

Revised: January 21, 2019

Product Information

Goat Anti-Rabbit IgG (H+L), Whole Antibody

CF® Dye	Ex/Em	Catalog No.	Size
CF®350		20141-1	50 uL
	347/448 nm	20141	500 uL
		20141-1mg	1 mg (lyophilized)
CF®405S		20082-1	50 uL
	404/431 nm	20082	500 uL
		20082-1mg	1 mg (lyophilized)
CF®405M	408/452 nm	20181-1	50 uL
		20181	500 uL
		20181-1mg	1 mg (lyophilized)
CF®405L	395/545 nm	20409-50uL	50 uL
		20409-500uL	500 uL
		20409-1mg	1 mg (lyophilized)
CF®488A	490/515 nm	20012-1	50 uL
		20012	500 uL
		20012-1mg	1 mg (lyophilized)
		20387-50uL	50 uL
CF®514	516/548 nm	20387-500uL	500 uL
		20387-1mg	1 mg (lyophilized)
		20366-50uL	50 uL
CF®532	527/558 nm	20366-500uL	500 uL
		20366-1mg	1 mg (lyophilized)
	541/560 nm	20309-1	50 uL
CF®543		20309	500 uL
		20309-1mg	1 mg (lyophilized)
	555/565 nm	20033-1	50 uL
CF®555		20033	500 uL
		20033-1mg	1 mg (lyophilized)
	562/583 nm	20102-1	50 uL
CF®568		20102	500 uL
		20102-1mg	1 mg (lyophilized)
	593/614 nm	20112-1	50 uL
CF®594		20112	500 uL
		20112-1mg	1 mg (lyophilized)
CF®633	630/650 nm	20122-1	50 uL
		20122	500 uL
		20122-1mg	1 mg (lyophilized)
	642/662 nm	20202-1	50 uL
CF®640R		20202	500 uL
		20202-1mg	1 mg (lyophilized)
CF®647	650/665 nm	20043-1	50 uL
		20043	500 uL
		20043-1mg	1 mg (lyophilized)
CF®660C	667/685 nm	20053-1	50 uL
		20053	500 uL
		20053-1mg	1 mg (lyophilized)

CF® Dye	Ex/Em	Catalog No.	Size
		20055-1	50 uL
CF®660R	663/682 nm	20055	500 uL
		20055-1mg	1 mg (lyophilized)
CF®750	755/777 nm	20073-1	50 uL
		20073	500 uL
		20073-1mg	1 mg (lyophilized)
CF®790	784/806 nm	20379-50uL	50 uL
	N/A	20185-1	50 uL
Biotin		20185	500 uL
		20185-1mg	1 mg (lyophilized)
APC	633,640/660	20412-100uL	100 uL
AFC	nm	20412-500uL	500 uL
AP	N/A	20465-100uL	100 uL
AF		20465-1mL	1 mL
HRP	N/A	20402-1mL	1 mL
		20402-100uL	100 uL
		20402-1mg	1 mg (lyophilized)
R-PE	496, 546,	20353-1mL	1 mL
R-PE	565/578 nm	20353-200uL	200 uL

Concentration

Liquid format:

CFdye and Biotin: 2 mg/mL in pH \sim 7.4 PBS containing 50% glycerol, 2 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

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

APC: 0.5 mg/mL in pH \sim 7.4 PBS containing 50% glycerol, 2 mg/ml bovine serum albumin (lgG-free and protease-free) and 0.05% sodium azide.

AP: 1 mg/mL in pH \sim 7.4 PBS containing 40% glycerol, 5 mg/ml bovine serum albumin (lgG-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 (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 dH $_2$ O 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 dH $_2$ O 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. Caution: 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 is an affinity purified goat antibody that recognizes rabbit IgG. This antibody is not highly-cross adsorbed against other species. Visit www.biotium.com to see our highly cross-adsorbed secondary antibodies for multiple labeling and isotype-specific secondary antibodies.

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.

Related Products

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.