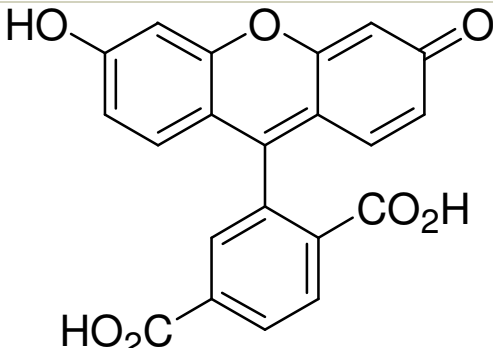


6-Carboxyfluorescein (single isomer)

Instruction Manual

Catalog Number	PK-CA707-51020
Description	Free acid form of 6-carboxyfluorescein single isomer. 6-FAM is one of the most popular green fluorescent reagents used for in situ labeling peptides, proteins and nucleotides. It has also been used to prepare various small fluorescent molecules.
Quantity	100 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (pH 9.0) = 495/519 nm; Extinction coefficient ϵ = 82,000 (pH 9.0)
Molecular Structure	
Molecular Weight / Molecular Formula	376.3 Da; C ₂₁ H ₁₂ O ₇
Purity	>98% (as determined by HPLC)
Appearance / Formulation / Solubility	Yellow-orange solid; soluble in DMSO, DMF and water (pH > 6.0). pKa = 6.5
Storage & Stability	Store at 4°C. Protect from light, especially when in solution).
Applications	Fluorescent pH indicator
References	Nucleic Acids Res 27, 1912-8(1999) Appl Environ Microbiol 62, 1347-53 (1996)
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.