

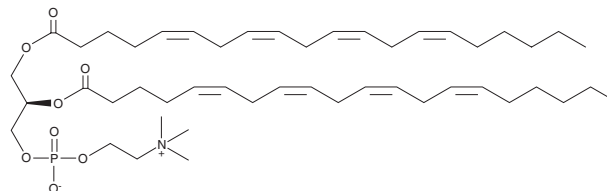
# PRODUCT INFORMATION



## 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-tetraenoyloxy)propyl (2-(trimethylammonio)ethyl) phosphate  
**Synonyms:** DAPC, 1,2-Diarachidonoyl-*sn*-glycero-3-Phosphatidylcholine, 20:4 (*cis*) PC  
**MF:** C<sub>48</sub>H<sub>80</sub>NO<sub>8</sub>P  
**FW:** 830.1  
**Purity:** ≥95%  
**Supplied as:** A neat oil  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

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

### 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.<sup>1,2</sup>

### References

- 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).
- 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, 03/17/2021

#### 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