

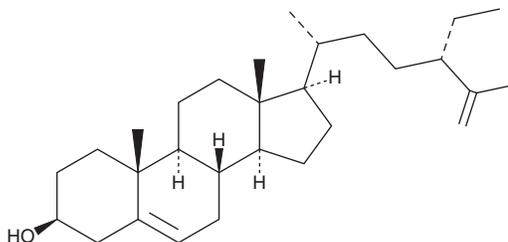
PRODUCT INFORMATION



Clerosterol

Item No. 35862

CAS Registry No.: 2364-23-0
Formal Name: (3 β ,24S)-stigmasta-5,25-dien-3-ol
Synonyms: (-)-Clerosterol, NSC 144945,
 $\Delta^{5,25}$ -Stigmastadienol
MF: C₂₉H₄₈O
FW: 412.7
Purity: \geq 95%
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

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

Description

Clerosterol is a phytosterol that has been found in *C. fragile* and has diverse biological activities.^{1,2} It is active against *T. b. brucei* (IC₅₀ = 53.6 μ g/ml).¹ Clerosterol inhibits the proliferation of A2058 melanoma cells (IC₅₀ = 150 μ M).² It induces apoptosis, cell cycle arrest at the sub-G₁ phase, and mitochondrial membrane depolarization in A2058 cells when used at a concentration of 150 μ M.

References

1. Hoet, S., Pieters, L., Muccioli, G.G., *et al.* Antitrypanosomal activity of triterpenoids and sterols from the leaves of *Strychnos spinosa* and related compounds. *J. Nat. Prod.* **70(8)**, 1360-1363 (2007).
2. Kim, A.D., Lee, Y., Kang, S.-H., *et al.* Cytotoxic effect of clerosterol isolated from *Codium fragile* on A2058 human melanoma cells. *Mar. Drugs* **11(2)**, 418-430 (2013).

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, 06/03/2024

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