

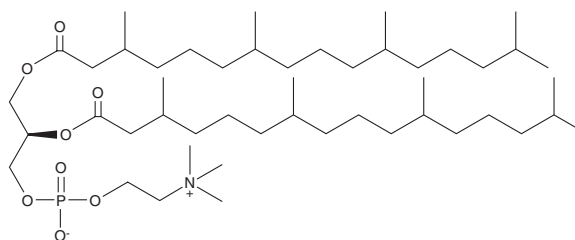
PRODUCT INFORMATION



1,2-Diphytanoyl-*sn*-glycero-3-PC

Item No. 25598

CAS Registry No.: 207131-40-6
Formal Name: (7R)-4-hydroxy-N,N,N,12,16,20,24-heptamethyl-10-oxo-7-[(3,7,11,15-tetramethyl-1-oxohexadecyl)oxy]-3,5,9-trioxa-4-phosphapentacosan-1-aminium, 4-oxide, inner salt
Synonyms: Diphytanoyl Phosphatidylcholine, 1,2-DPhPC, 4ME 16:0 PC
MF: C₄₈H₉₆NO₈P
FW: 846.3
Purity: ≥98%
Supplied as: A solution in chloroform
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

1,2-Diphytanoyl-*sn*-glycero-3-PC is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice.

Description

1,2-Diphytanoyl-*sn*-glycero-3-PC is a phospholipid containing the tetramethylated long-chain (16:0) diphytanic acid at the *sn*-1 and *sn*-2 positions. It has been used in the creation of lipid bilayer membranes.^{1,2} 1,2-Diphytanoyl-*sn*-glycero-3-PC, in combination with LPS, has been used to form the outer leaflet of a tethered bilayer lipid membrane as a model system to examine the effect of antibiotics on Gram-negative bacteria membranes.²

References

1. Schuster, B., Pum, D., Braha, O., *et al.* Self-assembled α -hemolysin pores in an S-layer-supported lipid bilayer. *Biochim. Biophys. Acta.* **1370**(2), 280-288 (1998).
2. Andersson, J., Fuller, M.A., Wood, K., *et al.* A tethered bilayer lipid membrane that mimics microbial membranes. *Phys. Chem. Chem. Phys.* **20**(18), 12958-12969 (2018).

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, 03/28/2022

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