

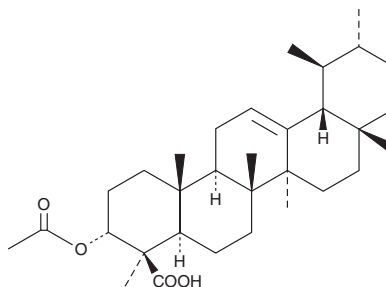
PRODUCT INFORMATION



β -acetyl-Boswellic Acid

Item No. 11674

CAS Registry No.: 5968-70-7
Formal Name: (3 α ,4 β)-3-(acetyloxy)-urs-12-en-23-oic acid
Synonyms: ABA, 3-O-acetyl- β -Boswellic acid, β -Boswellic acid acetate
MF: C₃₂H₅₀O₄
FW: 498.7
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years
Item Origin: Plant/*Boswellia serrata*



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

Laboratory Procedures

β -acetyl-Boswellic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the β -acetyl-boswellic acid in the solvent of choice. β -acetyl-Boswellic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of β -acetyl-boswellic acid in ethanol is approximately 5 mg/ml and approximately 25 mg/ml in DMSO and DMF.

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

Description

Boswellic acids are pentacyclic triterpenes first isolated from the resin of *Boswellia* plants. They have anti-inflammatory and anti-arthritis actions in mice.^{1,2} β -acetyl-Boswellic acid is a non-reducing inhibitor of 5-lipoxygenase (5-LO) that shows selectivity for 5-LO over 12-LO or COX.^{3,4} When studied as a 1:1 mixture with α -acetyl-boswellic acid, it inhibits topoisomerase II activity and arrests cell cycling or induces apoptosis in cancer cells.^{5,6}

References

1. Wang, Q., Pan, X., Wong, H.H., et al. *Osteoarthritis Cartilage* **22(1)**, 128-132 (2014).
2. Mostafa, D.M., Ammar, N.M., Basha, M., et al. *Drug Deliv.* **22(6)**, 748-756 (2015).
3. Safayhi, H., Mack, T., Sabieraj, J., et al. *J. Pharmacol. Exp. Ther.* **261**, 1143-1146 (1992).
4. Ammon, H.P., Safayhi, H., Mack, T., et al. *J. Ethnopharmacol.* **38(2-3)**, 113-119 (1993).
5. Zhao, W., Entschladen, F., Liu, H., et al. *Cancer Detect. Prev.* **27(1)**, 67-75 (2003).
6. Xia, L., Chen, D., Han, R., et al. *Mol. Cancer Ther.* **4(3)**, 381-388 (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, 06/18/2024

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