

Donkey Anti-Mouse IgG (H+L), Highly Cross-Adsorbed (Min X Rat)

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



Product Description

This is a highly cross-adsorbed donkey anti-mouse IgG (H L) secondary antibody labeled with our bright and photostable CF® Dyes and other labels. To minimize cross-reactivity, the antibody has been adsorbed against bovine, chicken, goat, guinea pig, Syrian hamster, horse, human, rabbit, rat and sheep serum proteins.

- Highly cross-adsorbed for specific staining with minimal background
- Available in 17 bright and photostable CF® Dyes
- Biotin, AP, and HRP conjugates also available
- Suitable for western, immunofluorescence, and immunohistology

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	Donkey
Antibody reactivity (target)	Mouse IgG
Species reactivity	Mouse
Cross adsorption	Bovine, Chicken, Goat, Guinea pig, Horse, Human, Rabbit, Rat, Sheep, Syrian hamster
Concentration	2 mg/mL, 1 mg/mL (HRP, AP conjugates)
Antibody/conjugate formulation	AP conjugates: 1 mg/mL in PBS/40% glycerol/5 mg/mL BSA/0.05% azide, 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.

Donkey Anti-Mouse IgG (H+L), Highly Cross-Adsorbed (Min X Rat)

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®350	347/448 nm	50 uL (100 ug)	20350-1	CF®350 Features
0.5 mL (1 mg)	20350			
1 mg (lyophilized)	20350-1mg			
CF®430	426/498 nm	50 uL (100 ug)	20461-50uL	CF®430 Features
0.5 mL (1 mg)	20461-500uL			
1 mg (lyophilized)	20461-1mg			
CF®488A	490/515 nm	50 uL (100 ug)	20014-1	CF®488A Features
0.5 mL (1 mg)	20014			
1 mg (lyophilized)	20014-1mg			
CF®514	516/548 nm	50 uL (100 ug)	20483-50uL	CF®514 Features
0.5 mL (1 mg)	20483-500uL			
1 mg (lyophilized)	20483-1mg			
CF®543	541/560 nm	50 uL (100 ug)	20305-1	CF®543 Features
0.5 mL (1 mg)	20305			
1 mg (lyophilized)	20305-1mg			
CF®555	555/565 nm	50 uL (100 ug)	20037-1	CF®555 Features
0.5 mL (1 mg)	20037			
1 mg (lyophilized)	20037-1mg			
CF®568	562/583 nm	50 uL (100 ug)	20105-1	CF®568 Features
0.5 mL (1 mg)	20105			
1 mg (lyophilized)	20105-1mg			
CF®583R	585/609 nm	50 uL (100 ug)	20894-50uL	CF®583R Features
0.5 mL (1 mg)	20894-500uL			
CF®594	593/614 nm	50 uL (100 ug)	20115-1	CF®594 Features
0.5 mL (1 mg)	20115			
1 mg (lyophilized)	20115-1mg			
CF®633	630/650 nm	50 uL (100 ug)	20124-1	CF®633 Features
0.5 mL (1 mg)	20124			
1 mg (lyophilized)	20124-1mg			
CF®640R	642/662 nm	50 uL (100 ug)	20177-1	CF®640R Features
0.5 mL (1 mg)	20177			
1 mg (lyophilized)	20177-1mg			
CF®647	650/665 nm	50 uL (100 ug)	20046-1	CF®647 Features
0.5 mL (1 mg)	20046			
1 mg (lyophilized)	20046-1mg			
CF®660R	663/682 nm	50 uL (100 ug)	20388-50uL	CF®660R Features
0.5 mL (1 mg)	20388-500uL			
1 mg (lyophilized)	20388-1mg			
CF®680R	680/701 nm	50 uL (100 ug)	20194-1	CF®680R Features
0.25 mL (500 ug)	20194			

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®725 0.25 mL (500 ug)	729/750 nm 20504-250uL	50 uL (100 ug)	20504-50uL	CF®725 Features
CF®740 0.25 mL (500 ug)	742/767 nm 20991-250uL	50 uL (100 ug)	20991-50uL	CF®740 Features
CF®790	784/806 nm	50 uL (100 ug)	20363-50uL	CF®790 Features
CF®800	797/816 nm	50 uL (100 ug)	20835-50uL	CF®800 Features
Biotin 0.5 mL (1 mg) 1 mg (lyophilized)	N/A 20190 20190-1mg	50 uL (100 ug)	20190-1	
HRP 1 mL (1 mg) 1 mg (lyophilized)	N/A 20404-1mL 20404-1mg	100 uL (100 ug)	20404-100uL	
AP 1 mL (1 mg)	N/A 20466-1mL	100 uL (100 ug)	20466-100uL	

View our full selection of bright and specific [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 curated [CF® Dye references](#).

This datasheet was generated on June 3, 2026 at 04:05:37 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/donkey-anti-mouse-igg-hl-highly-cross-adsorbed/>