

Intercept T20 (TBS) Antibody Diluent

P/N	Description
927-65001	500 mL
927-65003	3 x 500 mL
927-65010	10 x 500 mL

Intercept T20 (PBS) Antibody Diluent

P/N	Description
927-75001	500 mL
927-75003	3 x 500 mL
927-75010	10 x 500 mL

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) 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 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.

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).

Related Products

See 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) 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