

PRODUCT INFORMATION



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

Item No. 25660

CAS Registry No.: 70812-59-8

Formal Name: (5Z,8Z,11Z,14Z)-5,8,11,14-eicosatetraenoic acid, (1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]methyl]-2-[(1-oxohexadecyl)oxy]ethyl ester

Synonyms: 1-Hexadecanoyl-2-Eicosatetraenoyl-*sn*-glycero-3-Phosphatidylethanolamine, 1-Hexadecanoyl-2-Eicosatetraenoyl-*sn*-glycero-3-Phosphoethanolamine, 1-Palmitoyl-2-arachidonoyl PE, PAPE, PE(16:0/20:4), 16:0/20:4-PE

MF: C₄₁H₇₄NO₈P

FW: 740.0

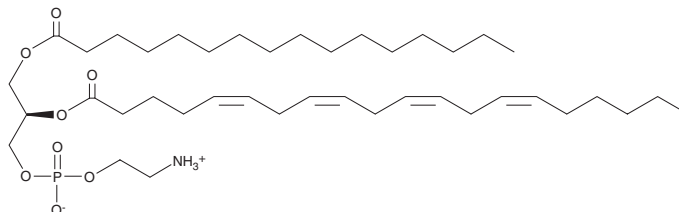
Purity: ≥95%

Supplied as: A solution in methanol:chloroform (7:3) from 100% chloroform

Storage: -80°C

Stability: ≥2years

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



Laboratory Procedures

1-Palmitoyl-2-arachidonoyl-*sn*-glycero-3-PE is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice.

Description

1-Palmitoyl-2-arachidonoyl-*sn*-glycero-3-PE is a phospholipid that contains palmitic acid (Item No. 10006627) and arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) at the *sn*-1 and *sn*-2 positions, respectively. It has been found in egg yolk and isolated rat retina.^{1,2}

References

1. Sun, N., Chen, J., Bao, Z., *et al.* Egg yolk phosphatidylethanolamine: Extraction optimization, antioxidative activity, and molecular structure profiling. *J. Food Sci.* **84**(5), 1002-1011 (2019).
2. Gugiu, B.G., Mesaros, C.A., Sun, M., *et al.* Identification of oxidatively truncated ethanolamine phospholipids in retina and their generation from polyunsaturated phosphatidylethanolamines. *Chem. Res. Toxicol.* **19**(2), 262-71 (2006).

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, 02/21/2024

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