the Multiply result by a large number of pixels, so that the background area is expanded to cover the object area.



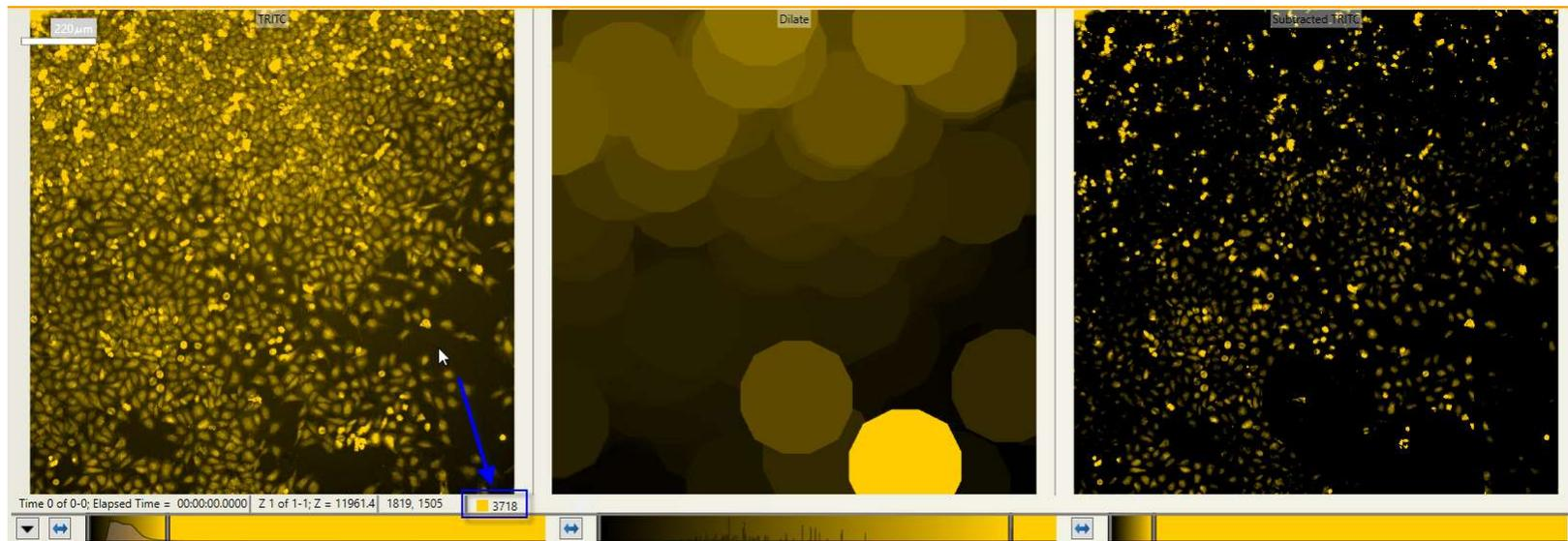
# Step 7: Modify Image > Subtract

- Subtract the Dilate result from the original image.
- The resulting image will have a background at or very close too zero.



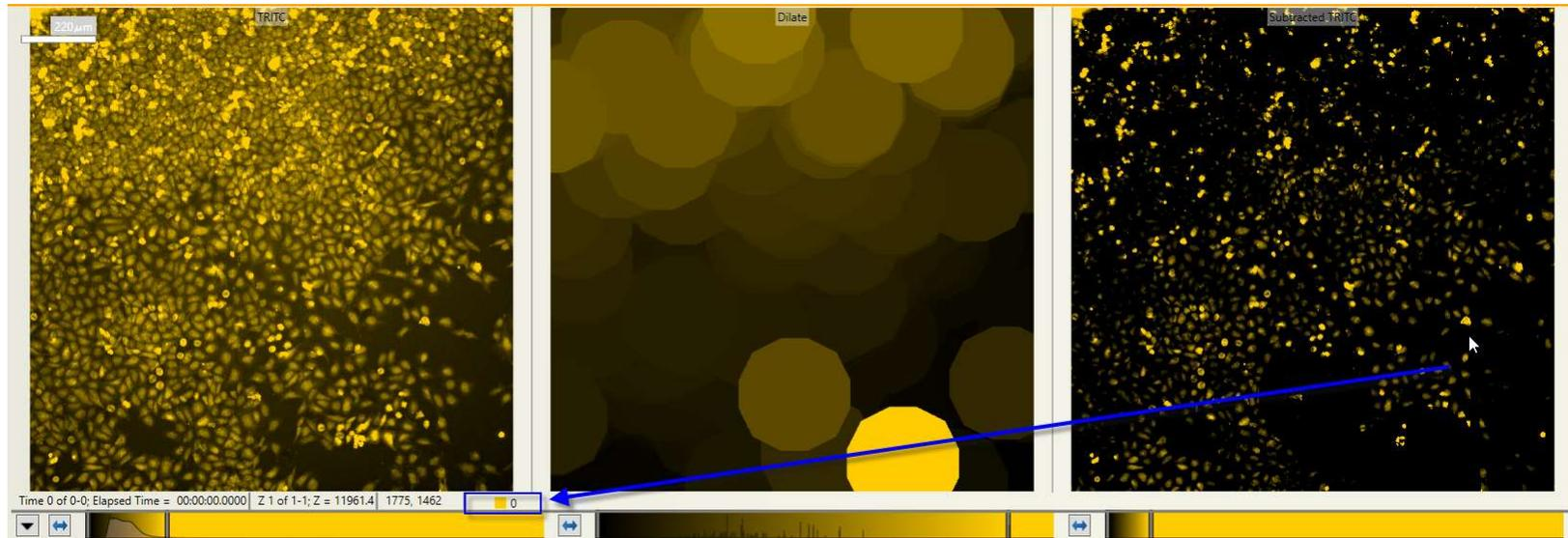
# Step 7: Modify Image > Subtract

- Subtract the Dilate result from the original image.
- The resulting image will have a background at or very close too zero.



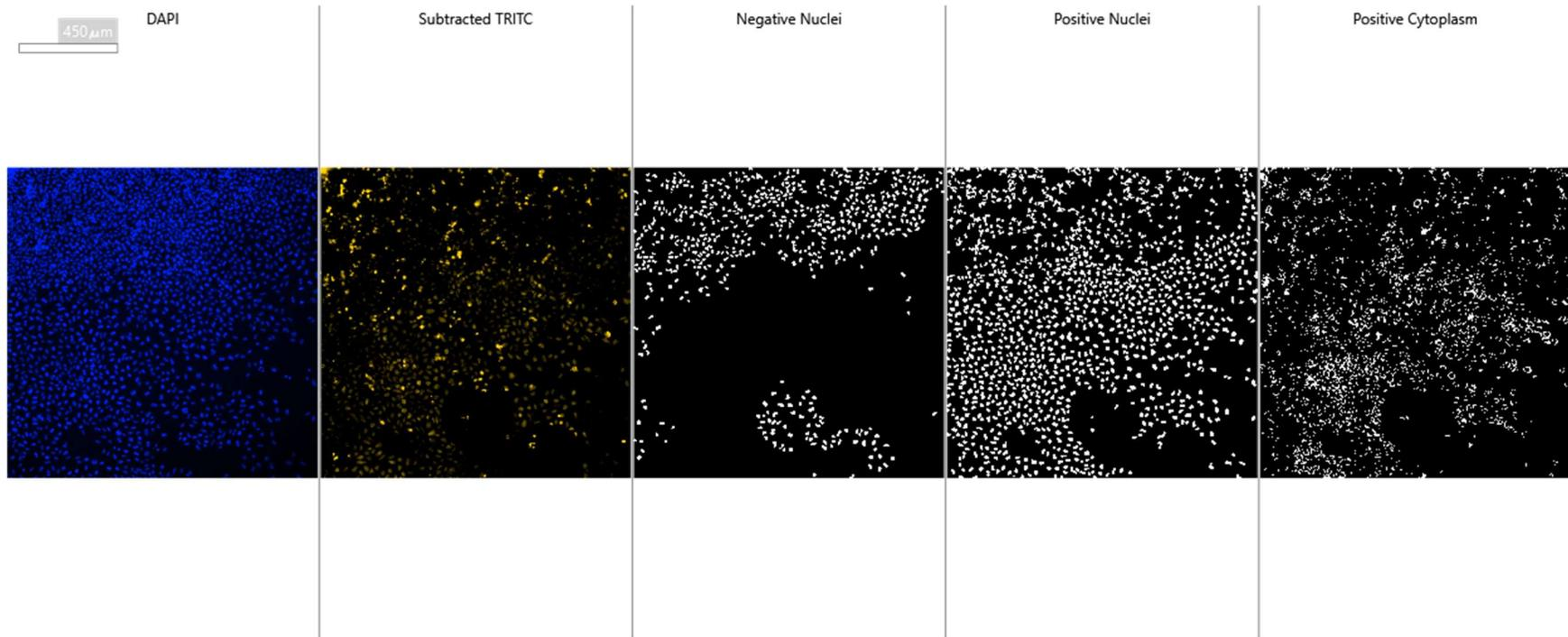
# Step 7: Modify Image > Subtract

- Subtract the Dilate result from the original image.
- The resulting image will have a background at or very close too zero.



# Step 8: Find Objects

- Add segmentation routine of interest.



# Step 9: Measure

- Set up measurements, making sure to select the subtract result image for intensity measurements

11 Measure Mask

Measurement Inputs

Standard Area Value

Create Object Overlay

Objects to Measure

Mask of Objects:

Image to Measure:  ... X

+ Add feature group.

Description:

Objects and features used for measurements.

Apply



# Step 9: Measure

- Set up measurements, making sure to select the subtract result image for intensity measurements

220 μm

