

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



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

### Product Description

This is a highly cross-adsorbed goat anti-rat 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, human, and rabbit serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 14 bright and photostable CF® Dyes
- APC, HRP, and R-PE conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

**Note:** Conjugates of blue fluorescent dyes like CF@350 and CF@405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

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

#### Product attributes

<b>Antibody type</b>	Secondary
<b>Clonality</b>	Polyclonal
<b>Host species</b>	Goat
<b>Antibody reactivity (target)</b>	Rat IgG
<b>Species reactivity</b>	Rat
<b>Cross adsorption</b>	Bovine, Horse, Human, Rabbit
<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, R-PE conjugates: PBS/2 mg/mL BSA/0.05% azide, 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.

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	0.5 mL (1 mg)	<a href="#">20147</a>	<a href="#">CF@350 Features</a>
1 mg	<a href="#">20147-1mg</a>			
50 uL (100 ug)	<a href="#">20147-1</a>			
CF@405M	408/452 nm	0.5 mL (1 mg)	<a href="#">20374-500uL</a>	<a href="#">CF@405M Features</a>
1 mg	<a href="#">20374-1mg</a>			
50 uL (100 ug)	<a href="#">20374-50uL</a>			
CF@488A	490/515 nm	0.5 mL (1 mg)	<a href="#">20023</a>	<a href="#">CF@488A Features</a>
1 mg	<a href="#">20023-1mg</a>			
50 uL (100 ug)	<a href="#">20023-1</a>			
CF@543	541/560 nm	0.5 mL (1 mg)	<a href="#">20321</a>	<a href="#">CF@543 Features</a>
1 mg	<a href="#">20321-1mg</a>			
50 uL (100 ug)	<a href="#">20321-1</a>			
CF@555	555/565 nm	0.5 mL (1 mg)	<a href="#">20233</a>	<a href="#">CF@555 Features</a>
1 mg	<a href="#">20233-1mg</a>			
50 uL (100 ug)	<a href="#">20233-1</a>			
CF@568	562/583 nm	0.5 mL (1 mg)	<a href="#">20096</a>	<a href="#">CF@568 Features</a>
1 mg	<a href="#">20096-1mg</a>			
50 uL (100 ug)	<a href="#">20096-1</a>			
CF@583R	585/609 nm	50 uL (100 ug)	<a href="#">20906-50uL</a>	<a href="#">CF@583R Features</a>
0.5 mL (1 mg)	<a href="#">20906-500uL</a>			
CF@594	593/614 nm	0.5 mL (1 mg)	<a href="#">20155</a>	<a href="#">CF@594 Features</a>
1 mg	<a href="#">20155-1mg</a>			
50 uL (100 ug)	<a href="#">20155-1</a>			
CF@633	630/650 nm	0.5 mL (1 mg)	<a href="#">20133</a>	<a href="#">CF@633 Features</a>
1 mg	<a href="#">20133-1mg</a>			
50 uL (100 ug)	<a href="#">20133-1</a>			
CF@640R	642/662 nm	0.5 mL (1 mg)	<a href="#">20088</a>	<a href="#">CF@640R Features</a>
1 mg	<a href="#">20088-1mg</a>			
50 uL (100 ug)	<a href="#">20088-1</a>			
CF@647	650/665 nm	0.5 mL (1 mg)	<a href="#">20283</a>	<a href="#">CF@647 Features</a>
1 mg	<a href="#">20283-1mg</a>			
50 uL (100 ug)	<a href="#">20283-1</a>			
CF@660C	667/685 nm	0.5 mL (1 mg)	<a href="#">20370-500uL</a>	<a href="#">CF@660C Features</a>
1 mg	<a href="#">20370-1mg</a>			
50 uL (100 ug)	<a href="#">20370-50uL</a>			
CF@680	681/698 nm	0.25 mL	<a href="#">20069</a>	<a href="#">CF@680 Features</a>
50 uL (100 ug)	<a href="#">20069-1</a>			
CF@770	770/797 nm	0.25 mL	<a href="#">20383-250uL</a>	<a href="#">CF@770 Features</a>
50 uL (100 ug)	<a href="#">20383-50uL</a>			
R-PE	496, 546, 565/578 nm	200 uL (100 ug)	<a href="#">20354-200uL</a>	
1 mL (500 ug)	<a href="#">20354-1mL</a>			

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
APC 0.5 mL (250 ug)	650/660 nm <a href="#">20413-500uL</a>	100 uL (50 ug)	<a href="#">20413-100uL</a>	
HRP 1 mL (1 mg) 1 mg	N/A <a href="#">20406-1mL</a> <a href="#">20406-1mg</a>	100 uL (100 ug)	<a href="#">20406-100uL</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 June 3, 2026 at 03:15:13 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-rat-igghl-highly-cross-adsorbed/>