

Imports Cellomics DIB Files into a Local, Network, or Database Location

Overview

Use the Import Cellomics Data command to convert sets of Cellomics DIB files into data sets in the database that can be accessed using either the Plate Data Utilities (DB) command or the Review Plate Data (DB) command. The command creates MetaXpress HTD files based on the data in the DIB files and converts the images to the TIFF format. It then moves the data into your MetaXpress screening database or a selected local or network folder.

Note: Cellomics DIB files do not include image calibration data. Unless you enter a specific calibration when you import the data, a default calibration of 1 μm = 1 pixel will be applied to the imported images.

Instructions

Use the following procedure to import Cellomics data into MetaXpress:

1. From the Apps Menu, select Plate Data Utilities. The Plate Data Utilities dialog box opens.
2. Click *Import Images*. The Import Images dialog box opens.
3. Click *Import Cellomics Data*. The Import Cellomics Data dialog box opens.
4. Click *Select Directory*. The Browse for Folder dialog box opens.
5. Navigate to the local or network folder containing the Cellomics data to import using the Browse for Folder dialog box, then click *OK*. The path is displayed next to the *Select Directory* button and the Cellomics files are displayed in the *Cellomics Data Files* field, as shown in Figure 1:

Document ID
D50027

Product
MetaXpress

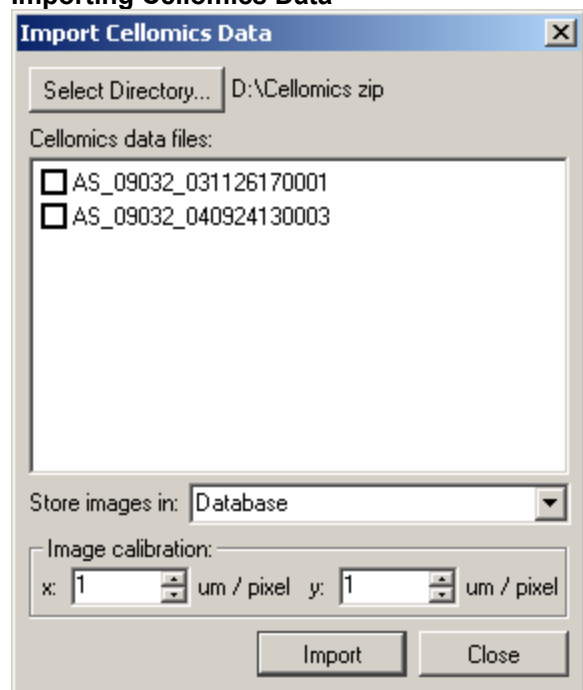
Created
16-December-2004

Last Reviewed
16-December-2004

Keyword: metaxpress cellomics database import

Issue Type: dbs database

Figure 1
Importing Cellomics Data



6. Select the check boxes next to the Cellomics file sets to import in the *Cellomics Data Files* field.
7. Select a location (either the database or an available local or network folder) to import the Cellomics data to from the *Store images in* drop-down box.
8. Enter a calibration ratio (ums per pixel) for the x and y values of the imported images in the *Image calibration* fields.
9. Click *Import*. The files are imported and will be available for review in the Review Plate Data (DB) dialog box.
10. Click *Close* to close the dialog box.