

PRODUCT INFORMATION



1-Palmitoyl-2-oleoyl-*sn*-glycero-3-phosphoserine

Item No. 46036

CAS Registry No.: 40290-44-6

Formal Name: 2S-amino-5-hydroxy-10-oxo-8R-[[[(1-oxohexadecyl)oxy]methyl]-4,6,9-trioxa-5-phosphaheptacos-18Z-enoic acid, 5-oxide

Synonyms: 1-Hexadecanoyl-2-(9Z-octadecenoyl)-*sn*-glycero-3-phosphoserine, 1-Palmitoyl-2-oleoyl Phosphatidylserine, 1,2-POPS, PS(16:0/18:1), PS(16:0/18:1(9Z)), PtdSer-(1-palmitoyl, 2-oleoyl)

MF: $C_{40}H_{76}NO_{10}P$

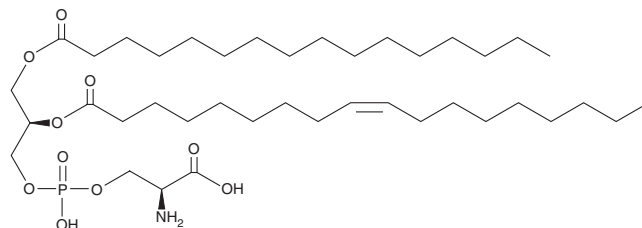
FW: 762.0

Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°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-phosphoserine is supplied as a solid. A stock solution may be made by dissolving the 1-palmitoyl-2-oleoyl-*sn*-glycero-3-phosphoserine in the solvent of choice, which should be purged with an inert gas. 1-Palmitoyl-2-oleoyl-*sn*-glycero-3-phosphoserine is sparingly soluble (1-10 mg/ml) in methanol.

Description

1-Palmitoyl-2-oleoyl-*sn*-glycero-3-phosphoserine is an anionic phospholipid that contains palmitic acid (Item Nos. 10006627 | 43561) and oleic acid (Item Nos. 90260 | 24659) at the *sn*-1 and *sn*-2 positions, respectively. It has been used in the synthesis of lipid membranes and synthetic high-density lipoproteins (HDLs).^{1,2}

References

1. Silin, V.I. and Hoogerheide, D.P. pH dependent electrical properties of the inner- and outer- leaflets of biomimetic cell membranes. *J. Colloid Interface Sci.* **594**, 279-289 (2021).
2. Yu, M., Dorsey, K.H., Halseth, T., *et al.* Enhancement of anti-inflammatory effects of synthetic high-density lipoproteins by incorporation of anionic lipids. *Mol. Pharm.* **20(11)**, 5454-5462 (2023).

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, 06/17/2026

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