www.biotium.com





Product Information

Goat Anti-Rabbit IgG (H+L), Whole Antibody, Highly Cross-Adsorbed (min x Hu, Ms, Rt)

CF® Dye	Ex/Em	Catalog No.	Size
CF®350	347/448 nm	20144-1	50 uL
		20144	500 uL
		20144-1mg	1 mg (lyophilized)
	408/452 nm	20373-50uL	50 uL
CF®405M		20373-500uL	500 uL
		20373-1mg	1 mg (lyophilized)
	426/498 nm	20460-50uL	50 uL
CF®430		20460-500uL	500 uL
		20460-1mg	1 mg (lyophilized)
	490/515 nm	20019-1	50 uL
CF®488A		20019	500 uL
		20019-1mg	1 mg (lyophilized)
	527/558 nm	20469-50uL	50 uL
CF®532		20469-500uL	500 uL
		20469-1mg	1 mg (lyophilized)
		20300-1	50 uL
CF®543	541/560 nm	20300	500 uL
		20300-1mg	1 mg (lyophilized)
	555/565 nm	20232-1	50 uL
CF®555		20232	500 uL
		20232-1mg	1 mg (lyophilized)
	562/583 nm	20103-1	50 uL
CF®568		20103	500 uL
		20103-1mg	1 mg (lyophilized)
	593/614 nm	20113-1	50 uL
CF®594		20113	500 uL
		20113-1mg	1 mg (lyophilized)
	630/650 nm	20123-1	50 uL
CF®633		20123	500 uL
		20123-1mg	1 mg (lyophilized)
	642/662 nm	20176-1	50 uL
CF®640R		20176	500 uL
		20176-1mg	1 mg (lyophilized)
CF®647	650/665 nm	20282-1	50 uL
		20282	500 uL
		20282-1mg	1 mg (lyophilized)
CF®660C	667/685 nm	20369-50uL	50 uL
		20369-500uL	500 uL
		20369-1mg	1 mg (lyophilized)

CF® Dye	Ex/Em	Catalog No.	Size
CF®680	681/698 nm	20067-1	50 uL
		20067	250 uL
CF®680R	680/701 nm	20193-1	50 uL
		20193	250 uL
CF®770	770/797 nm	20078-1	50 uL
		20078	250 uL
CF®790	784/806 nm	20343-50uL	50 uL
CF®800	797/816 nm	20832-50uL	50 uL
	N/A	20186-1	50 uL
Biotin		20186	500 uL
		20186-1mg	1 mg (lyophilized)
HRP	N/A	20403-100uL	100 uL
		20403-1mL	1 mL
		20403-1mg	1 mg (lyophilized)

Hu: human; Ms: mouse; Rt: rat

Concentration

Liquid format:

CFdye and Biotin: 2 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.

HRP: 1 mg/mL in pH \sim 7.4 PBS containing 50% glycerol and 15 mg/ml bovine serum albumin (IgG-free and protease-free).

Lyophilized format (after reconstitution):

CFdye and Biotin: 2 mg/mL in pH \sim 7.4 PBS containing 15 mg/mL bovine serum albumin (IgG-free and protease-free) and 20 mg/mL trehalose.

HRP: 1 mg/mL in pH ~7.4 PBS containing and 10 mg/mL trehalose.

Storage and Handling

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.

CF® Dye and biotin conjugates: add 0.5 mL dH2O HRP conjugate: add 1 mL dH2O

Add the indicated volume of water directly to the vial containing the lyophilized antibody and mix gently to dissolve. Store at -20°C and protect fluorescent conjugates 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.

Product Description

Goat anti-rabbit IgG (H+L), whole antibody, highly cross-adsorbed is an affinity purified goat antibody that recognizes rabbit IgG. This antibody is highly-cross adsorbed for multiple labeling, and has minimal cross-reactivity with human, mouse, or rat serum 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 Flow cytometry: 1 ug/10⁶ cells Near-infrared western detection: 50-100 ng/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.

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 ^{н⊤} Pap Pen 2.5 mm tip, ~400 uses	
22006	Super ^{н⊤} 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 pending US and international patents. Alexa Fluor is a registered trademark of Molecular Probes; CYDYE is a registered trademark of GE Healthcare; Draq7 is a registered trademark of Biostatus Ltd; DyLight is a registered trademark of Pierce Biotechnology; TWEEN is a registered trademark of Uniqema Americas LLC.