

PRODUCT INFORMATION

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

Item No. 39190

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

Synonyms: 1,2-Dicaproyl-*sn*-glycero-3-phosphatidylserine, 1,2-Dicaproyl-*sn*-glycero-3-phosphoserine, 1,2-Dihexanoyl-*sn*-glycero-3-phosphatidylserine, 1,2-Dihexanoyl-*sn*-glycero-3-phospho-L-serine, 1,2-Dihexanoyl-*sn*-glycero-3-phosphoserine, PS(6:0/6:0), 6:0/6:0-PS

MF: C₁₈H₃₃NO₁₀P • XNa

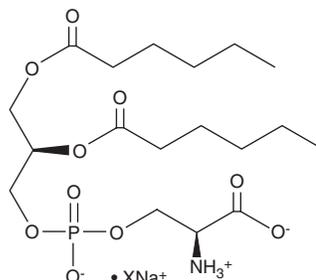
FW: 454.4

Purity: ≥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

1,2-Dihexanoyl-*sn*-glycero-3-PS (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 1,2-dihexanoyl-*sn*-glycero-3-PS (sodium salt) in the solvent of choice, which should be purged with an inert gas. 1,2-Dihexanoyl-*sn*-glycero-3-PS (sodium salt) is soluble in methanol.

Description

1,2-Dihexanoyl-*sn*-glycero-3-PS is a phospholipid containing hexanoic acid at the *sn*-1 and *sn*-2 positions.

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, 03/18/2025

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