

PRODUCT INFORMATION



1-Palmitoyl-2-Oleoyl-*sn*-glycero-3-PC

Item No. 15102

CAS Registry No.: 26853-31-6

Formal Name: 1-palmitoyl-2-oleoyl-*sn*-glycero-3-phosphatidylcholine

Synonyms: 1,2-Hexadecanoyl-2-Octadecenoyl-*sn*-glycero-3-Phosphatidylcholine,
1,2-Hexadecanoyl-2-Octadecenoyl-*sn*-glycero-3-Phosphocholine,
PC(16:0/18:1), 16:0/18:1-PC, 1,2-POPC

MF: C₄₂H₈₂NO₈P

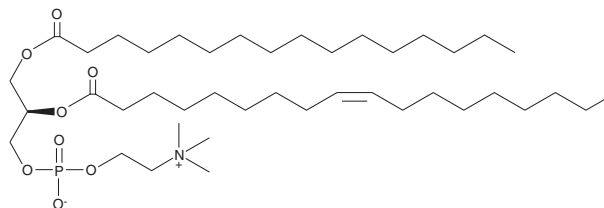
FW: 760.1

Purity: ≥95%

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

Description

1,2-POPC is a phospholipid containing 16:0 and 18:1 fatty acids at the *sn*-1 and *sn*-2 positions, respectively. It belongs to a class of phospholipids that are a major component of biological membranes.^{1,2} This compound can be used for liposome production in order to study the properties of lipid bilayers.

References

1. Moreno, M.J., Estronca, L.M.B.B., and Vaz, W.L.C. Translocation of phospholipids and dithionite permeability in liquid-ordered and liquid-disordered membranes. *Biophys. J.* **91**(3), 873-881 (2006).
2. Heberle, F.A. and Feigenson, G.W. Phase separation in lipid membranes. *Cold Spring Harb. Perspect. Biol.* **3**(4), 1-13 (2011).

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, 03/08/2024

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