

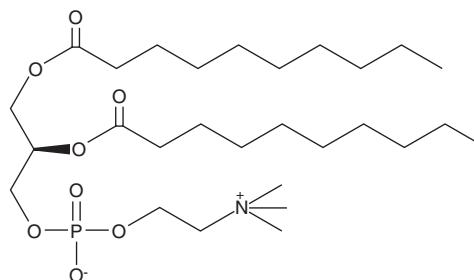
PRODUCT INFORMATION



1,2-Didecanoyl PC

Catalog No. 10009879

CAS Registry No.: 3436-44-0
Formal Name: 1,2-bis(O-decanoyl)-sn-glyceryl-3-phosphorylcholine
Synonyms: 1,2-Didecanoyl Phosphatidylcholine, Didecanoyl Lecithin, 1,2-DPC
MF: C₂₈H₅₆NO₈P
FW: 565.7
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

1,2-Didecanoyl PC (DPC) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of DPC in ethanol is approximately 30 mg/ml and approximately 20 mg/ml in DMSO and DMF.

DPC is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of DPC should be diluted with the aqueous buffer of choice. DPC has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Phosphatidylcholine species are a common class of phospholipids that comprise the mammalian cell membrane. DPC is a synthetic, less hydrophobic phospholipid that has been found to be useful for enhancing the absorption of peptide drugs and hormones such as insulin. Thus, the bioavailability of intranasally applied human growth hormone is enhanced when coadministered with the absorption enhancer DPC.¹

Reference

1. Agerholm, C., Bastholm, L., Johansen, P.B., *et al.* Epithelial transport and bioavailability of intranasally administered human growth hormone formulated with the absorption enhancers didecanoyl-L- α -phosphatidylcholine and α -cyclodextrin in rabbits. *J. Pharm. Sci.* **83(12)**, 1706-1711 (1994).

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, 03/25/2022

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