

Product Information

Goat Anti-Rabbit IgG (H+L), Whole Antibody, Highly Cross-Adsorbed (min x Hu, Ms, Rt) CF® Dye Single Label Conjugates for (d)-STORM

CF® Dye	Ex/Em (nm)	Cat. No.	Unit Size
CF®505	505/519	20877-50uL	50 uL
		20877-500uL	500 uL
CF®535ST	535/569	20822-50uL	50 uL
		20822-500uL	500 uL
CF®568	562/584	20801-50uL	50 uL
		20801-500uL	500 uL
CF®583R	586/609	20793-50uL	50 uL
		20793-500uL	500 uL
CF®597R	597/619	20797-50uL	50 uL
		20797-500uL	500 uL
CF®647	652/668	20809-50uL	50 uL
		20809-500uL	500 uL
CF®660C	667/685	20813-50uL	50 uL
		20813-500uL	500 uL
CF®680	681/698	20818-50uL	50 uL
		20818-500uL	500 uL
CF®750	755/779	20826-50uL	50 uL
		20826-500uL	500 uL

Hu: human; Ms: mouse; Rt: rat

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-rabbit 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 human, mouse, and rat 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.

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.

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
20792... 20876	Goat Anti-Mouse IgG (H+L), Highly Cross-Adsorbed, CF® Dye Conjugates, Single Label for STORM
20794... 20878	Donkey Anti-Mouse IgG (H+L), Highly Cross-Adsorbed, CF® Dye Single Label Conjugates 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 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.