

PRODUCT INFORMATION



DSPE-mPEG(2000) (sodium salt)

Item No. 37148

CAS Registry No.: 247925-28-6

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

Synonyms: 1,2-Dioctadecanoyl-*sn*-glycero-3-Phosphoethanolamine-N-mPEG(2000), 1,2-Distearoyl Phosphatidylethanolamine-N-mPEG(2000), 1,2-Distearoyl-*sn*-glycero-3-Phosphoethanolamine-N-methoxy-Polyethylene Glycol-2000, 1,2-DSPE-mPEG(2000), PE(18:0/18:0)-mPEG(2000), PtdEth-(1,2-Distearoyl)-N-mPEG(2000)

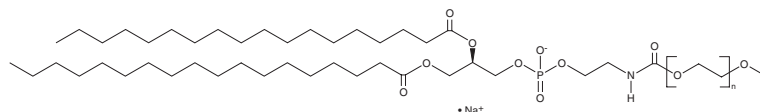
MF: $(C_{2}H_{4}O)_n C_{43}H_{83}NO_{10}P \cdot Na$

Purity: $\geq 95\%$

Supplied as: A solid

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

Stability: ≥ 2 years



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

Laboratory Procedures

DSPE-mPEG(2000) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the DSPE-mPEG(2000) (sodium salt) in the solvent of choice, which should be purged with an inert gas. DSPE-mPEG(2000) (sodium salt) is slightly soluble in dimethyl formamide and water.

Description

DSPE-mPEG(2000) is a PEGylated form of 1,2-distearoyl-*sn*-glycero-3-PE (DSPE; Item No. 15095).¹ It has been used in the generation of micelles and in combination with other lipids in the formation of liposomes. Liposomes containing DSPE-mPEG(2000) and encapsulating the DNA topoisomerase inhibitor doxorubicin (Item No. 15007) and the flavonoid quercetin (Item No. 10005169) reduce tumor growth in an MCF-7/adr multidrug-resistant breast cancer mouse xenograft model.

Reference

1. Zhang, J., Luo, Y., Zhao, X., et al. Co-delivery of doxorubicin and the traditional Chinese medicine quercetin using biotin-PEG₂₀₀₀-DSPE modified liposomes for the treatment of multidrug resistant breast cancer. *RSC Adv.* **6**(114), 113173–113184 (2016).

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, 04/11/2025

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