

CF@DI Aminoxy

CF@DI Aminoxy derivatives are useful for fluorescently labeling aldehyde/ketone groups on biomolecules such as polysaccharides, glycoproteins or antibodies to generate size- and charge-matched conjugates.



Product attributes

Chemical reactivity (reacts with)	Aldehydes/ketones
Colors	Green, Orange-red, Far-red
Functional group	Aminoxy (hydroxylamine)
Storage Conditions	Store at -10 to -35 °C, Protect from light

Product Description

CF@DI aminoxy dyes react with aldehyde/ketone groups on proteins or other biomolecules to form a stable oxime bond. The CF@DI dyes (CF@488DI, CF@555DI and CF@647DI) are designed to match each other for molecular weight and charge. The molecular weights of the same protein conjugated to different CF@DI dye colors are close to each other and the pI of the labeled proteins remains the same.

- React with carbonyl moieties of aldehyde or ketone groups on proteins and other biomolecules.
- Size- and charge-matched fluorescent aminoxy dyes.
- Available in 3 colors for multiplex analysis.

We also offer our original [CF@ Dye Aminoxy](#) with a wide selection of colors. CF@ Dyes are next-generation fluorescent dyes with advantages in brightness, photostability, and water solubility compared to other commercially available fluorescent dyes. For more information, download the [CF@ Dye Brochure](#).

Product	Ex/Em	MW (g/mol)	Size	Catalog No.
CF@488DI Aminoxy	483/508 nm	~633	1 mg	92177
CF@555DI Aminoxy	547/572 nm	~639	1 mg	92178
CF@647DI Aminoxy	639/668 nm	~621	1 mg	92179

References

1. Organic Letters (2017) 19, 12: 3179-3182. [DOI:10.1021/acs.orglett.7b01198](https://doi.org/10.1021/acs.orglett.7b01198)

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

This datasheet was generated on November 21, 2024 at 03:14:22 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/cfdi-dye-aminoxy/>