

# PRODUCT INFORMATION



## DSPE-mPEG(5000) (sodium salt)

Item No. 34652

CAS Registry No.: 459428-35-4

Formal Name:  $\alpha$ -[6-hydroxy-6-oxido-1,12-dioxo-9-[(1-oxooctadecyl)oxy]-5,7,11-trioxa-2-aza-6-phosphanonacos-1-yl]- $\omega$ -methoxy-poly(oxy-1,2-ethanediyl), monosodium salt

Synonyms: 1,2-Dioctadecanoyl-*rac*-glycero-3-Phosphoethanolamine-N-mPEG(5000), 1,2-Distearoyl Phosphatidylethanolamine-N-mPEG(5000), 1,2-Distearoyl-*rac*-glycero-3-Phosphoethanolamine-N-methoxy-Polyethylene Glycol-5000, 1,2-DSPE-mPEG(5000), mPEG-5000-DSPE, PE(18:0/18:0)-mPEG(5000), PtdEth-(1,2-Distearoyl)-N-mPEG(5000)

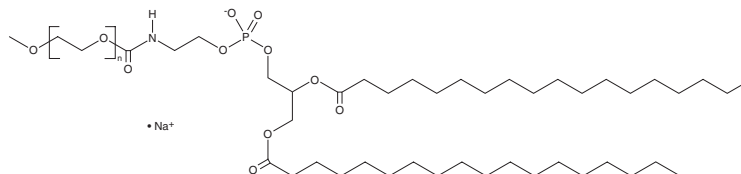
MF:  $(C_2H_4O)_n C_{43}H_{83}NO_{10}P \cdot Na$

Purity:  $\geq 85\%$

Supplied as: A solid

Storage:  $-20^\circ\text{C}$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

DSPE-mPEG(5000) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the DSPE-mPEG(5000) (sodium salt) in the solvent of choice, which should be purged with an inert gas. DSPE-mPEG(5000) (sodium salt) is soluble in the organic solvent ethanol at a concentration of approximately 3 mg/ml.

### Description

DSPE-mPEG(5000) is a PEGylated form of 1,2-distearoyl-*rac*-glycero-3-PE (1,2-DSPE).

#### 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, 04/10/2025

#### 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