

How to convert a multi-line or traced line region into a trace region

Problem

While trying to outline an object the user chose to use the multi-line

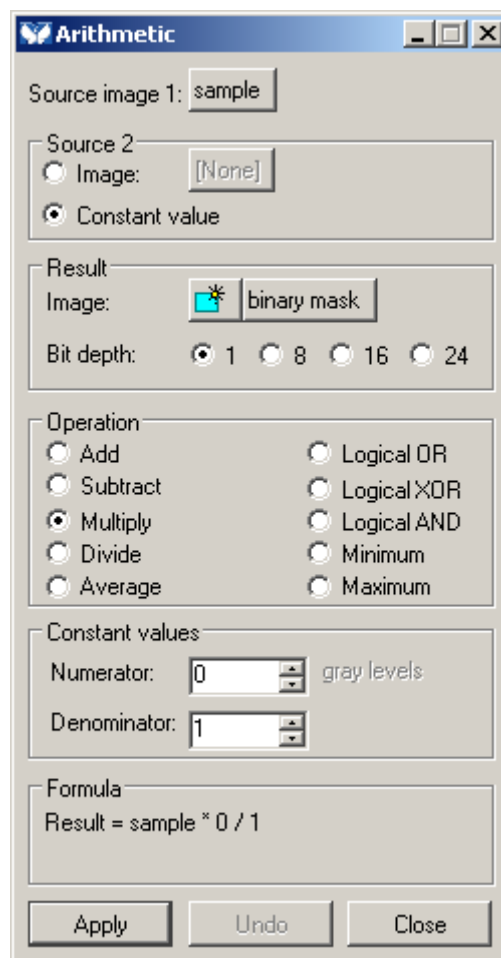
region  or traced line region  instead of a trace region



. The result was that the region selected the pixels underneath the line region, not the area encompassed by the region.

Solution

1. Deselect the region of interest.
2. Create a binary copy of the image by multiplying the image by 0 (Arithmetic dialog) and selecting a binary result image.



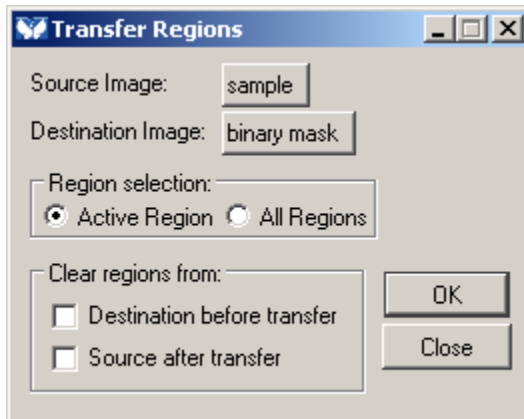
ARTICLE #
#20073

PRODUCTS
MetaMorph®
MetaVue®

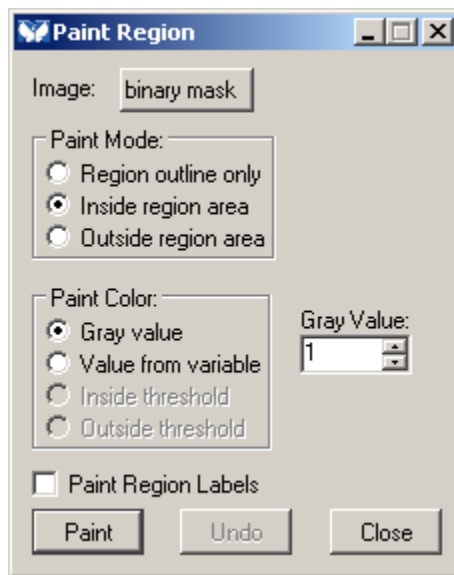
CREATED
5-May-2008

LAST UPDATED
5-May-2008

- Transfer the region of interest from the original image to the binary mask image using Transfer Regions (Region menu).



- On the binary image paint the region grey value 1 using Paint Region (Display menu, Graphics sub-menu).



- Clear the region from the binary image using Clear All Regions (Region menu)
- Choose Create Regions Around Objects (Region menu) to create a Traced region.
- Transfer the region to whichever image window you wish to use it on.