

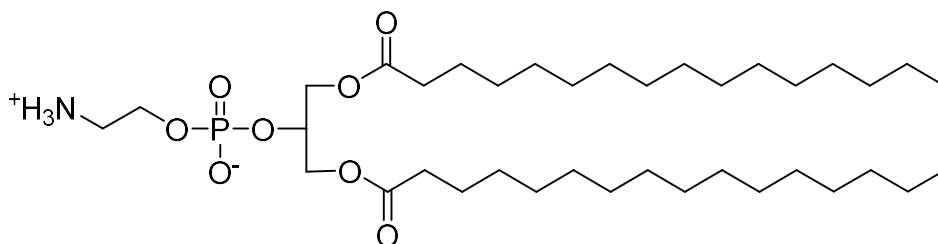
1,3-Dipalmitoyl-*sn*-glycero-2-phosphoethanolamine (1,3-DPPE)

Catalog number: L-2216

Molecular Formula: C₃₇H₇₄NO₈P

MW: 691.96

CAS: 67303-93-9



Alternate Name: 1,3-Dipalmitoyl phosphatidylethanolamine

Solubility: Chloroform > 10 mg/mL

Storage and Handling: Store dry at 4 °C. Stock solutions should be stored frozen (-20 °C or below).

Background: Phosphatidylethanolamine (PE) is the second most abundant phospholipid in mammalian cell membranes. It is formed by two pathways, the decarboxylation of phosphatidylserine or the reaction of CDP-ethanolamine and diacylglycerol. PE is one of the major components of cell membranes in bacteria and in nervous system.

References: 1) Vance, J.E., G. Tassavera. (2013) "Formation and function of phosphatidylserine and phosphatidylethanolamine in mammalian cells" *Biochim. Biophys. Acta* 1831(3):543-554.
2) Rockefeller, P, et al. (2015) "Phosphatidylethanolamine positively regulates autophagy and longevity" *Cell Death and Differentiation* 22: 499-508.
3) Vance, J.E. (2008) "Phosphatidylserine and phosphatidylethanolamine in mammalian cells: two metabolically related aminophospholipids." *J. Lipid Res.* 49(7): 1377-1387.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

