

PRODUCT INFORMATION



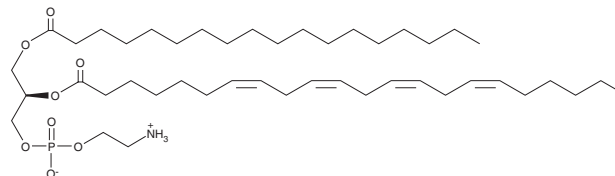
1-Stearoyl-2-Adrenoyl-*sn*-glycero-3-PE

Item No. 33661

CAS Registry No.: 131350-53-3

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

Synonyms: 1-Octadecanoyl-2-Docosatetraenoyl-
sn-glycero-3-Phosphatidylethanolamine,
1-Octadecanoyl-2-Docosatetraenoyl-*sn*-glycero-
3-Phosphoethanolamine, 1-Octadecanoyl-2-
Docosatetraenoyl-*sn*-glycero-3-PE, 1-Stearoyl-2-
Docosatetraenoyl-*sn*-glycero-3-PE, PE(18:0/22:4)



MF: $C_{45}H_{82}NO_8P$

FW: 796.1

Purity: $\geq 95\%$

Supplied as: A 50 mg/ml solution in chloroform

Storage: -20°C

Stability: ≥ 2 years

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

Description

1-Stearoyl-2-adrenoyl-*sn*-glycero-3-PE is a phospholipid that contains stearic acid (Item No. 10011298) and adrenic acid (Item No. 90300) at the *sn*-1 and *sn*-2 positions, respectively. 1-Stearoyl-2-adrenoyl-*sn*-glycero-3-PE levels are inversely correlated with subject age in mitochondria isolated from human post-mortem hippocampus.¹

Reference

1. Hancock, S.E., Friedrich, M.G., Mitchell, T.W., *et al.* Decreases in phospholipids containing adrenic and arachidonic acids occur in the human hippocampus over the adult lifespan. *Lipids* **50**(9), 861-872 (2015).

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, 11/20/2025

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