



Glowing products for science

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.



Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody type	Secondary
Host species	Goat
Antibody reactivity (target)	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
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@405S	404/431 nm	50 uL (100 ug)	20488-50uL	CF@405S Features
0.5 mL (1 mg)	20488-500uL			
1 mg (lyophilized)	20488-1mg			
CF@405M	408/452 nm	50 uL (100 ug)	20487-50uL	CF@405M Features
0.5 mL (1 mg)	20487-500uL			
1 mg (lyophilized)	20487-1mg			
CF@488A	490/515 nm	50 uL (100 ug)	20489-50uL	CF@488A Features
0.5 mL (1 mg)	20489-500uL			
1 mg (lyophilized)	20489-1mg			
CF@543	541/560 nm	50 uL (100 ug)	20490-50uL	CF@543 Features
0.5 mL (1 mg)	20490-500uL			
1 mg (lyophilized)	20490-1mg			
CF@555	555/565 nm	50 uL (100 ug)	20491-50uL	CF@555 Features
0.5 mL (1 mg)	20491-500uL			
1 mg (lyophilized)	20491-1mg			
CF@568	562/583 nm	50 uL (100 ug)	20492-50uL	CF@568 Features
0.5 mL (1 mg)	20492-500uL			
1 mg (lyophilized)	20492-1mg			
CF@583R	585/609 nm	50 uL (100 ug)	20900-50uL	CF@583R Features
0.5 mL (1 mg)	20900-500uL			
CF@594	593/614 nm	50 uL (100 ug)	20493-50uL	CF@594 Features
0.5 mL (1 mg)	20493-500uL			
1 mg (lyophilized)	20493-1mg			
CF@640R	642/662 nm	50 uL (100 ug)	20494-50uL	CF@640R Features
0.5 mL (1 mg)	20494-500uL			
1 mg (lyophilized)	20494-1mg			
CF@647	650/665 nm	50 uL (100 ug)	20495-50uL	CF@647 Features
0.5 mL (1 mg)	20495-500uL			
1 mg (lyophilized)	20495-1mg			
CF@660R	663/682 nm	50 uL (100 ug)	20496-50uL	CF@660R Features
0.5 mL (1 mg)	20496-500uL			
1 mg (lyophilized)	20496-1mg			
CF@660C	667/685 nm	50 uL (100 ug)	20497-50uL	CF@660C Features
0.5 mL (1 mg)	20497-500uL			
1 mg	20497-1mg			
CF@680	681/698 nm	50 uL (100 ug)	20499-50uL	CF@680 Features
0.5 mL (1 mg)	20499-500uL			
CF@680R	680/701 nm	50 uL (100 ug)	20498-50uL	CF@680R Features
0.5 mL (1 mg)	20498-500uL			
CF@770	770/797 nm	50 uL (100 ug)	20500-50uL	CF@770 Features
0.5 mL (1 mg)	20500-500uL			

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](#).

