

Product Information

Bovine Anti-Goat IgG (H+L), Whole Antibody, CF® Dye Conjugates

CF® Dye	Ex/Em	Catalog No.	Size
CF@488A	490/515 nm	20293-1	50 uL
		20293	500 uL
		20293-1mg	1 mg (lyophilized)
CF@543	541/560 nm	20313-1	50 uL
		20313	500 uL
		20313-1mg	1 mg (lyophilized)
CF@568	562/583 nm	20294-1	50 uL
		20294	500 uL
		20294-1mg	1 mg (lyophilized)
CF@594	593/614 nm	20295-1	50 uL
		20295	500 uL
		20295-1mg	1 mg (lyophilized)
CF@633	630/650 nm	20296-1	50 uL
		20296	500 uL
		20296-1mg	1 mg (lyophilized)
CF@640R	642/662 nm	20297-1	50 uL
		20297	500 uL
		20297-1mg	1 mg (lyophilized)

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

Bovine anti-goat IgG (H+L), whole antibody is an affinity purified bovine antibody that recognizes goat IgG. To minimize cross-reactivity, the antibody is highly cross-adsorbed against bovine, chicken, guinea pig, horse, human, mouse, rabbit, rat, and Syrian hamster 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

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.