

Revised: November 6, 2019



Product Information

BCN (Bicyclo[6.1.0]nonyne) Conjugates

Unit Size: 0.5 mg for CF® dye BCN

1.0 mg for Biotin BCN and MTS-BCN

Technical Summary

Cat. No.	BCN conjugate	Abs/Em (nm)	Membrane Permeability	Extinction coefficient	MW
92113	CF®405S	404/431	Impermeant	33,000	~788
92114	CF®405M	408/452	Impermeant	41,000	~720
96070	CF®440	440/515	Permeant	40,000	~698
92075	CF®488A	490/515	Impermeant	70,000	~1132
96026	CF®500	500/510	Permeant	70,000	~510
92076	CF®568	562/583	Impermeant	100,000	~932
92077	CF®594	593/614	Impermeant	115,000	~947
92078	CF®640R	642/662	Impermeant	105,000	~1050
96059	CF®647	650/665	Impermeant	240,000	~1203
96027	CF®650	650/670	Permeant	100,000	~717
96058	CF®680	681/698	Impermeant	210,000	~3371
92079	CF®680R	680/701	Impermeant	140,000	~1130
92169	Biotin				~463
96023	MTS				~331

Storage and Handling

Store BCN conjugates at - 20° C, protected from light. Product is guaranteed for at least 12 months from date of receipt when stored as recommended. Stock solution may be prepared in DMSO or dH₂O and can be stored at ≤ -20°C for at least 12 months.

Product Description

BCN (Bicyclo[6.1.0]nonyne) conjugates react with azide to form 1,2,3-triazole by copper-free 1,3-dipolar Huisgen cycloaddition. This copper-free bioorthogonal reaction can be used with live cells or cell extracts when there are concerns about native protein function loss with copper.

CF® dyes are Biotium's next-generation fluorescent dyes with exceptional brightness and photostability. CF®440 BCN, CF®500 BCN, and CF®650 BCN are membrane-permeant for intracellular copper-free reaction with azide.

Related Products

Catalog number	Product
92092	CF®405M Azide
92080	CF®488A Azide
92081	CF®555 Azide
92082	CF®568 Azide
92083	CF®594 Azide
92084	CF®647 Azide
92085	CF®640R Azide
92094	CF®660C Azide
92086	CF®488A Alkyne
92168	Biotin 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
92167	Biotin Azide
22004	Ultrafiltration Vial, 10K MWCO (5 per pack)
22018	Ultrafiltration Vial, 3K MWCO (5 per pack)
90082	DMSO, Anhydrous
22013	Bovine Serum Albumin, Fraction V
22014	Bovine Serum Albumin, 30% Solution
22020	10X Phosphate Buffered Saline
41024-4L	Water, Ultrapure Molecular Biology Grade

Please visit www.biotium.com to view our full selection of innovative products for life science research.

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.