

PRODUCT INFORMATION



DSPE-PEG(5000)-tetrazine

Item No. 46431

Formal Name: (2R)-3-((((14-(4-(1,2,4,5-tetrazin-3-yl)phenyl)-4,13-dioxo-6,9-dioxo-3,12-diazatetradecyl)oxy)(hydroxy)phosphoryl)oxy)propane-1,2-diyl distearate

Synonyms: DSPE-PEG_{5k}-Tre, DSPE-PEG(5000)-Tz

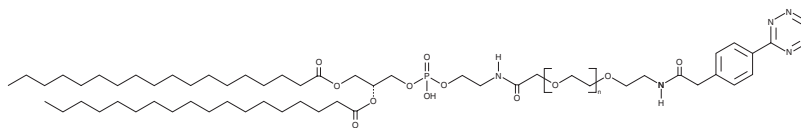
MF: (C₂H₄O)_nC₅₅H₉₅N₆O₁₁P

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

DSPE-PEG(5000)-tetrazine is supplied as a solid. A stock solution may be made by dissolving the DSPE-PEG(5000)-tetrazine in the solvent of choice, which should be purged with an inert gas. DSPE-PEG(5000)-tetrazine is soluble (≥10 mg/ml) in DMSO.

Description

DSPE-PEG(5000)-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 as a reactive anchor lipid by direct incorporation into the plasma membranes of CD8⁺ T cells.¹ DSPE-PEG(5000)-tetrazine, via membrane incorporation, has been used to conjugate T cells to liposomes containing DSPE-bicyclo[6.1.0]nonyne (DSPE-BCN) and encapsulating avasimibe (Item No. 18129) for tumor immunotherapy in mice.

Reference

1. Hao, M., Hou, S., Li, W., *et al.* Combination of metabolic intervention and T cell therapy enhances solid tumor immunotherapy. *Sci. Transl. Med.* **12**(571), eaaz6667 (2020).

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, 08/07/2026

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