

PRODUCT INFORMATION



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

Item No. 20953

CAS Registry No.: 17688-29-8

Formal Name: (R)-2,3-bis(((5Z,8Z,11Z,14Z)-icosa-5,8,11,14-tetraenoyl)oxy)propyl (2-(trimethylammonio)ethyl) phosphate
Synonyms: DAPC, 1,2-Diarachidonoyl-*sn*-glycero-3-Phosphatidylcholine, 20:4 (cis) PC

MF: $C_{48}H_{80}NO_8P$

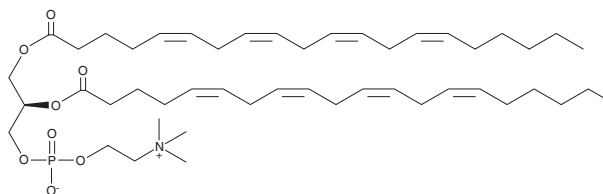
FW: 830.1

Purity: $\geq 90\%$

Supplied as: A 10 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥ 1 year



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

Description

1,2-Diarachidonoyl-*sn*-glycero-3-PC is a phospholipid containing the polyunsaturated long-chain (20:4) arachidonic 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.^{1,2}

References

1. Schwieger, C., Achilles, A., Scholz, S., *et al.* Binding of amphiphilic and triphilic block copolymers to lipid model membranes: The role of perfluorinated moieties. *Soft Matter* **10**(33), 6111-6362 (2014).
2. de Sa, M.M., Sresht, V., Rangel-Yagui, C.O., *et al.* Understanding miltefosine-membrane interactions using molecular dynamics simulations. *Langmuir* **31**(15), 4503-4512 (2015).

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, 11/12/2025

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