

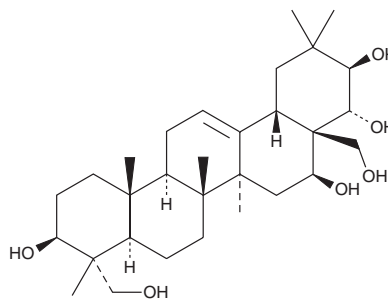
PRODUCT INFORMATION



Gymnemagenin

Item No. 11713

CAS Registry No.: 22467-07-8
Formal Name: 4 α -olean-12-ene-3 β ,16 β ,21 β ,22 α ,23,28-hexol
MF: C₃₀H₅₀O₆
FW: 506.7
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

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

Gymnemagenin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, gymnemagenin should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Gymnemagenin has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Gymnemagenin is the aglycone core of gymnemic acids A and B, triterpenoid sweetness inhibitors derived from *G. sylvestre*.¹ It is used as a biomarker for the quantification of gymnemic acids in medicinal plant extracts.²

References

1. Suttisri, R., Lee, I.S., and Kinghorn, A.D. Plant-derived triterpenoid sweetness inhibitors. *J. Ethnopharmacol.* **47(1)**, 9-26 (1995).
2. Mandal, V., Dewanjee, S., and Mandal, S.C. Microwave-assisted extraction of total bioactive saponin fraction from *Gymnema sylvestre* with reference to gymnemagenin: A potential biomarker. *Phytochem Anal.* **20(6)**, 491-497 (2009).

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, 10/26/2022

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