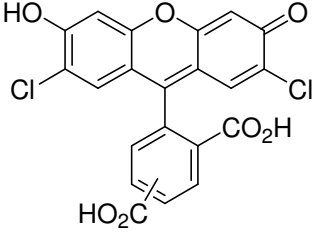


(Mixed isomer)

## Instruction Manual

Catalog Number	PK-CA707-51015
Description	5-(and-6)-carboxy-2',7'-dichlorofluorescein is similar to carboxyfluorescein (PK-CA707-51013) but has lower pKa (4.8). Thus, this water-soluble tracer is useful for detecting pH in more acidic environment. Please see also Cat.No. PK-CA707-51016.
Quantity	200 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em} = 504/529$ nm Extinction coefficient $\epsilon = 90,000$
Molecular Structure	
Molecular Weight / Molecular Formula	445 Da; C <sub>21</sub> H <sub>10</sub> Cl <sub>2</sub> O <sub>7</sub>
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Yellow solid; soluble in DMSO, DMF or basic water (pH >7).
Storage & Stability	Store at 4°C. Protect from light (especially when in solution).
Applications	Fluorescent pH indicator
References	<ol style="list-style-type: none"> <li>1. Bramhall, J., et al. Biochemistry 26, 6330(1987)</li> <li>2. Knight, C.G., et al. Biochim. Biophys. Acta 985, 75(1989)</li> <li>3. Nedergaard, M. et al. Anal. Biochem. 187, 109(1990)</li> </ol>
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.