



Glowing products for science

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

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

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. To minimize cross-reactivity, the antibody has been adsorbed against bovine, horse, human, rabbit, and swine serum. This antibody is not cross-adsorbed against rat. For staining rat tissue, or co-staining with rat antibodies, we recommend using [Goat Anti-Mouse IgG \(H L\), Highly Cross-Adsorbed \(Min x Rat\)](#).

- Highly-cross adsorbed for specific staining with minimal background
- Available with 19 bright and photostable CF® Dyes
- Biotin and HRP conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Note: Conjugates of blue fluorescent dyes like CF®350 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 [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.

References

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



Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Mouse IgG
Species reactivity	Mouse
Cross adsorption	Bovine, Horse, Human, Rabbit, Swine
Concentration	2 mg/mL, 1 mg/mL (HRP, AP conjugates)
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, HRP conjugates: PBS/50% glycerol/15 mg/mL BSA, HRP conjugates (lyophilized): PBS/10 mg/mL trehalose after reconstitution
Secondary/tag antibody applications	ELISA, Flow cytometry, IHC, IF (cells or tissue sections), Western blot

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20143-1 20143 20143-1mg	CF@350 Features
CF@405M	408/452 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20182-1 20182 20182-1mg	CF@405M Features
CF@430	426/498 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20459-50uL 20459-500uL 20459-1mg	CF@430 Features
CF@488A	490/515 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20018-1 20018 20018-1mg	CF@488A Features
CF@532	527/558 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20468-50uL 20468-500uL 20468-1mg	CF@532 Features
CF@543	541/560 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20299-1 20299 20299-1mg	CF@543 Features
CF@555	555/565 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20231-1 20231 20231-1mg	CF@555 Features
CF@568	562/583 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20101-1 20101 20101-1mg	CF@568 Features
CF@583R	585/609 nm	50 uL (100 ug) 0.5 mL (1 mg)	20904-50uL 20904-500uL	CF@583R Features
CF@594	593/614 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20111-1 20111 20111-1mg	CF@594 Features
CF@633	630/650 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20121-1 20121 20121-1mg	CF@633 Features
CF@640R	642/662 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20175-1 20175 20175-1mg	CF@640R Features
CF@647	650/665 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20281-1 20281 20281-1mg	CF@647 Features
CF@660C	667/685 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20052-1 20052 20052-1mg	CF@660C Features
CF@680R	680/701 nm	50 uL (100 ug) 0.25 mL (500 ug)	20192-1 20192	CF@680 Features
CF@680	681/698 nm	50 uL (100 ug) 0.25 mL (500 ug)	20065-1 20065	CF@680R Features
CF@750	755/777 nm	50 uL (100 ug) 250 uL (100 ug)	20463-50uL 20463-250uL	CF@750 Features
CF@770	770/797 nm	50 uL (100 ug) 0.25 mL (500 ug)	20077-1 20077	CF@770 Features
CF@790	784/806 nm	50 uL (100 ug)	20342-50uL	CF@790 Features
Biotin	N/A	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20184-1 20184 20184-1mg	
HRP	N/A	100 uL (100 ug) 1 mL (1 mg) 1 mg (lyophilized)	20401-100uL 20401-1mL 20401-1mg	

This datasheet was generated on April 27, 2024 at 07:39:46 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-mouse-igg-hi-highly-cross-absorbed/>