

PRODUCT INFORMATION



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

Item No. 41382

CAS Registry No.: 96998-01-5

Formal Name: 4Z,7Z,10Z,13Z,16Z,19Z-docosahexaenoic acid, (1R)-1-[[[(2-aminoethoxy)hydroxyphosphinyl]oxy]methyl]-2-[(1-oxooctadecyl)oxy]ethyl ester

Synonyms:

18:0/22:6-PE, PE(18:0/22:6), 1-Stearoyl-2-Docosahexaenoyl-*sn*-glycero-3-Phosphoethanolamine, 1-Stearoyl-2-Docosahexaenoyl-*sn*-glycero-3-Phosphatidylethanolamine

MF: $C_{45}H_{78}NO_8P$

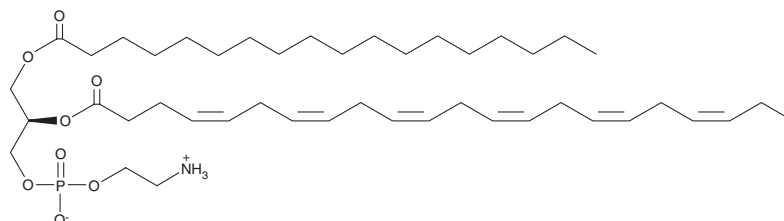
FW: 792.1

Purity: $\geq 95\%$

Supplied as: A solution in methanol:chloroform (30:70) and 0.1% amylenes

Storage: -20°C

Stability: ≥ 1 year



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

Laboratory Procedures

1-Stearoyl-2-docosahexaenoyl-*sn*-glycero-3-PE is supplied as a solution in methanol:chloroform (30:70) and 0.1% amylenes. To change the solvent, simply evaporate the methanol:chloroform (30:70) and 0.1% amylenes under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. 1-Stearoyl-2-docosahexaenoyl-*sn*-glycero-3-PE is sparingly soluble (1-10 mg/ml) in methanol and slightly soluble (0.1-1 mg/ml) in choloform.

Description

1-Stearoyl-2-docosahexaenoyl-*sn*-glyero-3-PE is a phospholipid that contains stearic acid (Item No. 10011298) and docosahexaenoic acid (DHA; Item No. 90310) at the *sn*-1 and *sn*-2 positions, respectively.

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, 09/24/2024

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