

Product Information

Donkey Anti-Rat IgG (H+L), Whole Antibody, Highly Cross-Adsorbed (min x Bv, Ch, Gt, GP, Hs, Hu, Ms, Rb, Sh, SHm)

CF® Dye	Ex/Em	Catalog No.	Size
CF@350	347/448 nm	20361-50uL	50 uL
		20361-500uL	500 uL
		20361-1mg	1 mg (lyophilized)
CF@405S	404/431 nm	20419-50uL	50 uL
		20419-500uL	500 uL
		20419-1mg	1 mg (lyophilized)
CF@488A	490/515 nm	20027-1	50 uL
		20027	500 uL
		20027-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20320-1	50 uL
		20320	500 uL
		20320-1mg	1 mg (lyophilized)
CF@568	562/583 nm	20092-1	50 uL
		20092	500 uL
		20092-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20159-1	50 uL
		20159	500 uL
		20159-1mg	1 mg (lyophilized)
CF@633	630/650 nm	20137-1	50 uL
		20137	500 uL
		20137-1mg	1 mg (lyophilized)
CF@640R	642/662 nm	20199-1	50 uL
		20199	500 uL
		20199-1mg	1 mg (lyophilized)
CF@647	650/665 nm	20843-50uL	50 uL
		20843-500uL	500 uL
		20843-1mg	1 mg (lyophilized)
CF@660R	663/682 nm	20390-50uL	50 uL
		20390-500uL	500 uL
		20390-1mg	1 mg (lyophilized)
CF@680	681/698 nm	20417-50uL	50 uL
		20417-250uL	250 uL
CF@750	755/777 nm	20857-50uL	50 uL
		20857-250uL	250uL

Bv: bovine; Ch: chicken; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; 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

Donkey anti-rat IgG (H+L), whole antibody, highly cross-adsorbed is an affinity purified donkey antibody that recognizes rat IgG. This antibody is highly-cross adsorbed for multiple labeling, and has minimal cross-reactivity with bovine, chicken, guinea pig, Syrian hamster, horse, human, mouse, rabbit, and sheep 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, download our Protocols for Antibody-Based Detection from the secondary antibody product webpage.

Accessory products for immunofluorescence

Cat. No.	Product Name	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 or 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	
22003	Mini-Cell Scrapers	<ul style="list-style-type: none"> • Collect cells from 96-well to 24-well plates • Polypropylene, sterile, disposable
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
23006	Flow Cytometry Fixation/Permeabilization Kit	<ul style="list-style-type: none"> • Ready-to-use fix/perm buffers for intracellular staining
22020	10X Phosphate Buffered Saline (PBS) 4L Cubitainer®	<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	

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, 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; TWEEN is a registered trademark of Uniqema Americas LLC.