

Goat Anti-Mouse IgG (H+L), Highly Cross-Adsorbed (Min X Rat)



Highly cross-adsorbed goat anti-mouse IgG (H L) secondary antibody labeled with our superior CF® Dyes.

Product Description

This is a highly cross-adsorbed goat anti-mouse IgG (H L) secondary antibody labeled with our bright and photostable CF® Dyes and other labels. To minimize cross-reactivity, the antibody has been adsorbed against bovine, chicken, goat, guinea pig, Syrian hamster, horse, human, rabbit, rat and sheep serum proteins.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 10 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology

Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. View our full selection of bright and specific [Secondary Antibodies](#), or search our catalog using our [Antibody Finder](#). Alternatively, you can view our [secondary antibody product listings](#) with catalog numbers. CF® Dyes offer exceptional brightness and photostability. For more information see our [CF® Dye technology page](#).

Storage and Handling Liquid format: Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Liquid format antibodies contain 50% glycerol and will not freeze at -20°C. Lyophilized format: Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Reconstitute antibodies in water using the indicated volumes below: CF® Dye and biotin conjugates: add 0.5 mL dH₂O HRP or DNP conjugates: add 1 mL dH₂O Add the indicated volume of water directly to the vial containing the lyophilized antibody and mix gently to dissolve. Store reconstituted antibody at -20°C and protect from light. Aliquot to avoid repeated freeze/thaw cycles. Alternatively, an equal volume of glycerol can be mixed with the reconstituted antibody so that it will remain liquid at -20°C. Optional: A preservative such as 0.05% sodium azide (final concentration) can be added to CF® Dye and biotin conjugates. Do not add sodium azide to HRP conjugates. **Note:** Storage of the antibody for more than a day at final working dilution is not recommended. CF is a registered trademark of Biotium, Inc.

Product attributes

| | |
|-------------------------------------|---|
| Clonality | Polyclonal |
| Antibody type | Secondary |
| Concentration | 2 mg/mL |
| Antibody/conjugate formulation | Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution |
| Species reactivity | Mouse |
| Secondary/tag antibody applications | Flow cytometry, IHC, IF (cells or tissue sections), Western blot |
| Host species | Goat |
| Antibody reactivity (target) | Mouse IgG |
| Cross adsorption | Bovine, Chicken, Goat, Guinea pig, Horse, Human, Rabbit, Rat, Sheep, Syrian hamster |

References

Download a list of [CF® dye references](#).

| Conjugation | Ex/Em | Size | Catalog No. | Dye Features |
|-------------|------------|--------------------|-----------------------------|----------------------------------|
| CF®405S | 404/431 nm | 50 uL (100 ug) | 20830-50uL | CF®405S Features |
| | | 0.5 mL (1 mg) | 20830-500uL | |
| | | 1 mg (lyophilized) | 20830-1mg | |
| CF®405M | 408/452 nm | 50 uL (100 ug) | 20340-50uL | CF®405M Features |
| | | 0.5 mL (1 mg) | 20340-500uL | |
| | | 1 mg (lyophilized) | 20340-1mg | |
| CF®488A | 490/515 nm | 50 uL (100 ug) | 20302-1 | CF®488A Features |
| | | 0.5 mL (1 mg) | 20302 | |
| | | 1 mg (lyophilized) | 20302-1mg | |
| CF®543 | 541/560 nm | 50 uL (100 ug) | 20328-1 | CF®543 Features |
| | | 0.5 mL (1 mg) | 20328 | |
| | | 1 mg (lyophilized) | 20328-1mg | |
| CF®568 | 562/583 nm | 50 uL (100 ug) | 20301-1 | CF®568 Features |
| | | 0.5 mL (1 mg) | 20301 | |
| | | 1 mg (lyophilized) | 20301-1mg | |
| CF®583R | 585/609 nm | 50 uL (100 ug) | 20903-50uL | CF®583R Features |
| | | 0.5 mL (1 mg) | 20903-500uL | |
| | | 1 mg (lyophilized) | 20903-1mg | |
| CF®594 | 593/614 nm | 50 uL (100 ug) | 20303-1 | CF®594 Features |
| | | 0.5 mL (1 mg) | 20303 | |
| | | 1 mg (lyophilized) | 20303-1mg | |
| CF®633 | 630/650 nm | 50 uL (100 ug) | 20341-50uL | CF®633 Features |
| | | 0.5 mL (1 mg) | 20341-500uL | |
| | | 1 mg (lyophilized) | 20341-1mg | |
| CF®640R | 642/662 nm | 50 uL (100 ug) | 20304-1 | CF®640R Features |
| | | 0.5 mL (1 mg) | 20304 | |
| | | 1 mg (lyophilized) | 20304-1mg | |
| CF®660C | 667/685 nm | 50 uL (100 ug) | 20368-50uL | CF®660C Features |
| | | 0.5 mL (1 mg) | 20368-500uL | |
| | | 1 mg (lyophilized) | 20368-1mg | |

To minimize crossreactivity, this antibody has been adsorbed against bovine, chicken, goat, guinea pig, hamster, horse, human, rabbit, rat and sheep serum.

Product link: <https://biotium.com/product/goat-anti-mouse-igg-hl-highly-cross-adsorbed-min-x-rat/>

