

PRODUCT INFORMATION



1-(1(Z)-Hexadecenyl)-2-linoleoyl-*sn*-glycero-3-phosphoethanolamine

Item No. 9004999

CAS Registry No.: 802981-98-2

Formal Name: 9Z,12Z-octadecadienoic acid, (1R)-2-[[[2-aminoethoxy]hydroxyphosphinyl]oxy]-1-[[[(1Z)-1-hexadecen-1-yloxy]methyl]ethyl ester

Synonyms: C16(plasm)-18:2-PE, 1-(1(Z)-Hexadecenyl)-2-(9(Z),12(Z)-Octadecadienoyl)-*sn*-glycero-3-phosphoethanolamine, 1-(1(Z)-Hexadecenyl)-2-linoleoyl Phosphatidylethanolamine, PE(P-16:0/18:2), PE(P-16:0/18:2(9Z,12Z)), PtdEtn-(1-(1(Z)hexadecenyl), 2-linoleoyl)

MF: $C_{39}H_{74}NO_7P$

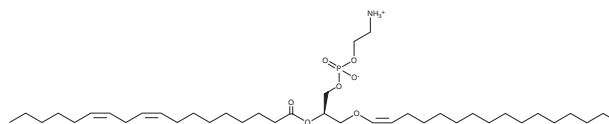
FW: 700.0

Purity: $\geq 95\%$ (9:1 mixture of *cis:trans*)

Supplied as: A 5 mg/ml solution in ethanol

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-(1(Z)-Hexadecenyl)-2-linoleoyl-*sn*-glycero-3-phosphoethanolamine is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvent such as chloroform purged with an inert gas can be used. 1-(1(Z)-Hexadecenyl)-2-linoleoyl-*sn*-glycero-3-phosphoethanolamine is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

1-(1(Z)-Hexadecenyl)-2-linoleoyl-*sn*-glycero-3-phosphoethanolamine is a plasmalogen that contains 1(Z)-hexadecenoic acid and linoleic acid (Item Nos. 90150 | 90150.1 | 21909) at the *sn*-1 and *sn*-2 positions, respectively.¹ It inhibits lead-induced decreases in viability and increases in reactive oxygen species (ROS) and oxidized phospholipids in HepG2 hepatocytes.

Reference

1. Wu, Y., Chen, Z., Darwish, W.S., *et al.* Choline and ethanolamine plasmalogens prevent lead-induced cytotoxicity and lipid oxidation in HepG2 cells. *J. Agric. Food Chem.* **67**(27), 7716-7725 (2019).

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, 04/29/2026

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