

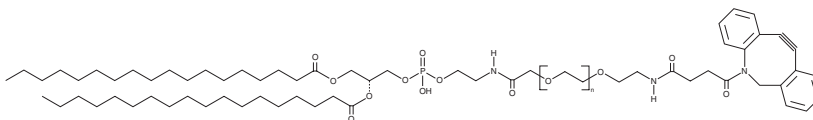
# PRODUCT INFORMATION



## DSPE-PEG(2000)-DBCO

Item No. 37798

CAS Registry No.: 3052817-89-4  
MF:  $(C_2H_4O)_n C_{64}H_{102}N_3O_{12}P$   
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^\circ C$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

DSPE-PEG(2000)-DBCO is supplied as a solid. A stock solution may be made by dissolving the DSPE-PEG(2000)-DBCO in the solvent of choice, which should be purged with an inert gas. DSPE-PEG(2000)-DBCO is soluble ( $\geq 10$  mg/ml) in ethanol.

### Description

DSPE-PEG(2000)-DBCO is a PEGylated and clickable form of 1,2-distearoyl-*sn*-glycero-3-PE (DSPE).

#### 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, 04/30/2025

#### 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