

Product Information

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

CF® Dye	Ex/Em	Catalog No.	Size
CF@350	347/448 nm	20142-1	50 uL
		20142	500 uL
		20142-1mg	1 mg (lyophilized)
CF@405S	404/431 nm	20416-50uL	50 uL
		20416-500uL	500 uL
		20416-1mg	1 mg (lyophilized)
CF@488A	490/515 nm	20016-1	50 uL
		20016	500 uL
		20016-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20314-1	50 uL
		20314	500 uL
		20314-1mg	1 mg (lyophilized)
CF@555	555/565 nm	20039-1	50 uL
		20039	500 uL
		20039-1mg	1 mg (lyophilized)
CF@568	562/583 nm	20106-1	50 uL
		20106	500 uL
		20106-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20116-1	50 uL
		20116	500 uL
		20116-1mg	1 mg (lyophilized)
CF@633	630/650 nm	20127-1	50 uL
		20127	500 uL
		20127-1mg	1 mg (lyophilized)
CF@640R	642/662 nm	20179-1	50 uL
		20179	500 uL
		20179-1mg	1 mg (lyophilized)
CF@647	650/665 nm	20048-1	50 uL
		20048	500 uL
		20048-1mg	1 mg (lyophilized)
CF@660C	667/685 nm	20051-1	50 uL
		20051	500 uL
		20051-1mg	1 mg (lyophilized)

CF® Dye	Ex/Em	Catalog No.	Size
CF@660R	663/682 nm	20391-50uL	50 uL
		20391-500ul	500 uL
		20391-1mg	1 mg (lyophilized)
CF@680	681/698 nm	20060-1	50 uL
		20060	250 uL
CF@680R	680/701 nm	20196-1	50 uL
		20196	500 uL
CF@750	755/777 nm	20362-50uL	50 uL
		20362-250uL	250 uL
CF@770	770/797 nm	20277-1	50 uL
		20277	500 uL
CF@790	784/806 nm	20345-50uL	50 uL
CF@800	797/816 nm	20834-50uL	50uL
Biotin	N/A	20189-1	50 uL
		20189	500 uL
		20189-1mg	1 mg (lyophilized)

Ch: chicken; GP: guinea pig; Hs: horse; Hu: human; Ms: mouse; Rb: rabbit; Rt: rat; SHm: Syrian hamster

Concentration

Liquid format:

CFdye 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.

Lyophilized format (after reconstitution):

CFdye 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.

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. 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

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

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
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.