



Glowing products for science

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

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

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



Product Description

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

- Highly-cross adsorbed for specific staining with minimal background
- Available in 15 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

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 [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

Antibody type	Secondary
Host species	Goat
Antibody reactivity (reacts)	Guinea Pig IgG
Cross adsorption	Bovine, Chicken, Goat, Horse, Human, Mouse, Rabbit, Rat, Sheep, Syrian hamster
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
Clonality	Polyclonal
Concentration	2 mg/mL
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot

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) 0.5 mL (1 mg) 1 mg (lyophilized)	20488-50uL 20488-500uL 20488-1mg	CF@405S Features
CF@405M	408/452 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20487-50uL 20487-500uL 20487-1mg	CF@405M Features
CF@488A	490/515 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20489-50uL 20489-500uL 20489-1mg	CF@488A Features
CF@543	541/560 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20490-50uL 20490-500uL 20490-1mg	CF@543 Features
CF@555	555/565 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20491-50uL 20491-500uL 20491-1mg	CF@555 Features
CF@568	562/583 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20492-50uL 20492-500uL 20492-1mg	CF@568 Features
CF@583R	585/609 nm	50 uL (100 ug) 0.5 mL (1 mg)	20900-50uL 20900-500uL	CF@583R Features
CF@594	593/614 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20493-50uL 20493-500uL 20493-1mg	CF@594 Features
CF@640R	642/662 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20494-50uL 20494-500uL 20494-1mg	CF@640R Features
CF@647	650/665 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20495-50uL 20495-500uL 20495-1mg	CF@647 Features
CF@660R	663/682 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20496-50uL 20496-500uL 20496-1mg	CF@660R Features
CF@660C	667/685 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg	20497-50uL 20497-500uL 20497-1mg	CF@660C Features
CF@680	681/698 nm	50 uL (100 ug) 0.5 mL (1 mg)	20499-50uL 20499-500uL	CF@680 Features
CF@680R	680/701 nm	50 uL (100 ug) 0.5 mL (1 mg)	20498-50uL 20498-500uL	CF@680R Features
CF@770	770/797 nm	50 uL (100 ug) 0.5 mL (1 mg)	20500-50uL 20500-500uL	CF@770 Features

This datasheet was generated on May 9, 2024 at 03:16:03 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/goat-anti-guinea-pig-igg-hl-highly-cross-adsorbed/>