



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.

References

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

Product attributes

Antibody type	Secondary
Host species	Goat
Antibody reactivity (species)	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

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 June 30, 2024 at 09:31:40 AM. 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/>