

# 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	Catalog No.	Size
CF@535ST	535/568 nm	20822-50uL	50 uL
		20822-500uL	500 uL
CF@568	562/583 nm	20801-50uL	50 uL
		20801-500uL	500 uL
CF@583R	586/609 nm	20793-50uL	50 uL
		20793-500uL	500 uL
CF@597R	597/619 nm	20797-50uL	50 uL
		20797-500uL	500 uL
CF@594ST	593/614 nm	20805-50uL	50 uL
		20805-500uL	500 uL
CF@647	650/665 nm	20809-50uL	50 uL
		20809-500uL	500 uL
CF@660C	667/685 nm	20813-50uL	50 uL
		20813-500uL	500 uL
CF@680	681/698 nm	20818-50uL	50 uL
		20818-500uL	500 uL
CF@750	755/777 nm	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 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](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.

### Related Products

Catalog Number	Product
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
23012	TrueBlack® IF Background Suppressor System (Permeabilizing)
23016	EverBrite™ Hardset Mounting Medium with NucSpot® 640
23017	EverBrite TrueBlack® Hardset Mounting Medium
23018	EverBrite TrueBlack® Hardset Mounting Medium with DAPI
23019	EverBrite TrueBlack® Hardset Mounting Medium with NucSpot® 640
40061	RedDot™2 Far Red Nuclear Counterstain
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ~400 uses
22006	Super <sup>HT</sup> 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
22030	AntiFix™ Universal Antigen Retrieval Buffer, 10X

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.

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.