

# PRODUCT INFORMATION

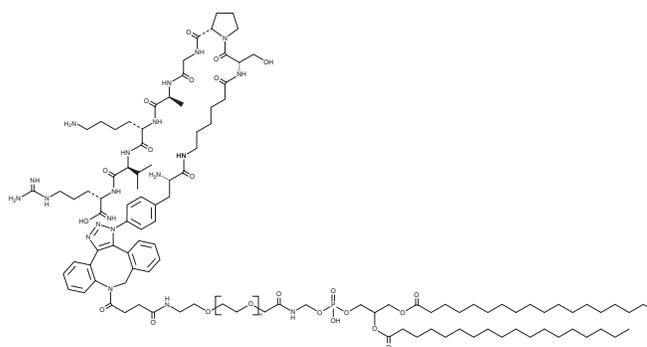


## DSPE-PEG(2000)-UNO Peptide

Item No. 41868

### Overview and Properties

<b>Contents:</b>	This vial contains 1 mg of DSPE-PEG(2000)-UNO click conjugate.
<b>Synonyms:</b>	CD206 Peptide Ligand, 1,2-Distearoyl Phosphatidylethanolamine-PEG(2000)-UNO, DSPE-PEG(2000)-SPGAKVR
<b>Peptide Sequence:</b>	SPGAKVR
<b>MF:</b>	$(C_2H_4O)_n C_{108}H_{175}N_{20}O_{22}P$
<b>Molecular Weight:</b>	3,958-4,224 (due to PEG congeners)
<b>Form:</b>	Solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	Dissolve conjugate in 0.2 ml ethanol prior to use.
<b>Applications:</b>	LNP formulations



### Description

DSPE-PEG(2000)-UNO is a peptide-ligand conjugate composed of DSPE-PEG(2000)-DBCO conjugated to azide-UNO peptide (Item No. 43998), a clickable form of a peptide that binds CD206 on tumor-associated macrophages (TAMs).<sup>1,2</sup>

### References

1. Scodeller, P., Simón-Gracia, L., Kopanchuk, S., *et al.* Precision targeting of tumor macrophages with a CD206 binding peptide. *Sci. Rep.* **7**(1), 14655 (2017).
2. Leland, 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, 01/28/2026

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