

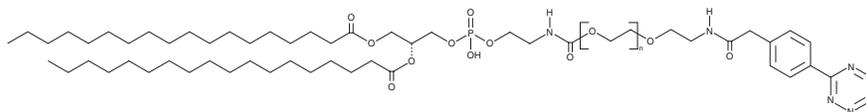
PRODUCT INFORMATION



DSPE-PEG(2000)-tetrazine

Item No. 45240

CAS Registry No.: 3101670-39-4
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 -[2-[[2-[4-(1,2,4,5-tetrazin-3-yl)phenyl]acetyl]amino]ethoxy]-poly(oxy-1,2-ethanediyl)]



Synonym: DSPE-PEG(2000)-Tz
MF: $(C_2H_4O)_n C_{54}H_{93}N_6O_{11}P$
Purity: $\geq 95\%$
Supplied as: A 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

DSPE-PEG(2000)-tetrazine is supplied as a solid. A stock solution may be made by dissolving the DSPE-PEG(2000)-tetrazine in the solvent of choice, which should be purged with an inert gas. DSPE-PEG(2000)-tetrazine is sparingly soluble (1-10 mg/ml) in ethanol.

Description

DSPE-PEG(2000)-tetrazine is a PEGylated and clickable form of 1,2-distearoyl-*sn*-glycero-3-PE (1,2-DSPE; Item No. 15095) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of small molecules *in vitro* and *in vivo*.¹ LNPs containing DSPE-PEG(2000)-tetrazine and encapsulating SN-38 (Item No. 15632) decrease the viability of HER2-overexpressing NIH3T6.7 cells pre-incubated with *trans*-cyclooctene-conjugated trastuzumab.

Reference

1. Yoo, J., Choi, S., Son, J., *et al.* Click chemistry-mediated tumor-targeting of SN38-loaded nanoparticles using trastuzumab. *Biochem. Biophys. Res. Commun.* **515**(1), 207-213 (2019).

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, 02/06/2026

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