

## **Intercept (PBS) Blocking Buffer and Antibody Diluent (927-76003)**

<b>P/N</b>	<b>Description</b>
<b>927-70001</b>	1 Intercept (PBS) Blocking Buffer, 500 mL
<b>927-75001</b>	2 Intercept T20 (PBS) Antibody Diluent , 500 mL

## **Intercept (TBS) Blocking Buffer and Antibody Diluent (927-66003)**

<b>P/N</b>	<b>Description</b>
<b>927-60001</b>	1 Intercept (TBS) Blocking Buffer, 500 mL
<b>927-65001</b>	2 Intercept T20 (TBS) Antibody Diluent , 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.

### **Specifications (Intercept T20 Antibody Diluent)**

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

- Shelf Life: See expiration date on bottle.
- Storage: 4 °C
- Contains a 0.05% concentration of Tween® 20.

**Note:** Because of the high purity and stability of the Tween 20 used in Intercept T20 Antibody Diluent, the concentration of 0.05% is sufficient. When using standard research grade Tween 20, use a concentration of 0.2%. See the NIR Western Blot Protocol ([licor.com/wbanalysis](http://licor.com/wbanalysis)) for more information.

- 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.

## **Description (Intercept T20 Antibody Diluent)**

Intercept T20 Antibody Diluents improve specificity of primary and secondary antibodies, reducing off-target effects. The diluent contains Intercept (PBS) Blocking Buffer or Intercept (TBS) Blocking Buffer preformulated with Tween® 20. When using a PVDF membrane, your secondary antibody diluent must contain a final concentration of 0.01 – 0.02% SDS.

Since there's no need to mix the diluent yourself, you can save time and reduce steps in your protocol that might lead to variability in your results.

## **Related Products**

See [licor.com/intercept](https://www.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) 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](http://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](mailto:biosales@licor.com)  
[licor.com/bio](http://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)