

Intercept (TBS) Blocking Buffer

P/N	Description
927-60000	125 mL
927-60001	500 mL
927-60003	3 x 500 mL
927-60010	10 x 500 mL
927-60025	25 x 500 mL
927-60050	50 x 500 mL

Intercept (PBS) Blocking Buffer

P/N	Description
927-70000	125 mL
927-70001	500 mL
927-70003	3 x 500 mL
927-70010	10 x 500 mL
927-70025	25 x 500 mL
927-70050	50 x 500 mL

Specifications (Intercept Blocking Buffer)

Note: *Shake well before each use.*

- Intercept Blocking Buffer is a ready-to-use formulation. For optimal performance, do not dilute.
- Shelf Life: See expiration date on bottle.
- Storage: 4 °C
- Suitable for a variety of applications, including: quantitative Western blots, chemiluminescent Western blots, In-Cell Western™ Assay, In-Gel Western, protein array, and glycoprotein detection.

Product Description (Intercept Blocking Buffer)

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. Intercept Blocking Buffer is available in ready-to-use formulations in tris-buffered saline or phosphate-buffered saline to provide optimal blocking conditions for antibodies requiring TBS-based or PBS-based buffer systems.

The blocker does not contain mammalian proteins.

The TBS blocking buffer is ideal for use with antibodies against phosphorylated protein targets.

Materials Required for Blocking Membrane

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

Guidelines

Use the same buffer system, PBS or TBS, for the entire Western blot process, including in wash solution.

Intercept T20 Antibody Diluents contain Tween® 20 and are available in TBS and PBS formulations, so you can use the appropriate buffer for your antibody diluent, without having to spend time measuring and adding Tween 20. When using a PVDF membrane, your secondary antibody diluent must contain a final concentration of 0.01 – 0.02% SDS.

Protocol

Note: Shake well before each use.

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.

Related Products

See licor.com/intercept for ordering information.

Blocking Buffer

- Intercept (PBS) Protein-Free Blocking Buffer
- Intercept (TBS) Protein-Free Blocking Buffer

Antibody Diluent

- Intercept T20 (PBS) Antibody Diluent
- Intercept T20 (TBS) Antibody Diluent

- Intercept T20 (PBS) Protein-Free Antibody Diluent
- Intercept T20 (TBS) Protein-Free Antibody Diluent

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 Intercept 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