

Components

P/N	Description
927-50100	125 mL
927-50000	500 mL
927-50003	3 x 500 mL
927-50010	10 x 500 mL

Specifications

- Odyssey Blocking Buffer (TBS) is a ready-to-use formulation, no dilution is required.
- Shelf Life: See expiration date on bottle.
- Storage: 4 °C

***Note:** Odyssey Blocking Buffer (TBS) contains 0.1% sodium azide and is not suitable for chemiluminescent detection.*

Product Description

Blocking buffers enhance the quality of Western blots by reducing background interference, increasing the signal-to-noise ratio, and promoting specific binding of the primary antibody while minimizing non-specific interactions. Odyssey Blocking Buffer (TBS) is a ready-to-use formulation in phosphate-buffered saline that provides optimal blocking conditions for antibodies requiring TBS-based buffer systems. The blocker does not contain mammalian proteins. It is suitable for a variety of applications, including In-Cell™ Western assays, protein arrays, ELISA, and transcription factor assays (EMSA).

Materials Required for Blocking Membrane

- Blotted PVDF or nitrocellulose membrane
- Western blot incubation box
- Orbital shaker for incubation at room temperature

Guidelines

If you're using Odyssey Blocking Buffer (TBS), use TBS-based buffers for the entire detection process and use TBS-based wash buffer.

Protocol

For a detailed Western blot protocol, see the *Near-Infrared (NIR) Western Blot Detection Protocol* (licor.com/wbanalysis). For troubleshooting ideas, see *Good Westerns Gone Bad: Tips to Make Your NIR Western Blot Great* (licor.com/GWGBIR).

Blocking Membrane

1. Place the PVDF or nitrocellulose membrane into the Western Blot Incubation Box and add undiluted Intercept Blocking Buffer to the membrane. Be sure to use sufficient blocking buffer to cover the membrane (a minimum of 0.4 mL/cm² is suggested).

Note: Do not use detergents like Tween® 20 or SDS during the blocking step, as this may generate a background signal. Detergents should only be added when diluting the primary and secondary antibodies.

2. Incubate for 1 hour at room temperature on an orbital shaker.
3. Proceed with Western blot detection protocol.

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, In-Cell Western, and Odyssey 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