

PRODUCT INFORMATION



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

Item No. 20960

CAS Registry No.: 326495-24-3

Formal Name: 9-octadecenoic acid (9Z)-(2R)-3-[[[(2,3-dihydroxypropoxy)hydroxyphosphinyl]oxy]-2-hydroxypropyl ester, monosodium salt

Synonyms: 18:1 Lyso-PG, 1-Oleoyl-2-hydroxy-*sn*-glycero-3-phospho-(1'-*rac*-glycerol), 1-Oleoyl-2-hydroxy-*sn*-glycero-3-phosphoglycerol

MF: $C_{24}H_{46}O_9P \cdot Na$

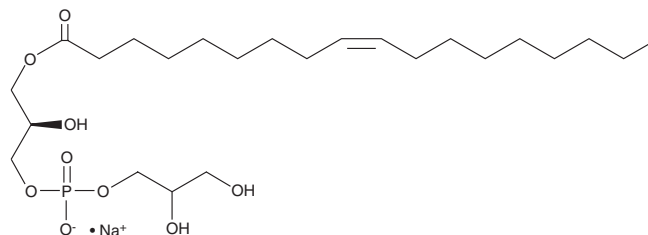
FW: 532.6

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

Supplied as: A 10 mg/ml solution in chloroform

Storage: $-80^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

1-Oleoyl-2-hydroxy-*sn*-glycero-3-PG (18:1 Lyso-PG) (sodium salt) is supplied as a solution in chloroform. To change the solvent, simply evaporate the 18:1 lyso-PG (sodium salt) under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 18:1 lyso-PG (sodium salt) in these solvents is approximately 16, 20, and 30 mg/ml.

Description

18:1 Lyso-PG is a lysophospholipid containing oleic acid (18:1) 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.^{1,2}

References

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).
2. Bulat, E. and Garrett, T.A. Putative N-acylphosphatidylethanolamine synthase from *A. thaliana* is a lysoglycerophospholipid acyltransferase. *J. Biol. Chem.* **286**(39), 33819-33831 (2011).

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, 01/07/2026

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