CF® Dye Hydrazide

Fluorescent CF® dye hydrazides are bright, extremely water-soluble, nontoxic molecules, useful as aldehyde-fixable cell tracers. They can also be used to label aldehyde or ketone groups of polysaccharides, carbohydrates or glycoproteins.



Product Description

Bright, extremely water-soluble and non-toxic CF® dye hydrazides are excellent fluorescent tracers for neurons or gap junction studies. The cell impermeant tracer dyes can be introduced into cells by microinjection. CF® dye hydrazides can also be used to fluorescently label aldehyde or ketone groups of carbohydrates or glycoproteins molecules after peroxidation with periodate.

- Fluorescently labeled low molecular weight, polar, non-toxic molecules; useful as aldehyde-fixable cell tracers.
- Membrane impermeant, can be easily loaded into cells by microinjection.
- Can be used to fluorescently label aldehyde or ketone groups of polysaccharides, carbohydrates, or glycoproteins.
- Bright, photostable and water-soluble CF® dyes are excellent options for fluorescent imaging.

For carbohydrates labeling applications, we recommend using <u>CF® dye aminooxy</u> forms, which are more reactive toward these groups than hydrazide forms. For fluorescence two-dimensional differential gel electrophoresis (2D-DIGE), we recommend <u>CFDI hydrazides</u>.

Superior CF® Dyes

CF® Dyes are Biotium's line of next-generation fluorescent dyes that have improved brightness, photostability and water solubility compared to other commercially available fluorescent dyes. Learn more about CF® Dyes. For more information download the CF® Dye Brochure.

CF Dye Hydrazides

Product	Ex/Em	Size	SKU	Purchase
CF®350 Hydrazide	347/448 nm	1 mg	92151	Purchase 92151
CF®405S Hydrazide	404/431 nm	1 mg	92183	Purchase 92183
CF®430 Hydrazide	426/498 nm	1 mg	96063	Purchase 96063
CF®440 Hydrazide	440/515 nm	1 mg	96064	Purchase 96064
CF®488A Hydrazide	490/515 nm	1 mg	92152	Purchase 92152
CF®555 Hydrazide	555/565 nm	1 mg	92153	Purchase 92153
CF®568 Hydrazide	562/583 nm	1 mg	92154	Purchase 92154
CF®594 Hydrazide	593/614 nm	1 mg	92158	Purchase 92158
CF®633 Hydrazide	630/650 nm	1 mg	92156	Purchase 92156
CF®640R Hydrazide	642/662 nm	1 mg	92157	Purchase 92157
CF®647 Hydrazide	650/665 nm	1 mg	92136	Purchase 92136
CF®660R Hydrazide	663/682 nm	1 mg	96024	Purchase 96024
CF®680R Hydrazide	680/701 nm	1 mg	96025	Purchase 96025

References

- 1. Cellular and Molecular Bioengineering (2019)12:121-130. DOI: 10.1007/s12195-018-0555-6
- 2. Journal of Controlled Release (2017) 263: 46-56. DOI: 10.1016/j.jconrel.2017.02.025
- 3. Int Forum Allergy Rhinol. (2017) 7:177- 184. DOI: 10.1002/alr.21865
- 4. Nature (2016) 538 (7623):96-98. DOI: 10.1038/nature19770

Download a list of CF® dye references.

This datasheet was generated on November 4, 2025 at 03:24:09 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cf-dye-hydrazide/

Product attributes

Call us: 800-304-5357

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Chemical reactivity (reacts with)	Aldehydes/ketones
Functional group	Hydrazide
Colors	Blue, Green, Red, Far-red, Near-infrared
Storage Conditions	Store at -10 to -35 °C, Protect from light

Email: btinfo@biotium.com