

# Product Information

## Goat Anti-Mouse IgG (H+L), Whole Antibody, Highly Cross-Adsorbed (min x Bv, Hs, Hu, Rb, Sw) CF® Dye Single Label Conjugates for (d)-STORM

CF® Dye	Ex/Em (nm)	Cat. No.	Unit Size
CF®505	505/519	20876-50uL	50 uL
		20876-500uL	500 uL
CF®535ST	535/569	20821-50uL	50 uL
		20821-500uL	500 uL
CF®568	562/584	20800-50uL	50 uL
		20800-500uL	500 uL
CF®583R	585/609	20792-50uL	50 uL
		20792-500uL	500 uL
CF®597R	597/619	20796-50uL	50 uL
		20796-500uL	500 uL
CF®647	652/668	20808-50uL	50 uL
		20808-500uL	500 uL
CF®660C	667/685	20812-50uL	50 uL
		20812-500uL	500 uL
CF®680	681/698	20817-50uL	50 uL
		20817-500uL	500 uL
CF®750	755/779	20825-50uL	50 uL
		20825-500uL	500 uL

Bv: bovine; Hs: horse; Hu: human; Rb: rabbit; Sw: swine

### Concentration

1 mg/mL in pH ~7.4 PBS containing 50% glycerol, 2 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

### Storage and Handling

Store at -20°C, protected from light. Product is stable for at least 12 months from date of receipt when stored as recommended. Antibodies contain 50% glycerol and will not freeze at -20°C.

**Note:** Storage of the antibody for more than a day at final working dilution is not recommended.

### Product Description

This product is prepared by single labeling (DOL=1) highly cross-adsorbed goat anti-mouse IgG (H+L) with CF® Dye for (d)-STORM super-resolution microscopy. This antibody is highly-cross adsorbed for multiple labeling, and has minimal cross-reactivity with bovine, horse, human, rabbit, and swine serum proteins.

CF® Dyes are a series of highly water-soluble fluorescent dyes spanning the visible and near-infrared spectrum for labeling biomolecules such as antibodies and nucleic acids. Developed in-house by our chemists, CF® Dyes offer superior brightness and signal-to-noise when compared to other commercially available fluorescent dyes. The CF® Dyes available with these secondary antibodies feature unique photoswitching properties that are ideal for STORM imaging. For more information about CF® Dyes, visit [www.biotium.com](http://www.biotium.com).

### Recommended Dilution Range

Fluorescence microscopy: 1-2 ug/mL

These concentrations are provided as a starting point for optimization; appropriate dilutions should be determined empirically. Generally IgG conjugates are used in the range of 1-10 ug/mL. For more detailed protocols for immunofluorescence staining for fluorescence microscopy, flow cytometry, or fluorescence-based western detection, please visit our website.

We do not currently offer detailed protocols for STORM and ask that customers look to the literature for guidance. For general protocols on immunofluorescence staining for fluorescence microscopy, flow cytometry, or fluorescence-based western detection, please visit our website.

### Related Products

Cat. No.	Product
20794... 20878	Donkey Anti-Mouse IgG (H+L), Highly Cross-Adsorbed, CF® Dye Single Label Conjugates for STORM
20793... 20877	Goat Anti-Rabbit IgG (H+L), Highly Cross-Adsorbed, CF® Dye Conjugates, Single Label for STORM
20795... 20879	Donkey Anti-Rabbit IgG (H+L), Highly Cross-Adsorbed, CF® Dye Conjugates, Single Label for STORM
20829... 20880	Donkey Anti-Goat IgG (H+L), Highly Cross-Adsorbed, CF® Dye Conjugates, Single Label for STORM
20837... 20881	Donkey Anti-Guinea Pig IgG (H+L), Highly Cross-Adsorbed, CF® Dye Conjugates, Single Label for STORM
00041, 00048	Phalloidin Conjugates (CF®647 and CF®680 for STORM)
30101-30104	MemBrite® Fix-ST membrane stains for STORM
30115-30118	ExoBrite™ STORM CTB EV Staining Kits
92500-92515	Mix-n-Stain™ Nanobody Labeling Kits
92549-92557	Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative
22030	AntiFix™ Universal Antigen Retrieval Buffer, 10X
22002	Tween®-20
23005	CoverGrip™ Coverslip Sealant
23007	TrueBlack® Lipofuscin Autofluorescence Quencher
23012	TrueBlack® IF Background Suppressor System (Permeabilizing)
29021...29073	Wheat Germ Agglutinin (WGA) Conjugates
40061	RedDot™2 Far Red Nuclear Counterstain
40083...41038	NucSpot® Nuclear Stains for dead or fixed cells

Please visit [www.biotium.com](http://www.biotium.com) to view our full selection of products featuring bright and photostable fluorescent CF® Dyes, including Mix-n-Stain™ antibody labeling kits, primary antibody conjugates, streptavidin, phalloidin, and other bioconjugates, as well as conjugates of biotin, HRP, AP, R-PE, APC and PerCP.

CF Dye technology is covered by US and international patents; TWEEN is a registered trademark of Uniqema Americas LLC. Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.