

(Single isomer)

Instruction Manual

Catalog Number	PK-CA707-51023
Description	The pKa of 5-carboxy-2',7'-dichlorosulfonefluorescein is near 4 and thus the dye is potentially an excellent pH indicator for acidic organelles. It is suitable for low pH detection.
Quantity	10 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (pH 9.0) = 509/532 nm; Extinction coefficient = 91,800 (pH 9.0)
Molecular Structure	
Molecular Weight / Molecular Formula	481.26 Da; C ₂₀ H ₁₀ Cl ₂ O ₈ S
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Yellow solid; soluble in DMSO, DMF or buffers with pH >6.0.
Storage & Stability	Store at 4°C. Protect from light (especially when in solution).
Applications	Fluorescent pH indicator
References	NA
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.