

PRODUCT INFORMATION



1,2-Dipalmitoyl-*rac*-glycero-3-PG (ammonium salt)

Item No. 30574

CAS Registry No.: 73548-70-6

Formal Name: hexadecanoic acid, 1-[[[(2,3-dihydroxypropoxy)hydroxyphosphinyl]oxy]methyl]-2-[(1-oxohexadecyl)oxy]ethyl ester, monoammonium salt

Synonyms: 1,2-Dipalmitoyl-*rac*-glycero-3-Phosphoglycerol, 1,2-DPPG, 16:0/16:0-PG, PG(16:0/16:0)

MF: $C_{38}H_{74}O_{10}P \cdot NH_4$

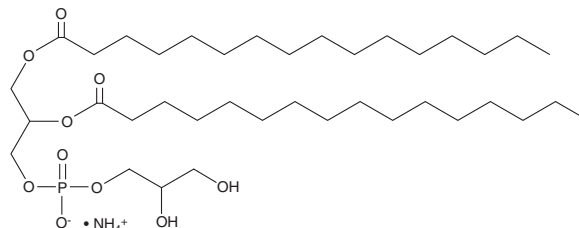
FW: 740.0

Purity: $\geq 95\%$

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-Dipalmitoyl-*rac*-glycero-3-PG (1,2-DPPG) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-DPPG in the solvent of choice, which should be purged with an inert gas. 1,2-DPPG is slightly soluble in ethanol and methanol.

Description

1,2-DPPG is a phosphoglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 and *sn*-2 positions. It has been used as a component of monolayers to study the membrane insertion ability of melittin (Item No. 17494) or apocytocrome c.¹

Reference

1. Han, X.H., Sui, S.F., and Yang, F.Y. A mini-trough for the study of membrane insertion ability of proteins. *Thin Solid Films* **284-285**, 789-792 (1996).

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, 11/03/2022

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