



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: [techsupport@biotium.com](mailto:techsupport@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 (Bos taurus), 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
<a href="#">CF@498</a>	498/519 nm	50 uL (50 ug)	<a href="#">20994-50uL</a>	
		0.5 mL (500 ug)	<a href="#">20994-500uL</a>	
<a href="#">CF@505</a>	505/519 nm	50 uL (50 ug)	<a href="#">20878-50uL</a>	
		0.5 mL (500 ug)	<a href="#">20878-500uL</a>	
<a href="#">CF@535ST</a>	535/568 nm	50 uL (50 ug)	<a href="#">20823-50uL</a>	<a href="#">CF@535ST Features</a>
		0.5 mL (500 ug)	<a href="#">20823-500uL</a>	
<a href="#">CF@568</a>	562/583 nm	50 uL (50 ug)	<a href="#">20802-50uL</a>	<a href="#">CF@568 Features</a>
		0.5 mL (500 ug)	<a href="#">20802-500uL</a>	
<a href="#">CF@583R</a>	586/609 nm	50 uL (50 ug)	<a href="#">20794-50uL</a>	<a href="#">CF@583R Features</a>
		0.5 mL (500 ug)	<a href="#">20794-500uL</a>	
<a href="#">CF@597R</a>	597/619 nm	50 uL (50 ug)	<a href="#">20798-50uL</a>	<a href="#">CF@597R Features</a>
		0.5 mL (500 ug)	<a href="#">20798-500uL</a>	
<a href="#">CF@647</a>	650/665 nm	50 uL (50 ug)	<a href="#">20810-50uL</a>	<a href="#">CF@647 Features</a>
		0.5 mL (500 ug)	<a href="#">20810-500uL</a>	
<a href="#">CF@660C</a>	667/685 nm	50 uL (50 ug)	<a href="#">20815-50uL</a>	<a href="#">CF@660C Features</a>
		0.5 mL (500 ug)	<a href="#">20815-500uL</a>	
<a href="#">CF@680</a>	681/698 nm	50 uL (50 ug)	<a href="#">20819-50uL</a>	<a href="#">CF@680 Features</a>
		0.5 mL (500 ug)	<a href="#">20819-500uL</a>	
<a href="#">CF@750</a>	755/777 nm	50 uL (50 ug)	<a href="#">20827-50uL</a>	<a href="#">CF@750 Features</a>
		0.5 mL (500 ug)	<a href="#">20827-500uL</a>	

## CF® Dye Secondary Antibodies, Single Label for STORM

Dye	<a href="#">Donkey Anti-Goat</a>	<a href="#">Donkey Anti-Guinea Pig</a>	<a href="#">Donkey Anti-Mouse</a>	<a href="#">Donkey Anti-Rabbit</a>	<a href="#">Goat Anti-Mouse</a>	<a href="#">Goat Anti-Rabbit</a>
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
<a href="#">CF@498</a>	<a href="#">20992</a>	<a href="#">20993</a>	<a href="#">20994</a>	<a href="#">20995</a>	<a href="#">20996</a>	<a href="#">20997</a>
<a href="#">CF@505</a>	<a href="#">20880</a>	<a href="#">20881</a>	<a href="#">20878</a>	<a href="#">20879</a>	<a href="#">20876</a>	<a href="#">20877</a>
<a href="#">CF@535ST</a>			<a href="#">20823</a>	<a href="#">20824</a>	<a href="#">20821</a>	<a href="#">20822</a>
<a href="#">CF@568</a>	<a href="#">20836</a>	<a href="#">20838</a>	<a href="#">20802</a>	<a href="#">20803</a>	<a href="#">20800</a>	<a href="#">20801</a>
<a href="#">CF@583R</a>			<a href="#">20794</a>	<a href="#">20795</a>	<a href="#">20792</a>	<a href="#">20793</a>
<a href="#">CF@597R</a>			<a href="#">20798</a>	<a href="#">20799</a>	<a href="#">20796</a>	<a href="#">20797</a>
<a href="#">CF@647</a>	<a href="#">20829</a>	<a href="#">20837</a>	<a href="#">20810</a>	<a href="#">20811</a>	<a href="#">20808</a>	<a href="#">20809</a>
<a href="#">CF@660C</a>			<a href="#">20815</a>	<a href="#">20816</a>	<a href="#">20812</a>	<a href="#">20813</a>
<a href="#">CF@680</a>			<a href="#">20819</a>	<a href="#">20820</a>	<a href="#">20817</a>	<a href="#">20818</a>
<a href="#">CF@750</a>			<a href="#">20827</a>	<a href="#">20828</a>	<a href="#">20825</a>	<a href="#">20826</a>

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 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](#).

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



