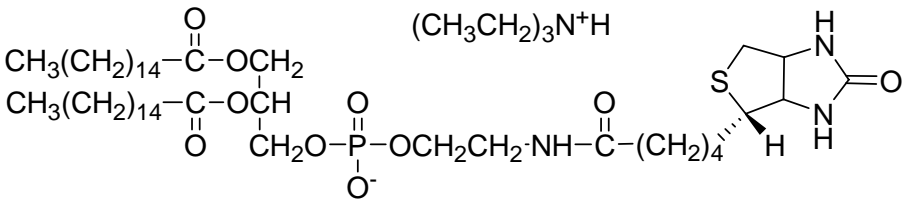


N-(biotinoyl)-1,2-dihexadecanoyl-*sn*-glycero-3-phosphoethanolamine, triethylammonium salt

Instruction Manual

Catalog Number	PK-CA707-60022
Description	Biotin-DHPE can be used to couple avidin or streptavidin to cell membranes or liposomes.
Quantity	10 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	1019.44 Da; C ₅₃ H ₁₀₃ N ₄ O ₁₀ PS
Purity	NA
Appearance / Formulation / Solubility	White solid; soluble in chloroform.
Storage & Stability	Store at -20C. 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.
Applications	See Description
References	1) Immunol. Meth. 158, 183(1993) 2) Anal. Biochem. 207, 341 (1992) 3) Biophys. J. 59, 387(1991)
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.