

PRODUCT INFORMATION

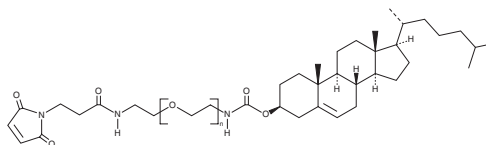


Cholesterol-PEG(2000)-Maleimide

Item No. 45804

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-(3-(2,5-dioxo-2,5-dihydro-1H-pyrrol-1-yl)propanamido)ethoxy)ethyl)carbamate
Synonym: Cholesterol-PEG₂₀₀₀-Maleimide
MF: (C₂H₄O)_nC₃₇H₅₇N₃O₅

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

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 cholesterol-PEG(2000)-maleimide can be prepared by directly dissolving the solid in aqueous buffers. Cholesterol-PEG(2000)-maleimide is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Cholesterol-PEG(2000)-maleimide is a PEGylated form of cholesterol (Item Nos. 9003100 | 39088 | 39448) that contains maleimide, which can react with the thiol group in cysteine residues to form a stable thioether bond.¹ It has been used to anchor an α₂-antiplasmin (α₂-AP) peptide to perfluorocarbon (PFC) nanoemulsions for use as thrombus-targeted contrast agents in a mouse model of pulmonary thromboembolism.

Reference

1. Temme, S., Grapentin, C., Quast, C., *et al.* Noninvasive imaging of early venous thrombosis by ¹⁹F magnetic resonance imaging with targeted perfluorocarbon nanoemulsions. *Circulation* **131**(16), 1405-1414 (2015).

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, 05/14x/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