

PRODUCT INFORMATION



1-Linoleoyl-d₁₁-2-Linoleoyl-*sn*-glycero-3-PC

Item No. 9004419

Formal Name: (7R,18Z,21Z)-4-hydroxy-N,N,N-trimethyl-10-oxo-7-[[[(9Z,12Z)-1-oxo-9,12-octadecadien-14,14,15,15,16,16,17,17,18,18,18-d₁₁-1-yl]oxy]-trioxa-4-phosphaheptacos-18,21-dien-1-aminium, inner salt, 4-oxide

Synonyms: 1,2-Dilinoleoyl-*sn*-glycero-3-Phosphocholine-d₁₁, DLPC-d₁₁, 18:2 (*cis*) PC-d₁₁

MF: C₄₄H₆₉D₁₁NO₈P

FW: 793.2

Chemical Purity: ≥95% (1,2-dilinoleoyl-*sn*-glycero-3-PC)

Deuterium

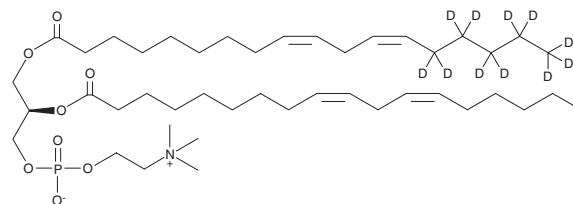
Incorporation: ≥99% deuterated forms (d₁-d₁₁); ≤1% d₀

UV/Vis.: λ_{max}: 231 nm

Supplied as: A 1 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.

Laboratory Procedures

1-Linoleoyl-d₁₁-2-Linoleoyl-*sn*-glycero-3-PC is intended for use as an internal standard for the quantification of 1,2-dilinoleoyl-*sn*-glycero-3-PC (Item No. 20954) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Description

1,2-Dilinoleoyl-*sn*-glycero-3-PC is a phospholipid containing the polyunsaturated long-chain (18:2) linoleic acid (Item Nos. 90150 | 90150.1 | 21909) inserted at the *sn*-1 and *sn*-2 positions. It has been 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, 01/13/2026

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