

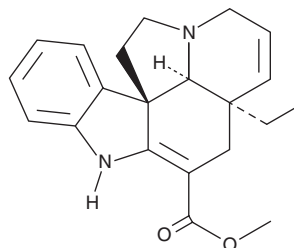
PRODUCT INFORMATION



Tabersonine

Item No. 26848

CAS Registry No.: 4429-63-4
Formal Name: (5 α ,12R,19 α)-2,3,6,7-tetradehydro-
aspidospermidine-3-carboxylic acid, methyl ester
Synonym: (-)-Tabersonine
MF: C₂₁H₂₄N₂O₂
FW: 336.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 226, 296, 329 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Malus* nut



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

Laboratory Procedures

Tabersonine is supplied as a crystalline solid. A stock solution may be made by dissolving the tabersonine in the solvent of choice, which should be purged with an inert gas. Tabersonine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of tabersonine in these solvents is approximately 2 mg/ml. Tabersonine is slightly soluble in ethanol.

Description

Tabersonine is an indole alkaloid that has been found in the beans of *Voacanga*.¹ It inhibits amyloid- β (1-42) (A β 42) fibril formation and disintegrates preformed A β 42 fibrils *in vitro* when used at a concentration of 10 μ M. Tabersonine is an intermediate in the biosynthesis of vindoline, a precursor in the synthesis of anticancer agents.²

References

1. Kai, T., Zhang, L., Wang, X., *et al.* Tabersonine inhibits amyloid fibril formation and cytotoxicity of A β (1-42). *ACS Chem. Neurosci.* **6**(6), 879-888 (2015).
2. Qu, Y., Easson, M.L.A.E., Froese, J., *et al.* Completion of the seven-step pathway from tabersonine to the anticancer drug precursor vindoline and its assembly in yeast. *Proc. Natl. Acad. Sci. U.S.A.* **112**(19), 6224-6229 (2015).

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, 11/16/2022

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