

# PRODUCT INFORMATION



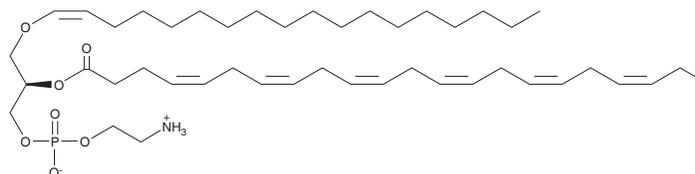
## 1-1(Z)-Octadecenyl-2-Docosahexaenoyl-*sn*-glycero-3-PE

Item No. 37138

**CAS Registry No.:** 206059-98-5  
**Formal Name:** 4Z,7Z,10Z,13Z,16Z,19Z-docosahexaenoic acid, (1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]methyl]-2-[(1Z)-1-octadecen-1-yloxy]ethyl ester

**Synonyms:** 1-1(Z)-Octadecenyl-2-Docosahexaenoyl-*sn*-glycero-3-Phosphatidylethanolamine, 1-1(Z)-Octadecenyl-2-Docosahexaenoyl-*sn*-glycero-3-Phosphoethanolamine, 18:0p/22:6-PE, C18(Plasm)-22:6-PE, PE(P-18:0/22:6)

**MF:** C<sub>45</sub>H<sub>78</sub>NO<sub>7</sub>P  
**FW:** 776.1  
**Purity:** ≥95%  
**Supplied as:** A 5 mg/ml solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Description

1-1(Z)-Octadecenyl-2-Docosahexaenoyl-*sn*-glycero-3-PE 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. Solvents such as chloroform purged with an inert gas can be used. The solubility of 1-1(Z)-Octadecenyl-2-Docosahexaenoyl-*sn*-glycero-3-PE in chloroform is approximately 1 mg/ml.

### Reference

1. Takahashi, R., Nakaya, M., Kotaniguchi, M., *et al.* Analysis of phosphatidylethanolamine, phosphatidylcholine, and plasmalogen molecular species in food lipids using an improved 2D high-performance liquid chromatography system. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **1077-1078**, 35-43 (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, 11/06/2025

### 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