

PRODUCT INFORMATION



Crassicauline A

Item No. 34674

CAS Registry No.: 79592-91-9
Formal Name: (1 α ,6 α ,14 α ,16 β)-20-ethyl-1,6,16-trimethoxy-4-(methoxymethyl)-aconitane-8,13,14-triol, 8-acetate 14-(4-methoxybenzoate)

Synonym: 3-Deoxyyunaconitine

MF: C₃₅H₄₉NO₁₀

FW: 643.8

Purity: \geq 98%

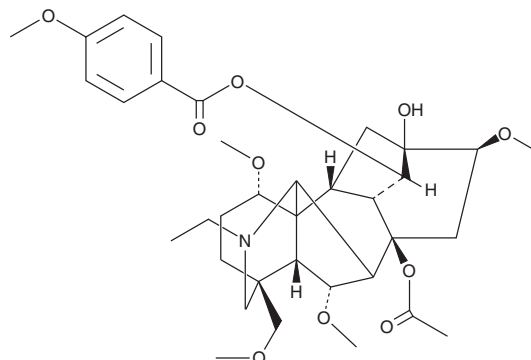
UV/Vis.: λ_{max} : 258 nm

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Aconitum transsectum*



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

Laboratory Procedures

Crassicauline A is supplied as a solid. A stock solution may be made by dissolving the crassicauline A in the solvent of choice, which should be purged with an inert gas. Crassicauline A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of crassicauline A in these solvents is approximately 5, 10, and 25 mg/ml, respectively.

Crassicauline A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, crassicauline A should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Crassicauline A has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Crassicauline A is a diterpene alkaloid originally isolated from *A. crassicaule* that has analgesic activity.^{1,2} It inhibits acetic acid-induced writhing in mice (ED_{50} = 0.048 mg/kg) and induces lethality in mice when administered at doses greater than 0.92 mg/kg.²

References

1. Fengpeng, W. and Qicheng, F. Alkaloids from roots of *Aconitum crassicaule*. *Planta Med.* **42(8)**, 375-379 (1981).
2. Wang, J.-L., Shen, X.-L., Chen, Q.-H., et al. Structure-analgesic activity relationship studies on the C₁₈- and C₁₉-diterpenoid alkaloids. *Chem. Pharmaceut. Bull.* **57(8)**, 801-807 (2009).

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, 11/22/2022

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