

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

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



Product Description

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

- Highly cross-adsorbed for specific staining with minimal background
- Available in 20 bright and photostable CF® Dyes
- Biotin and HRP conjugates also available
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

Note: Conjugates of blue fluorescent dyes like CF®350 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.

Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Goat
Antibody reactivity (target)	Rabbit IgG
Species reactivity	Rabbit
Cross adsorption	Human, Mouse, Rat
Concentration	2 mg/mL, 1 mg/mL (HRP, AP 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-Rabbit IgG (H+L), Highly Cross-Adsorbed

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®350	347/448 nm	50 uL (100 ug)	20144-1	CF®350 Features
		0.5 mL (1 mg)	20144	
		1 mg	20144-1mg	
CF®405M	408/452 nm	50 uL (100 ug)	20373-50uL	CF®405M Features
		0.5 mL (1 mg)	20373-500uL	
		1 mg	20373-1mg	
CF®430	426/498 nm	50 uL (100 ug)	20460-50uL	CF®430 Features
		1 mg	20460-1mg	
		0.5 mL (1 mg)	20460-500uL	
CF®488A	490/515 nm	50 uL (100 ug)	20019-1	CF®488A Features
		0.5 mL (1 mg)	20019	
		1 mg	20019-1mg	
CF®532	527/558 nm	50 uL (100 ug)	20469-50uL	CF®532 Features
		1 mg	20469-1mg	
		0.5 mL (1 mg)	20469-500uL	
CF®543	541/560 nm	50 uL (100 ug)	20300-1	CF®543 Features
		0.5 mL (1 mg)	20300	
		1 mg	20300-1mg	
CF®555	555/565 nm	50 uL (100 ug)	20232-1	CF®555 Features
		0.5 mL (1 mg)	20232	
		1 mg	20232-1mg	
CF®568	562/583 nm	50 uL (100 ug)	20103-1	CF®568 Features
		0.5 mL (1 mg)	20103	
		1 mg	20103-1mg	
CF®583R	585/609 nm	50 uL (100 ug)	20905-50uL	CF®583R Features
		0.5 mL (1 mg)	20905-500uL	
CF®594	593/614 nm	50 uL (100 ug)	20113-1	CF®594 Features
		0.5 mL (1 mg)	20113	
		1 mg	20113-1mg	
CF®633	630/650 nm	50 uL (100 ug)	20123-1	CF®633 Features
		0.5 mL (1 mg)	20123	
		1 mg	20123-1mg	
CF®640R	642/662 nm	50 uL (100 ug)	20176-1	CF®640R Features
		0.5 mL (1 mg)	20176	
		1 mg	20176-1mg	
CF®647	650/665 nm	50 uL (100 ug)	20282-1	CF®647 Features
		0.5 mL (1 mg)	20282	
		1 mg	20282-1mg	
CF®660C	667/685 nm	50 uL (100 ug)	20369-50uL	CF®660C Features
		0.5 mL (1 mg)	20369-500uL	
		1 mg	20369-1mg	

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®680	681/698 nm	50 uL (100 ug)	20067-1	CF®680 Features
		0.25 mL (500 ug)	20067	
CF®680R	680/701 nm	50 uL (100 ug)	20193-1	CF®680R Features
		0.25 mL (500 ug)	20193	
CF®740	742/767 nm	50 uL (100 ug)	20990-50uL	CF®740 Features
		250 uL (500 ug)	20990-250uL	
CF®770	770/797 nm	50 uL (100 ug)	20078-1	CF®770 Features
		0.25 mL (500 ug)	20078	
CF®790	784/806 nm	50 uL (100 ug)	20343-50uL	CF®790 Features
CF®800	797/816 nm	50 uL (100 ug)	20832-50uL	CF®800 Features
Biotin	N/A	50 uL (100 ug)	20186-1	
		0.5 mL (1 mg)	20186	
		1 mg	20186-1mg	
HRP	N/A	100 uL (100 ug)	20403-100uL	
		1 mL (1 mg)	20403-1mL	
		1 mg	20403-1mg	

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 [CF® dye references](#).

This datasheet was generated on December 18, 2025 at 01:28:50 AM. Visit product page to check for updated information before use.
 Product link: <https://legacy.biotium.com/product/goat-anti-rabbit-igg-hl-highly-cross-absorbed/>