

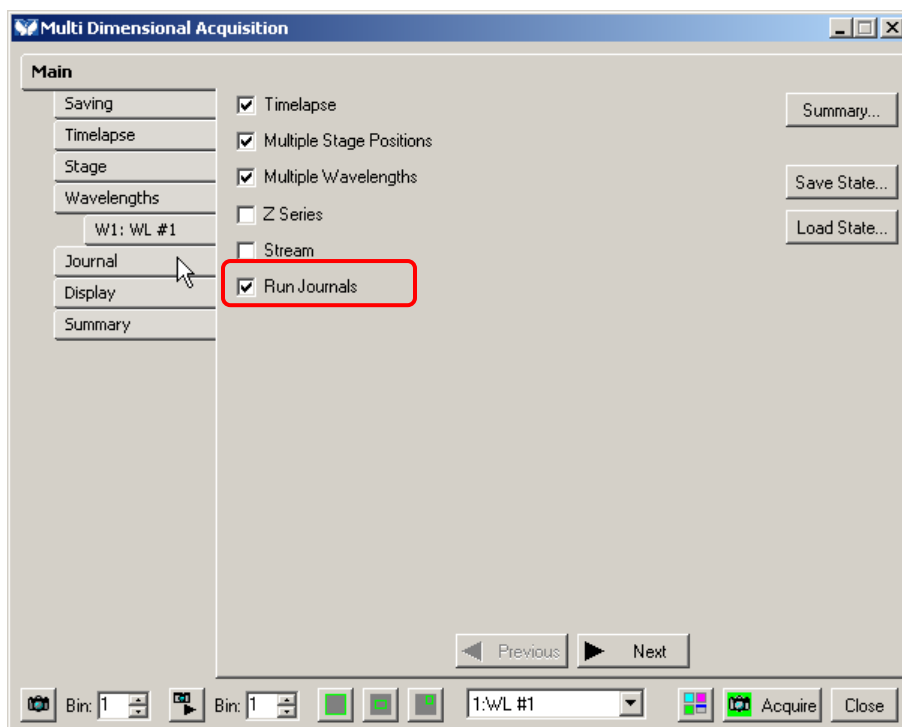
The use of journals during a Multi Dimensional Acquisition

Abstract

The Multi Dimensional Acquisition (MDA) dialog box provides tools and options for most multidimensional applications. However, there are a variety of applications and/or hardware configurations that are not easily handled by these tools and options alone. MetaMorph[®] software introduced the capability of using journals to further customize a multi dimensional experiment.

Instructions

To use journals in MDA, check the “Run Journals” box.



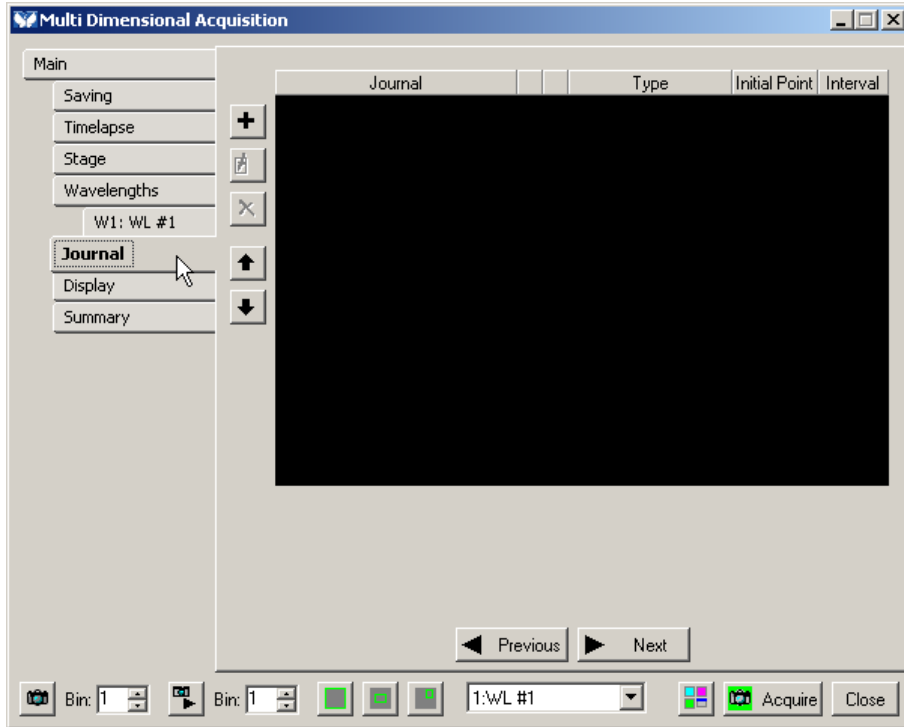
Click on the tab marked “Journal”. The following screen is what you should see if you have a version above 6.1.


ARTICLE #T20117

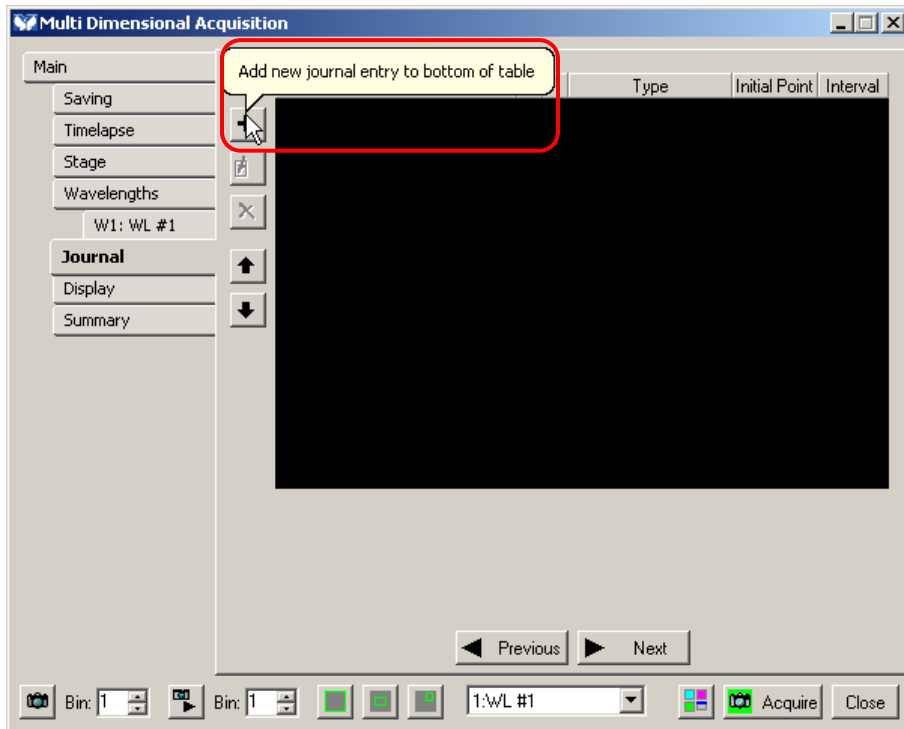
PRODUCTS
MetaMorph[®]
Version 6.1 or above

CREATED
23-June-2009

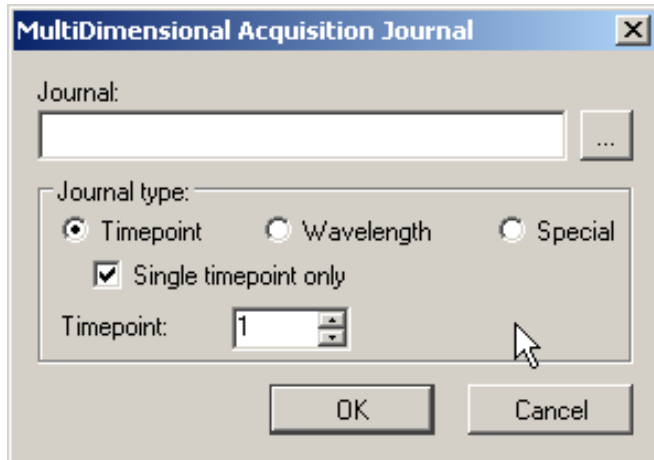
LAST UPDATED
23-June-2009



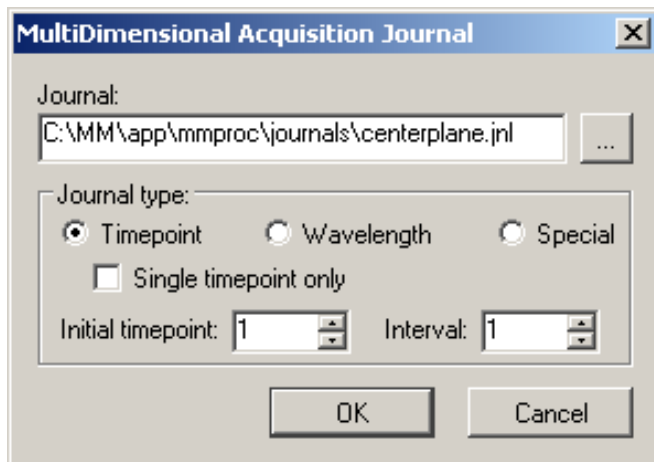
Click on the  sign to add a journal to run.

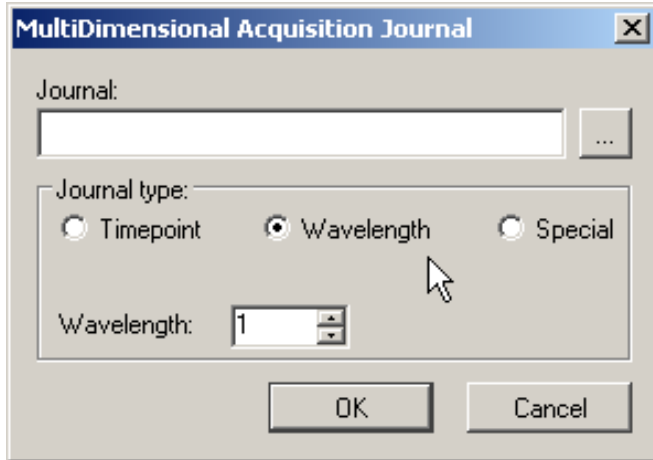


The journal can be run for certain timepoints, wavelengths or special circumstances, as illustrated below.

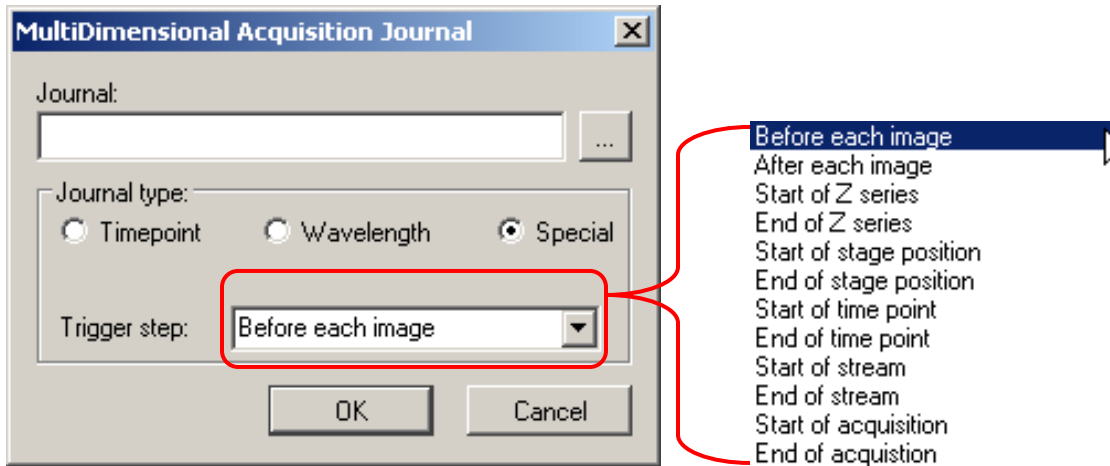


Select either a single timepoint for the journal to run on or select all timepoints by choosing an interval of one. By choosing an interval of two or more, the journal will run on the selected initial timepoint (not necessarily the first timepoint) then every second timepoint (interval of 2), every third timepoint (interval of 3), et cetera.




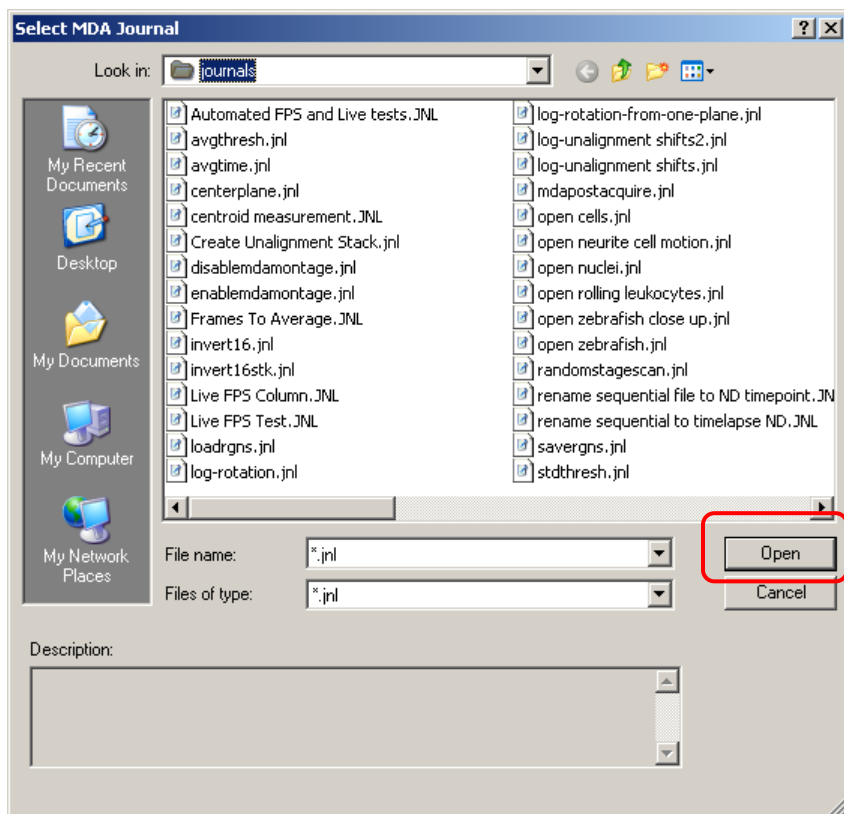


For the wavelength, only one wavelength can use the selected journal at one time. To have the journal run on multiple wavelengths, you must set it multiple times in the journal tab.

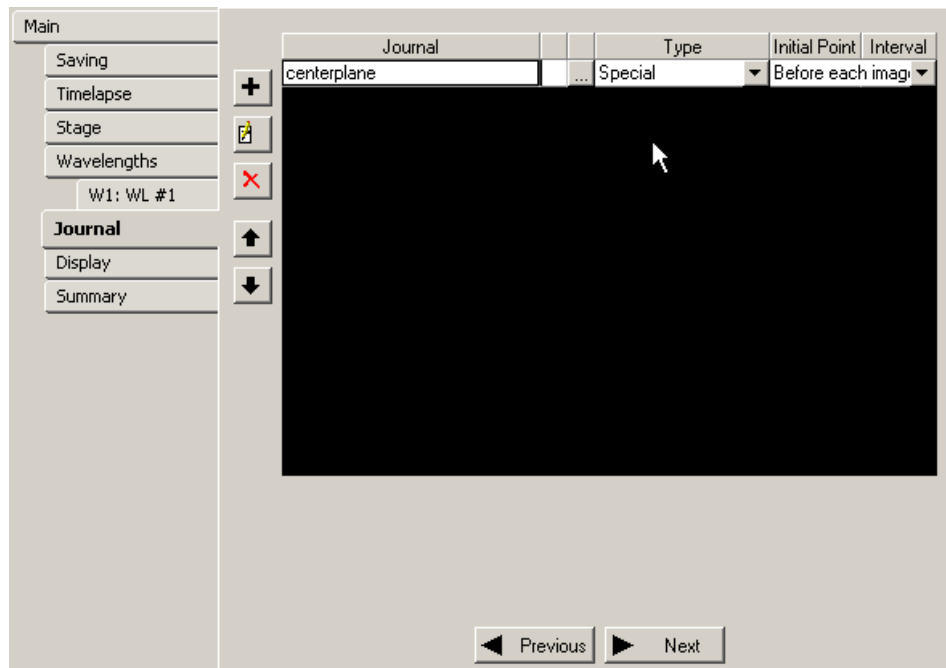


The special choice allows the user the greatest flexibility as to when a journal should run.

Push the button with 3 dots  to select the journal to run. Journals should be saved in <C:\MM\app\mmproc\journals> location on the computer running the device.



Once the journal is selected click on Open and the journal should appear in the Journal window of MDA.



This button allows the user to edit which journal is selected.



This button deletes the current journal selected.



These buttons move the selected journal up or down in the order.

Select the Summary tab and the summary should contain all information about the experiment as well as about the journal(s) selected to run during the MDA.

Multi Dimensional Acquisition Summary

MDA test

Directory: C:\MM\images\Testing

File Base Name: MDA test

Timelapse

6 time points every 50 ms, duration = 250 ms

Stage Positions

2 stage positions

Wavelengths

Wavelength 1

Illumination: WL #1

Collection: Every time point

Exposure: 25 ms

Gain: 0

Digitizer: 10 MHz (EM Gain)

Auto Focus: Off

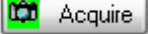
12 total images requiring 33.13 MB

Journals:

Before each image: centerplane

Description

Multi Dimensions Experiment

When all parameters of the Multi Dimensional Acquisition are correct, hit  to start the experiment.

For further assistance please contact Meta Imaging Series® support at:

800-635-5577 option 3-2-2 OR support.dtn@moldev.com.