

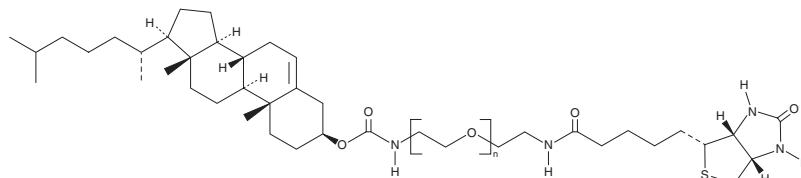
# PRODUCT INFORMATION



## 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



**Synonyms:** Biotin-PEG<sub>2000</sub>-Cholesterol, CholBio2k

**MF:** (C<sub>2</sub>H<sub>4</sub>O)<sub>n</sub>C<sub>40</sub>H<sub>66</sub>N<sub>4</sub>O<sub>4</sub>S

**FW:** 757.1

**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.<sup>1</sup> 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.<sup>2</sup>

### 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, 03/05/2024

#### 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