

PRODUCT INFORMATION



1,2-Didocosahexaenoyl-*sn*-glycero-3-PC

Item No. 10733

CAS Registry No.: 99296-81-8
Formal Name: (7R,13Z,16Z,19Z,22Z,25Z,28Z)-4-hydroxy-N,N,N-trimethyl-10-oxo-7-[[[(4Z,7Z,10Z,13Z,16Z,19Z)-1-oxo-4,7,10,13,16,19-docosahexaenyl]oxy]-3,5,9-trioxa-4-phosphahentriaconta-13,16,19,22,25,28-hexaen-1-aminium, inner salt, 4-oxide

Synonyms: 1,2-Didocosahexaenoyl PC, 1,2-Didocosahexaenoyl-*sn*-glycero-3-Phosphocholine, 1,2-Didocosahexaenoin-*sn*-glycero-3-Phosphatidylcholine, PC(22:6/22:6), 22:6/22:6-PC

MF: C₅₂H₈₀NO₈P

FW: 878.2

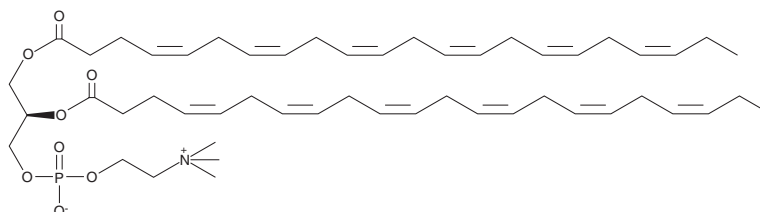
Purity: ≥95%

Supplied as: A solution in chloroform

Storage: -80°C

Stability: ≥2 years

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



Description

1,2-Didocosahexaenoyl-*sn*-glycero-3-PC is a phospholipid that contains docosahexaenoic acid (Item Nos. 90310 | 17950) at the *sn*-1 and *sn*-2 positions. It has been used in the formation of polyunsaturated endoplasmic reticulum-targeting liposomes (PERLs) that reduce cellular cholesterol levels in human peripheral blood mononuclear cells (PBMCs) and decrease the secretion of HIV-1 particles in PBMCs infected with HIV-1_{LAI}.¹ It has been used as a component of lipid monolayers and small unilamellar vesicles to study the effect of phospholipid composition on membrane fluidity.²

References

1. Pollock, S., Nichita, N.B., Böhmer, A., *et al.* Polyunsaturated liposomes are antiviral against hepatitis B and C viruses and HIV by decreasing cholesterol levels in infected cells. *Proc. Natl. Acad. Sci. USA.* **107**(40), 17176-17181 (2010).
2. Zerouga, M., Jenski, L.J., and Stillwell, W. Comparison of phosphatidylcholines containing one or two docosahexaenoic acyl chains on properties of phospholipid monolayers and bilayers. *Biochim Biophys Acta.* **1236**(2), 266-272 (1995).

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, 10/19/2023

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