

PRODUCT INFORMATION



1,2-Dioleoyl-*sn*-glycero-3-PG (sodium salt)

Item No. 20957

CAS Registry No.: 67254-28-8

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

Synonyms: 1,2-Dioleoyl-*sn*-glycero-3-phospho-
(1'-*rac*-glycerol), 1,2-Dioleoyl-*sn*-
glycero-3-phosphoglycerol,
1,2-DOPG, 18:1 (Δ^9 -*cis*) PG

MF: C₄₂H₇₈O₁₀P • Na

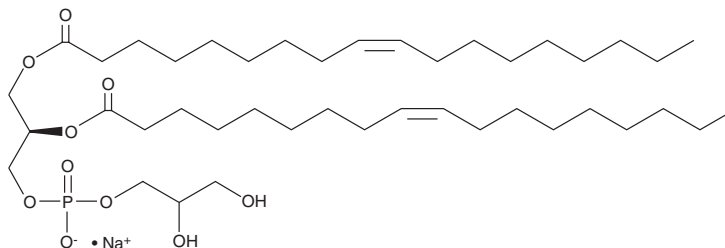
FW: 797.0

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -80°C

Stability: ≥2 years



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

Laboratory Procedures

1,2-Dioleoyl-*sn*-glycero-3-PG (DOPG) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the DOPG (sodium salt) in the solvent of choice, which should be purged with an inert gas. DOPG (sodium salt) is soluble in the organic solvent chloroform at a concentration of approximately 2 mg/ml.

Description

DOPG is a phospholipid containing the long-chain (18:1) fatty acid oleic acid (Item No. 90260) inserted at the *sn*-1 and *sn*-2 positions. It can be used in the generation of micelles, liposomes, and other artificial membranes.^{1,2}

References

1. Takahashi, T., Nomura, F., Yokoyama, Y., *et al.* Multiple membrane interactions and versatile vesicle deformations elicited by melittin. *Toxins (Basel)* **5**(4), 637-664 (2013).
2. Vequi-Suplicy, C.C., Riske, K.A., Knorr, R.L., *et al.* Vesicles with charged domains. *Biochim. Biophys. Acta* **1798**(7), 1338-1347 (2010).

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, 11/21/2022

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