

Chicken Anti-Mouse IgG (H+L)

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



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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

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.

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)	20208-1	CF®488A Features
0.5 mL (1 mg)	20208			
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)	20338-50uL	CF®568 Features
0.5 mL (1 mg)	20338-500uL			
1 mg (lyophilized)	20338-1mg			
CF®583R	585/609 nm	50 uL (100 ug)	20890-50uL	CF®583R Features
0.5 mL (1 mg)	20890-500uL			
CF®594	593/614 nm	50 uL (100 ug)	20221-1	CF®594 Features
0.5 mL (1 mg)	20221			
1 mg (lyophilized)	20221-1mg			
CF®633	630/650 nm	50 uL (100 ug)	20222-1	CF®633 Features
0.5 mL (1 mg)	20222			
1 mg (lyophilized)	20222-1mg			

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](#).