

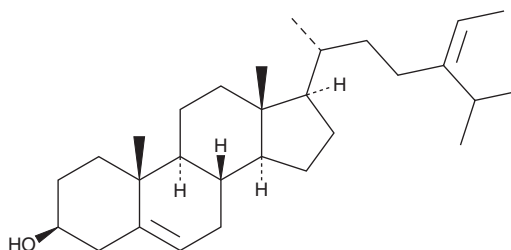
PRODUCT INFORMATION



Δ^5 -Avenasterol

Item No. 30154

CAS Registry No.: 18472-36-1
Formal Name: stigmasta-5,24(28)-dien-3 β -ol
Synonym: 24-ethylidene Cholesterol
MF: C₂₉H₄₈O
FW: 412.7
Purity: \geq 95% (mixture of isomers)
Supplied as: A solid
Storage: -20°C
Stability: \geq 2 years
Item Origin: Synthetic



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

Laboratory Procedures

Δ^5 -Avenasterol is supplied as a solid. A stock solution may be made by dissolving the Δ^5 -avenasterol in the solvent of choice, which should be purged with an inert gas. Δ^5 -Avenasterol is slightly soluble in chloroform and ethyl acetate.

Description

Δ^5 -Avenasterol is a phytosterol that has been found in various plant extracts, including *D. metel*, *D. innoxia*, *D. stramonium*, *D. tatula*, *H. muticus*, and *H. niger*.¹

Reference

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).

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, 03/30/2020

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