

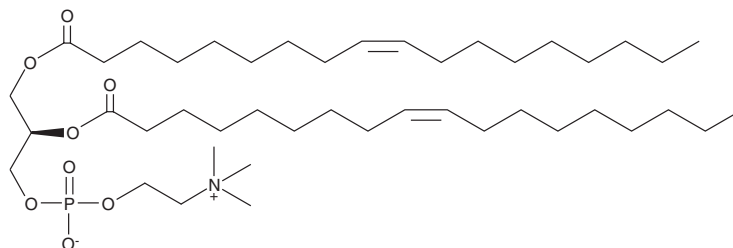
PRODUCT INFORMATION



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

Item No. 15098

CAS Registry No.: 4235-95-4
Formal Name: 1,2-dioleoyl-*sn*-glycero-3-phosphatidylcholine
Synonyms: *sn*-3-Dioleoyllecithin, 1,2-Dioleoyl-*sn*-glycero-3-Phosphocholine, DOPC, PDD 111
MF: C₄₄H₈₄NO₈P
FW: 786.1
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1,2-Dioleoyl-*sn*-glycero-3-PC (DOPC) is supplied as a crystalline solid. A stock solution may be made by dissolving the DOPC in the solvent of choice, which should be purged with an inert gas. DOPC is soluble in ethanol at a concentration of approximately 25 mg/ml.

Description

DOPC is a phospholipid containing the unsaturated long-chain (18:1) oleic acid inserted at the *sn*-1 and *sn*-2 positions. It is commonly used alone, or with other components, in the generation of micelles, liposomes, and other types of artificial membranes.¹⁻³

References

1. Konyakhina, T.M., Goh, S.L., Amazon, J., *et al.* Control of a nanoscopic-to-macroscopic transition: Modulated phases in four-component DSPC/DOPC/POPC/Chol giant unilamellar vesicles. *Biophys. J.* **101(2)**, L8-L10 (2011).
2. Rasch, M.R., Yu, Y., Bosoy, C., *et al.* Chloroform-enhanced incorporation of hydrophobic gold nanocrystals into dioleoylphosphatidylcholine (DOPC) vesicle membranes. *Langmuir* **28(36)**, 12971-12981 (2012).
3. Aguilar, L.F., Pino, J.A., Soto-Arriaza, M.A., *et al.* Differential dynamic and structural behavior of lipid-cholesterol domains in model membranes. *PLoS One* **7(6)**, e40254 (2012).

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, 01/10/2023

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