

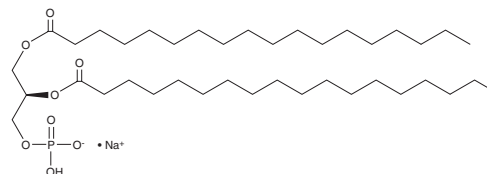
PRODUCT INFORMATION



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

Item No. 15083

CAS Registry No.: 108321-18-2
Formal Name: 1,2-distearoyl-*sn*-glycero-3-phosphate, monosodium salt
Synonyms: 1,2-Distearoyl-*sn*-glycero-3-phosphate, 1,2-Distearoyl-*sn*-glycero-3-phosphatidic Acid, 1,2-DSPA, 18:0/18:0-PA, PA (18:0/18:0)
MF: C₃₉H₇₆O₈P • Na
FW: 727.0
Purity: ≥98%
Supplied as: A crystalline 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-PA (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-distearoyl-*sn*-glycero-3-PA (sodium salt) in the solvent of choice, which should be purged with an inert gas. 1,2-Distearoyl-*sn*-glycero-3-PA (sodium salt) is soluble in the organic solvent chloroform at a concentration of approximately 1.6 mg/ml.

Description

1,2-Distearoyl-*sn*-glycero-3-PA (sodium salt) is a phospholipid containing a phosphatidic acid head group and 18:0 fatty acids at the *sn*-1 and *sn*-2 positions of the glycerol backbone. This negatively charged phospholipid has been used to study the dynamics of model lipid bilayers.^{1,2}

References

1. Li, S., Stein, A.J., Kruger, A., *et al.* Head groups of lipids govern the interaction and orientation between graphene oxide and lipids. *J. Phys. Chem. C* **117**(31), 16150-16158 (2013).
2. Bhagat, M. and Sofou, S. Membrane heterogeneities and fusogenicity in phosphatidylcholine-phosphatidic acid rigid vesicles as a function of pH and lipid chain mismatch. *Langmuir* **26**(3), 1666-1673 (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, 02/03/2025

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