

**926-32411**Highly Cross-Adsorbed Donkey Anti-Guinea Pig 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](http://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: Guinea Pig IgG, whole molecule

## Purity and Specificity

Isolation of specific antibodies was accomplished by immunoaffinity chromatography using antigens immobilized on agarose beads. Based on immunoelectrophoresis and/or ELISA, this antibody reacts with whole molecule guinea pig IgG as well as the light chains common to other guinea pig immunoglobulins. No reactivity was detected against non-immunoglobulin serum proteins. This antibody was tested by ELISA and/or solid phase adsorbed for minimal cross-reactivity with bovine, chicken, goat, Syrian hamster, horse, human, mouse, rabbit, rat, and sheep serum proteins, but may cross-react with immunoglobulins from other species. The conjugate has been specifically tested and qualified for Western blot and In-Cell Western applications.

## 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](http://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