

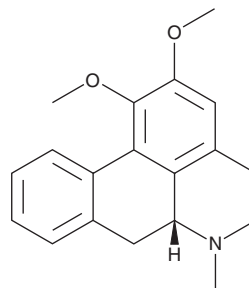
PRODUCT INFORMATION



Nuciferine

Item No. 34221

CAS Registry No.: 475-83-2
Formal Name: 5,6,6aR,7-tetrahydro-1,2-dimethoxy-6-methyl-4H-dibenzo[de,g]quinoline
Synonyms: (-)-Nuciferine, Sanjoinine E, VLT 049
MF: C₁₉H₂₁NO₂
FW: 295.4
Purity: ≥98%
UV/Vis.: λ_{max}: 272 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Nuciferine (Item No. 34221) is an analytical reference standard categorized as an alkaloid.¹ Nuciferine induces sedation, hypothermia, ptosis, and catalepsy in rodents. It has been found in refillable non-nicotinic vaping mixtures for electronic cigarettes.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 33839).

References

1. Bhattacharya, S.K., Bose, R., Ghosh, P., *et al.* Psychopharmacological studies on (-)-nuciferine and its Hofmann degradation product atherosperminine. *Psychopharmacology (Berl.)* **59(1)**, 29-33 (1978).
2. Holt, A.K., Poklis, J.L., and Peace, M.R. A retrospective analysis of chemical constituents in regulated and unregulated e-cigarette liquids. *Front. Chem.* **9**, 752342 (2021).

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, 06/28/2023

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