

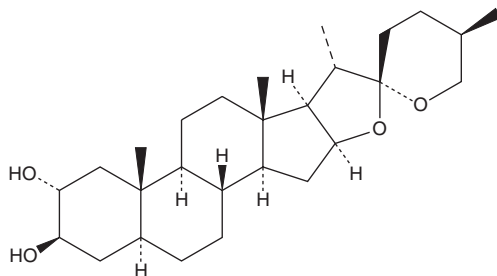
PRODUCT INFORMATION



Gitogenin

Item No. 34681

CAS Registry No.: 511-96-6
Formal Name: (2 α ,3 β ,5 α ,25R)-spirostan-2,3-diol
Synonyms: NSC 147752, 5 α -hydroxy Tigogenin
MF: C₂₇H₄₄O₄
FW: 432.6
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Tribulus terrestris* L.



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

Laboratory Procedures

Gitogenin is supplied as a solid. A stock solution may be made by dissolving the gitogenin in the solvent of choice, which should be purged with an inert gas. Gitogenin is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

Gitogenin is a sapogenin that has been found in *T. foenum-graecum* and has enzyme inhibitory and hormone-releasing activities.^{1,2} It is an inhibitor of UDP-glucuronosyltransferase (UGT) isoform UGT1A4 (IC₅₀ = 0.69 μ M).¹ Gitogenin (20 μ g/ml) stimulates the release of growth hormone in primary rat pituitary cells.²

References

1. Xu, M., Dong, P., Tian, X., *et al.* Drug interaction study of natural steroids from herbs specifically toward human UDP-glucuronosyltransferase (UGT) 1A4 and their quantitative structure activity relationship (QSAR) analysis for prediction. *Pharmacol. Res.* **110**, 139-150 (2016).
2. Shim, S.H., Lee, E.J., Kim, J.S., *et al.* Rat growth-hormone release stimulators from fenugreek seeds. *Chem. Biodivers.* **5**(9), 1753-1761 (2008).

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, 11/02/2022

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