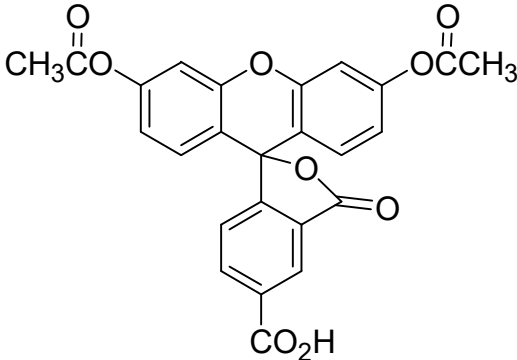


5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate (mixed isomer)

Instruction Manual

Catalog Number	PK-CA707-51016
Description	5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate is membrane-permeant and thus can be loaded into cells via incubation. The diacetate is readily hydrolyzed to 5-(and-6)-carboxy-2',7'-dichlorofluorescein (see Cat.No. PK-CA707-51015) by intracellular esterases. With a lower pKa (4.8), this water-soluble fluorescent tracer is useful for detecting pH in more acidic environment.
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em} = 495/529 \text{ nm}$
Molecular Structure	
Molecular Weight / Molecular Formula	529 Da; C ₂₅ H ₁₄ Cl ₂ O ₉
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Light-yellow solid; soluble in DMSO.
Storage & Stability	Store desiccated at RT or 4°C. Protect from light (especially when in solution).
Applications	NA
References	1) J. Cell Biol. 111, 877(1990) 2) Biochemistry 26, 6330(1987) 3) Biochim. Biophys. Acta 985, 75(1989) 4) Anal. Biochem. 187, 109(1990)
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.