

PRODUCT INFORMATION



1-Oleoyl-2-palmitoyl-*sn*-glycero-3-phosphoethanol (sodium salt)

Item No. 45069

CAS Registry No.: 2410279-88-6

Formal Name: 9Z-octadecenoic acid, (2R)-3-[(ethoxyhydroxyphosphinyl)oxy]2-[(1-oxohexadecyl)oxy]propyl ester, monosodium salt

Synonyms: 1-(9(Z)-Octadecenoyl)-2-hexadecanoyl-*sn*-glycero-3-phosphoethanol, 1-Oleoyl-2-Palmitoyl Phosphatidylethanol, PEth(18:1/16:0), PEth(18:1(9Z)/16:0), PtdEth-(1-oleoyl, 2-palmitoyl)

MF: $C_{39}H_{74}O_8P \cdot Na$

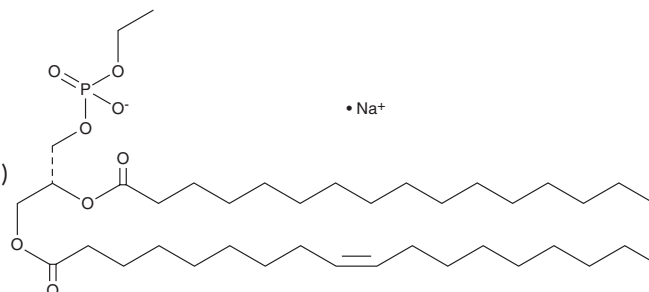
FW: 725.0

Purity: $\geq 99\%$

Supplied as: A 10 mg/ml solution in chloroform

Storage: $-20^{\circ}C$

Stability: ≥ 1 year



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

Description

1-Oleoyl-2-palmitoyl-*sn*-glycero-3-phosphoethanol is a glycerophospholipid and regioisomer of 1-palmitoyl-2-oleoyl-*sn*-glycero-3-phosphoethanol, which has been used as a biomarker of alcohol consumption.¹

Reference

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).

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