

Product Description

If your Odyssey Imager has an imaging drawer, place a sample you want to image on an Odyssey Imaging Tray and place the Tray in the imaging drawer. This pack (P/N 926-40020) has five trays.

Directions For Use

1. Use trays logo side up.
2. Lay sample flat on tray and within marks at each corner.
3. Place tray in instrument drawer. Notches in tray will align with tabs in drawer assembly.
4. Close drawer and proceed with imaging procedure.

Clean Tray

Trays may be discarded after use. Trays should be discarded if they are damaged or if an unacceptable level of residual signal remains after cleaning.

The following procedure may allow you to clean the tray sufficiently for reuse. If you choose to reuse the tray, it should be cleaned before and after each use.

1. Remove sample from tray.
Always handle membranes with forceps.
2. Clean tray with water and wipe with a lint-free wipe.
Warning: Do not use blue wipes.
3. Rinse with 70% ethanol or 70% isopropanol and lint-free wipes. Methanol may be used if necessary.
Warning: Do not use acetone or any other harsh solvents to clean the imaging tray.

Tray Disposal



Trays may be disposed with plastics, using "Other" classification.

LI-COR®

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

LI-COR Biosciences
4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com