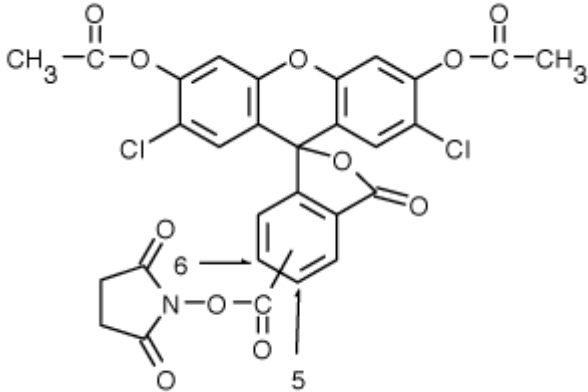


5-(and-6)-Carboxy-2',7'-dichloro-fluorescein diacetate, succinimidyl ester

Instruction Manual

Catalog Number	PK-CA707-90040
Description	5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate, succinimidyl ester is a useful fluorescent tracer that can passively diffuse into cells and covalently label intracellular proteins, resulting in long-term cell labeling. The reagent itself is colorless and nonfluorescent but becomes brightly green fluorescent once it is hydrolyzed by intracellular esterases. The fluorescence of the dye is relatively insensitive to pH change and the labeling has been shown to be stable to formaldehyde or glutaldehyde fixation condition.
Quantity	25 mg
Excitation / Emission Maxima	$\lambda_{exc}/\lambda_{em}$ (of hydrolysis product at pH 7) = 504/529 nm
Molecular Structure	
Molecular Weight / Molecular Formula	626 Da; C ₂₉ H ₁₇ Cl ₂ NO ₁₁
Purity	≥95% as determined by TLC.
Appearance / Formulation / Solubility	White solid; soluble in DMSO or DMF.
Storage & Stability	Store at -20°C. Protect from light, especially when in solution.
Applications	See Description
References	Development 116, 1087(1992)
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.