

Multi Dimensional Acquisition:

Alignment Dialog

Abstract

The Multi Dimensional Acquisition (MDA) dialog box provides tools and options for most multidimensional applications. This note will describe the Alignment dialog. For further information on Multi Dimensional Acquisition see T20117, T20121, T20122, T20123, T20124, T20125, T20127, T20128 and T20129.

The Alignment Dialog

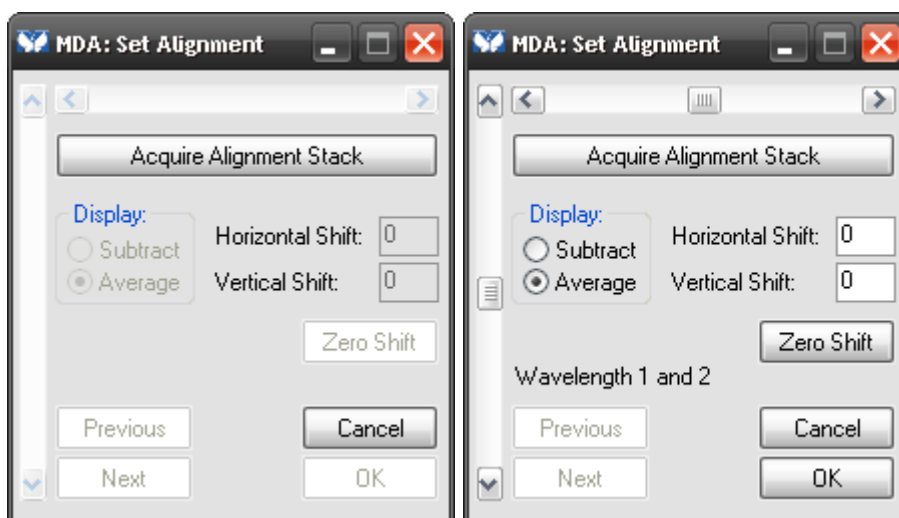
The Alignment dialog is where alignment is configured for the wavelengths in a multidimensional experiment. This dialog will only be displayed if the appropriate hardware is installed and detected by the MetaMorph® software.

ARTICLE #T20126

PRODUCTS
MetaMorph®
Version 7.0 or above

CREATED
27-August-2009

LAST UPDATED
27-August-2009



Acquire Alignment Stack

→ Acquires an image at each wavelength and builds a stack that can be used in adjusting alignment. Must be done first for the following options to become available.

Display

→ Selects the method to be used to display differences between the reference plane and the shifting plane:

Subtract

→ Uses subtraction to show the difference between the reference plane and the shifting plane. The planes will be aligned when there is nearly a uniform grayscale level throughout the entire image.

Average

- Uses averaging to display the offset between the two planes. The aligned plane should look like the original plane with as little blurring as possible.

Horizontal Shift (text box and slider)

- Adjusts the horizontal alignment of the plane in one-pixel increments

Vertical Shift (text box and slider)

- Adjusts the vertical alignment of the plane in one-pixel increments.

Zero Shift

- Resets the horizontal and vertical shift to zero.

Previous

- Places the previous plane in the alignment image window.

Next

- Places the next plane in the alignment image window.

Cancel

- Cancels your adjustments and closes the dialog box.

OK

- Applies the shift to all of the wavelengths.



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