

Product Information

Goat Anti-Mouse IgG1 F(ab'), Fc gamma specific

Conjugate	Ex/Em (nm)	Catalog no.	Unit size
CF@488A	490/515	20872-50uL	50 uL
CF@594	593/614	20873-50uL	50 uL

Concentration

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.

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. Antibodies contain 50% glycerol and will not freeze at -20°C.

Note: Storage of the antibody for more than a day at final working dilution is not recommended.

Product Description

This antibody is the F(ab') fragment of goat anti-mouse IgG1 Fc gamma-specific antibody labeled with CF@ dyes. Reacts with the Fc domain, but not the Fab portion, of mouse IgG1. May cross-react with other mouse IgG subclasses or immunoglobulins from other species. This antibody is not highly-cross adsorbed against other species. Visit www.biotium.com to see our highly cross-adsorbed secondary antibodies for multiple labeling.

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.

Accessory products for immunofluorescence

Cat. No.	Product	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	
40009, 40011, 40043	DAPI	<ul style="list-style-type: none"> • Popular UV (350 nm) excitable blue nuclear counterstains • Available in solid and liquid formats
40044, 40045,	Hoechst 33258	
40046, 40047	Hoechst 33342	
23005	CoverGrip™ Coverslip Sealant	<ul style="list-style-type: none"> • Replaces nail polish for coverslip sealing • Won't mix with aqueous mounting media
23014-T, 23014	TrueBlack® Plus Lipofuscin Autofluorescence Quencher	<ul style="list-style-type: none"> • Eliminates lipofuscin autofluorescence with less background than Sudan Black B • Reduces background from other sources • TrueBlack® Plus is the only quencher that works in aqueous buffer
23007	TrueBlack® Lipofuscin Autofluorescence Quencher	
23012	TrueBlack® IF Background Suppressor (Permeabilizing)	<ul style="list-style-type: none"> • Reduce background from non-specific antibody binding and charged fluorescent dyes • Block and permeabilize in just 10 minutes
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 • Convenient buffers and blocking agents for immunofluorescence or western
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	
22013	Bovine Serum Albumin Fraction V	
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.