

PRODUCT INFORMATION



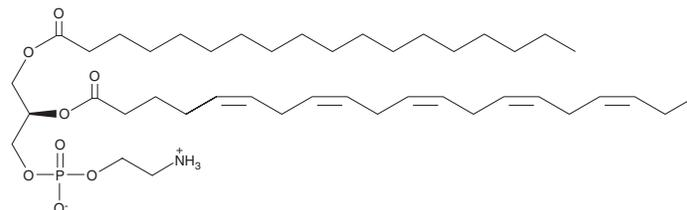
1-Stearoyl-2-Eicosapentaenoyl-*sn*-glycero-3-PE

Item No. 37702

CAS Registry No.: 96935-44-3
Formal Name: 5Z,8Z,11Z,14Z,17Z-eicosapentaenoic acid, (1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]methyl]-2-[(1-oxooctadecyl)oxy]ethyl ester

Synonyms: 1-Octadecanoyl-1-Eicosapentaenoyl-*sn*-glycero-3-Phosphatidylethanolamine, 1-Octadecanoyl-1-Eicosapentaenoyl-*sn*-glycero-3-Phosphoethanolamine, PE(18:0/20:5), 18:0/20:5-PE, 1-Stearoyl-2-Eicosapentaenoyl-*sn*-glycero-3-Phosphatidylethanolamine, 1-Stearoyl-2-Eicosapentaenoyl-*sn*-glycero-3-Phosphoethanolamine

MF: C₄₃H₇₆NO₈P
FW: 766.0
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-Stearoyl-2-eicosapentaenoyl-*sn*-glycero-3-PE is a phospholipid that contains stearic acid (Item No. 10011298) and eicosapentaenoic acid (EPA; Item Nos. 90110 | 90110.1 | 21908) at the *sn*-1 and *sn*-2 positions, respectively. It has been found in fish oils.¹

Reference

1. Yu, X., Wang, Q., Lu, W., *et al.* Fast and specific screening of EPA/DHA-enriched phospholipids in fish oil extracted from different species by HILIC-MS. *J. Agric. Food Chem.* **69(28)**, 7997-8007 (2021).

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