

PRODUCT INFORMATION



Ardisiacrispin A

Item No. 34615

CAS Registry No.: 23643-61-0
Formal Name: (3 β ,16 α ,20 β)-13,28-epoxy-3-[[O- β -D-glucopyranosyl-(1 \rightarrow 2)-O-[O- β -D-xylopyranosyl-(1 \rightarrow 2)- β -D-glucopyranosyl-(1 \rightarrow 4)]- α -L-arabinopyranosyl]oxy]-16-hydroxy-oleanan-29-al

Synonyms: Deglucocyclamin, LTS-4, Saxifragifolin B

MF: C₅₂H₈₄O₂₂

FW: 1,061.2

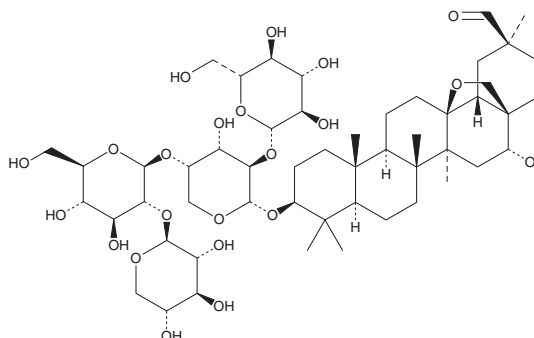
Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/Unspecified sp.



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

Laboratory Procedures

Ardisiacrispin A is supplied as a solid. A stock solution may be made by dissolving the ardisiacrispin A in the solvent of choice, which should be purged with an inert gas. Ardisiacrispin A is soluble in the organic solvent ethanol.

Description

Ardisiacrispin A is a triterpene saponin that has been found in *Myrsine australis* and has diverse biological activities.¹⁻³ It inhibits the cytopathic effects of poliovirus type I and herpes simplex virus 1 (HSV-1) in the filter paper disk assay when used at a concentration of 40 μ g/disk and is active against various *Candida* species (MICs = 80-160 μ g/ml).^{1,2} Ardisiacrispin A (10 and 13 μ M) induces apoptosis in HepG2 cells.³ It induces contractions in rat uterus preparations when used at a concentration of 7.5 μ M.²

References

1. Bloor, S.J. and Qi, L. Cytotoxic saponins from New Zealand *Myrsine* species. *J. Nat. Prod.* **57**(10), 1354-1360 (1994).
2. Çaliş, İ., Şatana, M.E., Yürüker, A., et al. Triterpene saponins from *Cyclamen mirabile* and their biological activities. *J. Nat. Prod.* **60**(3), 315-318 (1997).
3. Zhang, D.-M., Wang, Y., Tang, M.-K., et al. Saxifragifolin B from *Androsace umbellata* induced apoptosis on human hepatoma cells. *Biochem. Biophys. Res. Commun.* **362**(3), 759-765 (2007).

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, 10/06/2022

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