# **CF® Dye BCN**

CF® Dyes with a BCN group react with azides via a copper-free bioorthogonal reaction to label target molecules for surface staining of live cells or cell extracts. Membrane-permeant CF® Dye BCN conjugates for intracellular labeling reactions are also available.



#### **Product attributes**

Call us: 800-304-5357

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

Email: btinfo@biotium.com

### **Product Description**

CF® Dye BCN (bicyclo[6.1.0]nonyne) derivatives are alternatives to fluorescently labeled DIBO and DBCO. They react with azides to form 1,2,3-triazole by copper-free bioorthogonal 1,3-dipolar Huisgen cycloaddition. BCN dyes are useful for fluorescent labeling of live cell surface or cell extracts, particularly when there are concerns about native protein function loss with copper. Membrane-permeant CF® Dye BCN options are also available for intracellular labeling reactions.

- Fluorescent labeling of azide groups on target molecules via copper-free bioorthogonal chemistry.
- Alternative to fluorescently labeled DIBO and DBCO.
- Suitable for surface or intracellular labeling of live cells or cell extracts.
- Bright, photostable and water-soluble CF® Dyes are excellent options for fluorescent labeling.
- Available in several CF® Dye options including membrane-permeant forms for intracellular labeling.

Biotium also offers reactive <u>CF® Dye DBCQ</u> derivatives which react more rapidly with azide groups than other cyclooctynes through strain-promoted azide-alkyne cycloaddition (SPAAC).

#### Superior CF® Dyes

Biotium's next-generation CF® Dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about CF® Dyes. For more information download the CF® Dye Brochure.

## **CF® Dye BCN**

CF® Dye BCN	Ex/Em	Membrane Permeability	Size	Catalog No.	Dye Features
CF®405S	404/431 nm	Impermeant	0.5 mg	<u>92113</u>	CF®405S Features
CF®405M	408/452 nm	Impermeant	0.5 mg	<u>92114</u>	CF®405M Features
CF®440	440/515 nm	Permeant	0.5 mg	<u>96070</u>	CF®440 Features
CF®488A	490/515 nm	Impermeant	0.5 mg	<u>92075</u>	CF®488A Features
CF®500	502/511 nm	Permeant	0.5 mg	<u>96026</u>	
CF®568	562/583 nm	Impermeant	0.5 mg	<u>92076</u>	CF®568 Features
CF®594	593/614 nm	Impermeant	0.5 mg	<u>92077</u>	CF®594 Features
CF®640R	642/662 nm	Impermeant	0.5 mg	<u>92078</u>	CF®640R Features
CF®647	650/665 nm	Impermeant	0.5 mg	<u>96059</u>	CF®647 Features
CF®650	653/673 nm	Permeant	0.5 mg	<u>96027</u>	
CF®680	681/698 nm	Impermeant	0.5 mg	<u>96058</u>	CF®680 Features
CF®680R	680/701 nm	Impermeant	0.5 mg	<u>92079</u>	CF®680R Features

### References

- 1. Canadian Journal of Chemistry (2019),97(1):1-6. DOI: 10.1139/cjc-2018-0253.
- 2. eLife (2019),8: e50776. DOI: 10.7554/eLife.50776.001
- 3. Chem. Mater. (2019), 31:8035–8043. DOI: 10.1021/acs.chemmater.9b02485
- 4. Advanced functional materials, (2016),26(21):3612-3620. DOI: 10.1002/adfm.201505329

Download a list of CF® dye references.

This datasheet was generated on November 17, 2025 at 09:44:43 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cf-dye-bcn/">https://biotium.com/product/cf-dye-bcn/</a>