

PRODUCT INFORMATION



Vobtusine

Item No. 30057

CAS Registry No.: 19772-79-3
Formal Name: (2S,2aS,3'aR,4'aR,5aR,12bR,12'bR,12'cR,15aR,15'aS,17'aS)-2',3',4,4'a,5,6,8,12'c,13,13',14,14',17',17'a-tetradecahydro-12'c-hydroxy-9'-methoxy-spiro[1H,15aH-furo[2',3':7,8]indolizino[8,1-cd]carbazole-2(2aH),6'(7'H)-[4H,5H,15aH,16H]furo[2',3':7,8]indolizino[8,1-cd]pyrido[1,2,3-lm]carbazole]-7-carboxylic acid, methyl ester

Synonyms: NSC 180540, Vobtusin

MF: C₄₃H₅₀N₄O₆

FW: 718.9

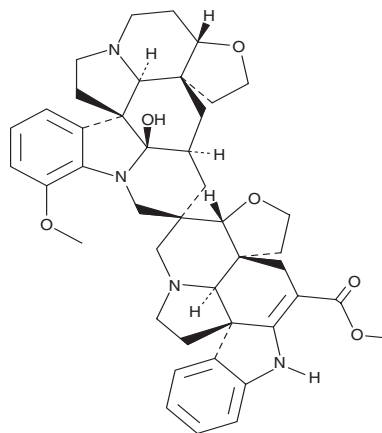
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Voacanga papuana*



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

Laboratory Procedures

Vobtusine is supplied as a solid. A stock solution may be made by dissolving the vobtusine in the solvent of choice, which should be purged with an inert gas. Vobtusine is soluble in ethanol, methanol, DMSO, and dimethylformamide.

Description

Vobtusine is an alkaloid that has been found in *V. africana* leaves.¹

Reference

1. Korocho, A.R., Juliani, H.R., Kulakowski, D., et al. *Voacanga africana*: Chemistry, quality and pharmacological activity. *African Natural Plant Products: New Discoveries and Challenges in Chemistry and Quality*. Juliani, H.R., Simon, J.E., and Ho, C.-T., editors, American Chemical Society (2009).

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, 01/04/2023

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