

PRODUCT INFORMATION



1-Palmitoyl-2-Oleoyl-*sn*-glycero-3-PA

Item No. 20963

CAS Registry No.: 62600-81-1

Formal Name: (1R)-1-[[[(1-oxohexadecyl)oxy]methyl]-2-(phosphonoxy)-9Z-octadecenoic acid ethyl ester

Synonyms: 16:0-18:1 PA, 16:0/18:1-PA, PA(16:0/18:1), 1-Hexadecanoyl-2-Octadecenoyl-*sn*-glycero-3-PA, 1-Hexadecanoyl-2-Octadecenoyl-*sn*-glycero-3-phosphate, 1-Hexadecanoyl-2-Octadecanoyl-*sn*-glycero-3-phosphatidic Acid, 1-Palmitoyl-2-Oleoyl-*sn*-glycero-3-phosphate, 1-Palmitoyl-2-Oleoyl-*sn*-glycero-3-Phosphatidic Acid, 1,2-POPA

MF: C₃₇H₇₁O₈P

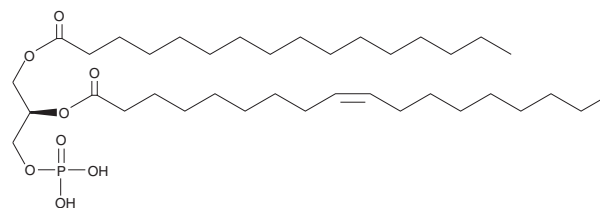
FW: 674.9

Purity: ≥95%

Supplied as: A 5 mg/ml solution in chloroform

Storage: -20°C

Stability: ≥4 years



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

Description

1,2-POPA is a phospholipid containing saturated palmitic acid (16:0) and monounsaturated oleic acid (18:1) inserted at the *sn*-1 and *sn*-2 positions, respectively. It can be used in the generation of micelles, liposomes, and other types of artificial membranes.¹

Reference

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).

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