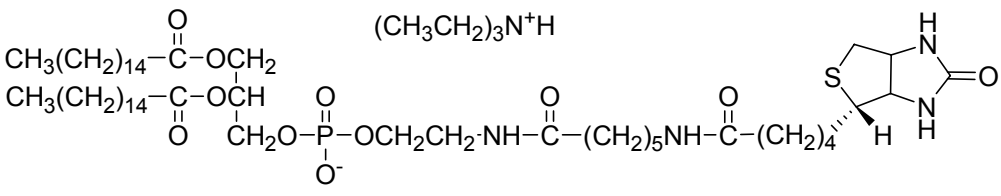


N-((6-(biotinoyl)amino)hexanoyl)-1,2-dihexadecanoyl-*sn*-glycero-3-phosphoethanolamine, triethylammonium salt

Instruction Manual

| | |
|---------------------------------------|---|
| Catalog Number | PK-CA707-60023 |
| Description | Biotin-X-DHPE can be used to couple avidin or streptavidin to cell membranes or liposomes. |
| Quantity | 5 mg |
| Excitation / Emission Maxima | $\lambda_{\text{ex}}/\lambda_{\text{em}}$ (in MeOH) = 362/433 nm |
| Molecular Structure |  |
| Molecular Weight / Molecular Formula | 1132.60 Da; C ₅₉ H ₁₁₄ N ₅ O ₁₁ PS |
| Purity | NA |
| Appearance / Formulation / Solubility | White 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. |
| 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.