

# Scan the Odyssey® Fc Imager Verification Plate (926-75002)

## Before You Begin

- » **Re-certify** the verification plate annually. Contact [biosupport@licor.com](mailto:biosupport@licor.com) to request the re-certification service (service number 926-75005).
- » **Handle** the verification plate with powder-free gloves.
- » **Store** the verification plate in the envelope it ships with.
- » **Only** touch the plate by the edges.
- » **Never** touch the glass or the stainless steel on the top of the plate.
- » **Never** attempt to clean the plate yourself. If the stainless steel face becomes contaminated, return the plate to LI-COR Biosciences for cleaning and re-certification.

## Position the Plate

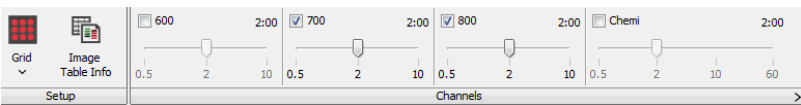
1. Ensure the instrument's imaging drawer is clean, dry, and free from any residue or dust.
2. Place the verification plate in the imaging drawer with the glass facing up.



## Begin Scan

**Note:** If you are using Image Studio 3 or earlier, please contact LI-COR for instructions.

1. Open Image Studio™ Software with a Work Area created specifically for verification plate images.
2. Connect to the instrument.
3. On the Acquire tab, ensure the following settings are chosen.



**Setup group:** Choose the Grid preset.


**Channels group:** Select the 700 and 800 channels and set the time to 2 minutes for both channels.

4. Click Start .

## Data Analysis

1. On the Analysis tab, ensure Grid is chosen in the Type group.
2. Ensure the following columns are in the Grid Spots table in the order shown.


Image Name	Channel	Spot Name	Signal	Total	Area	Bkgrnd.	Type	Channels	Instrument Name	Integration Times	Trim Mean
0000586_07	800	A01	0.372	0.383	256	0.0000434	Signal	700 800	OFC-0944	02:00 02:00	0.00150
0000586_07	800	A02	0.367	0.378	256	0.0000434	Signal	700 800	OFC-0944	02:00 02:00	0.00148
0000586_07	800	A03	0.365	0.376	256	0.0000434	Signal	700 800	OFC-0944	02:00 02:00	0.00147


Add any missing columns by clicking Columns  above and to the right of the tables.

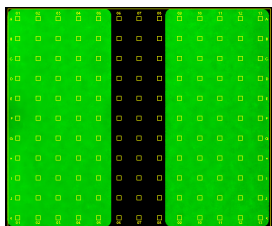
## Data Analysis (continued)





3. If you do not already have a Grid analysis template set up for analyzing verification plate images, set up the following template.

Rows	11	<input type="radio"/> Circle	↑	↓
Columns	13	<input checked="" type="radio"/> Square	Larger	Smaller
Grid	>	Spot		

- In the Grid group, set the number of Rows to 11 and the Columns to 13.
- In the Spot group, change the shape to Square.
- Also in the Spot group, click Smaller  until the Area value shown in the Grid Spots table is 256.
- Adjust the Grid Boundary until it is drawn closely around the two fluorescent regions.
  - The middle three columns of spots should be on the dark background of the plate.
  - All other spots should be inside the fluorescent regions.

**Note:** You may need to redraw the Grid Boundary to get the boundary and spots placed correctly. Click Redraw Boundary  and then click and drag a new boundary.



- Save the Grid Analysis settings as a template.
  - In the Grid group, click Custom  > Save Current Template .
  - Choose this template from the Template list the next time you need to analyze a verification plate. The settings in the Grid Analysis tab will be applied automatically.
- Assign background spots.
  - Click and drag to select the middle three columns of spots.
  - On the Grid Analysis tab, select User Defined from the background list.
  - Click OK to dismiss the User Defined Background dialog.
  - In the background group, click Assign Wells.
- Export data. Click Report  > Launch Spreadsheet  above and to the right of the tables.
- LI-COR provides an Excel® workbook for analyzing verification plate data. The workbook and instructions are available at [licor.com/ovp\\_fc](http://licor.com/ovp_fc).

© 2021 LI-COR, Inc. LI-COR, Odyssey, and Image Studio are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.



4647 Superior Street • Lincoln, NE 68504 • United States • Toll free: 800-645-4267  
Intl.: +1-402-467-0700 • [regulatory@licor.com](mailto:regulatory@licor.com)  
[licor.com/bio](http://licor.com/bio)

Doc #984-16285  
05/21