

Product Number
925-68031

Quantity: 0.1 mg
Concentration: 1 mg/mL

Lot Number:

Storage: 4 °C

December 2014

Updates available at:

<http://www.licor.com/packinserts>

Limitation of Liability and Limited Use Label License

LI-COR reagent products are offered for research purposes only and are not intended for human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the amount of product purchased and the components of the product in research conducted by the buyer (whether the buyer is a not-for-profit, academic or for-profit entity). The buyer shall not sell or otherwise transfer this product, its components, or materials made there from to any third party. Buyer shall not use this product or its components for commercial purposes. The term "commercial purposes" shall mean any activity by a party for consideration and may include, but is not limited to, use of the product or its components (i) in manufacturing, (ii) for therapeutic, diagnostic or prophylactic purposes, or (iii) for resale, whether or not such product or its components are resold for use in research. The use of this product by the buyer constitutes agreement with the terms of this limited use label license for LI-COR IRDye infrared dyes and reagent products. Inquiries regarding the licensing of one or more IRDye infrared dyes should be submitted by e-mail to busdev@licor.com.

LI-COR BIOSCIENCES DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR IRDYE INFRARED DYES AND REAGENT PRODUCTS. LI-COR BIOSCIENCES MAKES NO WARRANTIES OF ANY KIND, EITHER EXPRESS OR IMPLIED, AS TO ANY MATTER INCLUDING, BUT NOT LIMITED TO, WARRANTY OF FITNESS FOR PURPOSE, OR MERCHANTABILITY OR RESULTS OBTAINED FROM USE OF IRDYE INFRARED DYES. IN NO EVENT SHALL LI-COR BE LIABLE FOR LOST PROFITS, CONSEQUENTIAL, EXEMPLARY, SPECIAL, DIRECT, INCIDENTAL, OR PUNITIVE DAMAGES, OR ATTORNEY FEES, EVEN IF LI-COR HAD BEEN ADVISED OF, KNEW, OR SHOULD HAVE KNOWN, OF THE POSSIBILITIES THEREOF. NO EMPLOYEE, AGENT OR REPRESENTATIVE OF LI-COR HAS THE AUTHORITY TO BIND LI-COR TO ANY ORAL REPRESENTATION OR WARRANTY EXCEPT AS SPECIFICALLY SET FORTH HEREIN.

© 2014 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR, IRDye, and Odyssey are registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners. This product may be covered by US patents or patents pending. Visit www.licor.com/patents for details.

Doc # 988-14910

LI-COR

4647 Superior Street
Lincoln, Nebraska 68504 USA
North America: 800-645-4267
International: 402-467-0700
FAX: 402-467-0819

LI-COR GmbH Germany, Serving Europe, Middle East, and Africa: +49 (0) 6172 17 17 771
LI-COR UK Ltd. UK, Serving Denmark, Finland, Iceland, Ireland, Norway, UK, and Sweden: +44 (0) 1223 422104. All other countries, contact LI-COR Biosciences or a local LI-COR distributor: <http://www.licor.com/distributors>

www.licor.com/bio

IRDye® 680LT Streptavidin

Properties

Physical State: Liquid

Dye/Protein Ratio: moles of IRDye 680LT/mole Streptavidin

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 680LT Abs_{max} = 676 nm, Emission_{max} = 693 nm

IRDye 680LT Molecular Weight: 1402 g/mole

IRDye 680LT Extinction Coefficient = 250,000 M⁻¹ cm⁻¹ (in 1x PBS)

IRDye 680LT is not intended for In-Cell Western™ assays or *in vivo* applications.

Storage Conditions

Protect from light. Store vial at 4 °C. Expiration date is six months from date received.

Instructions

IRDye 680LT Streptavidin can be used as a secondary detection reagent for Western blotting and other applications using biotinylated conjugates. To reduce non-specific background staining due to possible aggregates in the solution, centrifuge briefly before use. A final concentration of 0.2 to 1.0 µg/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, 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.

Recommended Dilutions

Application	Suggested Range	Tween® 20* (v/v)	SDS* (v/v)
Odyssey® Western blot detection	1:1,000 – 1:5,000	0.1 – 0.2%	0.02 – 0.1%
Other	User optimized	User optimized	User optimized

* Added to reduce non-specific background staining.

Optimum dilutions will vary and should be determined empirically.