

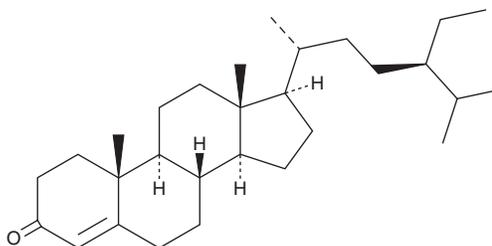
PRODUCT INFORMATION



β-Sitostenone

Item No. 36703

CAS Registry No.: 1058-61-3
Formal Name: stigmast-4-en-3-one
Synonym: NSC 49082
MF: C₂₉H₄₈O
FW: 412.7
Purity: ≥95%
UV/Vis.: λ_{max}: 241 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

β-Sitostenone is supplied as a solid. A stock solution may be made by dissolving the β-sitostenone in the solvent of choice, which should be purged with an inert gas. β-Sitostenone is soluble in organic solvents such as ethanol and dimethyl formamide (DMF). β-Sitostenone soluble in DMF at a concentration of approximately 0.33 mg/ml and is slightly soluble in ethanol.

Description

β-Sitostenone is a phytosterol that has been found in *A. occidentale* and has diverse biological activities.¹⁻⁴ It inhibits basal nitric oxide (NO) production in RAW 264.7 macrophages (IC₅₀ = 15.9 μM).¹ β-Sitostenone is active against *T. evansi* (IC₅₀ = 150 nM).² It reduces mean blood pressure in normotensive rats when administered at a dose of 2 mg/kg.³ β-Sitostenone (1.3 and 3 mg/kg) decreases blood glucose levels in dogs.⁴

References

1. Tewtrakul, S. and Subhadhirasakul, S. Effects of compounds from *Kaempferia parviflora* on nitric oxide, prostaglandin E₂ and tumor necrosis factor-alpha productions in RAW264.7 macrophage cells. *J. Ethnopharmacol.* **120(1)**, 81-84 (2008).
2. Nyunt, K.S., Elkhateeb, A., Tosa, Y., et al. Isolation of antitrypanosomal compounds from *Vitis repens*, a medicinal plant of Myanmar. *Nat. Prod. Commun.* **7(5)**, 609-610 (2012).
3. Barla, A., Birman, H., Kültür, Ş., et al. Secondary metabolites from *Euphorbia helioscopia* and their vasodepressor activity. *Turk. J. Chem.* **30(3)**, 325-332 (2006).
4. Alexander-Lindo, R.L., Morrison, E.Y.S.A., and Nair, M.G. Hypoglycaemic effect of stigmast-4-en-3-one and its corresponding alcohol from the bark of *Anacardium occidentale* (cashew). *Phytother. Res.* **18(5)**, 403-407 (2004).

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, 05/17/2023

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