

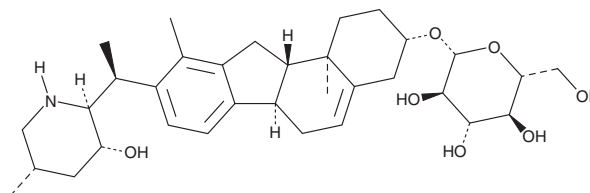
PRODUCT INFORMATION



Veratrosine

Item No. 34617

CAS Registry No.: 475-00-3
Formal Name: (3 β ,23 β)-14,15,16,17-tetradehydro-23-hydroxyveratraman-3-yl, β -D-glucopyranoside
Synonym: Veratramine 3-O-glucoside
MF: C₃₃H₄₉NO₇
FW: 571.7
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

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

Description

Veratrosine is a steroidal alkaloid that has been found in *Veratrum* and has diverse biological activities.^{1,2} It inhibits hedgehog (Hh) signaling in a reporter assay using NIH3T3 cells when used at a concentration of 8 μ M.¹ Veratrosine (1 mg/ml) inhibits aggregation of isolated rabbit platelets induced by arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607). It induces DNA damage in the cerebellum and cerebral cortex in mice when administered at doses of 0.25 and 2.5 μ mol/kg.² Veratrosine (1.5 and 4.5 μ mol/kg) decreases angiotensin II-induced increases in systolic blood pressure in mice.

References

1. Tang, J., Li, H.L., Shen, Y.H., *et al.* Antitumor and antiplatelet activity of alkaloids from *Veratrum dahuricum*. *Phytother. Res.* **24**(6), 821-826 (2010).
2. Cong, Y., Wu, Y., Shen, S., *et al.* A structure-activity relationship between the *Veratrum* alkaloids on the antihypertension and DNA damage activity in mice. *Chem. Biodivers.* **17**(2), e1900473 (2020).

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, 10/04/2022

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