

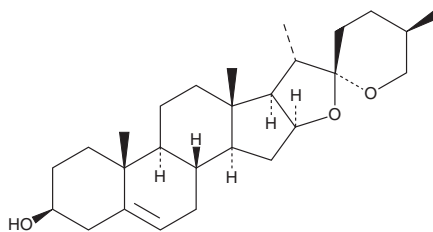
PRODUCT INFORMATION



Diosgenin

Item No. 19847

CAS Registry No.: 512-04-9
Formal Name: (3 β ,25R)-spirost-5-en-3-ol
Synonyms: Nitogenin, NSC 33396,
3 β -hydroxy-5-Spirostene
MF: C₂₇H₄₂O₃
FW: 414.6
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: Room temperature
Stability: \geq 4 years



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

Laboratory Procedures

Diosgenin is supplied as a crystalline solid. A stock solution may be made by dissolving the diosgenin in the solvent of choice. Diosgenin is soluble in organic solvents such as chloroform and dimethyl formamide, which should be purged with an inert gas. The solubility of diosgenin in these solvents is approximately 30 and 1 mg/ml, respectively.

Description

Diosgenin is a natural steroid sapogenin that has diverse biochemical and physiological actions. It has antioxidant and anti-inflammatory activities, is protective against metabolic and cardiovascular diseases, and can have immunomodulatory and estrogenic effects.¹⁻⁴

References

1. Lee, K. J., Oh, Y. C., Cho, W. K., *et al.* Antioxidant and anti-inflammatory activity determination of one hundred kinds of pure chemical compounds using offline and online screening HPLC assay. *Evid. Based Complement. Alternat. Med.* (2015).
2. Sung, B., Prasad, S., Yadav, V. R., *et al.* Cancer cell signaling pathways targeted by spice-derived nutraceuticals. *Nutr. Cancer* **64**(2), 173-197 (2012).
3. Chen, Y., Tang, Y. M., Yu, S. L., *et al.* Advances in the pharmacological activities and mechanisms of diosgenin. *Chin. J. Nat. Med.* **13**(8), 578-587 (2015).
4. Fuller, S., and Stephens, J. S. Diosgenin, 4-hydroxyisoleucine, and fiber from fenugreek: Mechanisms of actions and potential effects on metabolic syndrome. *Adv. Nutr.* **6**(2), 189-197 (2015).

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, 06/08/2023

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