

PRODUCT INFORMATION

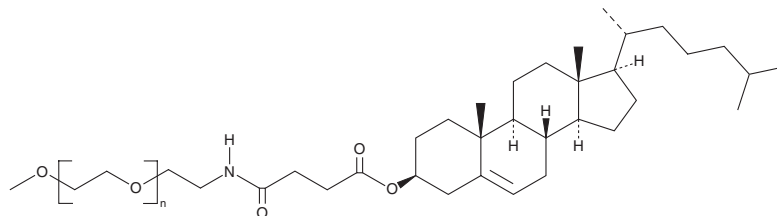


CLS-mPEG(2000)

Item No. 34647

Formal Name: (3S,8S,9S,10R,13R,14S,17R)-10,13-dimethyl-17-((R)-6-methylheptan-2-yl)-2,3,4,7,8,9,10,11,12,13,14,15,16,17-tetradecahydro-1H-cyclopenta[a]phenanthren-3-yl 4-((2-(2-methoxyethoxy)ethyl)amino)-4-oxobutanoate

Synonym: Cholesterol-mPEG
MF: $(C_2H_4O)_n C_{34}H_{57}NO_4$
Purity: $\geq 95\%$
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

CLS-mPEG(2000) is supplied as a solid. A stock solution may be made by dissolving the CLS-mPEG(2000) in the solvent of choice, which should be purged with an inert gas. CLS-mPEG(2000) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CLS-mPEG(2000) in these solvents is approximately 11 and 16 mg/ml, respectively. It is also slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CLS-mPEG(2000) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of CLS-mPEG(2000) in PBS (pH 7.2) is approximately 0.30 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

CLS-mPEG(2000) is a PEGylated form of cholesterol (Item No. 9003100). It has been used in the synthesis of liposomes for targeted drug delivery.¹

Reference

1. Zhao, X.B., Muthusamy, N., Byrd, J.C., *et al.* Cholesterol as a bilayer anchor for PEGylation and targeting ligand in folate-receptor-targeted liposomes. *J. Pharm. Sci.* **96(9)**, 2424-2435 (2007).

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, 04/11/2025

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