

PRODUCT INFORMATION



1-Palmitoyl-2-Docosahexaenoyl-*sn*-glycero-3-PE

Item No. 33440

CAS Registry No.: 96998-00-4
Formal Name: 4Z,7Z,10Z,13Z,16Z,19Z-docosahexaenoic acid, (1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]methyl]-2-[(1-oxohexadecyl)oxy]ethyl ester
Synonyms: 1-Hexadecanoyl-2-(Docosahexaenoyl)-*sn*-glycero-3-Phosphatidylethanolamine, 1-Hexadecanoyl-2-(Docosahexaenoyl)-*sn*-glycero-3-Phosphoethanolamine, 16:0/22:6-PE, PE(16:0/22:6)

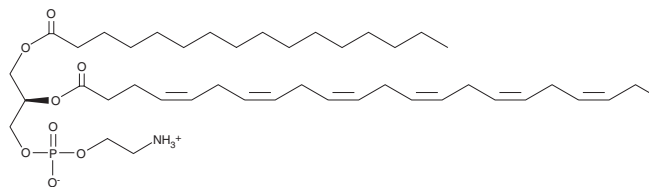
MF: C₄₃H₇₄NO₈P
FW: 764.0

Purity: ≥95%

Supplied as: A solution in chloroform

Storage: -80°C

Stability: ≥6 months



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

Description

1-Palmitoyl-2-docosahexaenoyl-*sn*-glycero-3-PE is a phospholipid that contains palmitic acid (Item No. 10006627) and docosahexaenoic acid (Item Nos. 90310 | 17950) at the *sn*-1 and *sn*-2 positions, respectively. It has been found in the nucleus, but not mitochondria, of mouse liver.¹ Levels of 1-palmitoyl-2-docosahexaenoyl-*sn*-glycero-3-PE are lower in pre-term infant plasma compared with maternal or fetal plasma.²

References

1. Aviram, R., Manella, G., Kopelman, N., *et al.* Lipidomics analyses reveal temporal and spatial lipid organization and uncover daily oscillations in intracellular organelles. *Mol. Cell* **62**(4), 636-648 (2016).
2. Bernhard, W., Raith, M., Koch, V., *et al.* Plasma phospholipids indicate impaired fatty acid homeostasis in preterm infants. *Eur. J. Nutr.* **53**(7), 1533-1547 (2014).

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, 04/01/2021

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