

## 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

<b>Antibody type</b>	Secondary
<b>Clonality</b>	Polyclonal
<b>Host species</b>	Chicken
<b>Antibody reactivity (target)</b>	Mouse IgG
<b>Species reactivity</b>	Mouse
<b>Cross adsorption</b>	Not cross-adsorbed
<b>Concentration</b>	2 mg/mL
<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
<b>Secondary/tag antibody applications</b>	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.

## Chicken Anti-Mouse IgG (H+L)

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

This datasheet was generated on November 21, 2024 at 03:17:13 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/chicken-anti-mouse-igg-hl/>