

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

Product attributes

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 (<i>Bos taurus</i>), 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

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	
CF@543	541/560 nm	50 uL (100 ug)	20445-50uL	CF@543 Features
		0.5 mL (1 mg)	20445-500uL	
		1 mg (lyophilized)	20445-1mg	
CF@594	593/614 nm	50 uL (100 ug)	20446-50uL	CF@594 Features
		0.5 mL (1 mg)	20446-500uL	
		1 mg (lyophilized)	20446-1mg	
CF@640R	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 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](#).

This datasheet was generated on November 21, 2024 at 03:42:59 AM. 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/>