

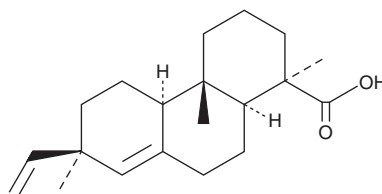
PRODUCT INFORMATION



Continentalic Acid

Item No. 35240

CAS Registry No.: 19889-23-7
Formal Name: (1R,4aS,4bR,7R,10aS)-7-ethenyl-1,2,3,4,4a,4b,5,6,7,9,10,10a-dodecahydro-1,4a,7-trimethyl-1-phenanthrenecarboxylic acid
Synonym: Pimaradienoic Acid
MF: C₂₀H₃₀O₂
FW: 302.5
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Aralia continentalis*



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

Laboratory Procedures

Continentalic acid is supplied as a solid. A stock solution may be made by dissolving the continentalic acid in the solvent of choice, which should be purged with an inert gas. Continentalic acid is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

Continentalic acid is a diterpene acid that has been found in *A. continentalis* and has diverse biological activities.¹⁻⁴ It is active against methicillin-susceptible *S. aureus* (MSSA) and clinical isolates of methicillin-resistant *S. aureus* (MRSA; MICs = 8 and 8-16 µg/ml, respectively).¹ Continentalic acid (200 µM) induces apoptosis in U-2932, OCI-LY1, and Ramos (RA 1) B cell lymphoma cells.² It also improves renal function and decreases renal malondialdehyde (MDA) and peroxidase levels, nitric oxide (NO) production, and neutrophil infiltration in a mouse model of *E. coli*- and LPS-induced renal injury.³

References

1. Jeong, S.-I., Han, W.-S., Yun, Y.-H., *et al.* Continentalic acid from *Aralia continentalis* shows activity against methicillin-resistant *Staphylococcus aureus*. *Phytother. Res.* **20(6)**, 511-514 (2006).
2. Jeon, B.-E., Kwon, C.-S., Lee, J.-E., *et al.* Anticancer activity of continentalic acid in B-cell lymphoma. *Molecules* **26(22)**, 6845 (2021).
3. Khan, A.M., Khan, A.U., Ali, H., *et al.* Continentalic acid exhibited nephroprotective activity against the LPS and *E. coli*-induced kidney injury through inhibition of the oxidative stress and inflammation. *Int. Immunopharmacol.* **80**, 106209 (2020).

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, 02/07/2024

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