

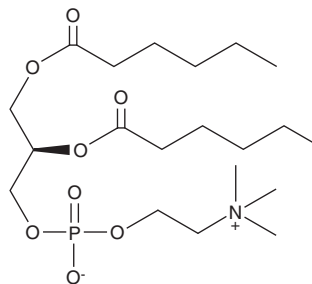
PRODUCT INFORMATION



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

Item No. 25585

CAS Registry No.: 34506-67-7
Formal Name: 4-hydroxy-N,N,N-trimethyl-10-oxo-7R-[(1-oxohexyl)oxy]-3,5,9-trioxa-4-phosphapentadecan-1-aminium, inner salt, 4-oxide
Synonyms: DHPC-C6, L-1,2-Dihexanoyllecithin, 1,2-Dihexanoyl-*sn*-glycero-3-Phosphocholine
MF: C₂₀H₄₀NO₈P
FW: 453.5
Purity: ≥95%
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-Dihexanoyl-*sn*-glycero-3-PC (DHPC-C6) is supplied as a crystalline solid. A stock solution may be made by dissolving the DHPC-C6 in the solvent of choice, which should be purged with an inert gas. DHPC-C6 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of DHPC-C6 in these solvents is approximately 30, 7, and 20 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of DHPC-C6 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of DHPC-C6 in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

DHPC-C6 is a synthetic phospholipid containing the short-chain (6:0) caproic acid inserted at the *sn*-1 and *sn*-2 positions. It is a substrate for phospholipase C isolated from *B. cereus* as well as phospholipase A₂ isolated from *A. halys blomhoffii* and *N. naja atra* snake venom.^{1,2} DHPC is commonly used in the generation of micelles, liposomes, and other types of artificial membranes.³

References

1. Little, C. Phospholipase C from *Bacillus cereus*. Action on some artificial lecithins. *Acta Chem. Scand. B* **31(4)**, 267-272 (1977).
2. Teshima, K., Kitagawa, Y., Samejima, Y., *et al.* Role of Ca²⁺ in the substrate binding and catalytic functions of snake venom phospholipases A₂. *J. Biochem.* **106(3)**, 518-527 (1989).
3. Batenjany, M.M., O'Leary, T.J., Levin, I.W., *et al.* Packing characteristics of two-component bilayers composed of ester- and ether-linked phospholipids. *Biophys. J.* **72(4)**, 1695-1700 (1997).

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, 11/08/2022

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