

PRODUCT INFORMATION



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

Item No. 15101

CAS Registry No.: 82765-47-7
Formal Name: (7R)-4-hydroxy-N,N,N-trimethyl-10-oxo-7-[(1-oxododecyl)oxy]-3,5,9-trioxa-4-phosphapentacosan-1-aminium, 4-oxide, inner salt

Synonyms: 1-Palmitoyl-2-Lauroyl-*sn*-glycero-3-Phosphatidylcholine, 1-Palmitoyl-2-Lauroyl-*sn*-glycero-3-Phosphocholine, 1,2-PLPC

MF: C₃₆H₇₂NO₈P

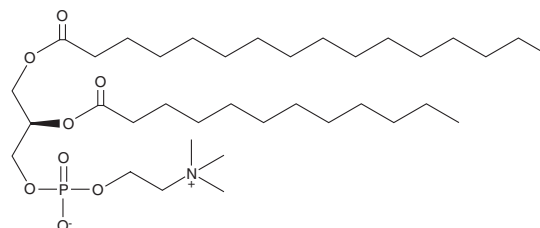
FW: 677.9

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

Description

1,2-PLPC is a phospholipid containing palmitoyl (16:0) and lauryl (12:0) acyl substituents at the *sn*-1 and *sn*-2 positions, respectively. This mixed-chain phosphatidylcholine has been used to study chain-chain contact interactions that contribute to the structural stability of lipid membrane bilayers.

Reference

1. Li, S., Wang, Z.-Q., Lin, H.-N., *et al.* Energy-minimized structures and packing states of a homologous series of mixed-chain phosphatidylcholines: A molecular mechanics study on the diglyceride moieties. *Biophys. J.* **65**(4), 1415-1428 (1993).

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, 11/01/2022

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