

# Verification Plate for the Odyssey® CLx Imager

## Before You Begin

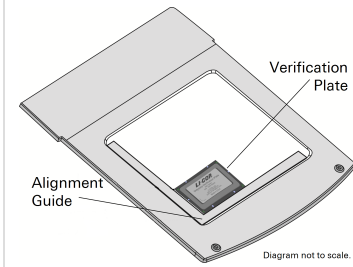
- » **Re-certify** the verification plate annually. Contact [biosupport@licor.com](mailto:biosupport@licor.com) to request the re-certification service (service number 926-75005).
- » **Handle** the verification plate with powder-free gloves.
- » **Store** the verification plate in the envelope it ships with.
- » **Only** touch the plate by the black outer casing.
- » **Never** touch the stainless steel part 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

Ensure the instrument's scan bed is clean, dry, and free of any residue or dust.

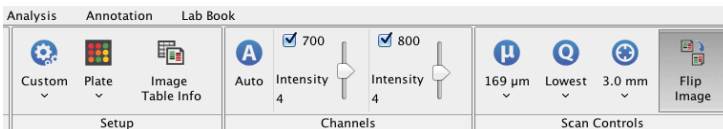
1. Place the alignment guide (PN 926-72200) in the front left corner of the instrument.
2. Place the verification plate with the stainless steel section down, the text facing up, and the left/bottom sides of the verification plate contacting the alignment guide.

Repeat the scan in the other three corners using the alignment guide to align the plate in each corner. Do not rotate the plate.



## Begin Scan

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 Plate preset.

**Channels group:** Deselect Auto and set intensity for both channels to 4.

**Scan Controls group:** Set the resolution to 169 µm, quality to Lowest, focus offset to 3.0mm, and enable Flip Image.

4. Click Start .

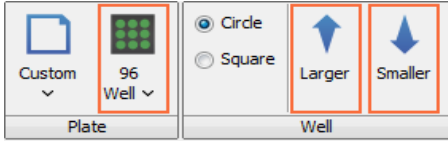
## Data Analysis

1. On the Analysis tab, ensure Plate is chosen in the Type group.
2. In the Plate Wells table, add the Trim Signal column.

**Image Studio 4.x or 5.x:** Click Columns  above and to the right of the tables.

**Image Studio 3.x or earlier:** On the Table tab, click Add/Remove Columns.

3. On the Plate Analysis tab, ensure the following are set up correctly.



**Plate group:** Choose the 96-Well Plate preset.

**Well group:** Click Larger/Smaller until the Area (in the Area column of the Plate Wells table) is 1649.

4. Assign background wells.
  - a. Click and drag to select the background wells (column 6).
  - b. On the Plate Analysis tab, select User Defined from the Background list.
  - c. Click OK to dismiss the User Defined Background dialog.
  - d. In the Background group, click Assign Wells.
5. Select images and choose data to export.
  - a. In the Images table, select the images to be analyzed. Click Filter > Selection.
  - b. In the Plate Wells table, clear the Display Current filter. Click Filter > Display Current.
6. Export Data.

Ensure columns in the Plate Wells table are in the order shown below.

Image Name	Channel	Signal	Total	Area	Type	Plate Name	Well Name	Focus	Instrument Name	Intensities	Trim Signal
0002466_01	700	402000	419000	1649	Signal	96 Well	H12	3.0	CLX-0012	4 4	407000
0002466_01	700	402000	419000	1649	Signal	96 Well	G12	3.0	CLX-0012	4 4	407000
0002466_01	700	401000	418000	1649	Signal	96 Well	G04	3.0	CLX-0012	4 4	407000
0002466_01	700	401000	418000	1649	Signal	96 Well	F01	3.0	CLX-0012	4 4	406000

**Image Studio 4.x or 5.x:** Click Report  > Launch Spreadsheet  above and to the right of the tables.

**Image Studio 3.x or earlier:** On the Table tab, click Launch Spreadsheet.

7. LI-COR provides an Excel® workbook for analyzing verification plate data. The workbook and instructions are available at [www.licor.com/ovp](http://www.licor.com/ovp).

© 2019 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 • Toll free: 800-645-4267 • Intl.: +1-402-467-0700 • [regulatory@licor.com](mailto:regulatory@licor.com)  
[licor.com/bio](http://licor.com/bio)