

Product Information

Goat Anti-Guinea Pig IgG (H+L), Whole Antibody, Highly Cross-Adsorbed (min x Bv, Ch, Gt, Hs, Hu, Ms, Rb, Rt, Sh, SHm)

CF® Dye	Ex/Em	Catalog No.	Size
CF@405M	408/452 nm	20487-50uL	50 uL
		20487-500uL	500 uL
		20487-1mg	1 mg (lyophilized)
CF@405S	404/431 nm	20488-50uL	50 uL
		20488-500uL	500 uL
		20488-1mg	1 mg (lyophilized)
CF@488A	490/515 nm	20489-50uL	50 uL
		20489-500uL	500 uL
		20489-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20490-50uL	50 uL
		20490-500uL	500 uL
		20490-1mg	1 mg (lyophilized)
CF@555	555/565 nm	20491-50uL	50 uL
		20491-500uL	500 uL
		20491-1mg	1 mg (lyophilized)
CF@568	562/583 nm	20492-50uL	50 uL
		20492-500uL	500 uL
		20492-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20493-50uL	50 uL
		20493-500uL	500 uL
		20493-1mg	1 mg (lyophilized)
CF@640R	642/662 nm	20494-50uL	50 uL
		20494-500uL	500 uL
		20494-1mg	1 mg (lyophilized)
CF@647	650/665 nm	20495-50uL	50 uL
		20495-500uL	500 uL
		20495-1mg	1 mg (lyophilized)
CF@660R	663/682 nm	20496-50uL	50 uL
		20496-500uL	500 uL
		20496-1mg	1 mg (lyophilized)
CF@660C	667/685 nm	20497-50uL	50 uL
		20497-500uL	500 uL
		20497-1mg	1 mg (lyophilized)
CF@680R	680/701 nm	20498-50uL	50 uL
		20498-500uL	500 uL
CF@680	681/698 nm	20499-50uL	50 uL
		20499-500uL	500 uL
CF@770	770/797 nm	20500-50uL	50 uL
		20500-500uL	500 uL

Bv: Bovine; Ch: chicken; Gt: goat; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Rt: rat; Sh: Sheep; SHm: Syrian hamster

Concentration

Liquid format:

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.

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, 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-guinea pig IgG (H+L), whole antibody, highly cross-adsorbed is an affinity purified goat antibody that recognizes guinea pig IgG. This antibody is highly-cross adsorbed for multiple labeling, and has minimal cross-reactivity with bovine, chicken, goat, Syrian hamster, horse, human, mouse, rabbit, rat and sheep serum proteins and non-immunoglobulin guinea pig 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.

Continued next page

Accessory products for immunofluorescence

Cat. No.	Product Name	Features
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 cyanine dyes like 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	
23005	CoverGrip™ Coverslip Sealant	<ul style="list-style-type: none"> • Replaces nail polish for coverslip sealing • Won't mix with aqueous mounting media
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
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, just 10 minutes for blocking & permeabilization • Complete system for blocking, permeabilizing, and antibody dilution • Non-mammalian blocking agents, for broad secondary antibody compatibility • Use together with TrueBlack@ Lipofuscin Autofluorescence Quencher to block antibody background and autofluorescence • For staining of cells or tissue sections
23013	TrueBlack@ WB Blocking Buffer Kit	<ul style="list-style-type: none"> • Blocks non-specific protein binding and background membrane fluorescence • Suppresses background caused by charged fluorescent dyes • Works as well or better than LI-COR's Odyssey@ Blocking Buffer, at a lower cost • Can be used with all types of fluorophores spanning the visible and NIR spectrum • Contains non-mammalian blocking agents, for broad secondary antibody compatibility • TBS based buffer suitable for phosphoprotein detection • Ready-to-use, no need to add detergents or optimize detergent concentration
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
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	
22003	Mini-Cell Scrapers	<ul style="list-style-type: none"> • Collect cells from 96-well to 24-well plates • Polypropylene, sterile, disposable
23006	Flow Cytometry Fixation/Permeabilization Kit	<ul style="list-style-type: none"> • Ready-to-use fix/perm buffers for intracellular staining for flow cytometry
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative	<ul style="list-style-type: none"> • EM-grade paraformaldehyde fixative, packaged under argon with no methanol preservative
22016	Permeabilization Buffer	<ul style="list-style-type: none"> • Convenient buffers and blocking agents for immunofluorescence or western
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)	

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.