

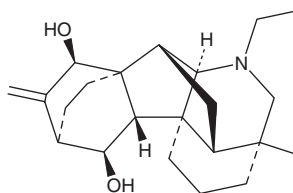
PRODUCT INFORMATION



Denudatine

Item No. 30904

CAS Registry No.: 26166-37-0
Formal Name: (3R,6aS,6bS,7S,8R,10R,10aS,11R,11aR,13R)-1-ethyl-dodecahydro-3-methyl-9-methylene-8,10a-ethano-11,3,6a-ethanylylidene-8H-indeno[2,1-b]azocine-7,10-diol



MF: C₂₂H₃₃NO₂
FW: 343.5
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥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

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

Description

Denudatine is an alkaloid that has been found in *A. jinyangense* and has antiarrhythmic activity.¹ *In vivo*, denudatine (50 and 100 mg/kg) prevents aconitine-induced arrhythmias in rats and increases survival in a rat model of calcium chloride-induced arrhythmias.

Reference

1. Jin, L.S., Zhou, Y.P., and Zeng, G.Y. Effects of denudatine on experimental arrhythmia and heart function. *Zhongguo Yao Li Xue Bao* **3(2)**, 104-108 (1982).

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, 10/24/2022

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