

Product Number

827-11081

Quantity: 25 µL

Storage: 4°C

Revised: January 2012

Updates available at:

<http://biosupport.licor.com>

Limitation of Liability and Limited Use Label License

LI-COR IRDye® Infrared Dyes and 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 therefrom, 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) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) 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.

© 2012 LI-COR, Inc. LI-COR is an ISO 9001 registered company. IRDye reagents, the Aeries and Odyssey Imaging systems are covered by U.S. and foreign patents and patents pending. LI-COR, IRDye, Odyssey, In-Cell Western, and Aeries are trademarks and registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

Doc #988-12729

LI-COR®

Experience Excellence

4647 Superior Street • P.O. Box 4000
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 UK, Ireland, and Scandinavia: +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 Goat (polyclonal) Anti-Rabbit IgG (H+L), Highly Cross Adsorbed – Sample Pack

Applications

IRDye 680LT secondary antibodies can be used for a variety of applications, including Western blotting. These antibodies are optimized for use with instruments within the Odyssey® and Aeries® family of imagers. Two-color detection can be achieved by multiplexing IRDye 680LT secondary antibodies with IRDye 800CW secondary antibodies. IRDye 680LT secondary antibodies are also suitable for immunofluorescent microscopy and other fluorescent imaging applications when using instrumentation with appropriate excitation and detection capabilities. Visit the LI-COR Biosciences website for more information. For detailed protocols, please reference the Application Protocols Manual included with all instrument installations, or download from <http://biosupport.licor.com>.
IRDye 680LT is not intended for *in vivo* applications.

Properties

Antibody Concentration: 1.0 mg/mL

Fluorophore: IRDye 680LT (MW 1402)

Excitation Wavelength: 679 nm (in PBS)

Emission Wavelength: 696 nm (in PBS)

Form of Antibody: IRDye 680LT purified immunoglobulin conjugate, in phosphate-buffered saline, pH 7.4. Contains 10 mg/mL BSA (free of IgG and protease) as stabilizer and 0.01% sodium azide as preservative.

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

Immunogen: Rabbit IgG

Purity and Specificity

Isolation of specific antibodies was accomplished by affinity chromatography using pooled IgG covalently linked to agarose. Based on ELISA and flow cytometry, this antibody reacts with the heavy and light chains of rabbit IgG, and with the light chains of rabbit IgM and IgA. This antibody was tested by Dot Blot and/or solid-phase adsorbed for minimal cross-reactivity with human, mouse, rat, sheep and chicken serum proteins, but may cross-react with immunoglobulins from other species. The conjugate has been specifically tested and qualified for Western blot.

Storage

Protect from light. Store at 4°C. This product is stable for up to 3 months at 4°C when stored as directed.

Recommended Dilutions

NOTE: When using PVDF membranes for Western blotting applications, SDS (final concentration of 0.01 - 0.02%) and Tween® 20 (final concentration of 0.1 - 0.2%) must be added during the detection incubation step to avoid non-specific background staining.

Application	Recommended
Odyssey Western blot detection	1:20,000
Other	User optimized

Optimum dilutions will vary and should be determined empirically.