

Part Numbers

927-40300 100 mL

927-40200 500 mL

Description

Casein blocking buffer is a blocking solution designed for Western Blot applications. It also serves as a diluent for antigens, antibodies, and other proteins.

Content

0.5% Casein (Hammarsten Grade) in 1X PBS.

Warning: Contains 0.1% sodium azide; do not use as HRP conjugated antibody diluent.

Appearance

Slightly yellow to colorless liquid.

Storage

Store refrigerated 2-8 °C.

Use

- **Preparation:** No dilution of blocking buffer required.
- **Incubation time:** Block according to standard laboratory procedures.

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 is a registered trademark 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