

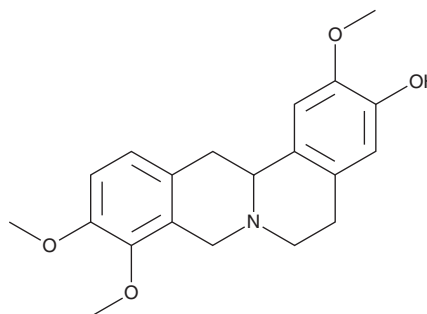
PRODUCT INFORMATION



Corypalmine

Item No. 42467

CAS Registry No.: 27313-86-6
Formal Name: 5,8,13,13a-tetrahydro-2,9,10-trimethoxy-6H-dibenzo[a,g]quinolizin-3-ol
Synonym: Tetrahydrojatrorrhizine
MF: $C_{20}H_{23}NO_4$
FW: 341.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Mahoniae Caulis*



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

Laboratory Procedures

Corypalmine is supplied as a solid. A stock solution may be made by dissolving the corypalmine in the solvent of choice, which should be purged with an inert gas. Corypalmine is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Corypalmine is an isoquinoline alkaloid that has been found in *F. recisa* and has receptor antagonist and antifungal activities.^{1,2} It is an antagonist of dopamine D_1 , D_2 , D_3 , and D_5 receptors (K_i s = 199, 42.4, 262, and 313 nM, respectively).¹ Corypalmine is active against *C. albicans*, *C. glabrata*, *C. krusei*, *C. parapsilosis*, and *C. neoformans* (MICs = 160, 160, 320, 160, and 80 $\mu\text{g/ml}$, respectively).²

References

1. Lee, D.Y.W., Liu, J., Zhang, S., et al. Asymmetric total synthesis of tetrahydroprotoberberine derivatives and evaluation of their binding affinities at dopamine receptors. *Bioorg. Med. Chem. Lett.* **27**(6), 1437-1440 (2017).
2. Rao, G.-X., Zhang, S., Wang, H.-M., et al. Antifungal alkaloids from the fresh rattan stem of *Fibraurea recisa* Pierre. *J. Ethnopharmacol.* **123**(1), 1-5 (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, 01/22/2025

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