



Product Information

CF™DI Dye Hydrazide

Unit Size: 1 mg

Technical Summary

Cat. No.	Dye	Ex/Em* (nm)	MW
92166	CF™488DI	483/508	574.0
92164	CF™555DI	547/572	564.7
92165	CF™647DI	639/668	562.7
92184	CF™660DI	661/703	627.0
92185	CF™680DI	676/711	676.9

^{*} in MeOH

Storage and Handling

Store CF™DI dye hydrazide at -20°C, protected from light. Stock solutions may be prepared in DMF or DMSO.

Product Description

CF™DI dye, hydrazide are CF™ dyes with a hydrazide reactive group. Hydrazide group react with molecules containing aldehyde or ketone groups to form a hydrazone bond.

The CF™DI dyes are designed to match for charge and molecular weight. When coupled to the carbonyl group of the proteins, the molecular weight of the labeled proteins are closed to each other and the pl of the labeled proteins remains the same.

Related Products

Biotium also offers a line of next-generation fluorescent CF™ dyes for labeling proteins and nucleic acids, with advantages in brightness, photostability, and water solubility compared to other fluorescent dyes. You may also be interested in the following related products from Biotium:

- CF™ dye succinimidyl (NHS) ester, hydrazide, maleimide, amine, aminooxy, alkyne, and azide derivatives
- CF™ dye protein labeling kits and Mix-n-Stain™ antibody labeling kits.
- CF™ dye secondary antibody conjugates and other bioconjugates
- · Please visit our website at www.biotium.com for details.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. CF dye and Mix-n-Stain are trademarks of Biotium.









