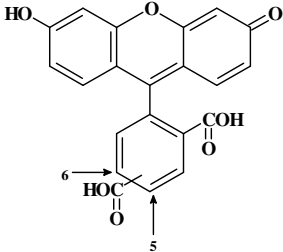


5-(and-6)-Carboxyfluorescein

Fluorescent pH indicator

Instruction Manual

Catalog Number	PK-CA707-51013
Description	Carboxyfluorescein has a pKa of 6.5 and can be used as a pH indicator. Its excitation spectrum and fluorescence response to pH are similar to those of BCECF. Carboxyfluorescein can also be used as a tracer dye. The dye is membrane-impermeant and can be loaded into cells by microinjection or scrape loading.
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (pH 9.0) = 492/514 nm; Extinction coefficient ϵ = 82,000
Molecular Structure	
Molecular Weight / Molecular Formula	376 Da; C ₂₁ H ₁₂ O ₇
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Yellow-orange solid; soluble in water with pH >6.0.
Storage & Stability	Store at 4°C. Protect from light, especially when in solution.
Applications	Fluorescent pH indicator
References	Weinstein, J.N., et al. Methods Enzymol. 128, 65(1986)
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.