

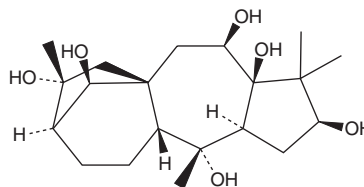
PRODUCT INFORMATION



Grayanotoxin III

Item No. 42645

CAS Registry No.: 4678-45-9
Formal Name: Grayanotoxane-3 β ,5,6 β ,10,14R,16-hexol
Synonyms: Andromedol, Deacetylandromedotoxin, Deacylasebotoxin I, GTX-III, NSC 323778
MF: C₂₀H₃₄O₆
FW: 370.5
Purity: $\geq$ 90%
Supplied as: A solution in water:acetonitrile
Storage: -20°C
Stability: $\geq$ 1 year



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

Description

Grayanotoxin III (GTX-III) is a diterpene toxin that has been found in *R. ponticum* nectar, a component of mad honey.¹ *In vivo*, GTX-III (0.1 and 0.25 mg/kg) induces salivation, emesis, and hind paw paralysis in mice.² It is lethal to mice (LD₅₀ = 0.84 mg/kg).

References

1. Kurtoglu, A.B., Yavuz, R., and Evrendilek, G.A. Characterisation and fate of grayanotoxins in mad honey produced from *Rhododendron ponticum* nectar. *Food Chem.* **161**, 47-52 (2014).
2. Ohgaki, T., Uchida, S., Meguri, H., *et al.* Preventive action of quisqualic acid against grayanotoxin-induced suppression of locomotor activity in mice. *Neuropharmacology* **27(10)**, 1045-1053 (1988).

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, 01/27/2025

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