

PRODUCT INFORMATION



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

Item No. 15096

CAS Registry No.: 61596-53-0
Formal Name: 1,2-diarachidoyl-*sn*-glycero-3-phosphatidylcholine

Synonyms: 1,2-DAPC, L- α -Diarachidonoyl lecithin, 1,2-Diarachidoyl-*sn*-glycero-3-Phosphocholine

MF: C₄₈H₉₆NO₈P

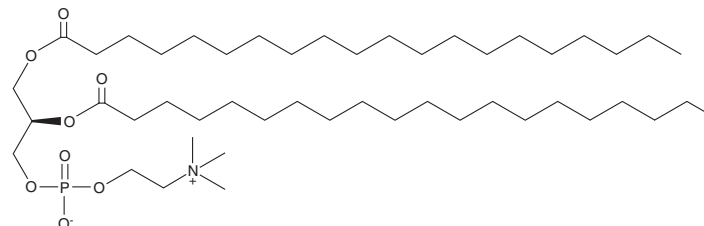
FW: 846.3

Purity: \geq 95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

1,2-Diarachidoyl-*sn*-glycero-3-PC (1,2-DAPC) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-DAPC in the solvent of choice. 1,2-DAPC is soluble in ethanol at a concentration of approximately 25 mg/ml.

Description

1,2-DAPC is a phospholipid containing the saturated long-chain (20:0) arachidic acid inserted at the *sn*-1 and *sn*-2 positions. It can be used in the generation of micelles, liposomes, and other types of artificial membranes.¹⁻³

References

1. Bagatolli, L.A. and Gratton, E. A correlation between lipid domain shape and binary phospholipid mixture composition in free standing bilayers: A two-photon fluorescence microscopy study. *Biophys. J.* **79**(1), 434-447 (2000).
2. Smith, E.A., van Gorkum, C.M., and Dea, P.K. Properties of phosphatidylcholine in the presence of its monofluorinated analogue. *Biophys. Chem.* **147**(1-2), 20-27 (2010).
3. Ritter, M., Schmidt, S., Jakab, M., et al. Evidence for the formation of symmetric and asymmetric DLPC-DAPC lipid bilayer domains. *Cell Physiol. Biochem.* **32**(1), 46-52 (2013).

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, 12/12/2022

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