

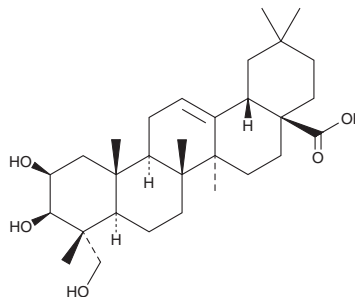
PRODUCT INFORMATION



Bayogenin

Item No. 35822

CAS Registry No.: 6989-24-8
Formal Name: (2 β ,3 β ,4 α)-2,3,23-trihydroxy-olean-12-en-28-oic acid
MF: C₃₀H₄₈O₅
FW: 488.7
Purity: \geq 90%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Bolbostemma paniculatum*



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

Laboratory Procedures

Bayogenin is supplied as a solid. A stock solution may be made by dissolving the bayogenin in the solvent of choice, which should be purged with an inert gas. Bayogenin is soluble in chloroform, DMSO, ethyl acetate, acetone, and dichloromethane.

Description

Bayogenin is a triterpene that has been found in *B. perennis* and has enzyme inhibitory activity.^{1,2} It inhibits rabbit muscle glycogen phosphorylase (IC₅₀ = 68 μ M).²

References

- Schöpke, T., Wray, V., Kunath, A., *et al.* Bayogenin and asterogenic acid glycosides from *Bellis perennis*. *Phytochemistry* **31(7)**, 2555-2557 (1992).
- Wen, X.-A., Liu, J., Zhang, L.-Y., *et al.* Synthesis and biological evaluation of arjunolic acid, bayogenin, hederagonic acid and 4-*epi*-hederagonic acid as glycogen phosphorylase inhibitors. *Chin. J. Nat. Med.* **8(6)**, 441-448 (2010).

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