

PRODUCT INFORMATION



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

Item No. 15097

CAS Registry No.: 18194-24-6

Formal Name: (7R)-4-hydroxy-N,N,N-trimethyl-10-oxo-7-[(1-oxotetradecyl)oxy]-3,5,9-trioxo-4-phosphatricosan-1-aminium, inner salt, 4-oxide

Synonyms: 1,2-Dimyristoyl-*sn*-glycero-3-Phosphocholine, 1,2-DMPC

MF: $C_{36}H_{72}NO_8P$

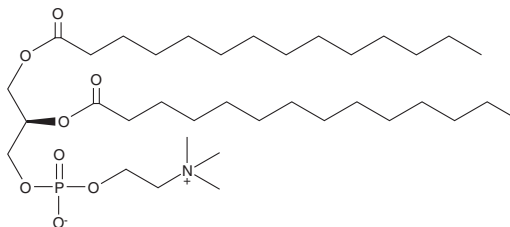
FW: 677.9

Purity: $\geq 98\%$

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

Description

1,2-DMPC is a phospholipid containing 14:0 fatty acids at the *sn*-1 and *sn*-2 positions. It is a synthetic phospholipid that has been used for liposome production in order to study the properties of lipid bilayer.¹⁻³

References

1. Uemura, A., Watarai, S., Iwasaki, T., *et al.* Induction of immune responses against glycosphingolipid antigens: Comparison of antibody responses in mice immunized with antigen associated with liposomes prepared from various phospholipids. *J. Vet. Med. Sci.* **67(12)**, 1197-1201 (2005).
2. Méléard, P., Gerbeaud, C., Pott, T., *et al.* Bending elasticities of model membranes: Influences of temperature and sterol content. *Biophys. J.* **72(6)**, 2616-2629 (1997).
3. Nagle, J.F. and Tristram-Nagle, S. Structure of lipid bilayers. *Biochim. Biophys. Acta.* **1469(3)**, 159-195 (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, 05/21/2024

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