

Pearl Imager Cart

This stainless steel cart for the Pearl Imager turns your optical imaging lab into a mobile workstation. The cart has a sturdy design with plenty of room to accommodate a wide variety of accessories, including:

- SmartFlow Anesthesia Suite
- Pearl Docking Station
- Computer
- Oxygen Tanks
- And more!



Install Rubber Shoes

The imaging cart includes new feet for your Pearl Imager.

1. Carefully remove each foot of the Pearl Imager by turning the foot counterclockwise (as you are looking at it).
2. Install the new feet on the bottom of the Pearl Imager by turning each foot clockwise (as you look at it).
3. Adjust feet appropriately to rest stably on the imaging cart.



© 2022 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