

PRODUCT INFORMATION



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

Item No. 25587

CAS Registry No.: 326495-21-0

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

Synonyms: 14:0 Lyso-PG, 1-Myristoyl-2-hydroxy-*sn*-glycero-3-phosphoglycerol, 1-Myristoyl-2-hydroxy-*sn*-glycero-3-phospho-(1'-*rac*-glycerol)

MF: $C_{20}H_{40}O_9P \cdot Na$

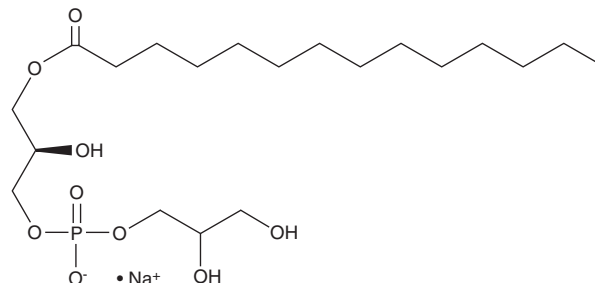
FW: 478.5

Purity: $\geq 95\%$ (9:1 mixture of 1-acyl and 2-acyl)

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-Myristoyl-2-hydroxy-*sn*-glycero-3-PG (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-myristoyl-2-hydroxy-*sn*-glycero-3-PG (sodium salt) in the solvent of choice. 1-Myristoyl-2-hydroxy-*sn*-glycero-3-PG (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 1-myristoyl-2-hydroxy-*sn*-glycero-3-PG (sodium salt) in these solvents is approximately 16, 30, and 20 mg/ml, respectively.

Description

1-Myristoyl-2-hydroxy-*sn*-glycero-3-PG is a lysophospholipid containing myristic acid (14:0) at the *sn*-1 position. It has been used in the generation of micelles, liposomes, and other types of artificial membranes, including lipid-based drug carrier systems.¹

Reference

1. Chadwick, A.C., Jensen, D.R., Peterson, F.C., *et al.* Expression, purification and reconstitution of the C-terminal transmembrane domain of scavenger receptor BI into detergent micelles for NMR analysis. *Protein Expr. Purif.* **107**, 35-42 (2015).

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, 07/10/2024

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