

# PRODUCT INFORMATION



## DSPE-PEG(2000)-Cyclo(RGDyK)

Item No. 41941

### Overview and Properties

**Contents:** This vial contains 200 µg of DSPE-PEG(2000)-cycloRGD conjugate.

**Synonyms:** 1,2-Distearoyl Phosphatidylethanolamine-PEG(2000)-cRGD Peptide, 1,2-DSPE-PEG(2000)-cRGDyK, DSPE-PEG(2000)-RGD, Integrin-Targeting Peptide Conjugate, PE(18:0/18:0)-PEG(2000)-CycloRGDyK

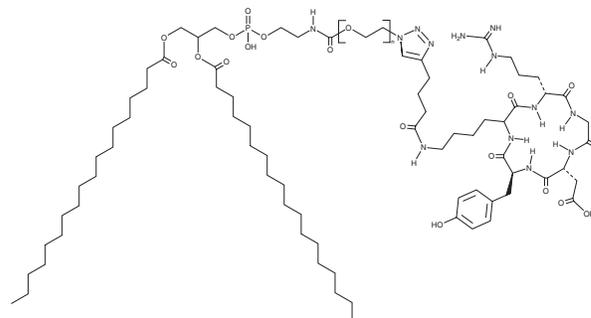
**MF:**  $(C_2H_4O)_n C_{75}H_{128}N_{13}O_{18}P$

**Form:** Solid

**Storage:** -20°C

**Stability:** ≥1 year

**Applications:** For dilution in ethanol and LNP formulations



### Description

DSPE-PEG(2000)-Cyclo(RGDyK) is a peptide-lipid conjugate. It was prepared from DSPE-PEG(2000)-azide (Item No. 40710) and alkyne-cyclo(RGDyK) (Item No. 41722), a clickable form of the peptide  $\alpha V\beta 3$  integrin ligand cyclo(RGDyK) (Item No. 35173), via copper click chemistry. DSPE-PEG(2000)-Cyclo(RGDyK) 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, 12/09/2025

**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