

PRODUCT INFORMATION



1,2-Dimyristoyl-*sn*-glycero-3-PS (sodium salt)

Item No. 20955

CAS Registry No.: 105405-50-3

Formal Name: (2S,8R)-5-oxide, 2-amino-5-hydroxy-11-oxo-8-[(1-oxotetradecyl)oxy]-4,6,10-trioxa-5-phosphatetrascanoic acid, monosodium salt

Synonyms: 1,2-Dimyristoyl-*sn*-glycero-3-phospho-L-serine, 1,2-Ditetradecanoyl-*sn*-glycero-phosphatidylserine, 1,2-Ditetradecanoyl-*sn*-glycero-phosphoserine, 1,2-DMPS, 14:0 PS, 14:0/14:0-PS, PS(14:0/14:0)

MF: $C_{34}H_{65}NO_{10}P \cdot Na$

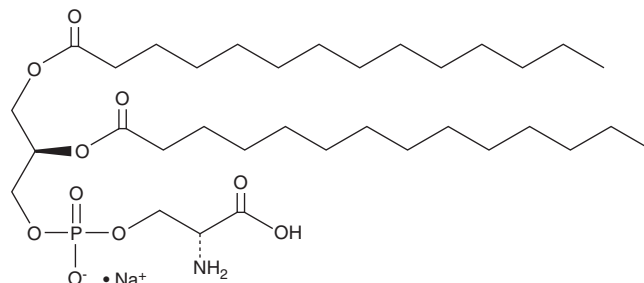
FW: 701.9

Purity: $\geq 95\%$

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

1,2-Dimyristoyl-*sn*-glycero-3-PS (1,2-DMPS) (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-DMPS (sodium salt) in the solvent of choice. 1,2-DMPS (sodium salt) is sparingly soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 0.5 mg/ml.

Description

1,2-DMPS is an anionic phospholipid containing myristic acid (14:0) inserted at the *sn*-1 and *sn*-2 positions. It can be used in the generation of micelles, liposomes, and other types of artificial membranes.^{1,2}

References

1. Mattai, J., Hauser, H., Demel, R.A., *et al.* Interactions of metal ions with phosphatidylserine bilayer membranes: Effect of hydrocarbon chain unsaturation. *Biochemistry* **28**(5), 2322-2330 (1989).
2. Tapia, M.J., Monteserin, M., Burrows, H.D., *et al.* Effect of the phospholipid chain length and head group on beta-phase formation of poly(9,9-dioctylfluorene) enclosed in liposomes. *Photochem. Photobiol.* **89**(6), 1471-1478 (2013).

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, 10/05/2022

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