

PRODUCT INFORMATION



Cholesterol-PEG(2000)-amine (hydrochloride)

Item No. 45631

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 (2-(2-aminoethoxy)ethyl) carbamate, monohydrochloride

MF: $(C_2H_4O)_n C_{30}H_{52}N_2O_2 \cdot HCl$

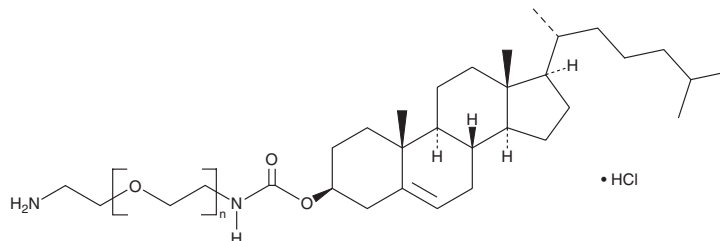
FW: 553.3

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

Cholesterol-PEG(2000)-amine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the cholesterol-PEG(2000)-amine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Cholesterol-PEG(2000)-amine (hydrochloride) is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

Cholesterol-PEG(2000)-amine is a PEGylated form of cholesterol (Item Nos. 9003100 | 39088). It has been used in the synthesis of conjugates containing cholesterol and pheophorbide a (Item No. 16072) for the formation of photosensitizing lipid nanoparticles (LNPs) with pro-oxidant and ferroptosis-inducing activity in cancer cells¹

Reference

1. Bae, G.H., Shin, S., Park, J.D., *et al.* Photosensitizing lipid nanoparticles for ferroptosis-enhanced photodynamic cancer therapy via GPX4 silencing. *Adv. Healthc. Mater.* **15**(7), e03748 (2026).

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, 04/21/2026

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