

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 16 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. 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](#).

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	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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@350	347/448 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20350-1 20350 20350-1mg	CF@350 Features
CF@430	426/498 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20461-50uL 20461-500uL 20461-1mg	CF@430 Features
CF@488A	490/515 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20014-1 20014 20014-1mg	CF@488A Features
CF@514	516/548 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20483-50uL 20483-500uL 20483-1mg	CF@514 Features
CF@543	541/560 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20305-1 20305 20305-1mg	CF@543 Features
CF@555	555/565 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20037-1 20037 20037-1mg	CF@555 Features
CF@568	562/583 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20105-1 20105 20105-1mg	CF@568 Features
CF@583R	585/609 nm	50 uL (100 ug) 0.5 mL (1 mg)	20894-50uL 20894-500uL	CF@583R Features
CF@594	593/614 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20115-1 20115 20115-1mg	CF@594 Features
CF@633	630/650 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20124-1 20124 20124-1mg	CF@633 Features
CF@640R	642/662 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20177-1 20177 20177-1mg	CF@640R Features
CF@647	650/665 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20046-1 20046 20046-1mg	CF@647 Features
CF@660R	663/682 nm	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20388-50uL 20388-500uL 20388-1mg	CF@660R Features
CF@680R	680/701 nm	50 uL (100 ug) 0.25 mL (500 ug)	20194-1 20194	CF@680R 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	N/A	50 uL (100 ug) 0.5 mL (1 mg) 1 mg (lyophilized)	20190-1 20190 20190-1mg	
HRP	N/A	100 uL (100 ug) 1 mL (1 mg) 1 mg (lyophilized)	20404-100uL 20404-1mL 20404-1mg	
AP	N/A	100 uL (100 ug) 1 mL (1 mg)	20466-100uL 20466-1mL	

This datasheet was generated on July 5, 2024 at 11:31:48 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/donkey-anti-mouse-igg-hl-highly-cross-adsorbed/>