

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

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



Product Description

This is a highly cross-adsorbed rabbit anti-mouse 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 11 bright and photostable CF® Dyes
- 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. 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.

Product attributes

Antibody type	Secondary
Clonality	Polyclonal
Host species	Rabbit
Antibody reactivity (target)	Mouse IgG
Species reactivity	Mouse
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

References

Download a list of [CF® dye references](#).

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®350	347/448 nm	50 uL (100 ug)	20149-1	CF®350 Features
		0.5 mL (1 mg)	20149	
		1 mg	20149-1mg	
CF®488A	490/515 nm	50 uL (100 ug)	20026-1	CF®488A Features
		0.5 mL (1 mg)	20026	
		1 mg	20026-1mg	
CF®543	541/560 nm	50 uL (100 ug)	20307-1	CF®543 Features
		0.5 mL (1 mg)	20307	
		1 mg	20307-1mg	
CF®555	555/565 nm	50 uL (100 ug)	20235-1	CF®555 Features
		0.5 mL (1 mg)	20235	
		1 mg	20235-1mg	
CF®568	562/583 nm	50 uL (100 ug)	20093-1	CF®568 Features
		0.5 mL (1 mg)	20093	
		1 mg	20093-1mg	
CF®583R	585/609 nm	50 uL (100 ug)	20911-50uL	CF®583R Features
		0.5 mL (1 mg)	20911-500uL	
		1 mg	20911-500uL	
CF®594	593/614 nm	50 uL (100 ug)	20158-1	CF®594 Features
		0.5 mL (1 mg)	20158	
		1 mg	20158-1mg	
CF®633	630/650 nm	50 uL (100 ug)	20136-1	CF®633 Features
		0.5 mL (1 mg)	20136	
		1 mg	20136-1mg	
CF®640R	642/662 nm	50 uL (100 ug)	20200-1	CF®640R Features
		0.5 mL (1 mg)	20200	
		1 mg	20200-1mg	
CF®647	650/665 nm	50 uL (100 ug)	20285-1	CF®647 Features
		0.5 mL (1 mg)	20285	
		1 mg	20285-1mg	
CF®680	681/698 nm	50 uL (100 ug)	20061-1	CF®680 Features
		0.25 mL (500 ug)	20061	

This datasheet was generated on July 17, 2024 at 08:30:27 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rabbit-anti-mouse-igg-hl-highly-cross-adsorbed/>

