

LI-COR® IRDye®

Infrared Imaging Reagents

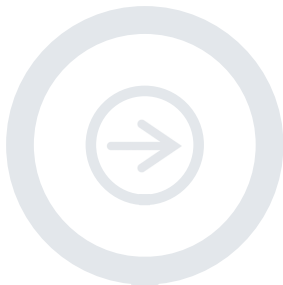
**IRDye 680 Conjugated Goat (polyclonal)
Anti-Mouse IgG, Highly Cross Adsorbed**

Product Number: 926-32220

Lot Number:

Quantity: 0.5 mg

Shipping: RT



LI-COR®

Biosciences

Product: IRDye 680 Conjugated Goat Anti-Mouse IgG (H+L), Cross-adsorbed

Applications

IRDye 680 antibodies can be used for a variety of applications, including Western blotting, In-Cell Western assays, In-Gel Westerns, and many others. Visit the LI-COR® website for more information. For detailed protocols, please reference the Odyssey® and/or Aeries® protocol manuals included with all instrument installations or download them from the LI-COR support site at: <http://biosupport.licor.com>

Two-color detection can be achieved by multiplexing IRDye 680 conjugates with IRDye 800CW or IRDye 800 antibody conjugates. IRDye™ 680 conjugates are also suitable for immunofluorescent microscopy and other fluorescent imaging applications, when using instrumentation with appropriate excitation and detection capabilities.

Antibody Concentration: 1.0 mg/ml when reconstituted as directed

Fluorophore: IRDye 680 (MW 950)

Excitation Wavelength: 683 nm (in PBS)

Emission Wavelength: 710 nm (in PBS)

Fluorophore/Protein Ratio:

_____ moles IRDye 680: 1 mole IgG

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

(Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)

Immunogen: Mouse IgG, whole molecule

Purity and Specificity: Isolation of specific antibodies was accomplished by immunoaffinity chromatography using antigens immobilized on agarose beads. Based on immunoelectrophoresis, this antibody reacts with the heavy chains of mouse IgG as well as the light chains common to most mouse immunoglobulins. No reactivity was detected against non-immunoglobulin serum proteins. This antibody was tested by ELISA and/or solid phase adsorbed for minimal cross-reactivity with human, bovine, horse, rabbit and swine serum proteins, but may cross-react with immunoglobulins from other species. The conjugate has been specifically tested and qualified for Western blot and In Cell Western applications.

Reconstitution and Storage: *Protect from light.* Store at 4°C prior to reconstitution. Reconstitute contents of vial with 0.5 mL sterile distilled water. Mix gently by inverting, and allow to rehydrate for at least 30 minutes before use. Centrifuge product if solution is not completely clear after standing at room temperature. Dilute only immediately prior to use. This product is stable for up to 3 months at 4°C when stored undiluted as directed. For extended storage, aliquot and freeze at -20°C; avoid repeated freeze-thaw cycles.

Continued

Recommended Dilutions:

Odyssey Western blot detection	1:15,000 Recommended (Suggested dilution range: 1:5,000 – 1:25,000)
In-Cell Western detection	1:200 Recommended (Suggested dilution range: 1:200 – 1:800)
Other applications	User optimized

Optimum dilutions will vary and should be determined empirically.

These products are offered for research and development purposes only. They cannot be used to manufacture derivatives or products for re-sale.

IRDye 680 and IRDye Infrared Dyes are covered by U.S. and foreign patents and patents pending. These products are distributed and sold for research purposes only and are not intended for diagnostic or therapeutic use. Purchase does not include or carry any right or license to use, develop or otherwise exploit products for commercial purposes. Inquiries regarding the licensing of one or more IRDye Infrared Dyes should be submitted by e-mail to busdev@licor.com

LIMITED WARRANTY AND LIMITATION OF DAMAGES

LI-COR IRDYE® INFRARED DYES ARE OFFERED FOR RESEARCH PURPOSES ONLY AND THE USE THEREOF REMAINS THE RESPONSIBILITY OF PURCHASER. LI-COR DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR IRDYE™ INFRARED DYES. LI-COR WARRANTS ONLY THAT IRDYE™ INFRARED DYES WILL MEET THE SPECIFICATIONS PROVIDED BY LI-COR. THIS LIMITED WARRANTY PROVIDED IN THIS PARAGRAPH SHALL BE IN LIEU OF ALL OTHER WARRANTIES, EXPRESS OR IMPLIED, INCLUDING ANY IMPLIED WARRANTY OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, NON-INFRINGEMENT AND NON-INTERFERENCE WITH ENJOYMENT. LI-COR'S OBLIGATION TO USER FOR ALL BREACHES OF THIS AGREEMENT, INCLUDING WARRANTIES, WHENEVER ALLEGED OR RAISED, SHALL NOT EXCEED THE TOTAL AMOUNT OF PAYMENTS MADE TO LI-COR BY PURCHASER OF IRDYE® INFRARED DYES DURING THE FIRST YEAR OF THIS AGREEMENT. IN NO EVENT SHALL LI-COR BE LIABLE FOR LOST PROFITS, CONSEQUENTIAL, EXEMPLARY, SPECIAL, DIRECT, INCIDENTAL, OR PUNITIVE DAMAGES, OR ATTORNEY FEES UNDER ANY CIRCUMSTANCES, 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.

LI-COR, Odyssey and IRDye are trademarks or registered trademarks of LI-COR, Inc. All other trademarks belong to their respective owners.



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 and Africa: +49 (0) 6172 17 17 771
LI-COR UK Ltd. UK, Serving UK, Ireland, Scandinavia: +44 (0) 1223 422104
www.licor.com

Doc #988-08357
Rev 1 0106