

N-(Texas Red sulfonyl)-1,2-dihexadecanoyl-*sn*-glycero-3-phosphoethanolamine, triethylammonium salt*

Instruction Manual

Catalog Number	PK-CA707-60027
Description	Rhodamine-labeled glycerophosphoethanolamine lipids such as Rhodamine- DHPE (Cat.No. PK-CA707-60026), TRITC-DHPE (Cat.No. PK-CA707-60028) and Texas Red-DHPE (Cat.No. PK-CA707-60027) have been used as fluorescence energy acceptors in combination with NBDPE as the donor in membrane fusion assays using the principle of fluorescence energy transfer (FRET). These probes have also been used for following membrane trafficking during endocytosis.
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{\text{exc}}/\lambda_{\text{em}}$ (in MeOH) = 582/601 nm; Extinction Coefficient: 90,000
Molecular Structure	
Molecular Weight / Molecular Formula	1380.77 Da; C ₇₄ H ₁₁₇ N ₄ O ₁₄ PS ₂
Purity	>90% (as determined by TLC)
Appearance / Formulation / Solubility	Dark 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	<ol style="list-style-type: none"> 1) Biochemistry 24, 6390 (1985) 2) Biochemistry 21, 1720(1982) 3) Biochemistry 20, 4093(1981) 4) Biochim. Biophys. Acta 1103, 185(1992) 5) Eur. J. Cell Biol. 53, 173(1990)
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.

* Texas Red is a trademark of Molecular Probes.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.