

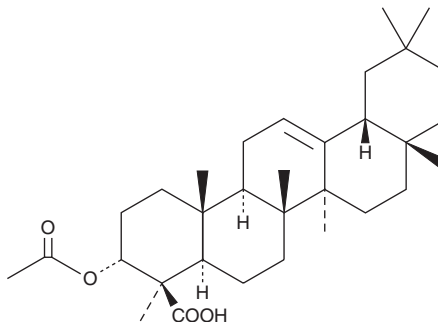
PRODUCT INFORMATION



α -acetyl Boswellic Acid

Item No. 11673

CAS Registry No.: 89913-60-0
Formal Name: (3 α ,4 β)-3-(acetyloxy)-olean-12-en-23-oic acid
Synonyms: ABA, 3-O-acetyl- α -Boswellic Acid, α -Boswellic Acid Acetate
MF: C₃₂H₅₀O₄
FW: 498.7
Purity: \geq 99%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Boswellia serrata*



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

Laboratory Procedures

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

Description

α -acetyl Boswellic acid is a pentacyclic triterpene and isomer of β -acetyl-boswellic acid (Item No. 11674) originally isolated from *B. serrata*.¹ It inhibits LPS-induced TNF- α mRNA expression in monocytes as well as NF- κ B activity in HEK293 cells when used at a concentration of 10 μ M.¹ α -acetyl Boswellic acid also inhibits phosphorylation of I κ B α mediated by I κ B kinases (IKKs) in monocytes.

Reference

1. Syrovets, T., Büchele, B., Krauss, C., *et al.* Acetyl-boswellic acids inhibit lipopolysaccharide-mediated TNF- α induction in monocytes by direct interaction with I κ B kinases. *J. Immunol.* **174**(1), 498-506 (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, 01/03/2023

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