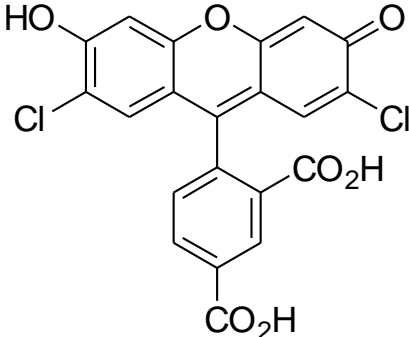


5-Carboxy-2',7'- dichlorofluorescein

(Single isomer)

Instruction Manual

Catalog Number	PK-CA707-51017
Description	This dye can be used as a low pH indicator similar to (PK-CA707-51015), and is also useful as a general green fluorescent label. Carboxy-2',7'-dichlorofluorescein is significantly more photo-stable than carboxyfluorescein.
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em} = 504/529$ nm (pH 8.0); Extinction coefficient = 90,000
Molecular Structure	
Molecular Weight / Molecular Formula	445 Da; C ₂₁ H ₁₀ Cl ₂ O ₇
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Orange solid; soluble in DMSO, DMF or basic water (pH >7).
Storage & Stability	Store at 4°C and protect from light.
Applications	Fluorescent pH indicator
References	1. Bramhall, J., et al. Biochemistry 26, 6330(1987) 2. Knight, C.G., et al. Biochim. Biophys. Acta 985, 75(1989) 3. Nedergaard, M. et al. 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.