

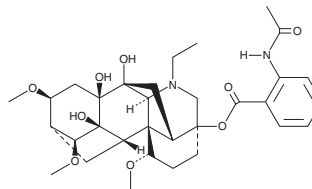
PRODUCT INFORMATION



Ranaconitine

Item No. 34774

CAS Registry No.: 1360-76-5
Formal Name: 20-ethyl-1 α ,14 α ,16 β -trimethoxy-aconitane-4,7,8,9-tetrol 4-[2-(acetylamino)benzoate]
Synonym: 7-hydroxy Lappaconitine
MF: C₃₂H₄₄N₂O₉
FW: 600.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 223, 252 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Aconitum carmichaelii*



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

Laboratory Procedures

Ranaconitine is supplied as a solid. A stock solution may be made by dissolving the ranaconitine in the solvent of choice, which should be purged with an inert gas. Ranaconitine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ranaconitine in these solvents is approximately 1, 10, and 30 mg/ml, respectively.

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

Description

Ranaconitine is an alkaloid that has been found in *Aconitum* and has analgesic and local anesthetic activities in mice (ED₅₀s = 6.99 and 0.167 μ mol/kg, respectively).¹

Reference

1. Bello-Ramírez, A.M. and Nava-Ocampo, A.A. A QSAR analysis of toxicity of *Aconitum* alkaloids. *Fundam. Clin. Pharmacol.* **18**(6), 699-704 (2004).

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, 12/08/2022

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