

CF® Dye DBCO

CF® Dye with dibenzocyclooctyne (i.e., DBCO, ADIBO) reacts with azide via copper-free click chemistry by strain-promoted azide-alkyne cycloaddition (SPAAC).



Product attributes

Storage Conditions	Store at -35°C to -5°C
Chemical reactivity (reacts with)	Azides/Picoly azides

Product Description

CF® Dye DBCO (Dibenzocyclooctyne) derivatives are highly efficient, fast-reacting alternatives to fluorescently labeled BCN and DIBO. These DBCO derivatives are specifically designed for copper-free click chemistry, reacting more rapidly with azide groups than other cyclooctynes through strain-promoted azide-alkyne cycloaddition (SPAAC). Unlike traditional click reactions, DBCO-azide reactions do not require toxic copper catalysts (CuAAC), making them more suitable for live-cell or *in vivo* experiments. CF® Dye DBCO is available in 14 derivatives.

- Fluorescent labeling of azide groups on target molecules
- Time-saving alternative to fluorescently labeled DIBO and BCN
- Suitable for labeling cell extracts or live cells without copper toxicity issues
- Wide selection of bright, photostable, and water-soluble CF® Dyes

DBCO dyes are useful for bioorthogonal reactions in combination with trans-cyclooctenes (TCO) and tetrazine chemistries, because DBCO does not interact with tetrazines unlike other cyclooctynes. DBCO conjugation reactions can be performed in aqueous buffers or in organic solvents, depending on the properties of the substrate molecules.

Superior CF® Dyes

Biotium offers a wide selection of CF® Dyes with the DBCO reactive group. Our outstanding series of CF® Dyes offer superior brightness, photostability, and signal-to-noise when compared to other fluorescent dyes. Learn more about [CF® Dyes](#).

CF® Dye DBCO

Product	Ex/Em	Size	Catalog No.	Dye Features
CF@405M	416/452 nm	1 mg	96110	CF@405M Features
CF@405S	411/431 nm	1 mg	96111	CF@405S Features
CF@488A	490/516 nm	1 mg	96112	CF@488A Features
CF@555	554/568 nm	1 mg	96113	CF@555 Features
CF@568	562/584 nm	1 mg	96114	CF@568 Features
CF@594	593/615 nm	1 mg	96115	CF@594 Features
CF@640R	642/663 nm	1 mg	96116	CF@640R Features
CF@647	652/668 nm	1 mg	96117	CF@647 Features
CF@660R	662/682 nm	1 mg	96118	CF@660R Features
CF@680	681/698 nm	1 mg	96119	CF@680 Features
CF@680R	680/701 nm	1 mg	96120	CF@680 Features
CF@700	696/721 nm	1 mg	96121	CF@700 Features
CF@740	742/767 nm	1 mg	96122	CF@740 Features
CF@750	755/779 nm	1 mg	96123	CF@750 Features

This datasheet was generated on December 21, 2024 at 05:37:04 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/cf-dye-dbc/>