

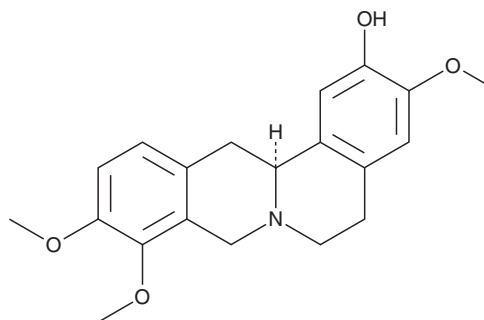
PRODUCT INFORMATION



(-)-Isocorypalmine

Item No. 29886

CAS Registry No.: 483-34-1
Formal Name: (13aS)-5,8,13,13a-tetrahydro-3,9,10-trimethoxy-6H-dibenzo[a,g]quinolizin-2-ol
Synonyms: L-Isocorypalmine, Isocorypalmine
MF: C₂₀H₂₃NO₄
FW: 341.4
Purity: ≥95%
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

(-)-Isocorypalmine is supplied as a solid. A stock solution may be made by dissolving the (-)-isocorypalmine in the solvent of choice, which should be purged with an inert gas. (-)-Isocorypalmine is soluble in methanol.

Description

(-)-Isocorypalmine is an isoquinoline alkaloid and a metabolite of L-tetrahydropalmitine (L-THP; Item No. 20535) that has been found in *Corydalis* and has diverse biological activities.¹⁻⁵ It binds to human dopamine D₁₋₅ receptors with K_i values of 6.2, 41.8, 37.3, 77.4, and 9.5 nM, respectively.² (-)-Isocorypalmine inhibits Epstein-Barr virus-early antigen (EBV-EA) activation induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in Raji cells (IC₅₀ = 300 mol ratio/32 pmol TPA), which is predictive for anti-tumor activity.³ It is cytotoxic to A549, SKOV3, SK-MEL-2, and HCT15 cancer cells (IC₅₀s = 67.32, 47.37, 47.66, and 67.32 μM, respectively).⁴ (-)-Isocorypalmine is active against clinical strains of *C. albicans*, *C. glabrata*s, *C. krusei*, *C. parapsilosis*, and *C. neoformans* (MICs = 40-320 μg/ml).⁵

References

1. Xiao, W., Zhuang, X., Shen, G., *et al.* Simultaneous quantification of L-tetrahydropalmitine and its urine metabolites by ultra high performance liquid chromatography with tandem mass spectrometry. *J. Sep. Sci.* **37(6)**, 696-703 (2014).
2. Lee, D.Y.W., Liu, J., Zhang, S., *et al.* Asymmetric total synthesis of tetrahydropyroberberine derivatives and evaluation of their binding affinities at dopamine receptors. *Bioorg. Med. Chem. Lett.* **27(6)**, 1437-1440 (2017).
3. Ito, C., Itoigawa, M., Tokuda, H., *et al.* Chemopreventive activity of isoquinoline alkaloids from *Corydalis* plants. *Planta Med.* **67(5)**, 473-475 (2001).
4. Kim, K.H., Piao, C., J., Choi, S.U., *et al.* New cytotoxic tetrahydropyroberberine-aporphine dimeric and aporphine alkaloids from *Corydalis turtchaninovii*. *Planta Med.* **76(15)**, 1732-1738 (2010).
5. 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, 12/09/2022

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