

Product Information

CF™ Dye Alkyne or Biotin Alkyne

Unit Size: 0.5 mg for CF™ dye alkyne
1.0 mg for Biotin alkyne

Technical Summary

Cat. No.	CF™ Dye	Abs _{max} (nm)	Em _{max} (nm)	Extinction coefficient	MW
92093	CF™405M	408	452	41,000	~539
92086	CF™488A	490	515	70,000	~951
92087	CF™555	555	565	150,000	~938
92088	CF™568	562	583	100,000	~751
92089	CF™594	593	614	115,000	~766
92091	CF™640R	642	662	105,000	~869
92090	CF™647	650	665	240,000	~908
92095	CF™660C	667	685	200,000	~3149
96004	CF™660R	663	682	100,000	~925
96005	CF™680	681	698	210,000	~3189
96006	CF™680R	680	701	140,000	~949
92168	Biotin	---	---	---	~458

Storage and Handling

Store CF™ dye alkyne or biotin alkyne 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 dH₂O and can be stored at ≤ -20°C for at least 12 months.

Product Description

For this set of products, we provide our CF™ dyes or biotin in the alkyne form. CF™ dyes are Biotium's line of next generation fluorescent dyes with combined advantages in brightness, photostability, and water solubility. CF™ dye alkyne or biotin alkyne can react with azide to form 1,2,3-triazole by copper(I) catalyzed 1,3-dipolar Huisgen cycloaddition. It can also function as monomer that allows copolymerization with other monomers to form a fluorescent polymer.

Other Related Products

You may also be interested in the following related products from Biotium:

- A full selection CF™ reactive dyes and CF™ dye conjugates
- CF™ dye-labeled azides

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.