

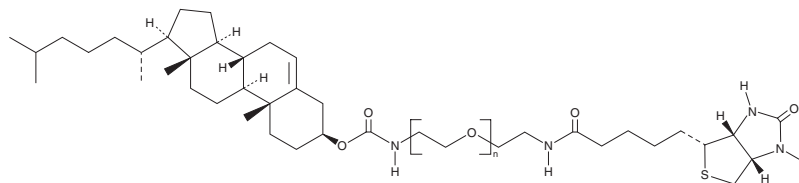
PRODUCT INFORMATION



Cholesterol-PEG(2000)-Biotin

Item No. 39036

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-(5-((3aR,4R,6aS)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanamido)ethoxy)ethyl)carbamate



Synonym: Biotin-PEG₂₀₀₀-Cholesterol

MF: (C₂H₄O)_nC₄₀H₆₆N₄O₄S

Purity: ≥90%

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)-biotin is supplied as a solid. A stock solution may be made by dissolving the cholesterol-PEG(2000)-biotin in the solvent of choice, which should be purged with an inert gas. Cholesterol-PEG(2000)-biotin is soluble in methanol.

Description

Cholesterol-PEG(2000)-biotin is a biotinylated and PEGylated form of cholesterol (Item Nos. 9003100 | 39088). It has been used as a lipid anchor to immobilize vesicles, liposomes, viral particles, bacteria, or yeast to microfluidic chips.¹ As a lipid anchor, cholesterol-PEG(2000)-biotin (10 μM) increases phycoerythrin-conjugated streptavidin density on the surface of mesenchymal stem cells greater than sulfo-NHS-LC-biotin.²

References

1. Kuhn, P., Eyer, K., Robinson, T., *et al.* A facile protocol for the immobilisation of vesicles, virus particles, bacteria, and yeast cells. *Integr. Biol. (Camb)*. **4**(12), 1550-1555 (2012).
2. Davis, K.A., Peng, H., Chelvarajan, L., *et al.* Increased yield of gelatin coated therapeutic cells through cholesterol insertion. *J. Biomed. Mater. Res. A*. **109**(3), 326-335 (2021).

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, 05/26/2026

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