

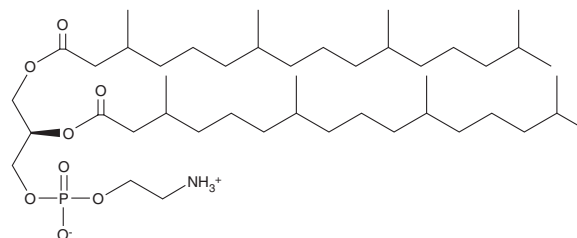
PRODUCT INFORMATION



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

Item No. 15094

CAS Registry No.: 201036-16-0
Formal Name: 1,2-diphytanoyl-*sn*-glycero-3-phosphatidylethanolamine
Synonyms: 1,2-Diphytanoyl-*sn*-glycero-3-Phosphoethanolamine, DiPPE, 1,2-DPyPE, 4ME 16:0 PE
MF: C₄₅H₉₀NO₈P
FW: 804.2
Purity: ≥95%
Supplied as: A 80:20 solution in chloroform:methanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

1,2-Diphytanoyl-*sn*-glycero-3-PE (1,2-DPyPE) is supplied as an 80:20 solution in chloroform:methanol. A stock solution may be made by dissolving the DPyPE in the solvent of choice, which should be purged with an inert gas. 1,2-DPyPE is soluble in the organic solvent chloroform at a concentration of approximately 3 mg/ml.

Description

Phosphatidylethanolamines (PEs) are phospholipids found in biological membranes that serve both structural and functional roles.^{1,2} Different types of PE are commonly used in the generation of micelles, liposomes, and other types of artificial membranes.^{3,4} 1,2-DPyPE is a phospholipid containing the tetramethylated long-chain (16:0) diphytanic acid inserted at the *sn*-1 and *sn*-2 positions.

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. Wellner, N., Diep, T.A., Janfelt, C., *et al.* N-acylation of phosphatidylethanolamine and its biological functions in mammals. *Biochim. Biophys. Acta*. **1831(3)**, 652-662 (2013).
3. Simões, S., Moreira, J.N., Fonseca, C., *et al.* On the formulation of pH-sensitive liposomes with long circulation times. *Adv. Drug Deliv. Rev.* **56(7)**, 947-965 (2004).
4. Fattal, E., Couvreur, P., and Dubernet, C. "Smart" delivery of antisense oligonucleotides by anionic pH-sensitive liposomes. *Adv. Drug Deliv. Rev.* **56(7)**, 931-946 (2004).

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, 01/16/2023

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