

PRODUCT INFORMATION



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

Item No. 25595

CAS Registry No.: 59403-52-0
Formal Name: (7R,12Z,15Z,18Z,21Z,24Z,27Z)-4-hydroxy-N,N,N-trimethyl-9-oxo-7-[[[(1-oxooctadecyl)oxy]methyl]-3,5,8-trioxa-4-phosphatriaconta-12,15,18,21,24,27-hexaen-1-aminium, 4-oxide, inner salt

Synonyms: DHA-PC, 18:0/22:6 PC, PC(18:0/22:6(4Z,7Z,10Z,13Z,16Z,19Z)), 1-Stearoyl-2-Docosahexaenoyl-*sn*-glycero-3-Phosphatidylcholine, 1-Stearoyl-2-Docosahexaenoyl-*sn*-glycero-3-Phosphocholine

MF: C₄₈H₈₄NO₈P

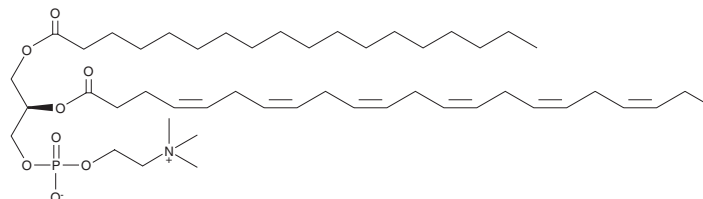
FW: 834.2

Purity: ≥95%

Supplied as: A solution in chloroform

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-docosahexaenoyl-*sn*-glycero-3-PC is a phospholipid containing stearic acid (Item No. 10011298) and docosahexaenoic acid (DHA; Item No. 90310) at the *sn*-1 and *sn*-2 positions, respectively. It has been used in the study of lipid membrane organization and dynamics.¹ The levels of 1-stearoyl-2-docosahexaenoyl-*sn*-glycero-3-PC are decreased in the L5 anterior horn in a superoxide dismutase 1 mutant (SOD1^{G93A}) transgenic mouse model of amyotrophic lateral sclerosis (ALS) at the terminal stage of disease.²

References

1. Leng, X., Kinnun, J.J., Cavazos, A.T., *et al.* All n-3 PUFA are not the same: MD simulations reveal differences in membrane organization for EPA, DHA and DPA. *Biochim. Biophys. Acta Biomembr.* **1860(5)**, 1125-1134 (2018).
2. Arima, H., Omura, T., Hayasaka, T., *et al.* Reductions of docosahexaenoic acid-containing phosphatidylcholine levels in the anterior horn of an ALS mouse model. *Neuroscience* **297**, 127-136 (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, 11/01/2022

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