

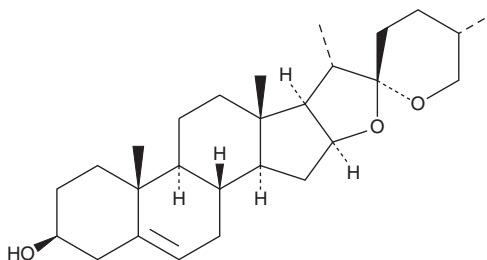
PRODUCT INFORMATION



Yamogenin

Item No. 36008

CAS Registry No.: 512-06-1
Formal Name: (3β,25S)-spirost-5-en-3-ol
Synonyms: Neodiosgenin, NSC 226132
MF: C₂₇H₄₂O₃
FW: 414.6
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Dioscorea collettii*



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

Laboratory Procedures

Yamogenin is supplied as a solid. A stock solution may be made by dissolving the yamogenin in the solvent of choice, which should be purged with an inert gas. Yamogenin is soluble in organic solvents such as ethanol and chloroform. The solubility of yamogenin in these solvents is approximately 1 and 10 mg/ml, respectively.

Description

Yamogenin is a steroidal sapogenin that has been found in fenugreek (*T. foenum graecum*) and has diverse biological activities.^{1,2} It inhibits activation of the liver X receptor (LXR) induced by the LXRα and LXRβ agonist T0901317 (Item No. 71810) in HepG2 cells in a reporter assay when used at concentrations of 5 or 10 μM.¹ Yamogenin (10 μM) decreases triglyceride levels in primary mouse hepatocytes. It is cytotoxic to HaCaT keratinocytes, as well as HeLa cervical and SKOV3 ovarian cancer cells (IC₅₀s = 16.4, 19.6, and 23.9 μg/ml, respectively).²

References

1. Moriwaki, S., Murakami, H., Takahashi, N., *et al.* Yamogenin in fenugreek inhibits lipid accumulation through the suppression of gene expression in fatty acid synthesis in hepatocytes. *Biosci. Biotechnol. Biochem.* **78(7)**, 1231-1236 (2014).
2. Stefanowicz-Hajduk, J., Kól-Kogus, B., Spzrak-Stefanowska, B., *et al.* Cytotoxic activity of standardized extracts, a fraction, and individual secondary metabolites from fenugreek seeds against SKOV-3, HeLa and MOLT-4 cell lines. *Pharm. Biol.* **59(1)**, 424-437 (2021).

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, 12/15/2022

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