

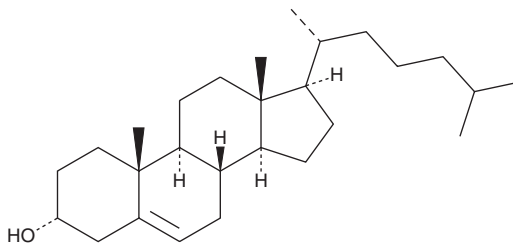
PRODUCT INFORMATION



Epicholesterol

Item No. 45822

CAS Registry No.: 474-77-1
Formal Name: (3 α)-cholest-5-en-3-ol
Synonyms: α -Cholesterol, 3 α -Cholesterol, 3-Epicholesterol
MF: C₂₇H₄₆O
FW: 386.7
Purity: $\geq 95\%$
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

Epicholesterol is supplied as a solid. A stock solution may be made by dissolving the epicholesterol in the solvent of choice, which should be purged with an inert gas. Epicholesterol is slightly soluble (0.1-1 mg/ml) in ethanol.

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.^{2,3} 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, 07/14/2026

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