

## 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

<b>Chemical reactivity (reacts with)</b>	Aldehydes/ketones
<b>Colors</b>	Green, Orange-red, Far-red
<b>Functional group</b>	Aminoxy (hydroxylamine)
<b>Storage Conditions</b>	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	<a href="#">92177</a>
CF@555DI Aminoxy	547/572 nm	~639	1 mg	<a href="#">92178</a>
CF@647DI Aminoxy	639/668 nm	~621	1 mg	<a href="#">92179</a>

## 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 July 31, 2025 at 03:28:00 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cfdi-dye-aminoxy/>