

Revised: October 26, 2017



Product Information

CF[™] Dye Picolyl Azide or Biotin Picolyl Azide

Unit Size: 0.5 mg

Technical Summary

Cat. No.	CF™ Dye	Abs _{max} (nm)	Em _{max} (nm)	Extinction coefficient	MW
92187	CF™488A	490	515	70,000	~1116
92188	CF™568	562	583	100,000	~916
92189	CF™594	593	614	115,000	~931
92190	CF™640R	642	662	105,000	~1034
92191	CF™647	650	665	240,000	~1187
96001	CF™660C	667	685	200,000	~3326
96002	CF™660R	663	682	100,000	~1090
96003	CF™680	681	698	210,000	~3355
96007	CF™680R	680	701	140,000	~1114
92186	Biotin				~694

Storage and Handling

Store $\overline{\mathsf{CF}}^{\mathsf{TM}}$ dye picolyl azide or biotin picolyl 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 dH₂O 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 picolyl azide form. CF[™] dyes are Biotium's line of next generation fluorescent dyes with combined advantages in brightness, photostability, and water solubility. CF[™] dye picolyl azide or biotin picolyl azide reacts with alkyne to form 1,2,3-triazole by 1,3-dipolar Huisgen cycloaddition through the use of much lower copper(I) concentration without sacrificing reaction efficiency.

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 alkynes

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.