

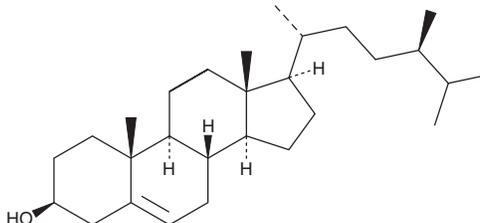
PRODUCT INFORMATION



Campesterol

Item No. 41695

CAS Registry No.: 474-62-4
Formal Name: (3 β ,24R)-ergost-5-en-3-ol
Synonyms: Campesterin, NSC 224330
MF: C₂₈H₄₈O
FW: 400.7
Purity: \geq 98% (mixture of epimers)
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/Soybean



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

Laboratory Procedures

Campesterol is supplied as a solid. A stock solution may be made by dissolving the campesterol in the solvent of choice, which should be purged with an inert gas. Campesterol is sparingly soluble (1-10 mg/ml) in chloroform and methanol.

Description

Campesterol is a phytosterol that has been found in *D. innoxia* and has diverse biological activities.¹⁻⁴ It inhibits basic FGF-induced proliferation in, and vessel formation by, human umbilical vein endothelial cells (HUVECs) when used at concentrations of 1, 5, or 10 μ M.² Campesterol (15 μ M) decreases the proliferation of MCF-7 and ZR-75-1 breast cancer cells.³ It reduces the levels of estrogen receptor α (ER α) in, and the diameter of, patient-derived breast cancer organoids when used at a concentration of 10 μ M. Campesterol has been used as a marker of cholesterol intestinal absorption in humans.⁴

References

1. Ramadan, M.F., Zayed, R., and El-Shamy, H. Screening of bioactive lipids and radical scavenging potential of some solanaceae plants. *Food Chem.* **103(3)**, 885-890 (2007).
2. Choi, J.M., Lee, E.O., Lee, H.J., et al. Identification of campesterol from *Chrysanthemum coronarium* L. and its antiangiogenic activities. *Phytother. Res.* **21(10)**, 954-959 (2007).
3. Majumder, R., Banerjee, S., Mandal, M., et al. A virtual drug discovery screening illuminates campesterol as a potent estrogen receptor alpha inhibitor in breast cancer. *J. Med. Chem.* **67(12)**, 10321-10335 (2024).
4. Wu, A.H., Ruan, W., Todd, J., et al. Biological variation of β -sitosterol, campesterol, and lathosterol as cholesterol absorption and synthesis biomarkers. *Clin. Chim. Acta* **430**, 43-47 (2014).

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, 12/18/2025

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