

PRODUCT INFORMATION



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

Item No. 20954

CAS Registry No.: 998-06-1

Formal Name: (7R,18Z,21Z)-4-hydroxy-N,N,N-trimethyl-10-oxo-7-
[[[(9Z,12Z)-1-oxo-9,12-octadecadien-1-yl]oxy]-trioxa-4-
phosphaheptacos-18,21-dien-1-aminium, inner salt, 4-oxide

Synonyms: 1,2-Dilinoleoyl-*sn*-glycero-3-Phosphatidylcholine,
DLPC, 18:2 (*cis*) PC

MF: C₄₄H₈₀NO₈P

FW: 782.1

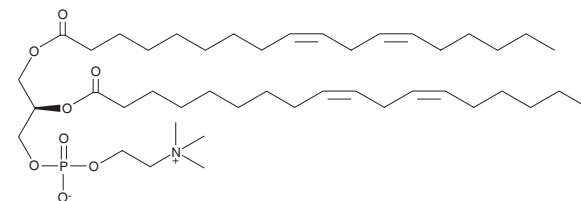
Purity: ≥95%

UV/Vis.: λ_{max}: 231 nm

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.

Laboratory Procedures

1,2-Dilinoleoyl-*sn*-glycero-3-PC is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice. 1,2-Dilinoleoyl-*sn*-glycero-3-PC is soluble in the organic solvent ethanol at a concentration of approximately 25 mg/ml.

Description

1,2-Dilinoleoyl-*sn*-glycero-3-PC is a phospholipid containing the polyunsaturated long-chain (18:2) linoleic acid inserted at the *sn*-1 and *sn*-2 positions. It can be used to generate micelles, liposomes, and other types of artificial membranes.¹⁻³

References

1. Ritter, M., Schmidt, S., Jakab, M., *et al.* Evidence for the formation of symmetric and asymmetric DLPC-DAPC lipid bilayer domains. *Cell Physiol. Biochem.* **32**(1), 46-52 (2013).
2. Mazari, A., Iwamoto, S., and Yamauchi, R. Effects of linoleic acid position in phosphatidylcholines and cholesterol addition on their rates of peroxidation in unilamellar liposomes. *Biosci. Biotechnol. Biochem.* **74**(5), 1013-1017 (2010).
3. Bagatolli, L.A. and Gratton, E.A. Correlation between lipid domain shape and binary phospholipid mixture composition in free standing bilayers: A two-photon fluorescence microscopy study. *Biophys. J.* **79**(1), 434-447 (2000).

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, 02/19/2024

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