

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



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

### Product Description

This is a highly cross-adsorbed rabbit 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 human serum.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 6 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

### Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Rabbit
Antibody reactivity (target)	Rat IgG
Species reactivity	Rat
Cross adsorption	Human
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
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.

## Rabbit Anti-Rat 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="#">20025-1</a>	<a href="#">CF@488A Features</a>
		0.5 mL (1 mg)	<a href="#">20025</a>	
		1 mg	<a href="#">20025-1mg</a>	
CF@568	562/583 nm	50 uL (100 ug)	<a href="#">20094-1</a>	<a href="#">CF@568 Features</a>
		0.5 mL (1 mg)	<a href="#">20094</a>	
		1 mg	<a href="#">20094-1mg</a>	
CF@583R	585/609 nm	50 uL (100 ug)	<a href="#">20912-50uL</a>	<a href="#">CF@583R Features</a>
		0.5 mL (1 mg)	<a href="#">20912-500uL</a>	
CF@594	593/614 nm	50 uL (100 ug)	<a href="#">20157-1</a>	<a href="#">CF@594 Features</a>
		0.5 mL (1 mg)	<a href="#">20157</a>	
		1 mg	<a href="#">20157-1mg</a>	
CF@633	630/650 nm	50 uL (100 ug)	<a href="#">20135-1</a>	<a href="#">CF@633 Features</a>
		0.5 mL (1 mg)	<a href="#">20135</a>	
		1 mg	<a href="#">20135-1mg</a>	
CF@640R	642/662 nm	50 uL (100 ug)	<a href="#">20201-1</a>	<a href="#">CF@640R Features</a>
		0.5 mL (1 mg)	<a href="#">20201</a>	
		1 mg	<a href="#">20201-1mg</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 December 3, 2024 at 09:20:27 AM. Visit product page to check for updated information before use.

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