



Glowing products for science

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

Highly cross-adsorbed goat anti-human IgG (H L) secondary antibody labeled with our superior CF® Dyes.



Product Description

This is a highly cross-adsorbed goat anti-human IgG (H L) secondary antibody labeled with our bright and photostable CF® Dyes. To minimize cross-reactivity, the antibody has been adsorbed against bovine, horse, and mouse serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 11 bright and photostable CF® Dyes
- HRP and R-PE conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

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

Product attributes

Antibody type	Secondary, Anti-Human Immunoglobulin
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Human IgG
Species reactivity	Human
Cross adsorption	Bovine, Horse, Mouse
Concentration	2 mg/mL, 1 mg/mL (HRP, AP conjugates), 0.5 mg/mL (R-PE, APC conjugates)
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, HRP conjugates: PBS/50% glycerol/15 mg/mL BSA, HRP conjugates (lyophilized): PBS/10 mg/mL trehalose after reconstitution
Secondary/tag antibody applications	ELISA, 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-Human IgG (H+L), Highly Cross-Adsorbed

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@488A	490/515 nm	50 uL (100 ug)	20022-1	CF@488A Features
0.5 mL (1 mg)	20022			
1 mg	20022-1mg			
CF@543	541/560 nm	50 uL (100 ug)	20319-1	CF@543 Features
0.5 mL (1 mg)	20319			
1 mg	20319-1mg			
CF@555	555/565 nm	50 uL (100 ug)	20230-1	CF@555 Features
0.5 mL (1 mg)	20230			
1 mg	20230-1mg			
CF@568	562/583 nm	50 uL (100 ug)	20097-1	CF@568 Features
0.5 mL (1 mg)	20097			
1 mg	20097-1mg			
CF@583R	585/609 nm	50 uL (100 ug)	20901-50uL	CF@583R Features
0.5 mL (1 mg)	20901-500uL			
CF@594	593/614 nm	50 uL (100 ug)	20154-1	CF@594 Features
0.5 mL (1 mg)	20154			
1 mg	20154-1mg			
CF@633	630/650 nm	50 uL (100 ug)	20132-1	CF@633 Features
0.5 mL (1 mg)	20132			
1 mg	20132-1mg			
CF@640R	642/662 nm	50 uL (100 ug)	20081-1	CF@640R Features
0.5 mL (1 mg)	20081			
1 mg	20081-1mg			
CF@647	650/665 nm	50 uL (100 ug)	20280-1	CF@647 Features
0.5 mL (1 mg)	20280			
1 mg	20280-1mg			
CF@680	681/698 nm	50 uL (100 ug)	20287-1	CF@680 Features
0.25 mL (500 ug)	20287			
CF@770	770/797 nm	50 uL (100 ug)	20288-1	CF@770 Features
0.25 mL (500 ug)	20288			
R-PE	496, 546, 565/578 nm	200 uL (100 ug)	20355-200uL	
1 mL (500 ug)	20355-1mL			
HRP	N/A	100 uL (100 ug)	20470-100uL	
1 mL (1 mg)	20470-1mL			
1 mg	20470-1mg			

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 curated [CF® Dye references](#).

This datasheet was generated on June 3, 2026 at 04:07:01 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-human-igg-hl-highly-cross-adsorbed/>