

## Intercept (TBS) Protein-Free Blocking Buffer

P/N	Description
927-80000	125 mL
927-80001	500 mL
927-80003	3 x 500 mL
927-80010	10 x 500 mL

## Intercept (PBS) Protein-Free Blocking Buffer

P/N	Description
927-90000	125 mL
927-90001	500 mL
927-90003	3 x 500 mL
927-90010	10 x 500 mL

## Specifications



Intercept Protein-Free Blocking Buffers now use an eco-friendly ingredient. While this ingredient changes the appearance of the blocking buffer, the blocking buffer still performs the same. Check the certificate of analysis to see if your buffer uses this ingredient.

### **Note: Shake well before each use.**

- Intercept Protein-Free Blocking Buffers are ready-to-use. 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.
- Compatible with avidin-biotin detection systems. Unlike casein or milk-based blocking buffers, Intercept Protein-Free Blocking Buffer will not interfere with avidin-biotin detection systems.

## 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. Intercept Protein-Free Blocking Buffers are ready-to-use formulations in phosphate-buffered saline (PBS) or tris-buffered saline (TBS) to provide optimal blocking conditions for PBS or TBS systems.

Intercept Protein-Free Blocking Buffer is an excellent alternative when protein-based blocking buffers are contributing to cross-reactivity.

## 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 (PBS or TBS) throughout the entire Western blot process, including in blocking buffer, antibody diluent, and wash solutions.

Intercept T20 Protein-Free 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.

## 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](http://licor.com/wbanalysis)). For troubleshooting ideas, see *Good Westerns Gone Bad: Tips to Make Your NIR Western Blot Great* ([licor.com/GWGBIR](http://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/cm<sup>2</sup> 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.

## Creating Antibody Diluent

- Intercept T20 Protein-Free Antibody Diluent contains Tween® 20, so you don't have to spend time measuring and adding Tween 20. PBS and TBS formulations of Intercept T20 Protein-Free Antibody Diluent are available.
- If you're using a PVDF membrane, be sure to add SDS to a final concentration of 0.01 - 0.02% in Intercept Protein-Free Blocking Buffer or Intercept Protein-Free Antibody Diluent when creating secondary antibody diluent.

See the *Near-Infrared Western Blot Protocol* ([licor.com/wbanalysis](http://licor.com/wbanalysis)) for more detail.

## Related Products

See [licor.com/intercept](http://licor.com/intercept) for ordering information.

## **Blocking Buffer**

- Intercept (PBS) Blocking Buffer
- Intercept (TBS) 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](https://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.



© 2022 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](mailto:biosales@licor.com)  
[licor.com/bio](https://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](mailto: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](mailto:bio-eu@licor.com)