

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.

ntotlum

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 Email: techsupport@biotium.com

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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovany 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	490/515 nm	50 uL (100 ug)	20022-1	CF®488A Features
		0.5 mL (1 mg)	<u>20022</u>	
		1 mg	20022-1mg	
CF®543 541/560 nm	541/560 nm	50 uL (100 ug)	<u>20319-1</u>	CF®543 Features
		0.5 mL (1 mg)	<u>20319</u>	
		1 mg	20319-1mg	
CF®555 555/565 nn	555/565 nm	50 uL (100 ug)	<u>20230-1</u>	CF®555 Features
		0.5 mL (1 mg)	20230	
		1 mg	20230-1mg	
CF®568 562/583 r	562/583 nm	50 uL (100 ug)	<u>20097-1</u>	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®594 593/61	593/614 nm	50 uL (100 ug)	<u>20154-1</u>	CF®594 Features
		0.5 mL (1 mg)	<u>20154</u>	
		1 mg	20154-1mg	
CF®633 630/650 nm	630/650 nm	50 uL (100 ug)	<u>20132-1</u>	CF®633 Features
		0.5 mL (1 mg)	<u>20132</u>	
		1 mg	20132-1mg	
CF®640R 642/662 nm	642/662 nm	50 uL (100 ug)	<u>20081-1</u>	CF®640R Features
		0.5 mL (1 mg)	<u>20081</u>	
		1 mg	20081-1mg	
CF®647 650/6	650/665 nm	50 uL (100 ug)	<u>20280-1</u>	CF®647 Features
		0.5 mL (1 mg)	20280	
		1 mg	20280-1mg	
CF®680	681/698 nm	50 uL (100 ug)	<u>20287-1</u>	CF®680 Features
		0.25 mL (500 ug)	<u>20287</u>	
CF®770	770/797 nm	50 uL (100 ug)	<u>20288-1</u>	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	N/A	100 uL (100 ug)	20470-100uL	
		1 mL (1 mg)	20470-1mL	
		1 mg	20470-1mg	

View our full selection of <u>Secondary Antibodies</u>, or search our catalog using our <u>Antibody Finder</u>. Alternatively, you can view our <u>secondary antibody product listings</u> 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 October 27, 2025 at 03:18:01 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/