

CF® Dye Azide

CF® Dyes with an azide group. Can be used to label or detect alkyne, BCN or cyclooctyne-containing molecules by fluorescence spectroscopy, microscopy or flow cytometry following copper-catalyzed azide-alkyne cycloaddition.



Product attributes

Chemical reactivity (reacts with)	Alkynes, BCN
Functional group	Azide
Storage Conditions	Store at -10 to -35 °C, Protect from light

Product Description

CF® Dye azide can react with terminal alkynes to form 1,2,3-triazole via copper-catalyzed cycloaddition. They can also undergo Staudinger ligation with a phosphine containing compounds. The azide derivatives also react with cyclooctynes and BCN groups via copper-free bioorthogonal reaction.

- Fluorescent detection or labeling of alkyne, BCN or cyclooctyne groups on target molecules.
- 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](#).

We also offer [CF® Dye Picolyl Azide](#) derivatives for azide-alkyne cycloaddition reactions at lower copper concentrations, suitable for labeling live cells or other copper-sensitive systems.

CF® Dye Azide

CF® Dye Azide	Ex/Em	Size	Catalog No.	Dye Features
CF@405M	408/452 nm	0.5 mg	92092	CF@405M Features
CF@488A	490/515 nm	0.5 mg	92080	CF@488A Features
CF@532	527/558 nm	0.5 mg	92180	CF@532 Features
CF@543	541/560 nm	0.5 mg	92181	CF@543 Features
CF@555	555/565 nm	0.5 mg	92081	CF@555 Features
CF@568	562/583 nm	0.5 mg	92082	CF@568 Features
CF@594	593/614 nm	0.5 mg	92083	CF@594 Features
CF@640R	642/662 nm	0.5 mg	92085	CF@640R Features
CF@647	650/665 nm	0.5 mg	92084	CF@647 Features
CF@660R	663/682 nm	0.5 mg	92182	CF@660R Features
CF@660C	667/685 nm	0.5 mg	92094	CF@660C Features
CF@680	681/698 nm	0.5 mg	92119	CF@680 Features
CF@680R	680/701 nm	0.5 mg	96000	CF@680R Features

References

Nat Protocols (2017) 12(5), 1011-1028. [doi: 10.1038/nprot.2017.020](https://doi.org/10.1038/nprot.2017.020) (CF@405M azide, CF@568 azide, SIM)

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

This datasheet was generated on June 3, 2026 at 05:20:19 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf-dye-azide/>