

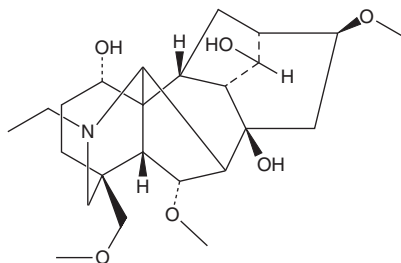
PRODUCT INFORMATION



Bullatine B

Item No. 36170

CAS Registry No.: 466-26-2
Formal Name: 20-ethyl-6,16-dimethoxy-4-(methoxymethyl)-(1 α ,6 α ,14 α ,16 β)-aconitane-1,8,14-triol
Synonym: Neoline
MF: C₂₄H₃₉NO₆
FW: 437.6
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/Aconitum



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

Laboratory Procedures

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

Description

Bullatine B is a diterpene alkaloid that has been found in *Aconitum* and has diverse biological activities.^{1,2} It inhibits voltage-gated sodium channel 1.7 (Na_v1.7) in HEK293 cells expressing the human channel (IC₅₀ = 25.8 nM) and scavenges DPPH (Item No. 14805) radicals by 65.3% at 1 mM. Bullatine B (10 mg/kg) increases the mechanical pain threshold in the von Frey test in streptozotocin-induced diabetic, but not non-diabetic, mice.²

References

1. Begum, S., Ali, A., Ahman, W., *et al.* Pharmacologically active C-19 diterpenoid alkaloids from the aerial parts of *Aconitum laeve* Royle. *Rec. Nat. Prod.* **8**(2), 83-92 (2014).
2. Nakatani, Y., Negoro, K., Yamauchi, M., *et al.* Neoline, an active ingredient of the processed aconite root in Goshajinkigan formulation, targets Na_v1.7 to ameliorate mechanical hyperalgesia in diabetic mice. *J. Ethnopharmacol.* **259**, 112963 (2020).

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, 12/19/2022

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