

925-32218

Highly Cross-Adsorbed Donkey Anti-Chicken IgG (H+L) 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: 0.1 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: Chicken IgY, whole molecule. IgY is the original designation for the IgG-like protein found in both serum and egg yolk

Purity and Specificity

The antibody was isolated from antisera by immunoaffinity chromatography using antigens coupled to agarose beads. Based on immunoelectrophoresis and/or ELISA, the antibody reacts with whole molecule chicken IgY. It also reacts with the light chains of other chicken 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 bovine, goat, guinea pig, Syrian hamster, horse, mouse, human, rabbit, rat, and sheep 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.

Reconstitute Secondary Antibody

1. Combine each 0.1 mg of secondary antibody with 0.1 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. Protect the solution from light.

Note: *If the solution is not completely dissolved 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.



© 2022 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