

# Goat Anti-Human IgG (H+L), Fc Gamma Specific, Highly Cross-Adsorbed

Highly cross-adsorbed goat anti-human IgG (H+L), Fc gamma specific secondary antibody labeled with our superior CF® dyes.

# **Product Description**

This is a highly cross-adsorbed goat anti-human IgG (H L), Fc gamma specific secondary antibody labeled with our superior CF® dyes. The antibody reacts specifically with the human IgG heavy chains and not with light chains or F(ab')2 fragment of human immunoglobulins. To minimize cross-reactivity, the antibodies are cross-adsorbed against bovine, horse, mouse and rabbit serum IgG prior to conjugation.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 5 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology

Clonality	Polyclonal		
Antibody type	Secondary, Anti-Human Immunoglobulin		
Concentration	2 mg/mL		
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		
Species reactivity	Human		
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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.		
Host species	Goat		
Antibody reactivity (target)	Human IgG Fc		
Cross adsorption	Bovine, Horse, Mouse, Rabbit		

Call us: 800-304-5357

Product attributes

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/515 nm	50 uL (100 ug)	20444-50uL	CF®488A Features
0.5 mL (1 mg)	20444-500uL			
1 mg (lyophilized)	20444-1mg			
<u>CF®543</u>	541/560 nm	50 uL (100 ug)	20445-50uL	CF®543 Features
0.5 mL (1 mg)	20445-500uL			
1 mg (lyophilized)	20445-1mg			
<u>CF®594</u>	593/614 nm	50 uL (100 ug)	20446-50uL	CF®594 Features
0.5 mL (1 mg)	20446-500uL			
1 mg (lyophilized)	20446-1mg			
<u>CF®640R</u>	642/662 nm	50 uL (100 ug)	20447-50uL	CF®640R Features
0.5 mL (1 mg)	20447-500uL			
1 mg (lyophilized)	20447-1mg			
CF®647	650/665 nm	50 uL (100 ug)	20448-50uL	CF®647 Features
0.5 mL (1 mg)	20448-500uL			
1 mg (lyophilized)	20448-1mg			

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 November 28, 2025 at 11:57:30 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/goat-anti-human-igg-hl-fc-gamma-specific-highly-cross-adsorbed/