

Pearl Imager® Compound Injection Clip

Components

- Package includes 25 Compound Injection Clips (P/N 9300-850)
- Package includes 100 Compound Injection Clips (P/N 9300-851)

Required Accessories (Not Included)

- Syringes with ODs: 0.22 – 0.3"
- Catheter with length at least 28 cm to allow clearance for the drawer to close

Description

- Compound Injection Clip accommodates syringes with ODs: 0.22 – 0.3"
- Compound Injection Clip attaches to the Pearl Imaging Bed for easy transfer from work area to instrument
- Compound Injection Clip attaches to the outside of the Pearl Imager drawer, allowing the drawer to be closed and imaging begun before administering the optical agent

Instructions

1. Insert the syringe in the Compound Injection Clip as shown (Figure 1).

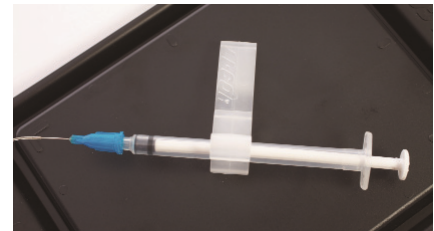


Figure 1.

2. Snap the Compound Injection Clip onto the outside edge of the Imaging Bed (Figure 2) while adjustments to the animal are made.

To protect the imaging bed from unwanted probe contamination, place a black drape (provided with Pearl Accessories) under the mouse's hindquarters, tail, and catheter/syringe assembly (Figure 2).



Figure 2.

3. Move the imaging bed to the Pearl Imager and secure the lever to initiate the flow of Isoflurane and heat to the Imaging Bed (Figure 3).



Figure 3.

4. Remove the Compound Injection Clip from the Imaging Bed and attach it to the front of the imaging drawer close to the catheterization point as shown (Figure 4).

Note: To avoid crimping the catheter when the drawer is closed, position 90% of the catheter's length inside the drawer front as shown (Figure 5).



Figure 4.

5. Start the desired image series using Image Studio™ Software. Once images are being captured, dispense the optical imaging agent quickly (within 5 seconds). If injection is too slow, the distribution effect will not be seen.

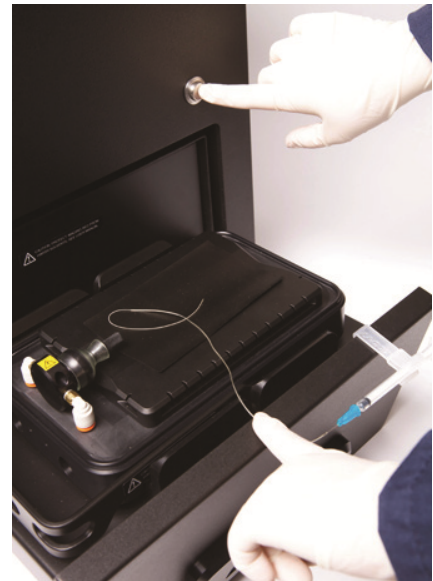


Figure 5.

LI-COR®

© 2020 LI-COR, Inc. LI-COR, Pearl, and Image Studio 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