

PRODUCT INFORMATION



DSPE-PEG(2000)-T7

Item No. 44433

Contents: This vial contains 200 µg or 1 mg of DSPE-PEG(2000)-T7 click conjugate as a lyophilized film.

Synonyms: 1,2-Distearoyl Phosphatidylethanolamine-PEG(2000)-T7, TfR Peptide Ligand, Transferrin Receptor Peptide Ligand

Peptide Sequence: FXGHAIYPRH (X = aminohexanoic acid)

MF: (C₂H₄O)_nC₁₂₁H₁₈₄N₂₃O₂₄P

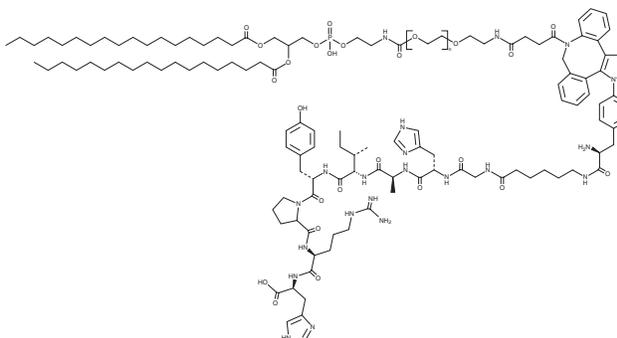
Molecular Weight: 3,839-5,222; 4,350.8 (Average due to PEG congeners)

Form: Solid

Storage: -20°C (as supplied)

Stability: ≥2 years

Applications: Suitable for LNP formulations, working concentration/dilution should be determined empirically.



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

Laboratory Procedures

DSPE-PEG(2000)-T7 is supplied as a solid. A stock solution may be made by dissolving the DSPE-PEG(2000)-T7 in the solvent of choice, which should be purged with an inert gas. DSPE-PEG(2000)-T7 is soluble in ethanol at a concentration of approximately 5 mg/ml.

Description

DSPE-PEG(2000)-T7 is a peptide-lipid conjugate that has been used for targeted delivery of liposomes *in vivo*.¹ It is composed of DSPE-PEG(2000)-DBCO conjugated to an azide-peptide containing the transferrin receptor-targeting peptide T7. Liposomes containing DSPE-PEG(2000)-T7 accumulate in tumors in an orthotopic C6 glioma mouse xenograft model. Liposomes containing DSPE-PEG(2000)-T7 and encapsulating doxorubicin increase survival in an orthotopic C6 glioma mouse xenograft model.

Reference

1. Zong, T., Mei, L., Gao, H., *et al.* Synergistic dual-ligand doxorubicin liposomes improve targeting and therapeutic efficacy of brain glioma in animals. *Mol. Pharm.* **11**(7), 2346-2357 (2014).

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, 01/27/2026

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