

CF®DI Hydrazide

CF®DI dye hydrazides react with carbonyl moieties of aldehyde or ketone groups. They are size- and charge-matched for detection of protein carbonylation by fluorescence two-dimensional differential gel electrophoresis (2D-DIGE).



Product attributes

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

Product Description

CF®DI dye hydrazides react with molecules containing aldehyde or ketone groups to form a hydrazone bond. The dyes are size- and charge-matched and do not significantly alter the molecular weight or pI of the labeled proteins. They are used for detection of protein carbonylation by fluorescence 2D-DIGE.

- React with carbonyl moieties of aldehyde or ketone groups.
- Size- and charge-matched, do not significantly alter the molecular weight or pI of the labeled protein.
- For detection of protein carbonylation by fluorescence 2D-DIGE.
- Available in several colors for detection of multiple samples on the same gel.

Also see our regular [CF® Dye Hydrazides](#), available in a wide selection of colors.

CFDI Hydrazide: Product & SKU Summary

Product	Conjugation	Ex/Em	Size	SKU	Purchase
CF®488DI Hydrazide	CF®488DI	483/508 nm	1 mg	92166	Purchase 92166
CF®555DI Hydrazide	CF®555DI	547/572 nm	1 mg	92164	Purchase 92164
CF®647DI Hydrazide	CF®647DI	639/668 nm	1 mg	92165	Purchase 92165
CF®660DI Hydrazide	CF®660DI	661/703 nm	1 mg	92184	Purchase 92184
CF®680DI Hydrazide	CF®680DI	676/711 nm	1 mg	92185	Purchase 92185

References

1. Free Radical Biology and Medicine (2020) 148:83-95. [DOI: 10.1016/j.freeradbiomed.2019.12.039](https://doi.org/10.1016/j.freeradbiomed.2019.12.039)
2. J Proteomics Computational Biol. (2016) 2(1):1-12. <http://fulir.irb.hr/id/eprint/3551>

This datasheet was generated on December 31, 2025 at 12:25:34 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cfdi-hydrazide/>