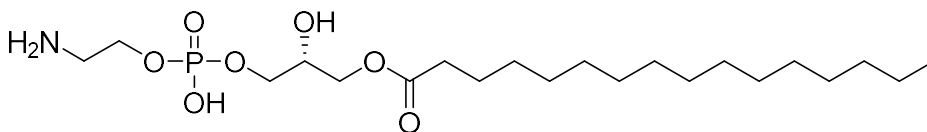


## 16:0-LysoPE (LPE)

**Catalog number:** L-2516



**Molecular Formula:** C<sub>21</sub>H<sub>43</sub>NNaO<sub>7</sub>P

**MW:** 453.55

**CAS#:** 53862-35-4

**Alternate Name:** 16:0 LPE, 1-Palmitoyl-2-hydroxy-*sn*-glycero-3-phospho-ethanolamine

**Solubility:** methanol, > 5 mg/mL

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

**Background:** Lysophospholipids are lipid mediators which elicit a diverse array of cellular responses through activation of phospholipid specific G protein-coupled receptors. Lysophosphatidylethanolamine (LPE) is a minor lipid found in low levels in plasma membranes and serum which is generated through PLA2. LPE has been shown to increase intracellular Ca<sup>2+</sup>, invasiveness, and migration in SKOV3 and MDA-MB-231 cells. In addition, LPE accelerates ripening and increase the shelf-life of some fruits and vegetables such as tomatoes, grapes, and cranberries.

**References:** 1). Kumiko Makide, et al "Novel lysophospholipid receptors: their structure and function" *Journal of Lipid Research*, **2014**, 55, 1986.

2) K.S. Park, et al, "Lysophosphatidylethanolamine stimulates chemotactic migration and cellular invasion in SK-OV3 human ovarian cancer cells: involvement of pertussis toxin-sensitive G-protein coupled receptor." *FEBS Lett.*, **2007**, 581, 4411-4416.

**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.

Technical Data Sheet, Rev 3, 6/3/2024 – **For research use only.** Not intended for diagnostic or therapeutic use.

