

PRODUCT INFORMATION



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

Item No. 37340

CAS Registry No.: 162440-05-3
Formal Name: 4,7R-dihydroxy-N,N,N-trimethyl-10-oxo-3,5,9-trioxa-4-phosphahentriaconta-13Z,16Z,19Z,22Z,25Z,28Z-hexaen-1-aminium, 4-oxide, inner salt

Synonyms: 1-Docosahexaenoyl-2-hydroxy-*sn*-glycero-3-Phosphocholine, LysoPC(22:6/0:0), PC(22:6/0:0), 22:6/0:0-PC

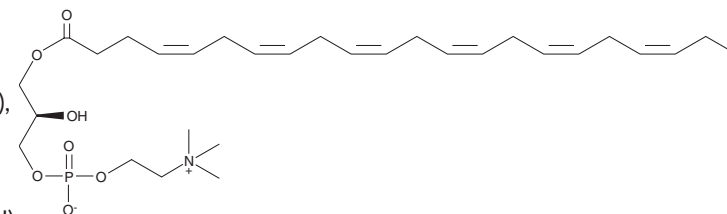
MF: C₃₀H₅₀NO₇P
FW: 567.7

Purity: ≥95% (as an 8:2 mixture of the 1-acyl and 2-acyl)

Supplied as: A solution in chloroform

Storage: -80°C

Stability: ≥1 year



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

Description

1-Docosahexaenoyl-2-hydroxy-*sn*-glycero-3-PC is a lysophospholipid containing docosahexaenoic acid (DHA; Item Nos. 90310 | 17950) at the *sn*-1 position that has been found in shark liver extract.¹ 1-Docosahexaenoyl-2-hydroxy-*sn*-glycero-3-PC decreases LPS-induced increases in TNF- α and IL-6 levels in RAW 264.7 cells in a concentration-dependent manner.² It decreases LPS-induced production of nitric oxide (NO) in RAW 264.7 cells when used at concentrations ranging from 12 to 200 μ M. 1-Docosahexaenoyl-2-hydroxy-*sn*-glycero-3-PC (15-500 μ g/kg) inhibits plasma leakage into the peritoneum in a mouse model of peritonitis induced by zymosan A (Item No. 21175).

References

1. Chen, S. and Li, K.W. Mass spectrometric identification of molecular species of phosphatidylcholine and lysophosphatidylcholine extracted from shark liver. *J. Agric. Food Chem.* **55**(23), 9670-9677 (2007).
2. Huang, L.S., Hung, N.D., Sok, D.-E., *et al.* Lysophosphatidylcholine containing docosahexaenoic acid at the *sn*-1 position is anti-inflammatory. *Lipids* **45**(3), 225-236 (2010).

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, 06/04/2024

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