

## 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	<a href="#">Purchase 92187</a>
CF@568 Picolyl Azide	CF@568	562/583 nm	0.5 mg	92188	<a href="#">Purchase 92188</a>
CF@594 Picolyl Azide	CF@594	593/614 nm	0.5 mg	92189	<a href="#">Purchase 92189</a>
CF@640R Picolyl Azide	CF@640R	642/662 nm	0.5 mg	92190	<a href="#">Purchase 92190</a>
CF@647 Picolyl Azide	CF@647	650/665 nm	0.5 mg	92191	<a href="#">Purchase 92191</a>
CF@660C Picolyl Azide	CF@660C	667/685 nm	0.5 mg	96001	<a href="#">Purchase 96001</a>
CF@660R Picolyl Azide	CF@660R	663/682 nm	0.5 mg	96002	<a href="#">Purchase 96002</a>
CF@680 Picolyl Azide	CF@680	681/698 nm	0.5 mg	96003	<a href="#">Purchase 96003</a>
CF@680R Picolyl Azide	CF@680R	680/701 nm	0.5 mg	96007	<a href="#">Purchase 96007</a>

## References

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

This datasheet was generated on December 7, 2025 at 10:42:58 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cf-dye-picolyl-azide/>