

# PRODUCT INFORMATION



## 1-Stearoyl-2-hydroxy-*sn*-glycero-3-PE

Item No. 25593

CAS Registry No.: 69747-55-3

Formal Name: octadecanoic acid, (2R)-3-[[[2-aminoethoxy]hydroxyphosphinyl]oxy]-2-hydroxypropyl ester

Synonyms: 18:0 LPE, 18:0 Lyso-PE, 1-Octadecanoyl-*sn*-glycero-3-phosphoethanolamine, 1-Stearoyl-2-hydroxy-*sn*-glycero-3-phosphatidylethanolamine, 1-Stearoyl LPE, 1-Stearoyl lyso-PE

MF: C<sub>23</sub>H<sub>48</sub>NO<sub>7</sub>P

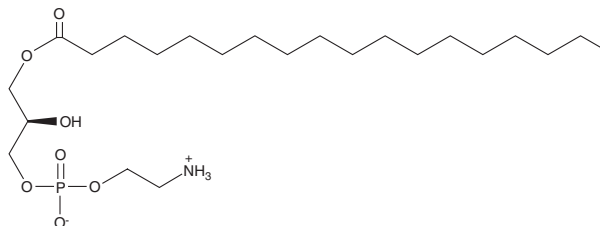
FW: 481.6

Purity: ≥95% (9:1 mixture of 1-acyl and 2-acyl)

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

1-Stearoyl-2-hydroxy-*sn*-glycero-3-PE (18:0 LPE) is supplied as a solid. A stock solution may be made by dissolving the 18:0 LPE in the solvent of choice, which should be purged with an inert gas. 18:0 LPE is slightly soluble in chloroform.

### Description

18:0 LPE is a lysophospholipid.<sup>1</sup> It induces transient increases in intracellular calcium in PC12 rat neuronal cells in a concentration-dependent manner, an effect that can be blocked by the lysophosphatidic acid receptor 1 (LPA<sub>1</sub>) receptor antagonist AM095 (Item No. 22141).

### Reference

1. Lee, J.M., Park, S.J., and Im, D.S. Lysophosphatidylethanolamine increases intracellular Ca<sup>2+</sup> through LPA<sub>1</sub> in PC-12 neuronal cells. *Biochem. Biophys. Res. Commun.* **461**(2), 378-382 (2015).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/16/2025

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM