

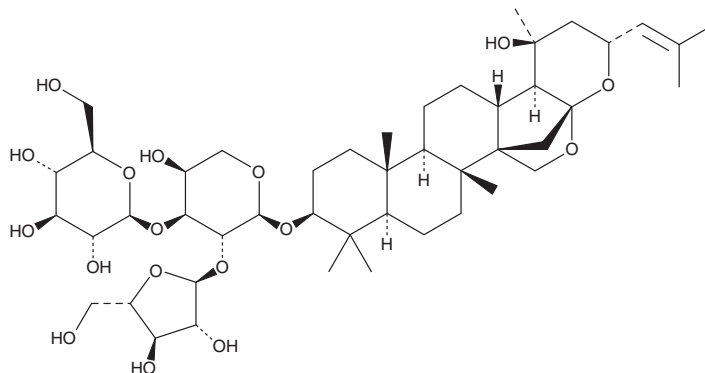
PRODUCT INFORMATION



Bacopaside X

Item No. 11851

CAS Registry No.: 94443-88-6
Formal Name: (3 β ,16 β ,23R)-16,23:16,30-diepoxy-20-hydroxydammar-24-en-3-yl O- α -L-arabinofuranosyl-(1 \rightarrow 2)-O- $[\beta$ -D-glucopyranosyl-(1 \rightarrow 3)]- α -L-arabinopyranoside
Synonyms: Bacopaside VII, Jujubogenin isomer of Bacopasaponin C
MF: C₄₆H₇₄O₁₇
FW: 899.1
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Bacopa monnieri*



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

Laboratory Procedures

Bacopaside X is supplied as a solid. A stock solution may be made by dissolving the bacopaside X in the solvent of choice, which should be purged with an inert gas. Bacopaside X is soluble in DMSO.

Description

Bacopaside X is a triterpenoid saponin that has been found in *B. monniera* and has antifungal and anticancer activities.¹⁻³ It inhibits the growth of the fungi *C. albicans*, *C. neoformans*, and *A. fumigatus* (MIC = 50 μ g/ml for all).² Bacopaside X decreases the proliferation of MDA-MB-231, SHG44, HCT8, A549, and PC3M cancer cell lines with IC₅₀ values of 14.3, 15.9, 9.8, 9.7, and 10.1 μ M, respectively.³ It reduces tumor growth by 84.13% in an S180 sarcoma mouse allograft model when administered at a dose of 50 μ mol/kg.

References

1. Zhou, Y., Shen, Y.-H., Zhang, C., et al. Triterpene saponins from *Bacopa monnieri* and their antidepressant effects in two mice models. *J. Nat. Prod.* **70**(4), 652-655 (2007).
2. Li, X.-C., ElSohly, H.N., Nimrod, A.C., et al. Antifungal jujubogenin saponins from *Colubrina retusa*. *J. Nat. Prod.* **62**(5), 674-677 (1999).
3. Peng, L., Zhou, Y., Kong, D.Y., et al. Antitumor activities of dammarane triterpene saponins from *Bacopa monniera*. *Phytother. Res.* **24**(6), 864-868 (2010).

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, 10/03/2022

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