

PRODUCT INFORMATION



Cimigenol-3-O- α -L-arabinoside

Item No. 42612

CAS Registry No.: 256925-92-5

Formal Name: (3 β ,15 α ,16 α ,23R,24S)-16,23:16,24-diepoxy-15,25-dihydroxy-9,19-cyclolanostan-3-yl α -L-arabinopyranoside

Synonyms: Cimicifugoside M, Cimiracemoside C, Cimigenol 3-O- α -L-arabinopyranoside

MF: C₃₅H₅₆O₉

FW: 620.8

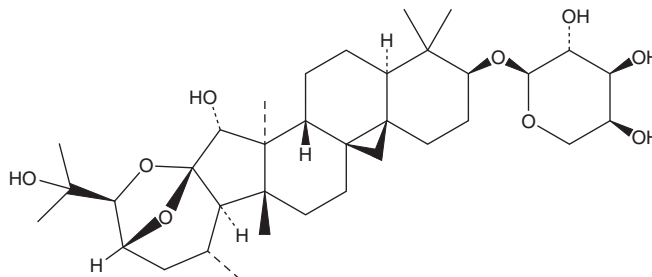
Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Cimicifugae rhizoma*



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

Laboratory Procedures

Cimigenol-3-O- α -L-arabinoside is supplied as a solid. A stock solution may be made by dissolving the cimigenol-3-O- α -L-arabinoside in the solvent of choice, which should be purged with an inert gas. Cimigenol-3-O- α -L-arabinoside is soluble in the organic solvent methanol.

Description

Cimigenol-3-O- α -L-arabinoside is a cycloartane glycoside that has been found in *C. racemosa*.¹

Reference

1. Watanabe, K., Mimaki, Y., Sakagami, H., *et al.* Cycloartane glycosides from the rhizomes of *Cimicifuga racemosa* and their cytotoxic activities. *Chem. Pharm. Bull. (Tokyo)* **50(1)**, 121-125 (2002).

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, 05/12/2026

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