

PRODUCT INFORMATION



Pulchinenoside A

Item No. 28545

CAS Registry No.: 129724-84-1
Formal Name: (3 β ,4 α)-3-[[2-O-(6-deoxy- α -L-mannopyranosyl)- α -L-arabinopyranosyl]oxy]-23-hydroxy-lup-20(29)-en-28-oic acid

Synonym: Anemoside A₃

MF: C₄₁H₆₆O₁₂

FW: 751.0

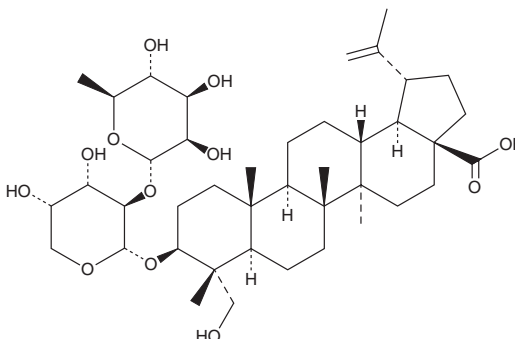
Purity: \geq 90%

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

Pulchinenoside A is supplied as a solid. A stock solution may be made by dissolving the pulchinenoside A in the solvent of choice, which should be purged with an inert gas. Pulchinenoside A is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of pulchinenoside A in these solvents is approximately 5 mg/ml.

Pulchinenoside A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pulchinenoside A should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Pulchinenoside A has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Pulchinenoside A is a triterpenoid saponin that has been found in *Pulsatilla chinensis* and has diverse biological activities.^{1,2} It inhibits NMDA-evoked currents in hippocampal neurons (IC₅₀s = 18.48 and 20.19 μ M at -50 and +50 mV, respectively) and inhibits NMDA-induced neuronal cell death when used at a concentration of 30 μ M.² Pulchinenoside A (30 and 100 mg/kg) increases the levels of serotonin (5-HT; Item No. 14332), 5-hydroxyindolacetic acid (5-HIAA; Item No. 22889), and norepinephrine (Item No. 16673) in the hippocampus and 5-HIAA and norepinephrine in the prefrontal cortex in mice. It reduces the time spent immobile in the forced swim test in mice when administered at doses of 30 and 100 mg/kg and increases the time spent in the target quadrant of the Morris water maze in mice at 100 mg/kg. Pulchinenoside A (12.6 and 25.2 mg/kg) reduces infarct volume in a middle cerebral artery occlusion (MCAO) rat model of ischemia. It also induces relaxation of precontracted isolated rat thoracic aorta and renal, mesenteric, and left coronary arteries in a concentration-dependent manner.¹

References

1. Zhang, D.-M., Lin, S.-M., Lau, C.-W., *et al.* Anemoside A₃-induced relaxation in rat renal arteries: Role of endothelium and Ca²⁺ channel inhibition. *Planta Med.* **76(16)**, 1814-1819 (2010).
2. Ip, F.C.F., Fu, W.-Y., Cheng, E.Y.L., *et al.* Anemoside A₃ enhances cognition through the regulation of synaptic function and neuroprotection. *Neuropsychopharmacology* **40(8)**, 1877-1887 (2015).

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, 12/05/2022

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