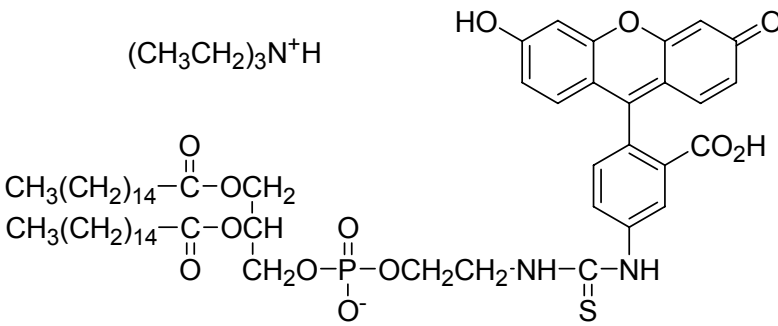


N-(fluorescein-5-thiocarbamoyl)-1,2-dihexadecanoyl-*sn*-glycero-3-phosphoethanolamine, triethylammonium salt

Instruction Manual

Catalog Number	PK-CA707-60024
Description	Fluorescein-DHPE is a useful surface pH indicator and has also been used for measuring lateral diffusion in membranes via fluorescence photobleaching recovery method.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em}$ (in MeOH) = 496/519 nm Extinction Coefficient $\geq 80,000$
Molecular Structure	 <p>(CH₃CH₂)₃N⁺H</p>
Molecular Weight / Molecular Formula	1182.53 Da; C ₆₄ H ₁₁₀ ON ₃ O ₁₃ PS
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	Orange solid; soluble in chloroform. Note: Chloroform may degrade the O-ring in the cap of the product vial. For long-term storage after reconstitution in chloroform, transfer the solution to a vial without an O-ring.
Storage & Stability	Store at -20°C. Protect from light.
Applications	See Description
References	1) Biochim. Biophys. Acta 939, 289(1988) 2) Biophys. J. 66, 25(1994) 3) J. Cell Biol. 103, 807(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.