

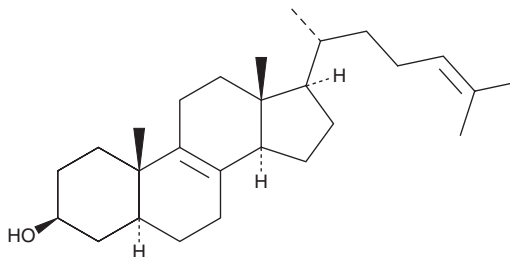
PRODUCT INFORMATION



Zymosterol

Item No. 45819

CAS Registry No.: 128-33-6
Formal Name: (5 α)-cholesta-8,24-dien-3 β -ol
MF: C₂₇H₄₄O
FW: 384.6
Purity: $\geq$ 95%
Supplied as: A solid
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

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

Description

Zymosterol is a precursor in the biosynthesis of cholesterol (Item Nos. 9003100 | 39088 | 39448).¹ It accumulates in the plasma membrane in fibroblasts and translocates to the endoplasmic reticulum (ER) during cholesterol synthesis.^{2,3}

References

1. Alexander, G.J. and Schwenk, E. Studies on biosynthesis of cholesterol. IX. Zymosterol as a precursor of cholesterol. *Arch. Biochem. and Biophys.* **66**(2), 381-387 (1957).
2. Lange, Y., Echevarria, F., and Steck, T.L. Movement of zymosterol, a precursor of cholesterol, among three membranes in human fibroblasts. *J. Biol. Chem.* **266**(32), 21439-21443 (1991).
3. Lange, Y. Cholesterol movement from plasma membrane to rough endoplasmic reticulum. Inhibition by progesterone. *J. Biol. Chem.* **269**(5), 3411-3414 (1994).

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, 06/10/2026

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