

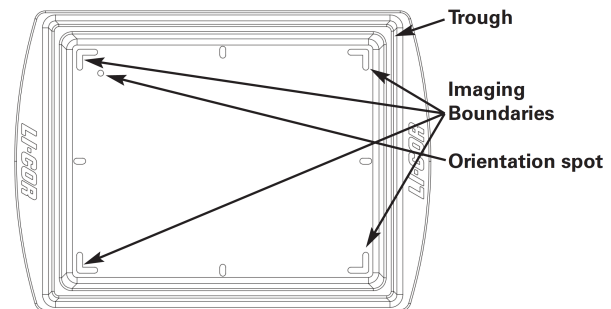
Organ Imaging Tray Description

5 pack (P/N 9300-836) or 25 pack (P/N 9300-837)

The Organ Imaging Tray is a flat polystyrene tray designed to be used on the Pearl Organ Tray Base (PN 9300-23). The Organ Imaging Tray provides a low near-infrared fluorescent background and a flat surface ideal for imaging excised organs. The trough will catch liquid and debris. Replace the Organ Imaging Tray frequently to ensure you have a low background imaging surface.

Instructions

1. Place the Organ Imaging Tray on the Organ Tray Base with the LI-COR logo facing up.
2. Place organ samples within the imaging boundary marks located in the four corners of the Organ Imaging Tray. The **imaging boundary marks** show the boundaries of the imaging area but are not present in the image. The **orientation spot** appears within the imaging area.
3. Place the Organ Imaging Tray and Base in the Pearl Imager according to instructions in the Pearl Imager Manual.
4. Close the drawer and acquire an image.



The top of the organ tray.

Cleaning

Clean the Organ Imaging Tray thoroughly using the Dye Decontamination Kit (P/N 9300-840) or discard and use a new Tray to prevent background contamination of the image from dye or dye-labeled debris.

LI-COR®

© 2019 LI-COR, Inc. LI-COR and Pearl 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