



1) Clean microplate () or membrane carrier ().

 Clean the bottom microplate surface with moist, lint-free paper.

 Clean membrane carrier with a mild laboratory detergent and rinse thoroughly with ultra-pure water. Do not use abrasive brushes or cleaning materials.

2) Press Eject. Load microplate or membrane carrier. Press Eject again.

 Orient **A1** on the microplate with **A1** on the microplate carrier, and lower the microplate between the guide pins until it rests on the microplate carrier.

 Place **Membranes** face down on the membrane carrier, avoiding air bubbles. Membranes (wet or dry) should be flat and within the boundaries of the carrier. For dry membranes, use a silicon mat (cleaned, rinsed, and dried) to flatten, if necessary. Place **gels** flat within the membrane carrier boundaries in any orientation, avoiding air bubbles. **To prevent spills, avoid excessive liquid in the membrane carrier.**

NOTE: Scanning (left to right) starts at the A1 location on the microplate and proceeds toward the front of the instrument one scan line at a time.



3) Open an existing Work Area or create a new one.

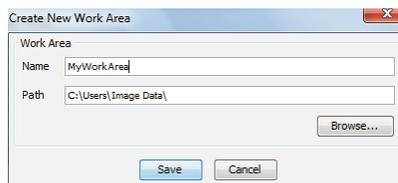
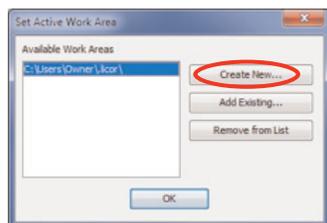
- a. Start Image Studio Software
- b. Choose Odyssey Sa as the acquisition system
- c. Open an existing Work Area or create a new one

1. To open an existing Work Area:

- a. Highlight available Work Area
- b. Click **OK**

2. To create a new Work Area:

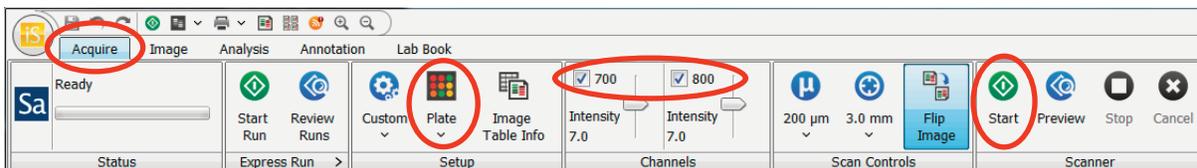
- a. Click **Create New**
- b. Browse to a folder on your hard drive to use as the Work Area
- c. Click on the **New Folder** icon and name it
- d. Click **Save**



- d. Enter User Name and Password. Click **OK** to continue.
- e. From the Setup Group, choose either:
 - i. Plate 200 μm
 1. Intensity: 7.0*
 2. Resolution: 200 μm
 3. Focus offset: 3.0 mm
 - ii. Membrane
 1. Intensity: 7.0*
 2. Resolution: 200 μm
 3. Focus offset: 3.0 mm
 - iii. Gels (Select **Membrane** and adjust Focus Offset as needed)
 1. Intensity: 7.0*
 2. Resolution: 200 μm
 3. Focus Offset: 3.0 mm + half the thickness of the gel (3.95 mm max)

* The recommended Intensity setting (in the Channels group) is 7.0. If a higher value is selected, the overall image will be brighter; however, if the Intensity is set too high, saturation may occur. Saturation is seen as a color change on the image.

- f. Select either:
 - i. **Plate, Western, or No Analysis** from the middle icon in the Setup group



- g. Click **Start** in the Scanner group
- h. Once scan has completed, press **Eject** and remove microplate or membrane carrier