

# Product Information

## CF® Dye Conjugated Monoclonal Mouse Anti-Biotin Antibodies

Catalog No.	Unit Size	Product Description
20203-1	50 uL	CF@405S Monoclonal Mouse Anti-Biotin
20203	0.25 mL	
20204-1	50 uL	CF@488A Monoclonal Mouse Anti-Biotin
20204	0.25 mL	
20205-1	50 uL	CF@594 Monoclonal Mouse Anti-Biotin
20205	0.25 mL	
20206-1	50 uL	CF@633 Monoclonal Mouse Anti-Biotin
20206	0.25 mL	
20207-1	50 uL	CF@640R Monoclonal Mouse Anti-Biotin
20207	0.25 mL	
20501-50uL	50 uL	CF@750 Monoclonal Mouse Anti-Biotin
20501-150uL	0.25 mL	
20367-50uL	50 uL	CF@770 Monoclonal Mouse Anti-Biotin
20367-250uL	0.25 mL	

**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.

**Clone:** 3D6.6

**Isotype:** Mouse IgG1, kappa

### Storage and Handling

Product is stable for at least 6 months at -20°C as an undiluted liquid. Storage of the antibody for more than a day at final working dilution is not recommended. Protect from light.

### Spectral Properties

$\lambda_{abs}/\lambda_{em}$  maxima in pH 7.4 PBS buffer

Conjugate	Abs (nm)	Em (nm)
CF@405S	404	431
CF@488A	490	515
CF@594	593	614
CF@633	630	650
CF@640R	642	662
CF@750	755	777
CF@770	770	797

### Product Description

Monoclonal mouse anti-biotin antibodies react with biotin conjugated to antibodies or other proteins. These monoclonal mouse anti-biotin antibodies are conjugated to Biotium's bright and photostable CF® dyes, and offer an alternative to streptavidin conjugates for detecting biotinylated probes.

### 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 antibody conjugates are used in the range of 1-10 ug/mL.

Monoclonal anti-biotin can be used as a primary antibody for staining of biotinylated proteins or other biotinylated targets in cells or tissue samples, or as a secondary antibody for detecting biotinylated primary antibodies.

For more detailed protocols for immunofluorescence staining for fluorescence microscopy, flow cytometry, or fluorescence-based western detection, please visit our website.

### Blocking of endogenous biotin

Cells or tissues contain endogenous biotin that can react with anti-biotin antibodies. We highly recommend blocking endogenous biotin to reduce background.

### Materials required but not provided

- Biotin blocking wash buffer: 1% (w/w) BSA and 0.05% (v/v) Tween® 20 in PBS
- Streptavidin, 0.1 mg/mL in biotin blocking wash buffer
- Biotin, 0.5 mg/mL in biotin blocking wash buffer

### Protocol

1. Fix, permeabilize, and perform immunofluorescence blocking of your cell or tissue samples following your preferred protocol.
2. Incubate samples with streptavidin solution for 15 minutes at room temperature.
3. Wash samples 3 x 5 minutes with biotin blocking wash buffer.
4. Incubate samples with biotin solution for 30 minutes at room temperature.
5. Wash samples 3 x 5 minutes with biotin blocking wash buffer.
6. Proceed to antibody incubation and washing according to your preferred immunofluorescence protocol.

## Related Products

Cat. No.	Product	Features
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, block &amp; permeabilize in just 10 minutes</li> <li>• Complete system for blocking, permeabilizing, and antibody dilution</li> <li>• Non-mammalian blocking agents, for broad secondary antibody compatibility</li> <li>• For immunofluorescence on cells or tissue sections</li> </ul>
23013	TrueBlack® WB Blocking Buffer Kit	<ul style="list-style-type: none"> <li>• Blocks as well or better than Odyssey® Blocking Buffer, at a lower price</li> <li>• Reduces non-specific protein bands and background over entire membrane</li> <li>• Suppresses background from charged dyes better than BSA, gelatin, or casein</li> <li>• Compatible with PVDF and nitrocellulose membranes</li> <li>• Contains no mammalian proteins, for broad antibody compatibility</li> <li>• For visible and near-IR fluorescent westerns</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>
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 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 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>
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>
40083	NucSpot® 470 Nuclear Stain,	<ul style="list-style-type: none"> <li>• Green fluorescent nuclear-specific counterstain for fixed cells or tissues</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	
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative	<ul style="list-style-type: none"> <li>• Ready-to-use fixation buffer</li> <li>• Methanol-free, prepared from EM grade paraformaldehyde</li> <li>• No glass ampoules to break, store in the original bottle</li> </ul>
22020	10X Phosphate Buffered Saline (PBS) 4L Cubitainer®	<ul style="list-style-type: none"> <li>• Convenient buffers and blocking agents for immunofluorescence or western</li> </ul>
22010	10X Fish Gelatin Blocking Agent	
22014	30% Bovine Serum Albumin Solution	

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.

CF dye technology is covered by pending US and international patents. Alexa Fluor, DyLight, and Image-iT are registered trademarks of Thermo Fisher Scientific; Cubitainer is a registered trademark of Hedwin Corporation. CY DYE is a registered trademark of GE Healthcare; Draq7 is a registered trademark of Biostatus Ltd; Odyssey is a registered trademark of LI-COR Biosciences.