

Product Information

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

Conjugate	Ex/Em	Catalog No.	Size
CF@350	347/448 nm	20140-1	50 uL
		20140	500 uL
		20140-1mg	1 mg (lyophilized)
CF@405S	404/431 nm	20080-1	50 uL
		20080	500 uL
		20080-1mg	1 mg (lyophilized)
CF@405M	408/452 nm	20180-1	50 uL
		20180	500 uL
		20180-1mg	1 mg (lyophilized)
CF@405L	395/545 nm	20408-50uL	50 uL
		20408-500uL	500 uL
		20408-1mg	1 mg (lyophilized)
CF@488A	490/515 nm	20010-1	50 uL
		20010	500 uL
		20010-1mg	1 mg (lyophilized)
CF@514	516/548 nm	20386-50uL	50 uL
		20386-500uL	500 uL
		20386-1mg	1 mg (lyophilized)
CF@532	527/558 nm	20365-50uL	50 uL
		20365-500uL	500 uL
		20365-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20306-1	50 uL
		20306	500 uL
		20306-1mg	1 mg (lyophilized)
CF@555	555/565 nm	20030-1	50 uL
		20030	500 uL
		20030-1mg	1 mg (lyophilized)
CF@568	562/583 nm	20100-1	50 uL
		20100	500 uL
		20100-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20110-1	50 uL
		20110	500 uL
		20110-1mg	1 mg (lyophilized)
CF@633	630/650 nm	20120-1	50 uL
		20120	500 uL
		20120-1mg	1 mg (lyophilized)
CF@640R	642/662 nm	20197-1	50 uL
		20197	500 uL
		20197-1mg	1 mg (lyophilized)
CF@647	650/665 nm	20040-1	50 uL
		20040	500 uL
		20040-1mg	1 mg (lyophilized)
CF@660C	667/685 nm	20050-1	50 uL
		20050	500 uL
		20050-1mg	1 mg (lyophilized)
CF@660R	663/682 nm	20054-1	50 uL
		20054	500 uL
		20054-1mg	1 mg (lyophilized)

Conjugate	Ex/Em	Catalog No.	Size
CF@750	755/777 nm	20070-1	50 uL
		20070	500 uL
		20070-1mg	1 mg (lyophilized)
CF@790	784/806 nm	20378-50uL	50 uL
Biotin	N/A	20183-1	50 uL
		20183	500 uL
		20183-1mg	1 mg (lyophilized)
APC	633,640/660 nm	20411-500uL	500 uL
		20411-100uL	100 uL
AP	N/A	20464-1mL	1 mL
		20464-100uL	100 uL
HRP	N/A	20400-1mL	1 mL
		20400-100uL	100 uL
		20400-1mg	1 mg (lyophilized)
R-PE	496, 546, 565/578 nm	20352-1mL	1 mL
		20352-200uL	200 uL

Concentration

Liquid format:

CF@ dye and biotin conjugates: 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 conjugates: 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 conjugates: 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.

AP conjugates: 1 mg/mL in pH ~7.4 PBS containing 40% glycerol, 5 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

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

Lyophilized format (after reconstitution):

CF@ dye and biotin conjugates: 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.

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

Storage and Handling

Store R-PE conjugates at 4°C, protected from light. Store all other conjugates 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 that contain glycerol 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. 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.

Continued next page

Product Description

Goat anti-mouse IgG (H+L), whole antibody is an affinity purified goat antibody that recognizes mouse 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.

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, DyLight, and Image-iT are registered trademarks of Thermo Fisher Scientific; Cubitainer is a registered trademark of Hedwin Corporation. CY DYE is a registered trademark of GE Healthcare; Draq7 is a registered trademark of Biostatus Ltd; Odyssey is a registered trademark of LI-COR Biosciences.

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	Features
23012	TrueBlack® IF Background Suppressor System (Permeabilizing)	<ul style="list-style-type: none"> • Suppresses background from non-specific antibody binding and charged fluorescent dyes • More efficient than Image-iT® FX, block & permeabilize in just 10 minutes • Complete system for blocking, permeabilizing, and antibody dilution • Non-mammalian blocking agents, for broad secondary antibody compatibility • For immunofluorescence on cells or tissue sections
23013	TrueBlack® WB Blocking Buffer Kit	<ul style="list-style-type: none"> • Blocks as well or better than Odyssey® Blocking Buffer, at a lower price • Reduces non-specific protein bands and background over entire membrane • Suppresses background from charged dyes better than BSA, gelatin, or casein • Compatible with PVDF and nitrocellulose membranes • Contains no mammalian proteins, for broad antibody compatibility • For visible and near-IR fluorescent westerns
23007	TrueBlack® Lipofuscin Autofluorescence Quencher	<ul style="list-style-type: none"> • Eliminates lipofuscin autofluorescence with less background than Sudan Black B • Reduces background from other sources • Can be used before or after IF staining
23001	EverBrite™ Mounting Medium	<ul style="list-style-type: none"> • Excellent protection from photobleaching • Compatible with a wide variety of dyes including cyanine dyes • Available in wet-set or hardset formulations • With or without DAPI
23002	EverBrite™ Mounting Medium with DAPI	
23003	EverBrite™ Hardset Mounting Medium	
23004	EverBrite™ Hardset Mounting Medium with DAPI	
23008	Drop-n-Stain EverBrite™ Mounting Medium	<ul style="list-style-type: none"> • Convenient dropper bottle for pipette-free mounting • Protects fluorescent dyes from rapid photobleaching • Compatible with Cy®3, Cy®5, and Alexa Fluor® 647, unlike VECTASHIELD® • Available with our without DAPI for nuclear counterstaining • Works as well as original EverBrite™, but less viscous for rapid DAPI staining
23009	Drop-n-Stain EverBrite™ Mounting Medium with DAPI	
23005	CoverGrip™ Coverslip Sealant	<ul style="list-style-type: none"> • Replaces nail polish for coverslip sealing • Won't mix with aqueous mounting media
40061	RedDot™2 Far Red Nuclear Counterstain	<ul style="list-style-type: none"> • Far-red nuclear dye for the Cy®5 channel • More specific than Draq®7
40083	NucSpot® 470 Nuclear Stain,	<ul style="list-style-type: none"> • Green fluorescent nuclear-specific counterstain for fixed cells or tissues
22005	Mini Super ^{HT} Pap Pen 2.5 mm tip, ~400 uses	<ul style="list-style-type: none"> • Create hydrophobic barriers around tissue sections • Heat-stable to 120°C • Insoluble in aqueous buffers, detergents, alcohol and acetone; can be removed with xylene
22006	Super ^{HT} Pap Pen 4 mm tip, ~800 uses	
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative	<ul style="list-style-type: none"> • Ready-to-use fixation buffer • Methanol-free, prepared from EM grade paraformaldehyde • No glass ampoules to break, store in the original bottle
22020	10X Phosphate Buffered Saline (PBS) 4L Cubitainer®	<ul style="list-style-type: none"> • Convenient buffers and blocking agents for immunofluorescence or western
22010	10X Fish Gelatin Blocking Agent	
22014	30% Bovine Serum Albumin Solution	

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.