

www.biotium.com

Revised: September 29, 2020

# **Product Information**

## CF® Dye Azide or Biotin Azide

### **Product List**

Catalog no.	Azide Conjugate	Unit size	Ex/Em (nm)	MW (free acid form)
92092	CF®405M	0.5 mg	408/452	~583
92080	CF®488A	0.5 mg	490/515	~996
92180	CF®532	0.5 mg	527/558	~768
92181	CF®543	0.5 mg	541/560	~969
92081	CF®555	0.5 mg	555/565	~983
92082	CF®568	0.5 mg	562/583	~796
92083	CF®594	0.5 mg	593/614	~811
92085	CF®640R	0.5 mg	642/662	~914
92084	CF®647	0.5 mg	650/665	~1067
92094	CF®660C	0.5 mg	667/685	~3194
92182	CF®660R	0.5 mg	663/682	~970
92119	CF®680	0.5 mg	681/698	~3,234
96000	CF®680R	0.5 mg	680/701	~994
92167	Biotin	1 mg		~445

#### Storage and Handling

Store CF® Dye Azide or Biotin Azide at -20°C, protected from light. Product is stable for at least 12 months from date of receipt if stored as recommended. Stock solution may be prepared in DMSO or dH<sub>2</sub>O and can be stored at  $\leq$  -20°C for at least 12 months.

#### **Product Description**

For this set of products, we provide our CF® dyes or biotin in the azide form. CF® dyes or biotin with an azide group can be used to label or detect alkyne, BCN, DBCO, or DIFO by fluorescence spectroscopy, microscopy, or flow cytometry. CF® Dye Azide or Biotin Azide can react with alkyne to form 1,2,3-triazole by copper(I) catalyzed 1,3-dipolar Huisgen cycloaddition. It can also label BCN, DBCO, and DIFO in a copper-free bioorthogonal 1,3-dipolar Huisgen cycloaddition or undergo Staudinger ligation with phosphine containing compounds.

CF® dyes are Biotium's line of next generation fluorescent dyes with combined advantages in brightness, photostability, and water solubility.

### **Related Products**

Catalog number	Product		
90082	DMSO, Anhydrous		
92187	CF®488A Picolyl Azide		
92188	CF®568 Picolyl Azide		
92189	CF®594 Picolyl Azide		
92190	CF®640R Picolyl Azide		
92191	CF®647 Picolyl Azide		
96001	CF®660C Picolyl Azide		
96002	CF®660R Picolyl Azide		
96003	CF®680 Picolyl Azide		
96007	CF®680R Picolyl Azide		
92086	CF®488A Alkyne		
92087	CF®555 Alkyne		
92088	CF®568 Alkyne		
92089	CF®594 Alkyne		
92090	CF®647 Alkyne		
92091	CF®640R Alkyne		
92093	CF®405M Alkyne		
92095	CF®660C Alkyne		
92113	CF®405S BCN		
92114	CF®405M BCN		
96070	CF®440 BCN		
92075	CF®488A BCN		
96026	CF®500 BCN		
92076	CF®568 BCN		
92077	CF®594 BCN		
92078	CF®640R BCN		
96059	CF®647 BCN		
96027	CF®650 BCN		
96058	CF®680 BCN		
92079	CF®680R BCN		
92168	Biotin Alkyne		
92169	Biotin BCN		
92186	Biotin Picolyl Azide		

Please visit our website at www.biotium.com for information on our life science research products, including a full selection of CF® reactive dyes and CF® dye conjugates including succinimidyl esters, maleimides, amines, and hydrazides. Biotium also offers convenient Mix-n-Stain<sup>™</sup> CF® Dye Antibody Labeling Kits for convenient and stable antibody labeling in only 30 minutes.

CF dye technology is covered by pending U.S. and international patents. Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.