

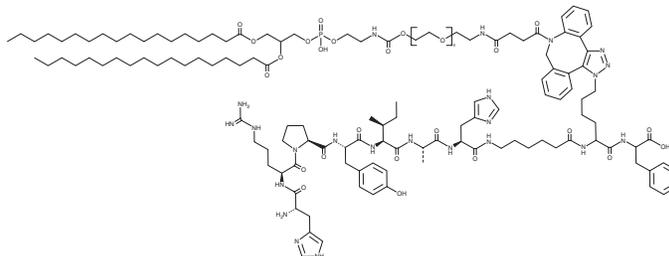
PRODUCT INFORMATION



DSPE-PEG(2000)-DT7

Item No. 42584

- Contents:** This vial contains 200 µg or 1 mg of DSPE-PEG(2000)-DT7 conjugate
- Synonyms:** 1,2-Distearoyl Phosphatidylethanolamine-PEG(2000)-DT7 Peptide,
1,2-DSPE-PEG(2000)-DT7,
PE(18:0/18:0)-PEG(2000)-DT7,
Transferrin Receptor-Targeting Peptide Conjugate
- Molecular Weight:** 4,014-5,070; 4,435 (Average due to PEG congeners)
- Form:** Solid
- Storage:** -20°C (as supplied)
- Stability:** ≥1 year
- Applications:** Suitable for LNP formulations, working concentration/dilution should be determined empirically.



Description

DSPE-PEG(2000)-DT7 is a peptide-lipid conjugate. It was prepared from DSPE-PEG(2000)-DBCO and a clickable form of DT7 peptide, a retro-inverso derivative of T7 peptide, a transferrin receptor (TfR) agonist. DSPE-PEG(2000)-DT7 is intended for use in the targeted delivery of lipid nanoparticles (LNPs) *in vitro* and *in vivo*.

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, 01/26/2026

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