

Product Information

Goat Anti-Rat IgG (H+L), Whole Antibody, Highly Cross-Adsorbed (min x Bv, Hs, Hu, Rb)

CF® Dye	Ex/Em	Catalog No.	Size
CF@350	347/448 nm	20147-1	50 uL
		20147	500 uL
		20147-1mg	1 mg (lyophilized)
CF@405M	408/452 nm	20374-50uL	50 uL
		20374-500uL	500 uL
		20374-1mg	1 mg (lyophilized)
CF@488A	490/515 nm	20023-1	50 uL
		20023	500 uL
		20023-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20321-1	50 uL
		20321	500 uL
		20321-1mg	1 mg (lyophilized)
CF@555	555/565 nm	20233-1	50 uL
		20233	500 uL
		20233-1mg	1 mg (lyophilized)
CF@568	562/583 nm	20096-1	50 uL
		20096	500 uL
		20096-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20155-1	50 uL
		20155	500 uL
		20155-1mg	1 mg (lyophilized)
CF@633	630/650 nm	20133-1	50 uL
		20133	500 uL
		20133-1mg	1 mg (lyophilized)
CF@640R	642/662 nm	20088-1	50 uL
		20088	500 uL
		20088-1mg	1 mg (lyophilized)
CF@647	650/665 nm	20283-1	50 uL
		20283	500 uL
		20283-1mg	1 mg (lyophilized)

CF® Dye	Ex/Em	Catalog No.	Size
CF@660C	667/685 nm	20370-50uL	50 uL
		20370-500uL	500 uL
		20370-1mg	1 mg (lyophilized)
CF@680	681/698 nm	20069-1	50 uL
		20069	250 uL
CF@770	770/797 nm	20383-50uL	50 uL
		20383-250uL	250 uL
APC	633,640/660 nm	20413-100uL	100 uL
		20413-500uL	500 uL
R-PE	496, 546, 565/578 nm	20354-200uL	200 uL
		20354-1mL	1 mL

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

Concentration

Liquid format:

CF dye: 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.

R-PE: 0.5mg/mL in pH ~7.4 PBS containing 2 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

APC: 0.5 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.

Lyophilized format (after reconstitution): 2 mg/mL in pH ~7.4 PBS containing 15 mg/mL bovine serum albumin (IgG-free and protease-free) and 20 mg/mL trehalose.

Storage and Handling

Store at -20°C (except for R-PE conjugates which is stored at 4°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.

Reconstitution (lyophilized format only): add 0.5 mL dH₂O and mix gently to dissolve. Store at -20°C, protected from light. Aliquot to avoid freeze-thaw cycles. Alternatively, glycerol can be added to the antibody so that it will not freeze at -20°C: add 0.25 mL dH₂O to the lyophilized antibody and mix gently to dissolve, then add 0.25 mL glycerol and mix well. Optional: a preservative may be added, such as 0.05% (final concentration) sodium azide.

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

Product Description

Goat anti-rat IgG (H+L), whole antibody, highly cross-adsorbed is an affinity purified goat antibody that recognizes rat IgG. This antibody is highly-cross adsorbed for multiple labeling, and has minimal cross-reactivity with bovine, horse, human, or rabbit 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 ^{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

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.