

Applications

VRDye 490 secondary antibodies can be used for a variety of applications, including Western blotting. Visit the LI-COR® website (licor.com/bio) for more information.

VRDye 490 Donkey anti-Goat IgG (H+L) (**P/N 926-49014**) is suitable for multiplex labeling experiments when combined with other secondary antibodies labeled with proper fluorescent dyes and when using instrumentation with appropriate excitation and detection capabilities. For example, two- or three-color detection can be achieved by multiplexing VRDye 490 secondary antibodies with IRDye® 680RD or IRDye 680LT and IRDye 800CW secondary antibodies for detection on the Odyssey® M Imager.

Specifications

- Quantity: 0.5 mg, lyophilized
- Antibody Concentration: 1.0 mg/mL when reconstituted as directed
- Fluorophore: VRDye 490 (1011 g/mole)
- Excitation Wavelength: 492 nm (in PBS)
- Emission Wavelength: 522 nm (in PBS)
- Immunogen: Goat IgG, whole molecule
- Form of Antibody: VRDye 490 conjugated purified immunoglobulin, lyophilized in phosphate buffered saline, pH 7.4. Contains 10 mg/mL Bovine serum albumin (BSA) (IgG and protease free) as stabilizer and 0.01% sodium azide as preservative after reconstitution.

Warning: *Contains sodium azide.*

Warning: *The BSA used as a stabilizer for this product is free of IgG. However, blockers and antibody diluents made with BSA and dry milk may contain IgG that reacts with anti-bovine IgG, anti-goat IgG, anti-horse IgG, and anti-sheep IgG antibodies. This can lead to a significant increase in background and/or reduction of secondary antibody titer for protein detection applications.*

Storage and Reconstitution

Lyophilized

- Storage temperature: 4 °C
- Storage conditions: Protect from light.
- Shelf life: This product is stable for six months when stored as recommended.

After Reconstitution

- Storage temperature: 4 °C
- Storage conditions: Protect from light.
- Shelf life: Expiration date is listed on the product's label.

When ready for use, prepare a 1.0 mg/mL stock solution by reconstituting contents of vial with 0.5 mL sterile distilled water. Mix gently by inverting. Allow to rehydrate for at least 30 minutes before use. Centrifuge product if solution is not completely transparent after standing at room temperature. Immediately dilute stock solution as recommended for each specific application prior to use. Do not re-use product.

Suggested Dilution Ranges

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

Application	Recommended	Suggested Range
Western blot detection	1:10,000	1:5,000 – 1:15,000
Other	User Optimized	

Purity and Specificity

The antibody was isolated from antisera by immunoaffinity chromatography using antigens coupled to agarose beads. Based on ELISA, this antibody reacts with heavy and light chains of goat IgG and sheep IgG. This antibody was tested by dot blot and/or solid phase adsorbed for minimal cross-reactivity with human, mouse, rabbit, rat, chicken, guinea pig, hamster, swine, and horse serum proteins but may cross-react with immunoglobulins from other species. The conjugate has been specifically tested and qualified for Western blot applications.

Limitation of Liability and Limited Use Label License

For Research Use Only. Not intended for human therapeutic or diagnostic use.

By using this product, you agree to the Limitation of Liability and the Limited Use Label License available online at licor.com/packinserts under the **Pack Inserts** heading (the "License"). The License is incorporated herein by this reference, and the License may be updated from time to time.



© 2021 LI-COR, Inc. LI-COR, Odyssey, IRDye, and VRDye are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com