

# Product Information

## Goat Anti-Mouse IgG (H+L) F(ab')<sub>2</sub> Fragment

CF® Dye	Ex/Em	Catalog No.	Size
CF@350	347/448 nm	20145-1	50 uL
		20145	0.25 mL
CF@488A	490/515 nm	20011-1	50 uL
		20011	0.25 mL
CF@543	541/560 nm	20329-1	50 uL
		20329	0.25 mL
CF@555	555/565 nm	20032-1	50 uL
		20032	0.25 mL
CF@568	562/583 nm	20109-1	50 uL
		20109	0.25 mL
CF@594	593/614 nm	20119-1	50 uL
		20119	0.25 mL
CF@633	630/650 nm	20130-1	50 uL
		20130	0.25 mL
CF@640R	642/662 nm	20086-1	50 uL
		20086	0.25 mL
CF@647	650/665 nm	20042-1	50 uL
		20042	0.25 mL
CF@680	681/698 nm	20063-1	50 uL
		20063	0.25 mL
Biotin	N/A	20188-1	50 uL
		20188	0.25 mL
APC	633, 640/660 nm	20415-100uL	100 uL
		20415-500uL	500 uL
R-PE	496, 546, 565/578 nm	20357-200uL	200 uL
		20357-500uL	500 uL

### Concentration

CF dyes 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.

R-PE: 0.5 mg/mL in pH ~7.4 PBS containing 2 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

APC: 0.5 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 (except for R-PE conjugates which are stored at 4°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')<sub>2</sub> fragment of affinity purified goat antibody that recognizes mouse IgG. This antibody is not highly-cross adsorbed against other species. Visit [www.biotium.com](http://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](http://www.biotium.com).

### Recommended Dilution Range

Fluorescence microscopy: 1-2 ug/mL

Flow cytometry: 1 ug/10<sup>6</sup> 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 Name	Features
23001	EverBrite™ Mounting Medium	<ul style="list-style-type: none"> <li>• Excellent protection from photobleaching</li> <li>• Compatible with a wide variety of dyes including cyanine dyes</li> <li>• Available in wet-set or hardset formulations</li> <li>• With or without DAPI</li> </ul>
23002	EverBrite™ Mounting Medium with DAPI	
23003	EverBrite™ Hardset Mounting Medium	
23004	EverBrite™ Hardset Mounting Medium with DAPI	
23005	CoverGrip™ Coverslip Sealant	<ul style="list-style-type: none"> <li>• Replaces nail polish for coverslip sealing</li> <li>• Won't mix with aqueous mounting media</li> </ul>
23007	TrueBlack™ Lipofuscin Autofluorescence Quencher	<ul style="list-style-type: none"> <li>• Eliminates lipofuscin autofluorescence with less background than Sudan Black B</li> <li>• Reduces background from other sources</li> <li>• Can be used before or after IF staining</li> </ul>
40061	RedDot™2 Far Red Nuclear Counterstain	<ul style="list-style-type: none"> <li>• Far-red nuclear dye for the Cy®5 channel</li> <li>• More specific than Draq®7</li> </ul>
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ~400 uses	<ul style="list-style-type: none"> <li>• Create hydrophobic barriers around tissue sections</li> <li>• Heat-stable to 120°C</li> <li>• Insoluble in aqueous buffers, detergents, alcohol and acetone; can be removed with xylene</li> </ul>
22006	Super <sup>HT</sup> Pap Pen 4 mm tip, ~800 uses	
22003	Mini-Cell Scrapers	<ul style="list-style-type: none"> <li>• Collect cells from 96-well to 24-well plates</li> <li>• Polypropylene, sterile, disposable</li> </ul>
23006	Flow Cytometry Fixation/Permeabilization Kit	<ul style="list-style-type: none"> <li>• Ready-to-use fix/perm buffers for intracellular staining</li> </ul>
22015	Fixation Buffer	<ul style="list-style-type: none"> <li>• Convenient buffers and blocking agents for immunofluorescence or western</li> </ul>
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](http://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.