

PRODUCT INFORMATION



Psychotridine

Item No. 31645

CAS Registry No.: 52617-25-1
Formal Name: 2,2',2'',2''',2''''',3,3',3'',3''',3''''',8,8',8'',8''',8''''',8a,8'a,8''a,8''''a,8''''''a-eicosahydro-1,1',1'',1''',1''''-pentamethyl-3a(1H),7':3'a,3''a(H,1''H):7'',3''''a(1''''H):7''''',3''''''a(1''''''H)-quinquepyrrolo[2,3-b]indole, stereoisomer

MF: C₅₅H₆₂N₁₀

FW: 863.2

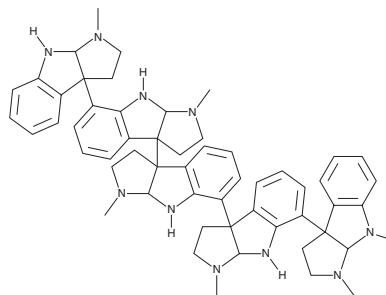
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Psychotria beccarioides*



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

Laboratory Procedures

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

Description

Psychotridine is an alkaloid that has been found in *P. forsteriana* and has diverse biological activities.¹⁻³ It inhibits ADP-, collagen-, or thrombin-induced aggregation of washed isolated human platelets with IC₅₀ values of 1.4, 1.4, and 3.9 μM, respectively.¹ Psychotridine (2.5 or 5 μM) is cytotoxic to HTC rat hepatocellular carcinoma cells.² It reduces paw licking induced by capsaicin (Item Nos. 92350 | 10010743) in mice when administered at doses of 0.5, 2.5, or 5 mg/kg.³

References

1. Beretz, A., Roth-Georger, A., Corre, G., *et al.* Polyindolinic alkaloids from *Psychotria forsteriana*. Potent inhibitors of the aggregation of human platelets. *Planta Med.* **51**(4), 300-303 (1985).
2. Roth, A., Kuballa, B., Bounthanh, C., *et al.* Cytotoxic activity of polyindoline alkaloids of *Psychotria forsteriana* (Rubiaceae) (1). *Planta Med.* **6**, 450-453 (1986).
3. Amador, T.A., Verotta, L., Nunes, D.S., *et al.* Involvement of NMDA receptors in the analgesic properties of psychotridine. *Phytomedicine* **8**(3), 202-206 (2001).

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, 12/08/2022

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