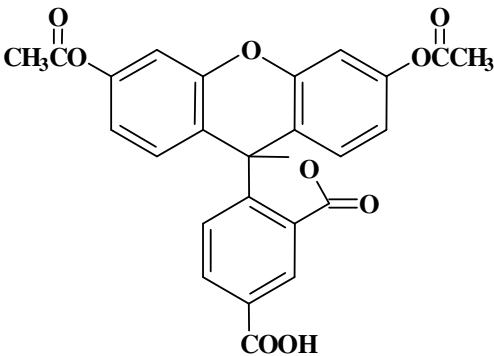


5-(and-6)-Carboxyfluorescein diacetate

Instruction Manual

Catalog Number	PK-CA707-51014
Description	5(6)-CFDA is membrane-permeant and thus can be loaded into cells via incubation. Once inside the cells, CFDA is hydrolyzed by intracellular esterases to carboxyfluorescein (Cat.No. PK-CA707-51013).
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{exc}/\lambda_{em} = 492/517$ nm (pH 9.0, after hydrolysis); Extinction coefficient = 78,000 (pH 9.0, after hydrolysis)
Molecular Structure	
Molecular Weight / Molecular Formula	460 Da; C ₂₅ H ₁₆ O ₉
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	Off-white solid; soluble in DMSO.
Storage & Stability	Stored desiccated at $\leq 4^{\circ}\text{C}$ upon receipt. Protect from light, especially when in solution.
Applications	Fluorescent pH indicator
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)
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.