

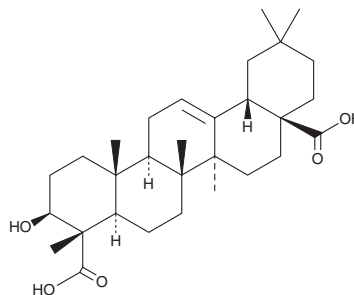
PRODUCT INFORMATION



Gypsogenic Acid

Item No. 36025

CAS Registry No.: 5143-05-5
Formal Name: (3 β ,4 α)-3-hydroxy-olean-12-ene-23,28-dioic acid
Synonym: Astrantiagenin J
MF: C₃₀H₄₆O₅
FW: 486.7
Purity: $\geq 95\%$
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

Gypsogenic acid is supplied as a solid. A stock solution may be made by dissolving the gypsogenic acid in the solvent of choice, which should be purged with an inert gas. Gypsogenic acid is slightly soluble in DMSO and methanol.

Description

Gypsogenic acid is a triterpene acid that has been found in *M. stenostachya* and has antibacterial and trypanocidal activities.^{1,2} It is active against the oral bacterial pathogens *E. faecalis*, *S. salivarius*, *S. sanguinis*, *S. mitis*, *S. mutans*, and *S. sobrinus* (MICs = 50-200 $\mu\text{g/ml}$).¹ Gypsogenic acid induces lysis of *T. cruzi* in isolated mouse blood (IC₅₀ = 56.6 μM).²

References

1. Cunha, L.C.S., e Silva, M.L.A., Furtado, N.A.J.C., et al. Antibacterial activity of triterpene acids and semi-synthetic derivatives against oral pathogens. *Z. Naturforsch. C. J. Biosci.* **62(9-10)**, 668-672 (2007).
2. Cunha, W.R., Martins, C.H., Ferreira, D.D.S., et al. *In vitro* trypanocidal activity of triterpenes from *Miconia* species. *Planta Med.* **69(5)**, 470-472 (2003).

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, 11/30/2022

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