PRODUCT INFORMATION



Tigogenin

Item No. 30137

CAS Registry No.:	77-60-1	
Formal Name:	(3β,5α,25R)-spirostan-3-ol	
Synonym:	NSC 93754	н /
MF:	C ₂₇ H ₄₄ O ₃	
FW:	416.6	
Purity:	≥90%	
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥4 years	но
Item Origin:	Natural/Source unknown	Н

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

Laboratory Procedures

Tigogenin is supplied as a solid. A stock solution may be made by dissolving the tigogenin in the solvent of choice, which should be purged with an inert gas. Tigogenin is soluble in the organic solvent dimethyl formamide at a concentration of approximately 2 mg/ml.

Description

Tigogenin is a steroidal sapogenin that has been found in A. sisalana and has diverse biological activities.¹⁻⁴ It increases the proliferation rate of, as well as mRNA levels of genes encoding the osteoblastic differentiation markers Cbfa1, collagen type I, and osteocalcin in, mouse bone marrow stromal cells (BMSCs) in a concentration-dependent manner.² It also increases matrix calcium deposition in BMSCs when used at concentrations of 30 and 90 μ M. Tigogenin is active against the fungus A. *fumigatus* (MIC₅₀ = 16 μ g/ml).³ It inhibits carrageenan-induced paw edema in rats when administered at a dose of 4.2 μ g/kg.

References

- 1. Cripps, A.L. and Blunden, G. A quantitative gas-liquid chromatographic method for the estimation of hecogenin and tigogenin in the leaves, juice and sapogenin concentrates of Agave sisalana. Steroids 31(5), 661-669 (1978).
- 2. Zhou, H., Yang, X., Wang, N., et al. Tigogenin inhibits adipocytic differentiation and induces osteoblastic differentiation in mouse bone marrow stromal cells. Mol. Cell Endocrinol. 270(1-2), 17-22 (2007).
- Upadhyay, S.K., Creech, C.C., Bowdy, K.L., et al. Synthesis and antifungal activity of functionalized 3 2,3-spirostane isomers. Bioorg. Med. Chem. Lett. 21(10), 2826-2831 (2011).
- 4. Peana, A.T., Moretti, M.D., Manconi, V., et al. Anti-inflammatory activity of aqueous extracts and steroidal sapogenins of Agave americana. Planta Med. 63(3), 199-202 (1997).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 12/12/2022

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