

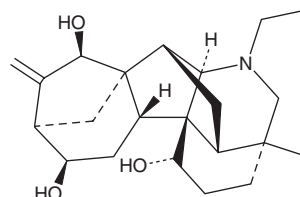
PRODUCT INFORMATION



12-Epinapelline

Item No. 36550

CAS Registry No.: 110064-71-6
Formal Name: (3R,6S,6aS,6bR,8R,9R,11R,11aR,12R,12aS,14R)-1-ethyldodecahydro-3-methyl-10-methylene-12,3,6a-ethanylylidene-9,11a-methanoazuleno[2,1-b]azocine-6,8,11-triol
Synonym: 12-*epi* Napelline
MF: C₂₂H₃₃NO₃
FW: 359.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Aconitum carmichaeli*



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

Laboratory Procedures

12-Epinapelline is supplied as a solid. A stock solution may be made by dissolving the 12-epinapelline in the solvent of choice, which should be purged with an inert gas. 12-Epinapelline is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

12-Epinapelline is a diterpenoid alkaloid that has been found in *Aconitum* and has antiarrhythmic activity.¹ It has antiarrhythmic activity in a rat model of aconitine-induced arrhythmia and antifibrillatory activity in a mouse model of aconitine-induced cardiac fibrillation (ED₅₀s = 8 and 20 mg/kg, respectively).

Reference

1. Dzhakhangirov, 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, 12/20/2022

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