



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

### Product attributes

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

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

This datasheet was generated on November 21, 2024 at 03:06:44 AM. Visit product page to check for updated information before use.

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