

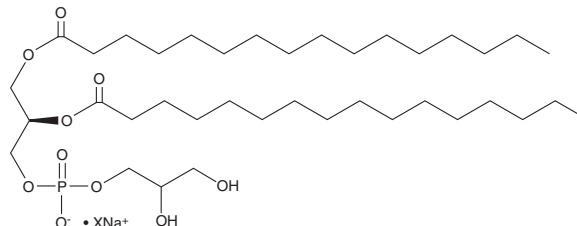
PRODUCT INFORMATION



1,2-Dipalmitoyl-*sn*-glycero-3-PG (sodium salt)

Item No. 15086

CAS Registry No.: 200880-41-7
Formal Name: 1,2-dipalmitoyl-*sn*-glycero-3-phospho-(1'-*rac*-glycerol), sodium salt
Synonyms: 1,2-Dihexadecanoyl-*sn*-glycero-3-phosphatidylglycerol, 1,2-Dipalmitoyl-*sn*-glycero-3-phospho-(1'-*rac*-glycerol), 1,2-Dihexadecanoyl-*sn*-glycero-3-phosphoglycerol, 1,2-DPPG, MG 6060LS, 16:0/16:0-PG, PG(16:0/16:0)
MF: C₃₈H₇₄O₁₀P • XNa
FW: 745.0
Purity: ≥98%
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-Dipalmitoyl-*sn*-glycero-3-PG (1,2-DPPG) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-DPPG (sodium salt) in the solvent of choice. 1,2-DPPG (sodium salt) is soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 2 mg/ml.

Description

1,2-DPPG is a phospholipid containing the long-chain (16:0) palmitic 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. Ege, C. and Lee, K.Y.C. Insertion of Alzheimer's Aβ40 peptide into lipid monolayers. *Biophys. J.* **87**, 1732-1740 (2004).
2. Mansour, H.M. and Zografi, G. The relationship between water vapor absorption and desorption by phospholipids and bilayer phase transitions. *J. Pharm. Sci.* **96**(2), 377-396 (2007).

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, 04/14/2025

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