

Product Information

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

For more detailed protocols for immunofluorescence staining for fluorescence microscopy, flow cytometry, or fluorescence-based western detection, please visit our website.

CF® Dye	Ex/Em	Catalog No.	Size
CF®568	562/583 nm	20836-50uL	50 uL
		20836-500uL	500 uL
CF®647	650/665 nm	20829-50uL	50 uL
		20829-500uL	500 uL

Ch: chicken; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Rt: rat; SHm: Syrian hamster

Concentration

Liquid format: 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. Liquid format 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, chicken, goat, guinea pig, horse, human, rabbit, sheep and syrian hamster proteins.

CF® dyes are superior to Alexa Fluor®, DyLight®, and Cy® dyes for antibody labeling by having combined advantages in brightness, photostability, specificity and novel features ideal for in vivo imaging. For more information about CF® dyes, download our CF® Dye Selection Guide at 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.

Accessory products for immunofluorescence

Cat. No.	Product Name
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
23007	TrueBlack™ Lipofuscin Autofluorescence Quencher
40061	RedDot™2 Far Red Nuclear Counterstain
22005	Mini Super ^{HT} Pap Pen 2.5 mm tip, ~400 uses
22006	Super ^{HT} Pap Pen 4 mm tip, ~800 uses
22003	Mini-Cell Scrapers
23006	Flow Cytometry Fixation/Permeabilization Kit
22015	Fixation Buffer
22016	Permeabilization Buffer
22017	Permeabilization and Blocking Buffer
22010	10X Fish Gelatin Blocking Agent
22011	Fish Gelatin Powder
22014	30% Bovine Serum Albumin Solution
22002	Tween®-20
22020	10X Phosphate-Buffered Saline (PBS)
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative

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.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

CF dye technology is covered by US and international patents. Alexa Fluor is a registered trademark of Molecular Probes; CYDYE is a registered trademark of GE Healthcare; DyLight is a registered trademark of Pierce Biotechnology; TWEEN is a registered trademark of Uniqema Americas LLC.