

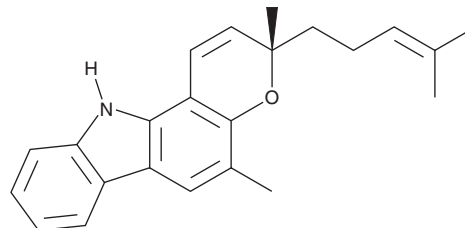
PRODUCT INFORMATION



Mahanimbine

Item No. 30897

CAS Registry No.: 21104-28-9
Formal Name: (3S)-3,11-dihydro-3,5-dimethyl-3-(4-methyl-3-penten-1-yl)-pyrano[3,2-a]carbazole
MF: C₂₃H₂₅NO
FW: 331.5
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 289 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Murraya koenigii*



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

Laboratory Procedures

Mahanimbine is supplied as a solid. A stock solution may be made by dissolving the mahanimbine in the solvent of choice, which should be purged with an inert gas. Mahanimbine is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

Mahanimbine is a carbazole alkaloid that has been found in *M. koenigii* and has diverse biological activities.¹⁻³ It is active against *S. aureus* and *S. pyogenes* (MIC₁₀₀ = 50 µg/ml for both).¹ Mahanimbine induces cell cycle arrest at the G₀/G₁ phase and apoptosis in Capan-2 and SW1190 cancer cells when used at a concentration of 7 µM, as well as inhibits proliferation in Capan-2, SW1190, BxPC-3, CFPAC-1, and HPAF-II cancer cells (IC₅₀s = 3.5, 3.5, 16, 64, and 32 µM, respectively).² It decreases body weight gain without affecting food intake and reduces increases in plasma total cholesterol and triglyceride levels in a rat model of high-fat diet-induced obesity when administered orally at a dose of 30 mg/kg.³

References

1. Ramsewak, R.S., Nair, M.G., Strasburg, G.M., *et al.* Biologically active carbazole alkaloids from *Murraya koenigii*. *J. Agric. Food Chem.* **47**(2), 444-447 (1999).
2. Pei, C., He, Q., Liang, S., *et al.* Mahanimbine exerts anticancer effects on human pancreatic cancer cells by triggering cell cycle arrest, apoptosis, and modulation of AKT/mammalian target of rapamycin (mTOR) and signal transducer and activator of transcription 3 (STAT3) signalling pathways. *Med. Sci. Monit.* **24**, 6975-6983 (2018).
3. Birari, R., Javia, V., and Bhutani, K.K. Antiobesity and lipid lowering effects of *Murraya koenigii* (L.) spreng leaves extracts and mahanimbine on high fat diet induced obese rats. *Fitoterapia* **81**(8), 1129-1133 (2010).

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, 08/25/2023

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