

PRODUCT INFORMATION



1,2-Distearoyl-*sn*-glycero-3-phosphoserine (sodium salt)

Item No. 38528

CAS Registry No.: 321595-13-5
Formal Name: (2S,8R)-4,6,10-trioxa-5-phosphaoctacosanoic acid, 2-amino-5-hydroxy-11-oxo-8-[(1-oxooctadecyl)oxy]-5-oxide, sodium salt
Synonyms: 1,2-Dioctadecanoyl-*sn*-glycero-3-phosphoserine, 1,2-Distearoyl Phosphatidylserine, 1,2-DSPS, PS(C18:0/C18:0), PtdSer-(1,2-distearoyl)

MF: C₄₂H₈₁NO₁₀P • xNa

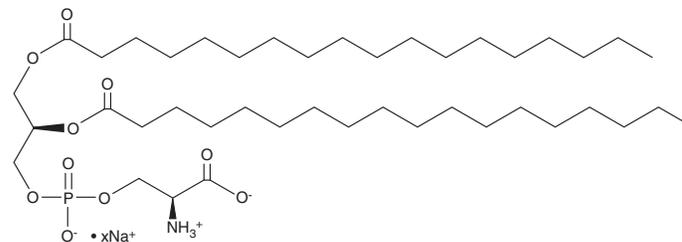
FW: 791.1

Purity: ≥90%

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

1,2-Distearoyl-*sn*-glycero-3-phosphoserine (1,2-DSPS) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 1,2-DSPS (sodium salt) in the solvent of choice, which should be purged with an inert gas. 1,2-DSPS (sodium salt) is slightly soluble (0.1-1 mg/ml) in chloroform:methanol.

Description

1,2-DSPS is a phospholipid containing stearic acid (Item No. 10011298) at the *sn*-1 and *sn*-2 positions. It has been used in the generation of liposomal nanoparticles for the delivery of siRNA *in vitro*.¹ Liposomal nanoparticles containing 1,2-DSPS and encapsulating siRNA targeting c-Rel decrease c-Rel levels in primary mouse macrophages.

Reference

1. Bressy, C., Zemani, A., Goyal, S., et al. Inhibition of c-Rel expression in myeloid and lymphoid cells with distearoyl -phosphatidylserine (DSPS) liposomal nanoparticles encapsulating therapeutic siRNA. *PLoS One* **17(12)**, e0276905 (2022).

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, 12/12/2025

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