

CF® Dye Alkyne

CF® Dye alkynes react with azide groups via copper-catalyzed bioorthogonal cycloaddition. Can be used to fluorescently detect or label azide groups on target molecules. CF® Dye alkynes can also be used as building blocks to form fluorescent polymers.



Product attributes

Chemical reactivity (reacts with)	Azides/Picolyl azides
Functional group	Alkyne
Storage Conditions	Store at -10 to -35 °C, Protect from light

Product Description

CF® Dye alkynes can react with azide groups to form 1,2,3-triazole via 1,3-dipolar Huisgen copper-catalyzed cycloaddition. They can also be used as monomeric building blocks to copolymerize with other monomers to form fluorescent polymers.

- Fluorescent detection or labeling azide groups on target molecules.
- Can be used as fluorescent monomers to form dye-labeled polymers.
- Bright, photostable and water-soluble CF® Dyes are excellent options for fluorescent labeling.

Superior CF® Dyes

CF® Dyes are Biotium's line of next-generation fluorescent dyes that have improved brightness, photostability and water solubility compared to Alexa Fluor®, DyLight®, and other fluorescent dyes.

Learn more about [CF® Dyes](#). For more information download the [CF® Dye Brochure](#).

CF® Dye Alkyne

CF® Dye Alkyne	Ex/Em	Size	Catalog No.	Dye Features
CF@405M	408/452 nm	0.5 mg	92093	CF@405M Features
CF@488A	490/515 nm	0.5 mg	92086	CF@488A Features
CF@555	555/565 nm	0.5 mg	92087	CF@555 Features
CF@568	562/583 nm	0.5 mg	92088	CF@568 Features
CF@594	593/614 nm	0.5 mg	92089	CF@594 Features
CF@640R	642/662 nm	0.5 mg	92091	CF@640R Features
CF@647	650/665 nm	0.5 mg	92090	CF@647 Features
CF@660R	663/682 nm	0.5 mg	96004	CF@660R Features
CF@660C	667/685 nm	0.5 mg	92095	CF@660C Features
CF@680	681/698 nm	0.5 mg	96005	CF@680 Features
CF@680R	680/701 nm	0.5 mg	96006	CF@680R Features

References

1. Dev Cell. (2019) 50(1):57-72.e6. [DOI: 10.1016/j.devcel.2019.04.035](#)
2. ACS Nano (2016) 10:8861–8870. [DOI: 10.1021/acs.nano.6b04748](#)

Download a list of [CF® dye references](#).

This datasheet was generated on January 17, 2026 at 02:49:49 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-alkyne/>