

Kit Components

P/N	Description
929-95100	Western Blot Storage Bags; 40 bags per pack

Specifications

- Storage: Room Temperature
- Impervious to light

Product Description

Western Blot Storage Bags are a convenient way to save, store, and organize Western blot membranes. The foil bags are lightproof and can accommodate standard size minigel blots (7 x 8cm) up to Criterion™ gel blots (8.5 x 13.5cm). Safeguarding blots in Western Blot Storage Bags prevents scratches and protects from photobleaching for long term storage.

Applications

Western Blot Storage Bags allow you to archive dry membranes.

- Dry fluorescent blots can be re-imaged.
- Fluorescent blots or chemiluminescent blots can be rehydrated and re-probed, when appropriate.

Note: See the Protocol section for specific instructions on how to archive blots you intend to re-probe.

Protocol

Warning: Blots that were stripped and re-probed cannot be used to generate quantitative results.

- To re-image chemiluminescent blots that were archived without first being stripped, you must add new HRP-labeled secondary antibody and new substrate. This process is likely to result in poor quality images.
- PVDF membranes will degrade over time, due to their chemical composition. This degradation will appear as background signal in the 700 nm channel around the edges of the blot.

Prepare Membrane for Archiving or Rescanning Only

1. Ensure that the membrane is completely dry.
2. Place the membrane between 2 pieces of standard, lab grade filter paper.
3. Place the membrane and filter paper in the bag and seal the bag, while removing air trapped in the bag.
4. Annotate the outside of the bag with appropriate experimental information using a suitable felt tip pen.
5. Store at room temperature in a dry place.

Prepare Membrane for Archiving and Re-probing

1. While the membrane is still wet, strip the membrane per the manufacturer's instructions.
2. Check for complete stripping.
3. Rinse the membrane with water or PBS/TBS.
4. Let the membrane dry completely.
5. Place the membrane between 2 pieces of standard, lab grade filter paper.
6. Place the membrane and filter paper in the bag and seal the bag, while removing air trapped in the bag.
7. Annotate the outside of the bag with appropriate experimental information using a suitable felt tip pen.
8. Store at room temperature in a dry place.

Re-probe Membrane

1. If using nitrocellulose, rehydrate the membrane in PBS/TBS for 5 minutes.
If using PVDF, briefly rehydrate in methanol before moving the membrane to PBS/TBS for 5 minutes.
2. Start the blocking step and continue through your standard Western blot procedure.

Limitation of Liability and Limited Use Label License

For Research Use Only. Not intended for human therapeutic or diagnostic use.

By using this product, you agree to the Limitation of Liability and the Limited Use Label License available online at licor.com/packinserts under the **Pack Inserts** heading (the "License"). The License is incorporated herein by this reference, and the License may be updated from time to time.



© 2020 LI-COR, Inc. LI-COR is a registered trademark 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