

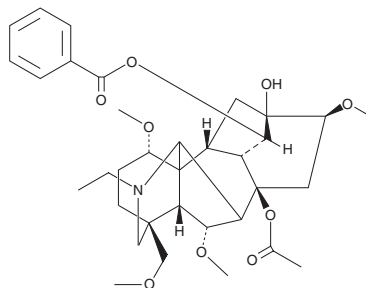
PRODUCT INFORMATION



3-Deoxyaconitine

Item No. 36151

CAS Registry No.: 3175-95-9
Formal Name: 20-ethyl-1 α ,6 α ,16 β -trimethoxy-4-(methoxymethyl)-aconitane-8,13,14 α ,15 α -tetrol, 8-acetate 14-benzoate
Synonym: Deoxyaconitine
MF: C₃₄H₄₇NO₁₀
FW: 629.7
Purity: $\geq$ 98%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years
Item Origin: Plant/*Aconitum carmichaelii* Debx



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

Laboratory Procedures

3-Deoxyaconitine is supplied as a solid. A stock solution may be made by dissolving the 3-deoxyaconitine in the solvent of choice, which should be purged with an inert gas. 3-Deoxyaconitine is soluble in DMSO.

Description

3-Deoxyaconitine is a diterpene alkaloid that has been found in *A. kusnezoffii*.¹⁻³ It induces spasms in isolated rabbit small intestine (EC₅₀ = 100 nM).² 3-Deoxyaconitine (0.375 mg/kg) also enhances lidocaine-induced sensory loss, weight-bearing capacity, and partial paralysis in rabbits.³

References

1. Wang, Y., Song, F., Xu, Q., *et al.* Characterization of aconitine-type alkaloids in the flowers of *Aconitum kusnezoffii* by electrospray ionization tandem mass spectrometry. *J. Mass. Spectrom.* **38**(9), 962-970 (2003).
2. Dzhakhangirov, F.N., Tursunkhodzhaeva, F.M., Sultankhodzhaev, M.N., *et al.* Spasmolytic activity of diterpenoid alkaloids and their derivatives. Structure-activity relationship. *Chem. Nat. Compd.* **49**(4), 702-706 (2013).
3. Komoda, Y., Nosaka, S., and Takenoshita, M. Enhancement of lidocaine-induced epidural anesthesia by deoxyaconitine in the rabbit. *J. Anesth.* **17**(4), 241-245 (2003).

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, 03/14/2023

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