

Revised: January 6, 2019

Product Information

CFDA (Carboxyfluorescein diacetate)

Catalog no.	Dye	Size
51018	5-CFDA (5-Carboxyfluorescein diacetate, single isomer)	100 mg
51021	6-CFDA (6-Carboxyfluorescein diacetate, single isomer)	100 mg
51014	5-(and 6)-Carboxyfluorescein diacetate (CFDA)	100 mg

Storage and Handling

Store at -20°C desiccated and protected from light. Product is stable for up to 12 months from date of receipt when stored as recommended as a solid.

Stock solutions may be prepared in anhydrous DMSO (catalog no. 90082). Solutions can be aliquoted and stored with desiccant and protected from light at -20°C, for 12 months.

Molecular Information: $C_{25}H_{16}O_{9}$ Molecular Weight: 460 Color and Form: Off-white solid Absorption/Emission: 492/517 nm (pH 9, after hydrolysis) Extinction Coefficient: 78,000 (pH 9, after hydrolysis)

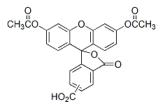


Figure 1. 5-(and 6)-Carboxyfluorescein diacetate (CFDA).

Product Description

Carboxyfluorescein diacetate (CFDA) is a membrane-permeant dye that fluoresces after being processed by intracellular esterases. Once processed, the carboxyfluorescein product has a pH-dependent spectral response similar to fluorescein, and requires cell membrane integrity to retain fluorescence within the cell. Therefore, CFDA can be used to monitor cell viability and intracellular pH.

References

1) Boitano, S., et al. *J. Cell Sci.* **98**, 343(1991). 2) Goodall, H., et al. *Nature* **295**, 524(1982). 3) Hansson, Y., et al. *J. Immunol. Meth.* **100**, 261(1987). 4) Bruning, J.W., et al. J. Immunol. Meth. **33**, 33(1980).

Related Products

Catalog number	Product
90082	DMSO, Anhydrous
90040	Carboxy-2',7'-dichlorofluorescein diacetate, succinimidyl ester, 5 and 6 isomers
90041	CFDA SE, 5 and 6 isomers
51015	5-(and-6)-Carboxy-2',7'-dichlorofluorescein
51016	5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate
51017	5-Carboxy-2',7'-dichlorofluorescein (single isomer)
51023	5-Carboxy-2',7'-dichlorosulfonefluorescein
51012	BCECF, AM Ester
51009	BCECF, AM Ester, 1 mg/mL in Anhydrous DMSO
51010	BCECF, Free Acid
30050	ViaFluor® CFSE Cell Proliferation Kit
30068	ViaFluor® 405 SE Cell Proliferation Kit
30086	ViaFluor® 488 SE Cell Proliferation Kit
32002- 32013	Live-or-Dye ™ Fixable Viability Staining Kits
30026	Calcein AM Cell Viability Assay

Please visit our website at www.biotium.com to view our full selection of cytoplasmic stains, including our amine-reactive ViaFluor® SE dyes, as well as our other viability and organelle stains for microscopy or flow cytometry. Also, see our full catalog of life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF® dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.