

PRODUCT INFORMATION



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

Item No. 15105

CAS Registry No.: 268550-95-4

Formal Name: 9Z-octadecenoic acid, (1R)-1-[[[(2,3-dihydroxypropoxy)hydroxyphosphinyl]oxy]methyl]-2-[(1-oxohexadecyl)oxy]ethyl ester, monosodium salt

Synonyms: 1-Palmitoyl-2-Oleoyl-*sn*-glycero-3-phosphoglycerol, 1-Palmitoyl-2-Oleoyl-*sn*-glycero-3-phospho-(1'-*rac*-glycerol), PG(16:0/18:1(9Z)), 1,2-POPG

MF: $C_{40}H_{76}O_{10}P \cdot Na$

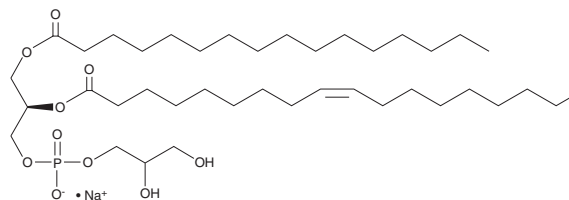
FW: 771.0

Purity: $\geq 98\%$

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-oleoyl-*sn*-glycero-3-PG (POPG) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the POPG (sodium salt) in the solvent of choice, which should be purged with an inert gas. POPG (sodium salt) is soluble in chloroform at a concentration of approximately 2 mg/ml.

Description

POPG is a phospholipid containing palmitoyl (16:0) and oleoyl (18:1) acyl substituents at the *sn*-1 and *sn*-2 positions, respectively. This mixed chain phosphatidylglycerol has been used to generate lipid membrane bilayers with controlled permeability and may be useful for various surfactant applications.¹⁻²

References

1. 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).
2. Dudia, A., Koëer, A., Subramaniam, V., et al. Biofunctionalized lipid-polymer hybrid nanocontainers with controlled permeability. *Nano Lett.* **8**(4), 1105-1110 (2008).

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, 11/01/2022

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