

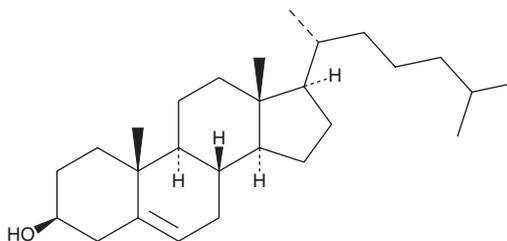
PRODUCT INFORMATION



Cholesterol

Item No. 9003100

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 is supplied as a solid. A stock solution may be made by dissolving the cholesterol in the solvent of choice, which should be purged with an inert gas. Cholesterol is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

Cholesterol is a major sterol produced in mammalian cells that is required for cell viability and proliferation.¹ It is a component of mammalian cell membranes that interacts with membrane phospholipids, sphingolipids, and proteins to influence their behavior. It is also a component of various lipid-based drug delivery (LBDD) systems, including liposomes and lipid nanoparticles (LNPs), where it has a role in membrane stability.² Cholesterol is a precursor of steroid hormones, bile acids, and the active form of vitamin D. Impaired cholesterol homeostasis is related to development of various diseases including fatty liver, diabetes, gallstones, dyslipidemia, atherosclerosis, heart attack, and stroke.³

References

1. Ohvo-Rekilä, H., Ramstedt, B., Leppimäki, P., *et al.* Cholesterol interactions with phospholipids in membranes. *Prog. Lipid Res.* **41(1)**, 66-97 (2002).
2. Tenchov, R., Bird, R., Curtze, A.E., *et al.* Lipid nanoparticles-from liposomes to mRNA vaccine delivery, a landscape of research diversity and advancement. *ACS Nano* **15(11)**, 16982-17015 (2021).
3. Yamanashi, Y., Takada, T., and Suzuki, H. Associations between lifestyle-related diseases and transporters involved in intestinal absorption and biliary excretion of cholesterol. *Biol. Pharm. Bull.* **41(1)**, 1-10 (2018).

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, 10/11/2022

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