



Glowing products for science

Donkey Anti-Mouse IgG (H+L), Highly Cross-Adsorbed, CF® Dye Conjugates, Single Label for STORM

Highly cross-adsorbed donkey anti-mouse IgG (H+L) secondary antibody with single CF® Dye label for STORM super-resolution microscopy.

Product Description

This is a highly cross-adsorbed donkey anti-mouse IgG (H L) secondary antibody that has a low degree of labeling (DOL) with one of our CF® Dyes validated for STORM.

- CF® Dye single label secondary antibody ideal for STORM imaging
- Highly cross-adsorbed for specific staining with minimal background
- Available in 10 CF® Dyes validated for STORM

Secondary antibodies with a low DOL, or number of dye molecules per antibody molecule, have been reported to be optimal for STORM ([Bittel et al. \(2015\) Proc. SPIE 9331](#)). This product is prepared by single labeling (DOL=1) of highly cross-adsorbed goat anti-mouse IgG (H L) with a selection of compatible CF® Dyes for **(d)-STORM super-resolution microscopy**. To minimize cross-reactivity, the antibody has been adsorbed against bovine, chicken, goat, guinea pig, horse, human, rabbit, rat, sheep, and Syrian hamster.

Learn more about [CF® Dyes for super-resolution microscopy](#).



Call us : [800-304-5357](tel:800-304-5357) Email: btinfo@biotium.com

Product attributes

Antibody reactivity (target)	Mouse IgG
Clonality	Polyclonal
Host species	Donkey
Cross adsorption	Bovine, Chicken, Goat, Guinea pig, Horse, Human, Rabbit, Rat, Sheep, Syrian hamster
Antibody type	Secondary
Concentration	1 mg/mL
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
Species reactivity	Mouse
Secondary/tag antibody applications	IF (cells or tissue sections), STORM
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, CF® Dye Conjugates, Single Label for STORM

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@498	498/519 nm	50 uL (50 ug)	20994-50uL	
		0.5 mL (500 ug)	20994-500uL	
CF@505	505/519 nm	50 uL (50 ug)	20878-50uL	
		0.5 mL (500 ug)	20878-500uL	
CF@535ST	535/568 nm	50 uL (50 ug)	20823-50uL	CF@535ST Features
		0.5 mL (500 ug)	20823-500uL	
CF@568	562/583 nm	50 uL (50 ug)	20802-50uL	CF@568 Features
		0.5 mL (500 ug)	20802-500uL	
CF@583R	586/609 nm	50 uL (50 ug)	20794-50uL	CF@583R Features
		0.5 mL (500 ug)	20794-500uL	
CF@597R	597/619 nm	50 uL (50 ug)	20798-50uL	CF@597R Features
		0.5 mL (500 ug)	20798-500uL	
CF@647	650/665 nm	50 uL (50 ug)	20810-50uL	CF@647 Features
		0.5 mL (500 ug)	20810-500uL	
CF@660C	667/685 nm	50 uL (50 ug)	20815-50uL	CF@660C Features
		0.5 mL (500 ug)	20815-500uL	
CF@680	681/698 nm	50 uL (50 ug)	20819-50uL	CF@680 Features
		0.5 mL (500 ug)	20819-500uL	
CF@750	755/777 nm	50 uL (50 ug)	20827-50uL	CF@750 Features
		0.5 mL (500 ug)	20827-500uL	

CF® Dye Secondary Antibodies, Single Label for STORM

Dye	Donkey Anti-Goat	Donkey Anti-Guinea Pig	Donkey Anti-Mouse	Donkey Anti-Rabbit	Goat Anti-Mouse	Goat Anti-Rabbit
Cross-adsorption	Ck, GP, Hs, Hu, Ms, Rb, Rt, SHm	Bv, Ck, Gt, Hs, Hu, Ms, Rb, Rt, Shp, SHm	Bv, Ck, Gt, GP, Hs, Hu, Rb, Rt, Shp, SHm	Bv, Ck, Gt, GP, Hs, Hu, Ms, Rt, Shp, SHm	Bv, Hs, Hu, Rb, Sw	Hu, Ms, Rt
CF@498	20992	20993	20994	20995	20996	20997
CF@505	20880	20881	20878	20879	20876	20877
CF@535ST			20823	20824	20821	20822
CF@568	20836	20838	20802	20803	20800	20801
CF@583R			20794	20795	20792	20793
CF@597R			20798	20799	20796	20797
CF@647	20829	20837	20810	20811	20808	20809
CF@660C			20815	20816	20812	20813
CF@680			20819	20820	20817	20818
CF@750			20827	20828	20825	20826

Bv: bovine; Ch: chicken; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Sh: sheep; SHm: Syrian hamster; Sw: swine; Rt: rat
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 [CF® Dye references](#).

1. Proc. SPIE 9331, (2015), DOI: [10.1117/12.2083209](#)

This datasheet was generated on March 31, 2025 at 09:39:15 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-cf-dye-storm/>

