

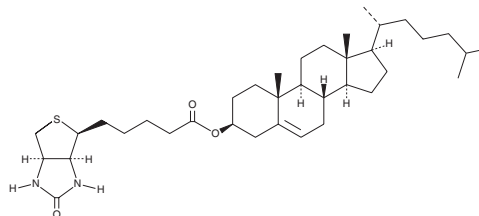
PRODUCT INFORMATION



Biotin-Cholesterol

Item No. 38852

CAS Registry No.: 2454045-15-7
Formal Name: (3 β)-cholest-5-en-3-ol, 3-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazole-4-pentanoate]
Synonym: Biotin-Provitamin D
MF: C₃₇H₆₀N₂O₃S
FW: 613.0
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

Biotin-cholesterol is supplied as a solid. A stock solution may be made by dissolving the biotin-cholesterol in the solvent of choice, which should be purged with an inert gas. Biotin-cholesterol is soluble in chloroform.

Description

Biotin-cholesterol is a biotinylated form of the major sterol cholesterol (Item No. 9003100). It has been used in the synthesis of ergosterol peroxide derivatives for cellular localization studies.¹ Biotin-cholesterol has also been used in the synthesis of biotin-conjugated liposomes and micelles for drug delivery to cancer cells *in vitro*.²

References

1. Ling, T., Lang, W.H., Martinez-Montemayor, M.M., *et al.* Development of ergosterol peroxide probes for cellular localisation studies. *Org. Biomol. Chem.* **17**(21), 5223-5229 (2019).
2. Li, W., Deng, L., Moosa, B., *et al.* "Nail" and "comb" effects of cholesterol modified NIPAm oligomers on cancer targeting liposomes. *Biomater. Sci.* **2**(4), 476-483 (2014).

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, 06/15/2023

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