# **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**

Call us: 800-304-5357

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

Email: btinfo@biotium.com

## **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	<u>92093</u>	CF®405M Features
<u>CF®488A</u>	490/515 nm	0.5 mg	<u>92086</u>	CF®488A Features
CF®555	555/565 nm	0.5 mg	<u>92087</u>	CF®555 Features
<u>CF®568</u>	562/583 nm	0.5 mg	<u>92088</u>	CF®568 Features
CF®594	593/614 nm	0.5 mg	<u>92089</u>	CF®594 Features
<u>CF®640R</u>	642/662 nm	0.5 mg	<u>92091</u>	CF®640R Features
CF®647	650/665 nm	0.5 mg	<u>92090</u>	CF®647 Features
CF®660R	663/682 nm	0.5 mg	<u>96004</u>	CF®660R Features
CF®660C	667/685 nm	0.5 mg	<u>92095</u>	CF®660C Features
<u>CF®680</u>	681/698 nm	0.5 mg	<u>96005</u>	CF®680 Features
CF®680R	680/701 nm	0.5 mg	<u>96006</u>	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/acsnano.6b04748

Download a list of CF® dye references.

This datasheet was generated on November 17, 2025 at 10:45:34 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cf-dye-alkyne/">https://biotium.com/product/cf-dye-alkyne/</a>