

PRODUCT INFORMATION



Pulsatilla Saponin D

Item No. 29455

CAS Registry No.: 68027-15-6

Formal Name: (3 β ,4 α)-3-[[O-6-deoxy- α -L-mannopyranosyl-(1 \rightarrow 2)-O- β -D-glucopyranosyl-(1 \rightarrow 4)]- α -L-arabinopyranosyl]oxy]-23-hydroxyolean-12-en-28-oic acid

Synonyms: Hederagenin 3-O- α -L-rhamnopyranosyl-(1 \rightarrow 2)- β -D-glucopyranosyl-(1 \rightarrow 4)]- α -L-arabinopyranoside, Hederacolchiside A, Pulchinoside B3

MF: C₄₇H₇₆O₁₇

FW: 913.1

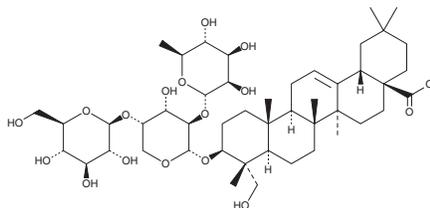
Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Pulsatilla chinensis*



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

Laboratory Procedures

Pulsatilla saponin D is supplied as a solid. A stock solution may be made by dissolving the pulsatilla saponin D in the solvent of choice, which should be purged with an inert gas. Pulsatilla saponin D is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pulsatilla saponin D in these solvents is approximately 10, 5, and 15 mg/ml, respectively.

Pulsatilla saponin D is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pulsatilla saponin D should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Pulsatilla saponin D has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Pulsatilla saponin D is an oleanane-type saponin that has been found in *P. koreana* and has anticancer activity.¹ It is cytotoxic to A549, SKOV3, SK-MEL-2, and HCT15 cells (EC₅₀s = 11.25, 13.17, 3.04, and 11.86 μ M, respectively). Pulsatilla saponin D (6 mg/kg per day) reduces tumor volume in a murine Lewis lung carcinoma model.

Reference

1. Bang, S.-C., Lee, J.-H., Song, G.-Y., *et al.* Antitumor activity of *Pulsatilla koreana* saponins and their structure-activity relationship. *Chem. Pharm. Bull. (Tokyo)*. **53(11)**, 1451-1454 (2005).

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, 11/04/2022

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