

PRODUCT INFORMATION



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

Item No. 24630

CAS Registry No.: 70897-27-7

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

Synonyms: 1,2-DHDPC, 1,2-Diheptadecanoyl-PC,
1,2-Diheptadecanoyl-*sn*-glycero-3-Phosphocholine,
PC(17:0/17:0)

MF: C₄₂H₈₄NO₈P

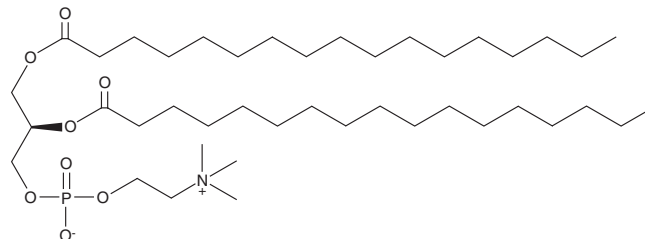
FW: 762.1

Purity: ≥98%

Supplied as: A 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-Diheptadecanoyl-*sn*-glycero-3-PC (1,2-DHDPC) is supplied as a solid. A stock solution may be made by dissolving the 1,2-DHDPC in the solvent of choice, which should be purged with an inert gas. 1,2-DHDPC is soluble in organic solvents such as methanol and dichloromethane.

Description

1,2-DHDPC is a synthetic phosphatidylcholine containing 17:0 fatty acids at the *sn*-1 and *sn*-2 positions. It has been used as an internal standard for the quantification of phosphatidylcholines by LC-MS.¹

Reference

1. Bird, S.S., Marur, V.R., Sniatynski, M.J., *et al.* Serum lipidomics profiling using LC-MS and high-energy collisional dissociation fragmentation: Focus on triglyceride detection and characterization. *Anal. Chem.* **83**(17), 6648-6657 (2011).

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, 12/19/2022

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