

**LI-COR® IRDye®**

---

Infrared Imaging Reagents

**IRDye® 680 Donkey Anti-Goat IgG (H+L)  
highly cross adsorbed**

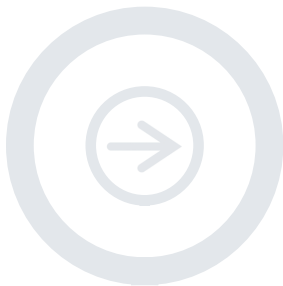
**Product Number:** 926-32224

**Lot Number:** B70820-02

**Quantity:** 0.5 mg

**Shipping:** RT

**Storage:** 4°C



Published September 2007.  
The most recent version of this  
pack insert is posted at  
<http://biosupport.licor.com/support>.

***LI-COR***®  
Biosciences

**Antibody Concentration:** 1.0 mg/ml when reconstituted as directed

**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:** Goat IgG, whole molecule

**Fluorophore:** IRDye 680

**Excitation Wavelength:** 683 nm (in PBS)

**Emission Wavelength:** 710 nm (in PBS)

**Fluorophore/Protein Ratio:** 2.4 moles IRDye 680:1 mole IgG

*Caution: Bovine serum albumin (BSA) and dry milk may contain IgG which reacts with anti-bovine IgG, anti-goat IgG, anti-horse IgG, and anti-sheep IgG antibodies. Therefore, use of BSA and/or dry milk to block or dilute these antibodies and/or your primary antibody may significantly increase background and/or reduce secondary antibody titer.*

**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® Imaging System 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 800CW conjugates with IRDye 680 antibody conjugates. IRDye 800CW conjugates are also suitable for immunofluorescent microscopy and other fluorescent imaging applications, when using instrumentation with appropriate excitation and detection capabilities.

**Purity and Specificity:** The antibody was isolated from antisera by immunoaffinity chromatography using antigens coupled to agarose beads. Based on immunoelectrophoresis, the antibody reacts with the heavy chains on goat IgG and with the light chains common to most goat immunoglobulins. No antibody was detected against non-immunoglobulin serum proteins. This antibody has been tested by ELISA and/or solid-phase adsorbed to ensure minimal cross-reaction with chicken, guinea pig, Syrian hamster, horse, human, mouse, rabbit, and rat serum proteins, but the antibody may cross-react with immunoglobulins from other species. The conjugate has been specifically tested and qualified for Western blot 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 or below; avoid repeated freeze-thaw cycles.

**Recommended Dilutions:** Odyssey Western blot detection: 1:15,000 recommended (suggested dilution range: 1:5,000 – 1:25,000). Other applications: User optimized. *Optimum dilutions will vary and should be determined empirically.*

**Note:** For *in vitro* research use only; not for diagnostic or therapeutic use. This product is not a medical device.

### **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. Buyer shall not determine the structure or otherwise reverse engineer this product. 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 DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR IRDYE® INFRARED DYES AND REAGENT PRODUCTS. LI-COR 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.

IRDye 800CW, IRDye 700DX, IRDye 800PA, IRDye 700PA and IRDye infrared dyes are covered by U.S. and foreign patents and patents pending. LI-COR, IRDye, Odyssey, Aeries, and In-Cell Western are trademarks or registered trademarks of LI-COR, Inc.

---

**LI-COR**®

Biosciences

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:\* +49 (0) 6172 17 17 771 • LI-COR UK Ltd.: +44 (0) 1223 422104

[www.licor.com](http://www.licor.com)

\*Serving Austria, Czech Republic, Germany, Hungary, Slovakia, and Switzerland

Doc # 988-09392

0907