

# Product Information

## Goat Anti-Mouse IgM (mu Chain), Whole Antibody

CF® Dye	Ex/Em	Catalog No.	Size
CF@488A	490/515 nm	20840-50uL	50 uL
		20840-250uL	250 uL
CF@555	555/565 nm	20485-50uL	50 uL
		20485-250uL	250 uL
CF@680	681/698 nm	20384-50uL	50 uL
		20384-250uL	250 uL
CF@680R	680/701 nm	20841-50uL	50 uL
		20841-250uL	250 uL
CF@770	770/797 nm	20385-50uL	50 uL
		20385-250uL	250 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. Liquid format 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

Goat Anti-Mouse IgM ( $\mu$  chain) CF® dye conjugates are prepared from affinity purified antibodies, and are suitable for immunofluorescence and other standard immunoassay applications. The antibodies react specifically with mouse IgM heavy chains ( $\mu$  chains) and not with immunoglobulin light chains.

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.

Continued next page

## 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	
23008	Drop-n-Stain EverBrite™ Mounting Medium	<ul style="list-style-type: none"> <li>• Convenient dropper bottle for pipette-free mounting</li> <li>• Protects fluorescent dyes from rapid photobleaching</li> <li>• Compatible with cyanine dyes like Cy@3, Cy@5, and Alexa Fluor@ 647, unlike VECTASHIELD@</li> <li>• Available with our without DAPI for nuclear counterstaining</li> <li>• Works as well as original EverBrite™, but less viscous for rapid DAPI staining</li> </ul>
23009	Drop-n-Stain EverBrite™ Mounting Medium	
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>
23012	TrueBlack@ IF Background Suppressor System (Permeabilizing)	<ul style="list-style-type: none"> <li>• Suppresses background from non-specific antibody binding and charged fluorescent dyes</li> <li>• More efficient than Image-iT@ FX, just 10 minutes for blocking &amp; permeabilization</li> <li>• Complete system for blocking, permeabilizing, and antibody dilution</li> <li>• Non-mammalian blocking agents, for broad secondary antibody compatibility</li> <li>• Use together with TrueBlack@ Lipofuscin Autofluorescence Quencher to block antibody background and autofluorescence</li> <li>• For staining of cells or tissue sections</li> </ul>
23013	TrueBlack@ WB Blocking Buffer Kit	<ul style="list-style-type: none"> <li>• Blocks non-specific protein binding and background membrane fluorescence</li> <li>• Suppresses background caused by charged fluorescent dyes</li> <li>• Works as well or better than LI-COR's Odyssey@ Blocking Buffer, at a lower cost</li> <li>• Can be used with all types of fluorophores spanning the visible and NIR spectrum</li> <li>• Contains non-mammalian blocking agents, for broad secondary antibody compatibility</li> <li>• TBS based buffer suitable for phosphoprotein detection</li> <li>• Ready-to-use, no need to add detergents or optimize detergent concentration</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 for flow cytometry</li> </ul>
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative	<ul style="list-style-type: none"> <li>• EM-grade paraformaldehyde fixative, packaged under argon with no methanol preservative</li> </ul>
22016	Permeabilization Buffer	<ul style="list-style-type: none"> <li>• Convenient buffers and blocking agents for immunofluorescence or western</li> </ul>
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)	

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.