

926-32350Highly Cross-Adsorbed Goat Anti-Mouse IgG₁ (γ₁ Chain Specific) Antibody
Conjugated to IRDye 800CW

Applications

IRDye 800CW secondary antibodies can be used for numerous applications, including: Western blotting, In-Cell Western™ Assay, In-Gel Western, and many others.

Two-color detection can be achieved by multiplexing IRDye 800CW secondary antibodies with IRDye 680RD or IRDye 680LT secondary antibodies. IRDye 800CW secondary antibodies are also suitable for immunofluorescent microscopy and other fluorescent imaging applications when using instrumentation with appropriate excitation and detection capabilities. Visit licor.com/support for protocols, example experiments, recommendations, and more technical information.

Storage

- Storage temperature: 4 °C
- Storage conditions: Protect from light and moisture.
- Shelf life: When stored as recommended, this product is stable for 3 months.

Specifications

- Fluorophore: IRDye 800CW (MW 1166)
- Size: 0.5 mg
- Antibody Concentration: 1.0 mg/mL (when reconstituted as directed)
- Form of Antibody: IRDye 800CW purified immunoglobulin conjugate, 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, after reconstitution.

Warning: *Contains sodium azide.*

- Excitation Wavelength: 778 nm (in PBS)
- Emission Wavelength: 795 nm (in PBS)
- Immunogen: Mouse IgG₁ paraproteins

Purity and Specificity

Isolation of specific antibodies was accomplished by affinity chromatography on mouse IgG₁ covalently linked to agarose. Based on ELISA and flow cytometry, this antibody reacts with the heavy chain of mouse IgG₁. This antibody was tested by Dot Blot and/or solid phase adsorbed for minimal cross-reactivity with mouse IgM, IgG_{2a}, IgG_{2b}, IgG₃ and IgA, pooled human sera, and purified human paraproteins. The conjugate has been specifically tested and qualified for Western blot and In-Cell Western applications. Results with your primary antibody may vary and specificity should be confirmed prior to performing two-color detection.

Reconstitute Secondary Antibody

1. Combine the contents of each vial of secondary antibody with 0.5 mL of sterile distilled water.
2. Gently mix by inverting.
3. Allow the solution to stand at room temperature for at least 30 minutes before use.

Note: If the solution is not completely transparent after standing at room temperature, centrifuge the solution.

When stored as recommended, this product is stable for 3 months.

Recommended Dilutions

Optimum dilutions will vary and should be determined empirically.

Avoid non-specific background staining on PVDF membranes: If you are making your own antibody diluent, add SDS (final concentration 0.01 – 0.02%) and Tween® 20 (final concentration 0.2%) to the diluent. If using Intercept® T20 Antibody Diluent, only add SDS (final concentration 0.01 – 0.02%).

Recommended Dilutions for Western blotting

Recommended	Suggested Range
1:20,000	1:5,000 – 1:25,000

Recommended Dilutions for the In-Cell Western™ Assay

Recommended	Suggested Range
1:800	1:200 – 1:1,200

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, In-Cell Western, Intercept, and IRDye 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
licor.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