

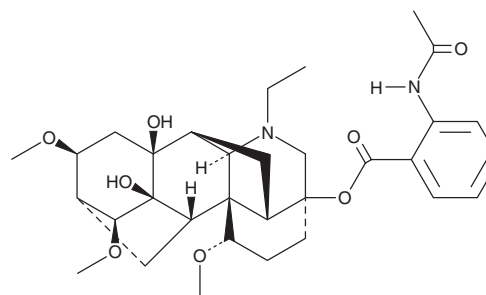
PRODUCT INFORMATION



Lappaconitine

Item No. 28747

CAS Registry No.: 32854-75-4
Formal Name: 20-ethyl-1 α ,14 α ,16 β -trimethoxy-aconitane-4,8,9-triol 4-[2-(acetylamino)benzoate]
MF: C₃₂H₄₄N₂O₈
FW: 584.7
Purity: \geq 90%
UV/Vis.: λ_{max} : 223, 253, 310 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/Aconitum sinomontanum



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

Laboratory Procedures

Lappaconitine is supplied as a solid. A stock solution may be made by dissolving the lappaconitine in the solvent of choice, which should be purged with an inert gas. Lappaconitine is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

Lappaconitine is a diterpene alkaloid that has been found in *A. sinomontanum* and has diverse biological activities.¹⁻³ It is an inhibitor of voltage-gated sodium channel 1.7 (Na_v1.7; IC₅₀ = 27.67 μ M for the human channel).² Lappaconitine has antiarrhythmic activity in a rat model of aconitine-induced arrhythmia with an ED₅₀ value of 0.05 μ g/kg.³ It increases the mechanical and thermal paw withdrawal thresholds in rats in a model of spinal nerve ligation-induced neuropathic pain (ED₅₀s = 1.1 and 1.6 mg/kg, respectively).¹

References

1. Sun, M.L., Ao, J.-P., Wang, Y.-R., *et al.* Lappaconitine, a C18-diterpenoid alkaloid, exhibits antihypersensitivity in chronic pain through stimulation of spinal dynorphin A expression. *Psychopharmacology (Berl.)* **235(9)**, 2559-2571 (2018).
2. Li, Y.-F., Zheng, Y.-M., Yu, Y., *et al.* Inhibitory effects of lappaconitine on the neuronal isoforms of voltage-gated sodium channels. *Acta Pharmacol. Sin.* **40(4)**, 451-459 (2019).
3. Dzhakhgairov, F.N., Sultankhodzhaev, M.N., Tashkhodzhaev, B., *et al.* Diterpenoid alkaloids as a new class of antiarrhythmic agents. Structure-activity relationship. *Chem. Nat. Compd.* **33(2)**, 190-202 (1997).

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, 10/10/2022

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