

MetaXpress[®] 6.5 Software Guide

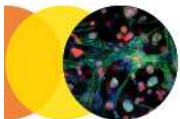
Search Plates

Date Revised 10/23/2018 Version A



Chapter Purpose

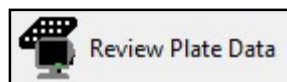
The purpose of this chapter is to explain the Search Plates feature added in MetaXpress 6.5.



Review Plate Data

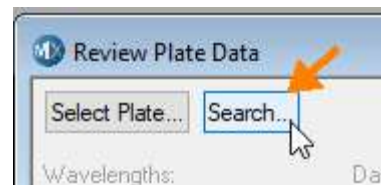
1. Open **Review Plate Data**

- In the main toolbar click on

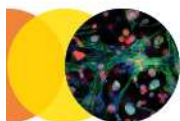


OR

- Under the **Screening** menu, select **Review Plate Data**



2. Click **Search**



Search Plates dialog

Review Plate Data - Search Plates

Search text: Start Date: End Date:

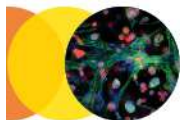
Search exact text Search current user only

Search Fields: Plate Name Plate Description Folder Name Barcode

Plate ID	Plate Name	Folder Name	Barcode	Plate Description	Creator...	Time Created

Double-click on search results to load a plate into Review Plate Data

- Enter the **Search text** and optional date range.
- Enable or disable Search options.
- Click **Search**.
- Optionally change default search options with **Save Reset State**.
- **NOTE:** If your database uses both 'Folder Name' and 'Experiment Set', the 'Folder Name' search will check both.



Search Results

The screenshot shows a software window titled "Review Plate Data - Search Plates". It features a search interface with the following elements:

- Search text:** A text box containing "transfluor".
- Start Date:** A date input field with the placeholder "yyyy-mm-dd".
- End Date:** A date input field with the placeholder "yyyy-mm-dd".
- Buttons:** "Search" and "Reset" buttons are located to the right of the date fields.
- Search Fields:** A section with checkboxes for "Plate Name", "Plate Description", "Folder Name", and "Barcode", all of which are checked.
- Search Options:** Checkboxes for "Search exact text" (unchecked) and "Search current user only" (checked).
- Save Reset State:** A button located to the right of the search fields.
- Table:** A table displaying search results with columns: Plate ID, Plate Name, Folder Name, Barcode, Plate Description, and a partially visible column on the right. The table contains three rows, with the second row (ID 11) highlighted in blue.

Plate ID	Plate Name	Folder Name	Barcode	Plate Description	Cn
10	TransfluorPitsBinned	Training		Transfluor Pits 40x 2x2 binning	Pa
11	TransfluorVesiclesBinned	Training		Transfluor Vesicles 20x 2x2 binning	Pa
155	EX2TransfluorVesicles	CME Training		EX2TransfluorVesicles	Pa

At the bottom of the window, there is a message: "Double-click on search results to load a plate into Review Plate Data" and a "Close" button.

- Double-click on a row to select it in the Review Plate Data
- Search results can remain open, in case you want to check multiple plates in the list



Support Resources

- F1 / HELP within MetaXpress® Software
- Support and Knowledge Base: <http://mdc.custhelp.com/>
- Request Support: <http://mdc.custhelp.com/app/ask>
- Find Contact information for your local team:
 - <https://www.moleculardevices.com/contact>
 - For North America, call 1 (800) 635-5577
 - Select options for Tech Support → Cellular Imaging Products → ImageXpress Instruments

