

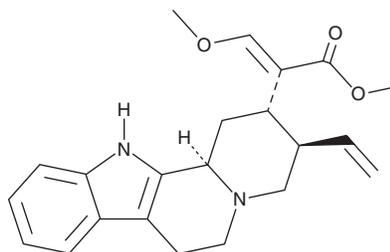
PRODUCT INFORMATION



Hirsuteine

Item No. 34939

CAS Registry No.: 35467-43-7
Formal Name: (αE,2S,3R,12bR)-3-ethenyl-1,2,3,4,6,7,12,12b-octahydro-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester
Synonym: 3-Epicorynantheine
MF: C₂₂H₂₆N₂O₃
FW: 366.5
Purity: ≥98%
UV/Vis.: λ_{max}: 226 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Uncaria rhynchophylla*



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

Laboratory Procedures

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

Description

Hirsuteine is an indole alkaloid that has been found in *U. sinensis* and has neuroprotective activities.^{1,2} It reduces glutamate-induced cytotoxicity in primary rat cerebellar granule neurons when used at concentrations ranging from 100 to 300 μM.¹ Hirsuteine (1-100 μM) inhibits nicotine-induced dopamine release in PC12 cells.²

References

1. Shimada, Y., Goto, H., Itoh, T., *et al.* Evaluation of the protective effects of alkaloids isolated from the hooks and stems of *Uncaria sinensis* on glutamate-induced neuronal death in cultured cerebellar granule cells from rats. *J. Pharm. Pharmacol.* **51(6)**, 715-722 (1999).
2. Watano, T., Nakazawa, K., Obama, T., *et al.* Non-competitive antagonism by hirsuteine of nicotinic receptor-mediated dopamine release from rat pheochromocytoma cells. *Jpn. J. Pharmacol.* **61(4)**, 351-356 (1993).

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, 10/14/2022

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