

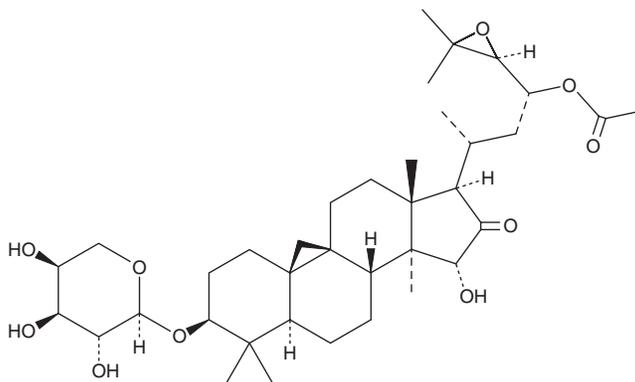
PRODUCT INFORMATION



Acetylcimigenol Arabinoside

Item No. 27596

CAS Registry No.: 402513-88-6
Formal Name: (3 β ,15 α ,23R,24S)-23-(acetyloxy)-3-(α -L-arabinopyranosyloxy)-24,25-epoxy-15-hydroxy-9,19-cyclolanostan-16-one
MF: C₃₇H₅₈O₁₀
FW: 662.9
Purity: \geq 98%
UV/Vis.: λ_{max} : 318 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Dioscorea opposita* Thunb.



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

Laboratory Procedures

Acetylcimigenol arabinoside is supplied as a solid. A stock solution may be made by dissolving the acetylcimigenol arabinoside in the solvent of choice, which should be purged with an inert gas. Acetylcimigenol arabinoside is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of acetylcimigenol arabinoside in these solvents is approximately 15 and 5 mg/ml, respectively.

Acetylcimigenol arabinoside is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, acetylcimigenol arabinoside should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Acetylcimigenol arabinoside 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

Acetylcimigenol arabinoside is a triterpene glycoside originally isolated from *Cimicifuga*.¹

Reference

1. Nian, Y., Zhang, X.-M., Li, Y., *et al.* Cycloartane triterpenoids from the aerial parts of *Cimicifuga foetida* Linnaeus. *Phytochemistry* **72**(11-12), 1473-1481 (2011).

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, 12/02/2022

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