

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



Highly cross-adsorbed donkey anti-rabbit 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-rabbit IgG (H L) secondary antibody that has a low degree of labeling (DOL) with one of our bright and photostable CF® Dyes.

- CF® Dye single label secondary antibody ideal for STORM imaging
- Highly cross-adsorbed for specific staining with minimal background
- Available in 9 bright and photostable CF® Dyes

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, mouse, rat, sheep, and Syrian hamster. Learn more about [CF® Dyes for super-resolution microscopy](#). 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.

Product attributes

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

References

Download a list of [CF® Dye references](#).

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@505	505/519 nm	50 uL (50 ug) 0.5 mL (500 ug)	20879-50uL 20879-500uL	
CF@535ST	535/568 nm	50 uL (50 ug) 0.5 mL (500 ug)	20824-50uL 20824-500uL	CF@535ST Features
CF@568	562/583 nm	50 uL (50 ug) 0.5 mL (500 ug)	20803-50uL 20803-500uL	CF@568 Features
CF@583R	586/609 nm	50 uL (50 ug) 0.5 mL (500 ug)	20795-50uL 20795-500uL	CF@583R Features
CF@597R	597/619 nm	50 uL (50 ug) 0.5 mL (500 ug)	20799-50uL 20799-500uL	CF@597R Features
CF@647	650/665 nm	50 uL (50 ug) 0.5 mL (500 ug)	20811-50uL 20811-500uL	CF@647 Features
CF@660C	667/685 nm	50 uL (50 ug) 0.5 mL (500 ug)	20816-50uL 20816-500uL	CF@660C Features
CF@680	681/698 nm	50 uL (50 ug) 0.5 mL (500 ug)	20820-50uL 20820-500uL	CF@680 Features
CF@750	755/777 nm	50 uL (50 ug) 0.5 mL (500 ug)	20828-50uL 20828-500uL	CF@750 Features