

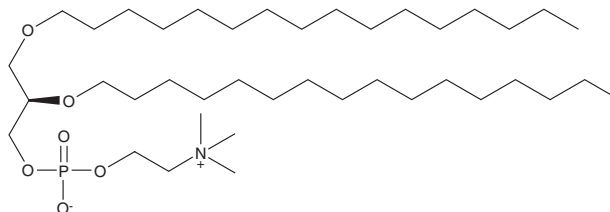
# PRODUCT INFORMATION



## 1,2-Dihexadecyl-*sn*-glycero-3-PC

Item No. 25588

**CAS Registry No.:** 36314-47-3  
**Formal Name:** (7R)-7-(hexadecyloxy)-4-hydroxy-N,N,N-trimethyl-3,5,9-trioxa-4-phosphapentacosan-1-aminium, 4-oxide, inner salt  
**Synonym:** 1,2-Dihexadecyl-*sn*-glycero-3-Phosphocholine  
**MF:** C<sub>40</sub>H<sub>84</sub>NO<sub>6</sub>P  
**FW:** 706.1  
**Purity:** ≥95%  
**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-Dihexadecyl-*sn*-glycero-3-PC is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dihexadecyl-*sn*-glycero-3-PC in the solvent of choice. 1,2-Dihexadecyl-*sn*-glycero-3-PC is soluble in the organic solvent ethanol, which should be purged with an inert gas, at a concentration of approximately 30 mg/ml.

### Description

1,2-Dihexadecyl-*sn*-glycero-3-PC is a synthetic ether-linked phospholipid containing hexadecyl groups at the *sn*-1 and *sn*-2 positions. It is commonly used in the generation of liposomes and artificial membranes to study membrane dynamics.<sup>1,2</sup>

### References

1. Fernández, M.S. and Juárez, J.A. Activity of phospholipase A<sub>2</sub> on a fluorescent substrate incorporated into non-hydrolyzable phospholipid liposomes. *Biochim. Biophys. Acta.* **1192(1)**, 132-142 (1994).
2. Mukherjee, S. and Chattopadhyay, A. Influence of ester and ether linkage in phospholipids on the environment and dynamics of the membrane interface: A wavelength-selective fluorescence approach. *Langmuir* **21(1)**, 287-293 (2005).

#### 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, 12/19/2022

#### 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