

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 attributes

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

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 [CF® dye aminoxy](#) forms, which are more reactive toward these groups than hydrazide forms. For fluorescence two-dimensional differential gel electrophoresis (2D-DIGE), we recommend [CFDI hydrazides](#).

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 June 16, 2026 at 09:30:38 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cf-dye-hydrazide/>