

Product Number

**926-54110**

Quantity: 25 µL

Storage: 4 °C

July 2013

Updates available at:

<http://biosupport.licor.com>

**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) 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 reagent products. LI-COR BIOSCIENCES DOES NOT PROVIDE RESEARCH ADVICE OR DETERMINE OR RECOMMEND ANY POTENTIAL USES FOR 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 REAGENT PRODUCTS. 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.

© 2013 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR and VRDye are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. For patent information, visit [www.licor.com/patents](http://www.licor.com/patents)

Doc #988-13598

**LI-COR**

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

# VRDye™ 549 Goat (polyclonal) anti-Mouse IgG (H+L), Highly Cross-Adsorbed – Sample Pack

## Applications

VRDye 549 secondary antibodies can be used for a variety of applications, including immunofluorescence microscopy, flow cytometry, and others. Visit the LI-COR® website (<http://www.licor.com/bio>) for more information.

VRDye 549 Goat anti-Mouse IgG (H+L) is suitable for multiplex labeling experiments when combined with other secondary antibodies labeled with proper fluorescent dyes, and using instrumentation with appropriate excitation and detection capabilities.

## Properties

**Antibody Concentration:** 1.0 mg/mL

**Fluorophore:** VRDye 549 (MW 1040)

**Excitation Wavelength:** 555 nm (in PBS)

**Emission Wavelength:** 569 nm (in PBS)

**Form of Antibody:** VRDye 549 purified immunoglobulin conjugate 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.

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

**Immunogen:** Mouse IgG paraproteins

## Purity and Specificity

Isolation of specific antibodies was accomplished by affinity chromatography using pooled mouse IgG covalently linked to agarose. Based on ELISA and flow cytometry, this antibody reacts with the heavy and light chains of mouse IgG<sub>1</sub>, IgG<sub>2a</sub>, IgG<sub>2b</sub>, and IgG<sub>3</sub>, and with the light chains of mouse IgM and IgA. This antibody was tested by dot blot and/or solid-phase adsorbed for minimal cross-reactivity with human, rabbit, goat, rat, and horse serum proteins, but may cross-react with immunoglobulins from other species. The conjugate has been specifically tested and qualified for immunofluorescence microscopy and flow cytometry.

## Storage

**Protect from light.** Store at 4 °C. This product is stable for up to 3 months at 4 °C.

## Recommended Dilutions

Application	Recommended	Suggested Range
Immunofluorescence Microscopy	1:400	1:100 - 1:1,000
Flow Cytometry	1:1,000	1:200 - 1:2,000
Other	User-optimized	

*Suggested working dilutions are given as a guide only. Optimum dilutions will vary and should be determined empirically.*