

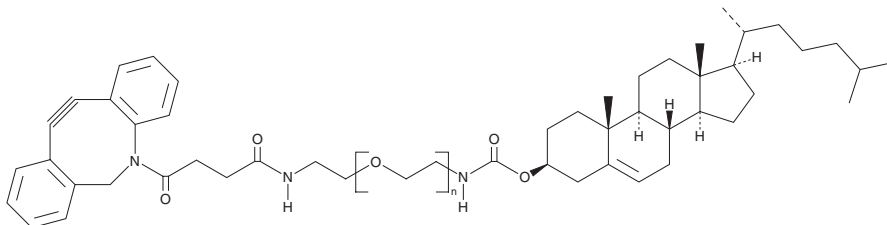
PRODUCT INFORMATION



Cholesterol-PEG(2000)-DBCO

Item No. 45803

Synonyms: Cholesterol-PEG₂₀₀₀-DBCO,
Cholesterol-PEG(2000)-
dibenzocyclooctyne
MF: (C₂H₄O)_nC₄₉H₆₅N₃O₄
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

Cholesterol-PEG(2000)-DBCO is a PEGylated and clickable form of cholesterol (Item Nos. 9003100 | 39088 | 39448). It has been used as a chemically reactive anchor lipid *in vitro* and is more efficiently incorporated into the cell membrane with less internalization than DSPE-PEG(2000)-DBCO (Item No. 44696).¹

Reference

1. Vabbilisetty, P., Boron, M., Nie, H., *et al.* Chemical reactive anchoring lipids with different performance for cell surface re-engineering application. *ACS Omega* **3**(2), 1589-1599 (2018).

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, 05/27/2026

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