

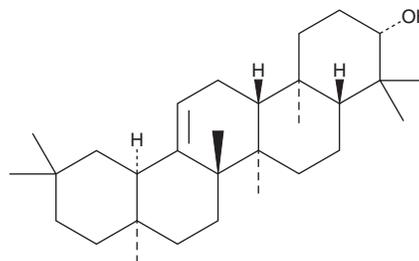
PRODUCT INFORMATION



β-Amyrin

Item No. 20949

CAS Registry No.: 559-70-6
Formal Name: (3β)-olean-12-en-3-ol
Synonyms: (+)-β-Amyrin, NSC 527971, Olean-12-en-3β-ol
MF: C₃₀H₅₀O
FW: 426.7
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

β-Amyrin is supplied as a crystalline solid. A stock solution may be made by dissolving the β-amyryn in the solvent of choice, which should be purged with an inert gas. β-Amyryn is soluble in organic solvents such as ethanol and dimethyl formamide (DMF). The solubility of β-amyryn in these solvents is approximately 1 and 3 mg/ml, respectively.

β-Amyryn is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, β-Amyryn should first be dissolved in DMF and then diluted with the aqueous buffer of choice. β-Amyryn has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

β-Amyryn is a natural pentacyclic triterpenoid commonly found in plants. It is a member of the oleanane group of secondary metabolites biosynthesized from squalene in the triterpenoid pathway. β-Amyryn inhibits collagen-induced platelet aggregation (IC₅₀ = 10.5 μM), inhibits acetylcholinesterase (IC₅₀ = 165 μM), and has antifungal and hepatoprotective effects.¹⁻⁴ Extracts of certain plants, used in herbal medicine to treat inflammation, have been shown to contain β-amyryn, suggesting that this compound contributes to anti-inflammatory actions.⁵

References

1. Ching, J., Chua, T.K., Chin, L.C., *et al.* Beta-amyryn from *Ardisia elliptica* Thunb. is more potent than aspirin in inhibiting collagen-induced platelet aggregation. *Indian J. Exp. Biol.* **48(3)**, 275-279 (2010).
2. Jabeen, K., Javid, A., Ahmad, E., *et al.* Antifungal compounds from *Melia azedarach* leaves for management of *Ascochyta rabiei*, the cause of chickpea blight. *Nat. Prod. Res.* **25(3)**, 264-276 (2011).
3. Oliveira, F.A., Chaves, M.H., Almeida, F.R., *et al.* Protective effect of alpha- and beta-amyryn, a triterpene mixture from *Protium heptaphyllum* (Aubl.) March. trunk wood resin, against acetaminophen-induced liver injury in mice. *J. Ethnopharmacol.* **98(1-2)**, 103-108 (2005).
4. Park, S.J., Ahn, Y.J., Oh, S.R., *et al.* Amyryn attenuates scopolamine-induced cognitive impairment in mice. *Biol. Pharm. Bull.* **37(7)**, 1207-1213 (2014).
5. Ferraz-Filha, Z.S., de Paula Michel Araújo, M.C., Ferrari, F.C., *et al.* *Tabebuia roseoalba*: In vivo hypouricemic and anti-inflammatory effects of its ethanolic extract and constituents. *Planta Med.* **82(16)**, 1395-1402 (2016).

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, 10/05/2022

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