

PRODUCT INFORMATION

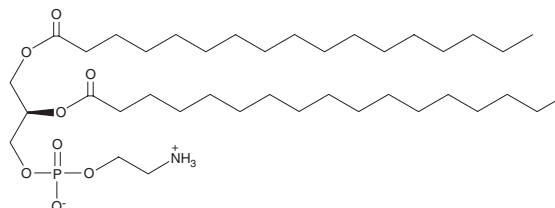


1,2-Diheptadecanoyl-*sn*-glycero-3-PE

Item No. 38319

CAS Registry No.: 140219-78-9
Formal Name: heptadecanoic acid, 1,1'-[(1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]methyl]-1,2-ethanediyl] ester
Synonyms: 17:0/17:0-PE, PE(17:0/17:0), 1,2-Heptadecyloyl-*sn*-glycero-3-Phosphoethanolamine, 1,2-Heptadecyloyl-*sn*-glycero-3-Phosphatidylethanolamine

MF: C₃₉H₇₈NO₈P
FW: 720.0
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1,2-Diheptadecanoyl-*sn*-glycero-3-PE is supplied as a solid. A stock solution may be made by dissolving the 1,2-diheptadecanoyl-*sn*-glycero-3-PE in the solvent of choice, which should be purged with an inert gas. 1,2-Diheptadecanoyl-*sn*-glycero-3-PE is lightly soluble (0.1-1 mg/ml) in chloroform and methanol.

Description

1,2-Diheptadecanoyl-*sn*-glycero-3-PE is a glycerophospholipid containing heptadecanoic acid (Item No. 19722) at the *sn*-1 and *sn*-2 positions. It has been used as an internal standard for the detection of phosphatidylethanolamines in human adipose tissue.¹

Reference

1. Al-Sari, N., Suvitaival, T., Mattila, I., *et al.* Lipidomics of human adipose tissue reveals diversity between body areas. *PLoS One* **15(6)**, e0228521 (2020).

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, 12/05/2025

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