

Organ Tray Base Description (P/N 9300-23)

The Organ Tray Base is installed on the Pearl Imager similarly to the Pearl Imaging Bed but does not provide heat or anesthetic gas. Organs are placed on a disposable Organ Imaging Tray (P/N 9300-836) that fits on the Organ Tray Base.

Note: *The Organ Tray Base should not be used for imaging without an Organ Imaging Tray due to optical focusing issues and potential dye contamination.*



Image showing the Organ Tray Base with Organ Tray.

Installation

See the Pearl Imager Manual.

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.



© 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