

PRODUCT INFORMATION



1,2-Dioleoyl-*sn*-glycero-3-phosphoethanol (sodium salt)

Item No. 45185

CAS Registry No.: 322647-53-0

Formal Name: 9-octadecenoic acid (9Z), (1R)-1-
[[[ethoxyhydroxyphosphinyl]oxy]
methyl]-1,2-ethanediyl ester,
monosodium salt

Synonyms: 1,2-Di-(9(Z)-Octadecenoyl)-
sn-glycero-3-phosphoethanol,
1,2-Dioleoyl Phosphatidylethanol,
PEth(18:1/18:1),
PEth(18:1(9Z)/18:1(9Z)),
PtdEth-(1,2-dioleoyl)

MF: C₄₁H₇₆O₈P • Na

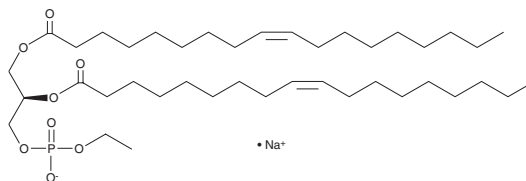
FW: 751.0

Purity: ≥98%

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

1,2-Dioleoyl-*sn*-glycero-3-phosphoethanol (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 1,2-dioleoyl-*sn*-glycero-3-phosphoethanol (sodium salt) in the solvent of choice, which should be purged with an inert gas. 1,2-Dioleoyl-*sn*-glycero-3-phosphoethanol (sodium salt) is sparingly soluble (1-10 mg/ml) in methanol.

Description

1,2-Dioleoyl-*sn*-glycero-3-phosphoethanol is a glycerophospholipid and homolog of 1-palmitoyl-2-oleoyl-*sn*-glycero-3-phosphoethanol, which has been used as a biomarker of alcohol consumption.¹ It has been found in postmortem human blood samples from individuals with moderate and high blood alcohol concentrations.²

References

1. Bantle, M., Long, J.O.T., Brydon, S.C., *et al.* Isomer composition assessment of synthetic phosphatidylethanol by collision-induced and ozone-induced dissociation mass spectrometry. *Drug Test. Anal.* **18**(4), 501-509 (2026).
2. Gnann, H., Weinmann, W., Engelmann, C., *et al.* Selective detection of phosphatidylethanol homologues in blood as biomarkers for alcohol consumption by LC-ESI-MS/MS. *J. Mass Spectrom.* **44**(9), 1293-1299 (2009).

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, 04/23/2026

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