



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) Email: [btinfo@biotium.com](mailto:btinfo@biotium.com)

### Product attributes

<b>Antibody type</b>	Secondary
<b>Host species</b>	Goat
<b>Antibody reactivity (target)</b>	Guinea Pig IgG
<b>Cross adsorption</b>	Bovine, Chicken, Goat, Horse, Human, Mouse, Rabbit, Rat, Sheep, Syrian hamster
<b>Antibody/conjugate formulation</b>	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution
<b>Clonality</b>	Polyclonal
<b>Concentration</b>	2 mg/mL
<b>Secondary/tag antibody applications</b>	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
<b>Product origin</b>	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
<a href="#">CF@405S</a>	404/431 nm	50 uL (100 ug)	<a href="#">20488-50uL</a>	<a href="#">CF@405S Features</a>
		0.5 mL (1 mg)	<a href="#">20488-500uL</a>	
		1 mg (lyophilized)	<a href="#">20488-1mg</a>	
<a href="#">CF@405M</a>	408/452 nm	50 uL (100 ug)	<a href="#">20487-50uL</a>	<a href="#">CF@405M Features</a>
		0.5 mL (1 mg)	<a href="#">20487-500uL</a>	
		1 mg (lyophilized)	<a href="#">20487-1mg</a>	
<a href="#">CF@488A</a>	490/515 nm	50 uL (100 ug)	<a href="#">20489-50uL</a>	<a href="#">CF@488A Features</a>
		0.5 mL (1 mg)	<a href="#">20489-500uL</a>	
		1 mg (lyophilized)	<a href="#">20489-1mg</a>	
<a href="#">CF@543</a>	541/560 nm	50 uL (100 ug)	<a href="#">20490-50uL</a>	<a href="#">CF@543 Features</a>
		0.5 mL (1 mg)	<a href="#">20490-500uL</a>	
		1 mg (lyophilized)	<a href="#">20490-1mg</a>	
<a href="#">CF@555</a>	555/565 nm	50 uL (100 ug)	<a href="#">20491-50uL</a>	<a href="#">CF@555 Features</a>
		0.5 mL (1 mg)	<a href="#">20491-500uL</a>	
		1 mg (lyophilized)	<a href="#">20491-1mg</a>	
<a href="#">CF@568</a>	562/583 nm	50 uL (100 ug)	<a href="#">20492-50uL</a>	<a href="#">CF@568 Features</a>
		0.5 mL (1 mg)	<a href="#">20492-500uL</a>	
		1 mg (lyophilized)	<a href="#">20492-1mg</a>	
<a href="#">CF@583R</a>	585/609 nm	50 uL (100 ug)	<a href="#">20900-50uL</a>	<a href="#">CF@583R Features</a>
		0.5 mL (1 mg)	<a href="#">20900-500uL</a>	
<a href="#">CF@594</a>	593/614 nm	50 uL (100 ug)	<a href="#">20493-50uL</a>	<a href="#">CF@594 Features</a>
		0.5 mL (1 mg)	<a href="#">20493-500uL</a>	
		1 mg (lyophilized)	<a href="#">20493-1mg</a>	
<a href="#">CF@640R</a>	642/662 nm	50 uL (100 ug)	<a href="#">20494-50uL</a>	<a href="#">CF@640R Features</a>
		0.5 mL (1 mg)	<a href="#">20494-500uL</a>	
		1 mg (lyophilized)	<a href="#">20494-1mg</a>	
<a href="#">CF@647</a>	650/665 nm	50 uL (100 ug)	<a href="#">20495-50uL</a>	<a href="#">CF@647 Features</a>
		0.5 mL (1 mg)	<a href="#">20495-500uL</a>	
		1 mg (lyophilized)	<a href="#">20495-1mg</a>	
<a href="#">CF@660R</a>	663/682 nm	50 uL (100 ug)	<a href="#">20496-50uL</a>	<a href="#">CF@660R Features</a>
		0.5 mL (1 mg)	<a href="#">20496-500uL</a>	
		1 mg (lyophilized)	<a href="#">20496-1mg</a>	
<a href="#">CF@660C</a>	667/685 nm	50 uL (100 ug)	<a href="#">20497-50uL</a>	<a href="#">CF@660C Features</a>
		0.5 mL (1 mg)	<a href="#">20497-500uL</a>	
		1 mg	<a href="#">20497-1mg</a>	
<a href="#">CF@680</a>	681/698 nm	50 uL (100 ug)	<a href="#">20499-50uL</a>	<a href="#">CF@680 Features</a>
		0.5 mL (1 mg)	<a href="#">20499-500uL</a>	
<a href="#">CF@680R</a>	680/701 nm	50 uL (100 ug)	<a href="#">20498-50uL</a>	<a href="#">CF@680R Features</a>
		0.5 mL (1 mg)	<a href="#">20498-500uL</a>	
<a href="#">CF@770</a>	770/797 nm	50 uL (100 ug)	<a href="#">20500-50uL</a>	<a href="#">CF@770 Features</a>
		0.5 mL (1 mg)	<a href="#">20500-500uL</a>	

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 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 March 29, 2025 at 01:37:33 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/>

