

CF® Dye Picolyl Azide

CF® dye picolyl azide derivatives fluorescently label alkyne groups on biomolecules via the standard azide-alkyne cycloaddition, but at low copper concentrations.



Product attributes

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

Product Description

CF® dye picolyl azide derivatives react with alkyne groups to form 1,2,3-triazole by 1,3-dipolar Huisgen cycloaddition. They offer benefits of the typical azide-alkyne cycloaddition but at a much lower copper (I) concentration without sacrificing reaction efficiency making it compatible with live cell labeling and copper-sensitive systems.

- Fluorescent detection or labeling of alkyne containing biomolecules in vitro or in vivo.
- Bioorthogonal azide-alkyne cycloaddition reaction at low copper (I) concentrations.
- 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 azide](#) derivatives for copper-catalyzed azide-alkyne cycloaddition reactions or copper-free bioorthogonal reactions with cyclooctynes and BCN group containing molecules.

CF Dye Picolyl Azide: Product & SKU Summary

Product	Conjugation	Ex/Em	Size	SKU	Purchase
CF@488A Picolyl Azide	CF@488A	490/515 nm	0.5 mg	92187	Purchase 92187
CF@568 Picolyl Azide	CF@568	562/583 nm	0.5 mg	92188	Purchase 92188
CF@594 Picolyl Azide	CF@594	593/614 nm	0.5 mg	92189	Purchase 92189
CF@640R Picolyl Azide	CF@640R	642/662 nm	0.5 mg	92190	Purchase 92190
CF@647 Picolyl Azide	CF@647	650/665 nm	0.5 mg	92191	Purchase 92191
CF@660C Picolyl Azide	CF@660C	667/685 nm	0.5 mg	96001	Purchase 96001
CF@660R Picolyl Azide	CF@660R	663/682 nm	0.5 mg	96002	Purchase 96002
CF@680 Picolyl Azide	CF@680	681/698 nm	0.5 mg	96003	Purchase 96003
CF@680R Picolyl Azide	CF@680R	680/701 nm	0.5 mg	96007	Purchase 96007

References

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

This datasheet was generated on June 9, 2026 at 05:50:49 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-picolyl-azide/>