

PRODUCT INFORMATION

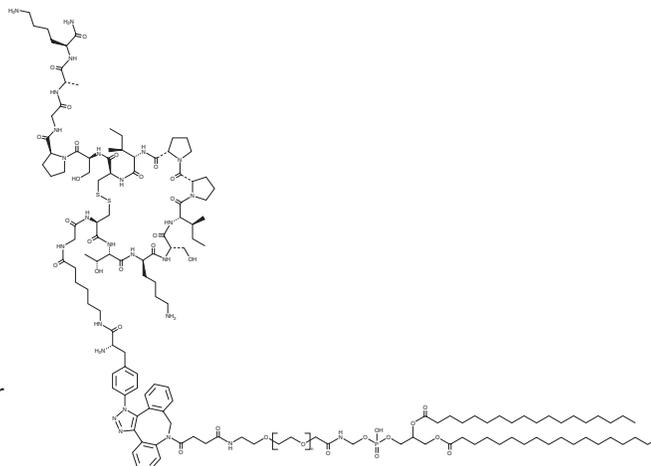


DSPE-PEG(2000)-MACTIDE

Item No. 43956

Overview and Properties

Contents:	This vial contains 200 µg or 1 mg of DSPE-PEG(2000)-MACTIDE click conjugate.
Synonyms:	1,2-Distearoyl Phosphatidylethanolamine-PEG(2000)-MACTIDE, 1,2-DSPE-PEG(2000)-CD206 Cyclic Peptide Ligand, DSPE-PEG(2000)-triazole-F-Ahx-G(CTKSIPPIC)SPGAK, DSPE-PEG(2000)-UNO-related Cyclic Peptide
Peptide Sequence:	Phe-Ahx-G[CTKSIPPIC]-SPGAK
MF:	$(C_2H_4O)_n C_{140}H_{225}N_{26}O_{32}PS_2$
Molecular Weight:	4,086-5,530 (due to PEG congeners)
Form:	Solid
Storage:	-20°C
Stability:	≥1 year
Storage Buffer:	Reconstitute lyophilized conjugate in 0.2 ml ethanol prior to use
Applications:	Suitable for LNP formulations



Description

DSPE-PEG(2000)-MACTIDE is a peptide-lipid conjugate composed of DSPE-PEG(2000)-DBCO conjugated to MACTIDE azide (Item No. 43997), a clickable form of a peptide that binds CD206 on tumor-associated macrophages (TAMs).¹ DSPE-PEG(2000)-MACTIDE is intended for use in the targeted delivery of lipid nanoparticles (LNPs) *in vitro* and *in vivo*.

Reference

1. Lepland, A., Peranzoni, E., Haljasorg, U., *et al.* Peptide-drug conjugate for therapeutic reprogramming of tumor-associated macrophages in breast cancer. *Adv. Sci. (Weinh)* **12**(10), e2410288 (2025).

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