

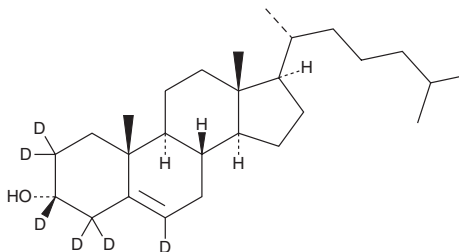
PRODUCT INFORMATION



Epicholesterol-d₆

Item No. 45739

CAS Registry No.: 92543-07-2
Formal Name: (3 α)-cholest-5-en-2,2,3,4,4,6-d₆-3-ol
Synonyms: α -Cholesterol-d₆, 3 α -Cholesterol-d₆, 3-Epicholesterol-d₆
MF: C₂₇H₄₀D₆O
FW: 392.7
Chemical Purity: \geq 98% (Epicholesterol)
Deuterium Incorporation: \geq 99% deuterated forms (d₁-d₆); \leq 1% d₀
Supplied as: A 1 mg/ml solution in ethanol
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Epicholesterol-d₆ is intended for use as an internal standard for the quantification of epicholesterol by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Description

Epicholesterol is an epimer of cholesterol (Item Nos. 9003100 | 39088 | 39448), where the C-3 hydroxyl group is in the α configuration.¹ It has been used to study the stereospecific effects of cholesterol on antibody recognition and receptor function.^{1,2} Epicholesterol has also been used to study the interactions of cholesterol-containing membranes with stereoisomeric ligands.¹

References

1. Tsuchiya, H. and Mizogami, M. Discrimination of stereoisomers by their enantioselective interactions with chiral cholesterol-containing membranes. *Molecules* **23**(1), 49 (2017).
2. Geva, M., Izhaky, D., Mickus, D.E., *et al.* Stereoselective recognition of monolayers of cholesterol, *ent*-cholesterol, and epicholesterol by an antibody. *Chembiochem.* **2**(4), 265-271 (2001).
3. Jafurulla, M., Rao, B.D., Sreedevi, S., *et al.* Stereospecific requirement of cholesterol in the function of the serotonin1A receptor. *Biochim. Biophys. Acta.* **1838**(1 Pt B), 158-163 (2013).

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