

Revised: April 4, 2024

# **Product Information**

# **CF® Dye Biotin**

## **Product List**

Cat. No.	Conjugate	Unit Size	Ex/Em (nm)	MW
80034	CF®488A	1 mg	490/516	~1182
80038	CF®532	1 mg	531/552	~954
80029	CF®568	1 mg	562/584	~981
80035	CF®594	1 mg	593/615	~997
80031	CF®633	1 mg	629/650	~1089
80032	CF®640R	1 mg	642/663	~1100
80036	CF®647	1 mg	652/668	~1254
80037	CF®680R	1 mg	680/701	~1180

### Storage and Handling

Store desiccated at 4°C. Protect from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

Stock solutions may be prepared in DMSO or DMF.

#### **Product Description**

Potential applications of Biotin-CF® Dyes include detection of biotin binding sites on avidin/streptavidin or in biological samples. In addition, Biotin-CF® Dyes can be used as a polar tracer by microinjection to study the morphology of cells. CF® Dyes are very photostable, which makes them suitable for studies where prolonged exposure to light may be necessary.

Browse our website to view our full selection of biotin and CF® Dye conjugates.

#### **Related Products**

Related Produ			
Cat. No.	Product		
9220892228	CF® Dye & Biotin SE Protein Labeling Kits		
00017	Biotin-XX-α-Bungarotoxin		
00020	Biotin-cAMP		
60022	Biotin-DHPE		
80019	Fluorescein Biotin		
80022	Biotin-Rhodamine 110		
90062	Biotin-4-Fluorescein		
92167	Biotin Azide		
92176	Biotin-XX Tyramide		
92168	Biotin Alkyne		
29030 29086	CF® Dye Streptavidin Conjugates		
29049	HRP Streptavidin		
29071	Alkaline Phosphatase Streptavidin		
29048	APC Streptavidin		
29044	R-PE Streptavidin		
40061	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO		
4008341040	NucSpot® Nuclear Stains		
23001	EverBrite™ Mounting Medium		
23002	EverBrite™ Mounting Medium with DAPI		
23003	EverBrite™ Hardset Mounting Medium		
23004	EverBrite™ Hardset Mounting Medium with DAPI		
23005	CoverGrip™ Coverslip Sealant		
22015	Fixation Buffer, 100 mL		
22016	Permeabilization Buffer, 100 mL		
22017	Permeabilization and Blocking Buffer, 100 mL		
22010	10% Fish Gelatin Blocking Buffer, 100 mL		
22011	Fish Gelatin Powder, 2 X 50 g		
22014	30% Bovine Serum Albumin Solution, 100 mL		
22002	Tween®-20, 50 mL		

Please visit our website at www.biotium.com for information on our life science research products, including fluorescent CF® Dye antibody conjugates and reactive dyes, biotinylation reagents with a full selection of reactive functional groups, as well as easy-to-use protein and antibody biotinylation kits, and kits for cell biology research.

CF dye technology is covered by pending US and international patents. TWEEN is a registered trademark of CRODA AMERICAS LLC.

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