

Product Information

CF® Dye Conjugated Monoclonal Mouse Anti-Biotin Antibodies

CF® Dye	Ex/Em (nm)	Cat. No.	Size
CF®405S	411/431	20203-1	50 uL
		20203	250 uL
CF®488A	490/516	20204-1	50 uL
		20204	250 uL
CF®568	562/584	20502-1	50 uL
		20502	250 uL
CF®594	593/615	20205-1	50 uL
		20205	250 uL
CF®633	629/650	20206-1	50 uL
		20206	250 uL
CF®640R	642/663	20207-1	50 uL
		20207	250 uL
CF®750	755/779	20501-50uL	50 uL
		20501-150uL	250 uL
CF®770	770/797	20367-50uL	50 uL
		20367-250uL	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.

Clone: 3D6.6

Isotype: Mouse IgG1, kappa

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.

Product Description

Monoclonal mouse anti-biotin antibodies bind to antibodies or other proteins conjugated to biotin. 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/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, 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. The only protocol supplied herein is for blocking endogenous biotin.

Protocol for 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
92244...92444	Mix-n-Stain™ Biotin Antibody Labeling Kit	<ul style="list-style-type: none"> Label your antibody with biotin in just 15 minutes without a purification step. Labeling tolerates many common buffer components including BSA and ascites.
29030...29086	Streptavidin Conjugates	<ul style="list-style-type: none"> High quality biotin-binding protein conjugated to Biotium's signature bright and photostable fluorescent CF® dyes, and a selection of other labels.
23012	TrueBlack® IF Background Suppressor System (Permeabilizing)	<ul style="list-style-type: none"> Suppresses background from non-specific antibody binding and charged fluorescent dyes More efficient than Image-iT® FX, block & permeabilize in just 10 minutes Complete system for blocking, permeabilizing, and antibody dilution Non-mammalian blocking agents, for broad secondary antibody compatibility For immunofluorescence on cells or tissue sections
23013	TrueBlack® WB Blocking Buffer Kit	<ul style="list-style-type: none"> Blocks as well or better than Odyssey® Blocking Buffer, at a lower price Reduces non-specific protein bands and background over entire membrane Suppresses background from charged dyes better than BSA, gelatin, or casein Compatible with PVDF and nitrocellulose membranes Contains no mammalian proteins, for broad antibody compatibility For visible and near-IR fluorescent westerns
23007	TrueBlack® Lipofuscin Autofluorescence Quencher	<ul style="list-style-type: none"> Eliminates lipofuscin autofluorescence with less background than Sudan Black B Reduces background from other sources Can be used before or after IF staining
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	
23008	Drop-n-Stain EverBrite™ Mounting Medium	<ul style="list-style-type: none"> Convenient dropper bottle for pipette-free mounting Protects fluorescent dyes from rapid photobleaching
23009	Drop-n-Stain EverBrite™ Mounting Medium with DAPI	<ul style="list-style-type: none"> Compatible with Cy®3, Cy®5, and Alexa Fluor® 647, unlike VECTASHIELD® Available with or without DAPI for nuclear counterstaining Works as well as original EverBrite™, but less viscous for rapid DAPI staining
23005	CoverGrip™ Coverslip Sealant	<ul style="list-style-type: none"> Replaces nail polish for coverslip sealing Won't leach into aqueous mounting media
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 Draq7®
40083...41038	NucSpot® Nuclear Stains	<ul style="list-style-type: none"> Fluorescent nuclear-specific counterstain for fixed cells or tissues, from green to far-red
23023	Mini Super ^{HT} Pap Pen 2.0	<ul style="list-style-type: none"> Create hydrophobic barriers around tissue sections, ~400 uses Heat-stable to 120°C Insoluble in aqueous buffers, detergents, alcohol and acetone; can be removed with xylene
23024	Super ^{HT} Pap Pen 2.0	
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative	<ul style="list-style-type: none"> Ready-to-use fixation buffer Methanol-free, prepared from EM grade paraformaldehyde No glass ampoules to break, store in the original bottle
22010	10X Fish Gelatin Blocking Agent	<ul style="list-style-type: none"> Convenient blocking agents for immunofluorescence or western
22014	30% Bovine Serum Albumin Solution	

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.

CF Dye technology is covered by US and international patents. Alexa Fluor, DyLight, and Image-iT are registered trademarks of Thermo Fisher Scientific; CY DYE is a registered trademark of Cytiva; Draq7 is a registered trademark of Biostatus Ltd; Odyssey is a registered trademark of LI-COR Biosciences; VECTASHIELD is a registered trademark of Vector Laboratories, Inc.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.