

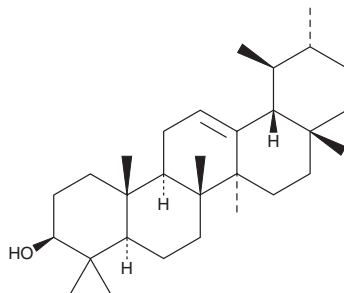
PRODUCT INFORMATION



α -Amyrin

Item No. 29831

CAS Registry No.: 638-95-9
Formal Name: (3 β)-urs-12-en-3-ol
Synonym: NSC 114787
MF: C₃₀H₅₀O
FW: 426.7
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/Burseraceae



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

Laboratory Procedures

α -Amyrin is supplied as a solid. A stock solution may be made by dissolving the α -amyrin in the solvent of choice, which should be purged with an inert gas. α -Amyrin is soluble in the organic solvent methanol.

Description

α -Amyrin is a triterpenoid that has been found in *B. copallifera* and has diverse biological activities.¹⁻⁴ It inhibits HIV-1 reverse transcriptase (IC₅₀ = 3.3 μ M).¹ α -Amyrin inhibits the proliferation of B16 2F2 melanoma cells and LPS-induced nitric oxide (NO) production in RAW 264.7 macrophages (IC₅₀s = 50 and 8.98 μ M, respectively).^{2,3} It reduces myeloperoxidase (MPO) activity, indicating a decrease in polymorphonuclear leukocyte (PMN) infiltration, and ear edema induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in mice (ID₅₀s = 0.45 and 0.31 mg/ear, respectively).⁴

References

1. Akihisa, T., Ogihara, J., Kato, J., *et al.* Inhibitory effects of triterpenoids and sterols on human immunodeficiency virus-1 reverse transcriptase. *Lipids* **36**(5), 507-512 (2001).
2. Hata, K., Hori, K., and Takahashi, S. Differentiation- and apoptosis-inducing activities by pentacyclic triterpenes on a mouse melanoma cell line. *J. Nat. Prod.* **65**(5), 645-648 (2002).
3. Romero-Estrada, A., Maldonado-Magaña, A., González-Christen, J., *et al.* Anti-inflammatory and antioxidative effects of six pentacyclic triterpenes isolated from the Mexican copal resin of *Bursera copallifera*. *BMC Complement. Altern. Med.* **16**(1), 422 (2016).
4. Otuki, M.F., Vieira-Lima, F., Malheiros, Â., *et al.* Topical antiinflammatory effects of the ether extract from *Protium kleinii* and α -amyrin pentacyclic triterpene. *Eur. J. Pharmacol.* **507**(1-3), 253-259 (2005).

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, 10/31/2023

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