

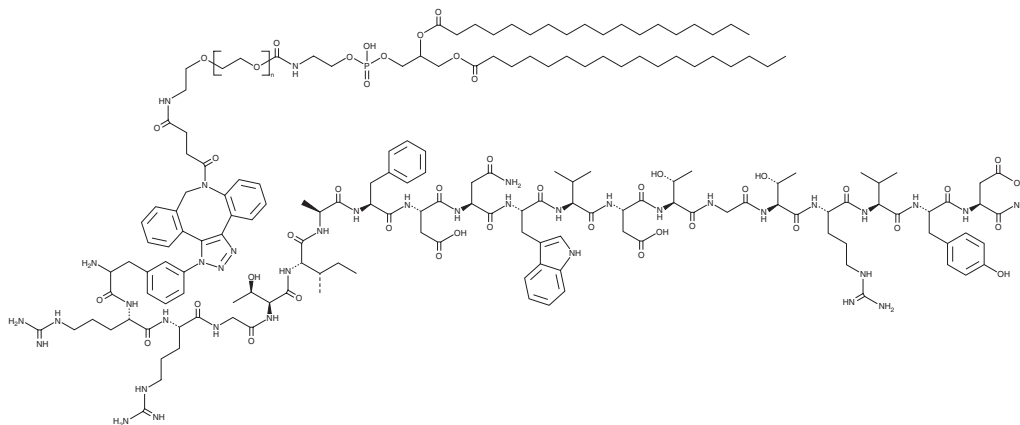
PRODUCT INFORMATION

DSPE-PEG(2000)-Pep2

Item No. 43421

Overview and Properties

Contents:	This vial contains 200 µg of DSPE-PEG(2000)-Pep2 click conjugate in 100 µl DMSO.
Synonyms:	1,2-Distearoyl Phosphatidylethanolamine-PEG(2000)-Pep2 Peptide, DSPE-PEG(2000)-RRGTIAFDNWVDTGTRVYDF, Scavenger Receptor B2/CD36 Targeting Peptide
Peptide Sequence:	RRGTIAFDNWVDTGTRVYDF
MF:	$(C_{24}H_{48}O)_n C_{170}H_{257}N_{38}O_{43}P$
Molecular Weight:	5,765 (Average)
Form:	Liquid
Storage:	-20°C
Stability:	≥1 year
Storage Buffer:	DMSO
Applications:	LNP formulations



Description

DSPE-PEG(2000)-Pep2 is a peptide-lipid conjugate that has been used for the targeted delivery of lipid nanoparticles (LNPs) *in vivo*.¹ It is composed of DSPE-PEG(2000)-DCBO conjugated to Pep2 peptide (Item No. 40258), a peptide ligand for the transmembrane glycoprotein CD36, also known as scavenger receptor B2, that selectively binds to CD36 over human serum albumin (HSA), IgG, and CD44.² LNPs containing DSPE-PEG(2000)-Pep2 and encapsulating the indoleamine 2,3-dioxygenase (IDO) inhibitor NLG-919 reduce the number of lung metastasis in a 4T1 murine mammary carcinoma model.¹

References

- Hu, C., Song, Y., Zhang, Y., *et al.* Sequential delivery of PD-1/PD-L1 blockade peptide and IDO inhibitor for immunosuppressive microenvironment remodeling via an MMP-2 responsive dual-targeting liposome. *Acta Pharm. Sin. B.* **13**(5), 2176-2187 (2023).
- Zheng, Y., Fang, X., Yang, Y., *et al.* Peptide-directed delivery of drug-loaded nanocarriers targeting CD36 overexpressing cells. *Colloid. Surface. A* **610**, 12590 (2021).

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