

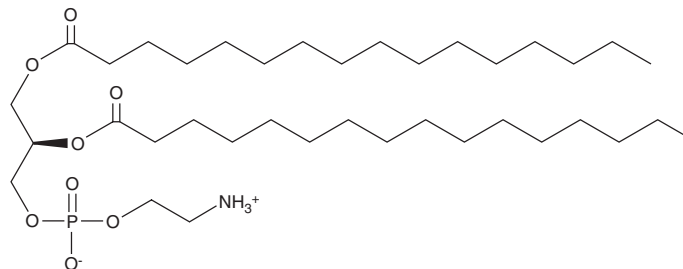
# PRODUCT INFORMATION



## 1,2-Dipalmitoyl-*sn*-glycero-3-PE

Item No. 15092

**CAS Registry No.:** 923-61-5  
**Formal Name:** hexadecanoic acid, 1,1'-[(1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]methyl]-1,2-ethanediyl] ester  
**Synonyms:** 1,2-Dipalmitoyl-*sn*-glycerol-3-Phosphoethanolamine, 1,2-DPPE  
**MF:** C<sub>37</sub>H<sub>74</sub>NO<sub>8</sub>P  
**FW:** 692.0  
**Purity:** ≥98%  
**Supplied as:** A crystalline 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

1,2-Dipalmitoyl-*sn*-glycero-3-PE is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dipalmitoyl-*sn*-glycero-3-PE in the solvent of choice, which should be purged with an inert gas. 1,2-Dipalmitoyl-*sn*-glycero-3-PE is soluble in the organic solvent chloroform at a concentration of approximately 3 mg/ml.

### Description

1,2-Dipalmitoyl-*sn*-glycero-3-PE (1,2-DPPE) is a naturally-occurring PE containing 16:0 fatty acids at the *sn*-1 and *sn*-2 positions. It belongs to a class of phospholipids that are the most abundant lipids in the inner leaflet of the plasma membrane.<sup>1</sup> 1,2-DPPE interacts with cholesterol to form a condensed lipid monolayer with tight hydrogen bonding of the 1,2-DPPE interheadgroups, resulting in a more fluid membrane that may aid in transport and signaling across the bilayer.<sup>2,3</sup>

### References

1. Vance, J.E. and Tasseva, G. Formation and function of phosphatidylserine and phosphatidylethanolamine in mammalian cells. *Biochim. Biophys. Acta* **1831(3)**, 543-554 (2013).
2. McQuaw, C.M., Sostarecz, A.G., Zheng, L., *et al.* Lateral heterogeneity of dipalmitoylphosphatidylethanolamine-cholesterol Langmuir-Blodgett films investigated with imaging time-of-flight secondary ion mass spectrometry and atomic force microscopy. *Langmuir* **21(3)**, 807-813 (2005).
3. Leekumjorn, S. and Sum, A.K. Molecular simulation study of structural and dynamic properties of mixed DPPC/DPPE bilayers. *Biophys. J.* **90(11)**, 3951-3965 (2006).

#### 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, 11/14/2022

#### 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