

Specifications

- Size: 0.5 mg
- Concentration: 1 mg/ml
- Physical State: Liquid
- Buffer: 10 mM Phosphate, 183 mM NaCl and, 2.7 mM KCl, pH 7.4
- Preservative: Sodium azide 0.005% (w/v)

Warning: *Sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.*

- IRDye 800CW Abs_{max} = 774 nm, Emission_{max} = 789 nm
- IRDye 800CW Molecular Weight: 1166 g/mole
- IRDye 800CW Extinction Coefficient = 240,000 M⁻¹ cm⁻¹ (in 1X PBS)
- Storage: Store vial at 4 °C **protected from light**. Expiration date is six months from date received.

Protocol

IRDye 800CW Streptavidin can be used as a secondary detection reagent for microscopy, protein arrays, In-Gel Westerns, and Western blotting. Centrifuge briefly before use to eliminate aggregates that may have formed in solution. This will reduce non-specific background staining.

Recommended Dilutions

A final concentration of 0.2 to 1.0 g/mL (1:1,000 to 1:5,000) is usually satisfactory for most applications. However, appropriate dilution may need to be determined empirically. For membrane-based applications and In-Gel Westerns, it is recommended to add SDS (0.02% to 0.1% final concentration) and Tween® 20 (0.1 to 0.2% final concentration) during the detection incubation step to reduce non-specific background staining.

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 and IRDye 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