

Product Information

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

Conjugate	Ex/Em	Catalog No.	Size
CF@350	347/448 nm	20350-1	50 uL
		20350	500 uL
		20350-1mg	1 mg (lyophilized)
CF@430	426/498 nm	20461-50uL	50 uL
		20461-500uL	500 uL
		20461-1mg	1 mg (lyophilized)
CF@488A	490/515 nm	20014-1	50 uL
		20014	500 uL
		20014-1mg	1 mg (lyophilized)
CF@514	516/548 nm	20483-50uL	50 uL
		20483-500uL	500 uL
		20483-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20305-1	50 uL
		20305	500 uL
		20305-1mg	1 mg (lyophilized)
CF@555	555/565 nm	20037-1	50 uL
		20037	500 uL
		20037-1mg	1 mg (lyophilized)
CF@568	562/583 nm	20105-1	50 uL
		20105	500 uL
		20105-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20115-1	50 uL
		20115	500 uL
		20115-1mg	1 mg (lyophilized)
CF@633	630/650 nm	20124-1	50 uL
		20124	500 uL
		20124-1mg	1 mg (lyophilized)
CF@640R	642/662 nm	20177-1	50 uL
		20177	500 uL
		20177-1mg	1 mg (lyophilized)
CF@647	650/665 nm	20046-1	50 uL
		20046	500 uL
		20046-1mg	1 mg (lyophilized)

Conjugate	Ex/Em	Catalog No.	Size
CF@660R	663/682 nm	20388-50uL	50 uL
		20388-500uL	500 uL
		20388-1mg	1 mg (lyophilized)
CF@680R	680/701 nm	20194-1	50 uL
		20194	500 uL
CF@790	784/806 nm	20363-50uL	50 uL
CF@800	797/816 nm	20835-50uL	50 uL
Biotin	N/A	20190-1	50 uL
		20190	500 uL
		20190-1mg	1 mg (lyophilized)
AP	N/A	20466-100uL	100 uL
		20466-1mL	1 mL
HRP	N/A	20404-100uL	100 uL
		20404-1mL	1 mL
		20404-1mg	1 mg (lyophilized)

Bv: bovine; Ch: chicken; Gt: goat; GP: guinea pig; Hs: horse; Hu: human; Rb: rabbit; Rt: rat; Sh: sheep; SHm: Syrian hamster

Concentration

Liquid format:

CF@ dye and Biotin: 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.

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: 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: 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: 1 mg/mL in pH ~7.4 PBS containing and 10 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):

CF® Dye and biotin conjugates: add 0.5 mL dH2O

HRP conjugate: add 1 mL dH2O

Add the indicated volume of water directly to the vial containing the lyophilized antibody and mix gently to dissolve. Store at -20°C and protect fluorescent conjugates from light. Aliquot to avoid repeated freeze/thaw cycles. Alternatively, an equal volume of glycerol can be mixed with the reconstituted antibody so that it will remain liquid at -20°C. Optional: a preservative such as 0.05% sodium azide (final concentration) can be added to CF® dye and biotin conjugates. 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

Donkey Anti-Mouse IgG (H+L) (Min x Rat), Whole Antibody, Highly Cross-Adsorbed is an affinity purified donkey antibody that recognizes mouse IgG. This antibody is highly-cross adsorbed for multiple labeling, and has minimal cross-reactivity with bovine, chicken, goat, guinea pig, Syrian hamster, horse, human, rabbit, rat, 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, please visit our website.

Related Products

Catalog Number	Product
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
22020	10X Phosphate-Buffered Saline (PBS)
22023	Paraformaldehyde, 4% in PBS, Ready-To-Use Fixative

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.