

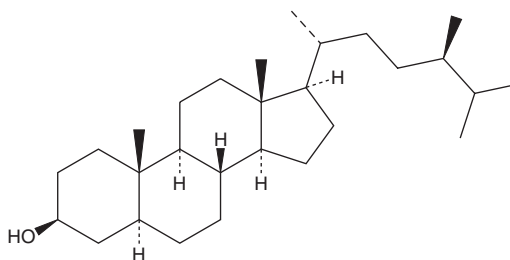
PRODUCT INFORMATION



Campestanol

Item No. 26538

CAS Registry No.: 474-60-2
Formal Name: (3 β ,5 α ,24R)-ergostan-3-ol
Synonym: 24-Methyl-5 α -cholestan-3 β -ol
MF: C₂₈H₅₀O
FW: 402.7
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Campestanol is supplied as a solid. A stock solution may be made by dissolving the campestanol in the solvent of choice, which should be purged with an inert gas. Campestanol is slightly soluble in chloroform.

Description

Campestanol is a phytosterol found in vegetables, fruits, nuts, and seeds.¹ Dietary administration of vegetable stanol esters (1% w/w) containing campestanol (30.1%) and sitostanol (65.7%; Item No. 26094) lower serum cholesterol, LDL, VLDL, and intermediate-density lipoprotein (IDL) cholesterol levels in an APOE*3-Leiden transgenic mouse model of atherosclerosis.² This combination of campestanol with sitostanol also reduces atherosclerotic lesion area by 91% and the proportion of lesions with extensive macrophage infiltration.

References

1. Gylling, H. and Miettinen, T.A. Cholesterol reduction by different plant stanol mixtures and with variable fat intake. *Metabolism* **48(5)**, 575-580 (1999).
2. Volger, O.L., Mensink, R.P., Plat, J., *et al.* Dietary vegetable oil and wood derived plant stanol esters reduce atherosclerotic lesion size and severity in apoE*3-Leiden transgenic mice. *Atherosclerosis* **157(2)**, 375-381 (2001).

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, 11/01/2022

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