

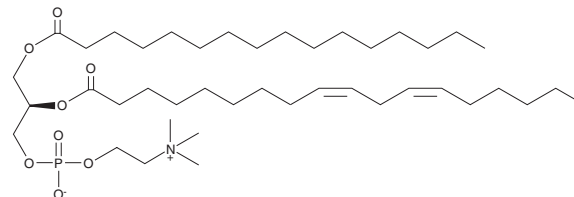
PRODUCT INFORMATION



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

Item No. 20962

CAS Registry No.: 17708-90-6
Formal Name: (7R,17Z,20Z)-4-hydroxy-N,N,N-trimethyl-9-oxo-7-[[[(1-oxohexadecyl)oxy]methyl]-3,5,8-trioxa-4-phosphahexacos-17,20-dien-1-aminium-4-oxide, inner salt
Synonyms: 1-Palmitoyl-2-Linoleoyl-*sn*-glycero-3-Phosphatidylcholine, 16:0-18:2 PC, PLPC
MF: C₄₂H₈₀NO₈P
FW: 758.1
Purity: ≥95%
Supplied as: A solution in chloroform
Storage: -80°C
Stability: ≥1 year



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

Laboratory Procedures

1-Palmitoyl-2-Linoleoyl-*sn*-glycero-3-PC (PLPC) 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. Solvents such as ethanol purged with an inert gas can be used. The solubility of PLPC in ethanol is approximately 50 mg/ml.

PLPC is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PLPC should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. PLPC has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PLPC is a phospholipid containing palmitic (16:0) and linoleic (18:2) acids inserted at the *sn*-1 and *sn*-2 positions, respectively. It can be used in the generation of micelles, liposomes, and other types of artificial membranes and has been particularly useful in the study of lipid peroxidation.¹⁻³

References

1. Conte, E., Megli, F.M., Khandelia, H., *et al.* Lipid peroxidation and water penetration in lipid bilayers: A W-band EPR study. *Biochimica et Biophysica Acta* **1828**, 510-517 (2013).
2. Milne, G.L., Seal, J.R., Havrilla, C.M., *et al.* Identification and analysis of products formed from phospholipids in the free radical oxidation of human low density lipoproteins. *J. Lipid Res.* **46**, 307-319 (2005).
3. Høyryup, P., Mouritsen, O.G., and Jørgensen, K. Phospholipase A₂ activity towards vesicles of DPPC and DMPC-DSPC containing small amounts of SMPC. *Biochim. Biophys. Acta* **1515(2)**, 133-143 (2001).

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, 11/02/2022

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