



Glowing products for science

## 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.

Call us : [800-304-5357](tel:800-304-5357) Email: [techsupport@biotium.com](mailto:techsupport@biotium.com)

### 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
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Host species	Goat
Antibody reactivity (target)	Mouse IgG
Cross adsorption	Bovine, Chicken, Goat, Guinea pig, Horse, Human, Rabbit, Rat, Sheep, Syrian hamster

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

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

To minimize crossreactivity, this antibody has been adsorbed against bovine, chicken, goat, guinea pig, hamster, horse, human, rabbit, rat and sheep serum. 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 dH<sub>2</sub>O

HRP or DNP conjugates: add 1 mL dH<sub>2</sub>O

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.

## References

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

This datasheet was generated on July 12, 2025 at 07:23:52 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/goat-anti-mouse-igg-hl-highly-cross-adsorbed-min-x-rat/>