Chicken Anti-Mouse IgG (H+L)

Chicken anti-mouse IgG (H L) secondary antibody labeled with our superior CF® Dyes.



Product Description

This is a chicken anti-mouse IgG (H L) secondary antibody labeled with our superior CF® Dyes.

- Available in 7 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.

Product attributes

Antibody type	Secondary		
Clonality	Polyclonal		
Host species	Chicken		
Antibody reactivity (target)	Mouse IgG		
Species reactivity	Mouse		
Cross adsorption	Not cross-adsorbed		
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 (Bos taurus), or recombinant BSA produced in Chipase hamster over years calls. Inquire for the specific let		

Call us: 800-304-5357 Email: btinfo@biotium.com

Chicken Anti-Mouse IgG (H+L)

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®350	347/448 nm	50 uL (100 ug)	20331-50uL	CF®350 Features
		0.5 mL (1 mg)	20331-500uL	
		1 mg (lyophilized)	20331-1mg	
CF®488A	490/515 nm	50 uL (100 ug)	<u>20208-1</u>	CF®488A Features
		0.5 mL (1 mg)	<u>20208</u>	
		1 mg (lyophilized)	20208-1mg	
CF®543	541/560 nm	50 uL (100 ug)	20334-50uL	CF®543 Features
		0.5 mL (1 mg)	20334-500uL	
		1 mg (lyophilized)	20334-1mg	
CF®568	562/583 nm	50 uL (100 ug)	<u>20338-50uL</u>	CF®568 Features
		0.5 mL (1 mg)	20338-500uL	
		1 mg (lyophilized)	20338-1mg	
CF®583R	585/609 nm	50 uL (100 ug)	<u>20890-50uL</u>	CF®583R Features
		0.5 mL (1 mg)	20890-500uL	
CF®594	593/614 nm	50 uL (100 ug)	<u>20221-1</u>	CF®594 Features
		0.5 mL (1 mg)	<u>20221</u>	
		1 mg (lyophilized)	20221-1mg	
CF®633	630/650 nm	50 uL (100 ug)	<u>20222-1</u>	CF®633 Features
		0.5 mL (1 mg)	<u>20222</u>	
		1 mg (lyophilized)	20222-1mg	

View our full selection of <u>Secondary Antibodies</u>, or search our catalog using our <u>Antibody Finder</u>. Alternatively, you can view our <u>secondary antibody product listings</u> 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 November 8, 2025 at 04:18:20 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/chicken-anti-mouse-igg-hl/