

PRODUCT INFORMATION



Hederacoside D

Item No. 34202

CAS Registry No.: 760961-03-3
Formal Name: (3 β ,4 α)-3-(α -L-arabinopyranosyloxy)-23-hydroxy-olean-12-en-28-oic acid, O-6-deoxy- α -L-mannopyranosyl-(1 \rightarrow 4)-O- β -D-glucopyranosyl-(1 \rightarrow 6)- β -D-glucopyranosyl ester

MF: C₅₃H₈₆O₂₂

FW: 1,075.2

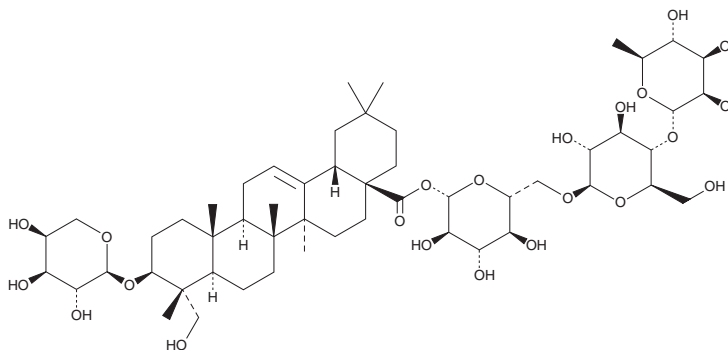
Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Hedera*



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

Laboratory Procedures

Hederacoside D is supplied as a solid. A stock solution may be made by dissolving the hederacoside D in the solvent of choice, which should be purged with an inert gas. Hederacoside D is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of hederacoside D in ethanol is approximately 1 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of hederacoside D can be prepared by directly dissolving the solid in aqueous buffers. The solubility of hederacoside D in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Hederacoside D is a triterpene saponin that has been found in *S. hexaphylla*.¹

Reference

1. Vinh, L.B., Jang, H.-J., Phong, N.V., *et al.* Isolation, structural elucidation, and insights into the anti-inflammatory effects of triterpene saponins from the leaves of *Stauntonia hexaphylla*. *Bioorg. Med. Chem. Lett.* **29**(8), 965-969 (2019).

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, 11/18/2022

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