

PRODUCT INFORMATION



1-Palmitoyl-2-hydroxy-*sn*-glycero-3-PG (sodium salt)

Item No. 20961

CAS Registry No.: 326495-22-1

Formal Name: hexadecanoic acid, (2R)-3-[[[2,3-dihydroxypropoxy]hydroxyphosphinyl]oxy]-2-hydroxypropyl ester, monosodium salt

Synonyms: 16:0 Lyso-PG, 1-Palmitoyl-2-hydroxy-*sn*-glycero-3-phospho-(1'-*rac*-glycerol), 1-Palmitoyl-2-hydroxy-*sn*-glycero-3-phosphoglycerol

MF: $C_{22}H_{44}O_9P \cdot Na$

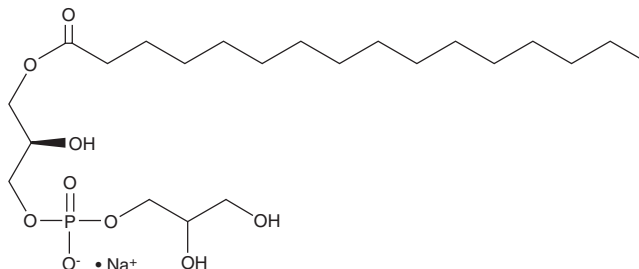
FW: 506.5

Purity: $\geq 95\%$

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

1-Palmitoyl-2-hydroxy-*sn*-glycero-3-PG (16:0 Lyso-PG) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 16:0 lyso-PG in the solvent of choice, which should be purged with an inert gas. 16:0 Lyso-PG is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 16:0 lyso-PG in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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 16:0 lyso-PG can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 16:0 lyso-PG in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

16:0 Lyso-PG is a lysophospholipid containing palmitic acid (16:0) (Item No. 10006627) at the *sn*-1 position. It can be used in the generation of micelles, liposomes, and other types of artificial membranes, including lipid-based drug carrier systems.¹

Reference

1. Jeschek, D., Lhota, G., Wallner, J., *et al.* A versatile, quantitative analytical method for pharmaceutical relevant lipids in drug delivery systems. *J. Pharm. Biomed. Anal.* **119**, 37-44 (2016).

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, 12/07/2022

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