

Multi Dimensional Acquisition:

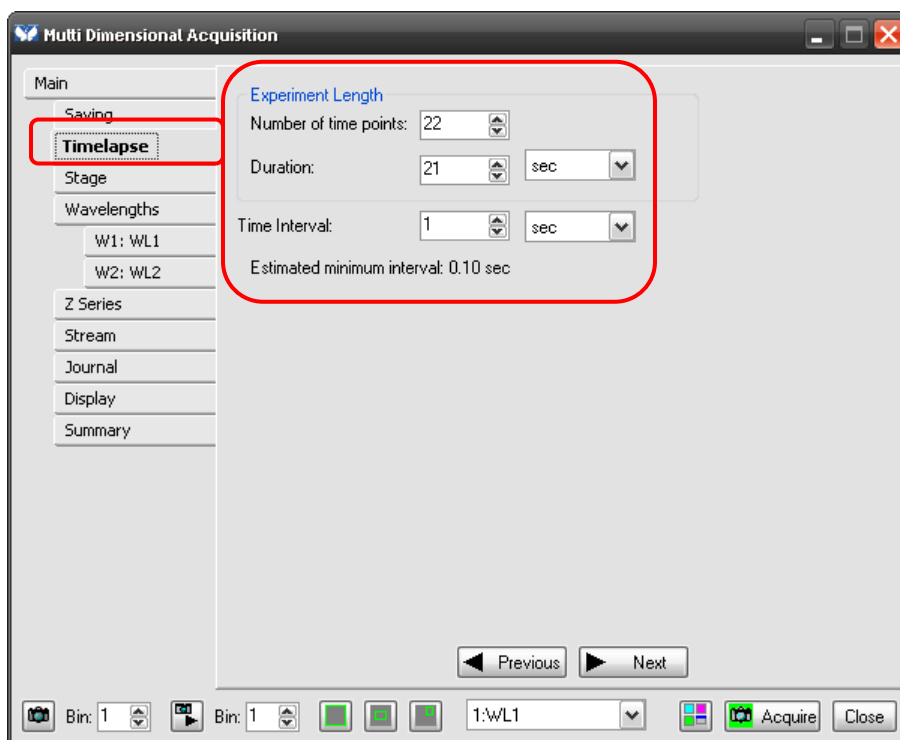
Timelapse Tab

Abstract

The Multi Dimensional Acquisition (MDA) dialog box provides tools and options for most multidimensional applications. This note will describe the Timelapse Tab. For further information on Multi Dimensional Acquisition see T20117, T20121, T20123, T20124, T20125, T20126, T20127, T20128 and T20129.

The Timelapse Tab

The Timelapse tab is where the timelapse experiment is configured.



Number of Time Points

- Specifies the total number of times you want an image to be acquired. Enter 0 if you want the experiment to run until halted.

Duration

- Specifies the time it will take to acquire the number of times points based on the interval. Changing this field updates the *Number of time points* field by calculating the number of time points from the time interval and the duration. Set the time units in the adjacent box. The available units are msec, seconds, minutes, and hours. This setting is not visible if either the *Number of time points* or the *Time Interval* is set to 0.

Time Interval

ARTICLE #T20122

PRODUCTS
MetaMorph®
Version 7.0 or above

CREATED
23-June-2009

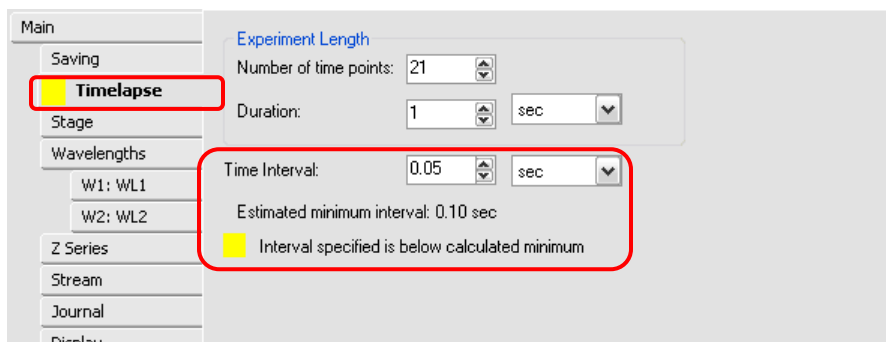
LAST UPDATED
25-August-2009

→ Specifies the amount of time to wait between acquisitions. In the box next to the Time Interval box, select the unit of time you want to use.

NOTE: If the acquisition cannot be accomplished in a time period shorter than the interval, acquisition will move to the next time point as soon as the current time point is completed.

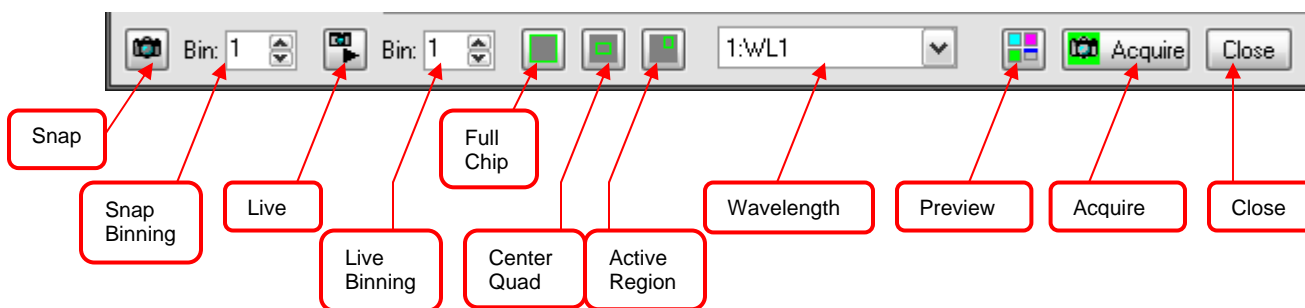
Estimate minimum interval

→ Indicates the estimated minimum time interval calculated from the sum of the exposure time for the wavelength multiplied by the number of images that need to be acquired for each loop.



→ If the specified interval is not sufficient, an appropriate warning message is displayed, such as *Interval specified is below the calculated minimum*.

Across the bottom of the dialog box are the functions for snapping an image, binning, live, live binning, chip dimensions, wavelength chooser, preview, acquire and close. These options are always available no matter which tab is selected.



If there are any further questions, please contact Technical support at 800-635-5577 x1820 or support.dtn@moldev.com.