

Intercept T20 (TBS) Protein-Free Antibody Diluent

P/N	Description
927-85001	500 mL
927-85003	3 x 500 mL
927-85010	10 x 500 mL

Intercept T20 (PBS) Protein-Free Antibody Diluent

P/N	Description
927-95001	500 mL
927-95003	3 x 500 mL
927-95010	10 x 500 mL

Specifications



The Intercept Protein-Free Blocking Buffer in the antibody diluent now uses an eco-friendly ingredient. While this ingredient changes the appearance of the diluent, the diluent still performs the same. Check the certificate of analysis to see if your diluent uses this ingredient.

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 Protein-Free 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) 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.

Description

Intercept T20 Antibody Diluents improve specificity of primary and secondary antibodies, reducing off-target effects. The diluent contains Intercept (PBS) Protein-Free Blocking Buffer or Intercept (TBS) Protein-Free Blocking Buffer preformulated with Tween® 20. This diluent can be used where animal-based products are prohibited.

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.

Intercept T20 Protein-Free Antibody Diluent is ideal for use with Intercept Protein-Free Blocking Buffer.

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](https://www.licor.com/wbanalysis)). For troubleshooting ideas, see *Good Westerns Gone Bad: Tips to Make Your NIR Western Blot Great* ([licor.com/GWGBIR](https://www.licor.com/GWGBIR)).

Preparing Secondary Antibody Diluent

If you're using a PVDF membrane, add SDS to a final concentration of 0.01 - 0.02% in Intercept Protein-Free Antibody Diluent when preparing the secondary antibody diluent.

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

Related Products

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

Blocking Buffer

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

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.



© 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
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