

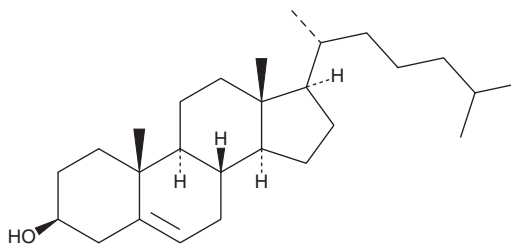
PRODUCT INFORMATION



Cholesterol (plant)

Item No. 39448

CAS Registry No.: 57-88-5
Formal Name: (3 β)-cholest-5-en-3-ol
Synonyms: NSC 8798, Provitamin D
MF: C₂₇H₄₆O
FW: 386.7
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

Cholesterol (plant) is supplied as a solid. A stock solution may be made by dissolving the cholesterol (plant) in the solvent of choice, which should be purged with an inert gas. Cholesterol (plant) is soluble in chloroform and methanol.

Description

Cholesterol is a minor sterol produced in plant cells and a component of plant surface lipids.¹⁻³ Cholesterol content in plants is generally several hundreds- to thousands-fold less than that of mammalian cells.¹ It is a precursor in the synthesis of various bioactive plant metabolites, including steroidal glycoalkaloids and phytoecdysteroids.²

References

1. Nakajima, N., Fujioka, S., Tanaka, T., *et al.* Biosynthesis of cholestanol in higher plants. *Phytochemistry* **60**(3), 275-279 (2002).
2. Sonawane, P.D., Pollier, J., Panda, S., *et al.* Plant cholesterol biosynthetic pathway overlaps with phytosterol metabolism. *Nat. Plants* **3**(1), 16205 (2016).
3. Behrman, E.J. and Gopalan, V. Cholesterol and plants. *J. Chem. Educ.* **82**(12), 1791-1793 (2005).

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, 10/24/2023

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