

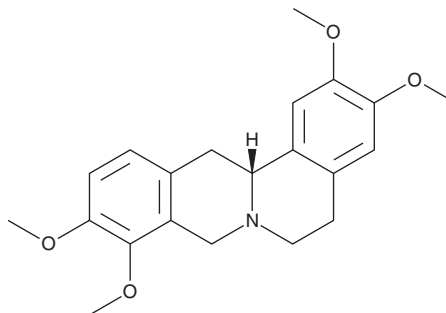
PRODUCT INFORMATION



D-Tetrahydropalmatine

Item No. 35577

CAS Registry No.: 3520-14-7
Formal Name: 5,8,13,13aR-tetrahydro-2,3,9,10-tetramethoxy-6H-dibenzo[a,g]quinolizine
Synonyms: (+)-Corydalis B, (+)-Tetrahydropalmatine, (R)-Tetrahydropalmatine, D-THP
MF: C₂₁H₂₅NO₄
FW: 355.4
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

D-Tetrahydropalmatine is supplied as a solid. A stock solution may be made by dissolving the D-tetrahydropalmatine in the solvent of choice, which should be purged with an inert gas. D-Tetrahydropalmatine is slightly soluble in chloroform, ethanol, and methanol.

Description

D-Tetrahydropalmatine is a tetrahydroberberine alkaloid that has been found in *Corydalis*.¹ *In vivo*, D-Tetrahydropalmatine (32 mg/kg) reverses apomorphine-induced inhibition of dopaminergic firing in the substantia nigra pars compacta (SNC) in paralyzed rats.²

References

1. Liao, J., Liang, W.-Z., and Tu, G.-S. Isolation and identification of eleven tertiary alkaloids in *Corydalis decumbens*. *J. Chin. Pharma. Sci.* **4**(2), 57-61 (1995).
2. Sun, B.C., Huang, K.X., and Jin, G.Z. Comparison of effects of tetrahydropalmatine enantiomers on firing activity of dopamine neurons in substantia nigra pars compacta. *Zhongguo Yao Li Xue Bao* **13**(4), 292-297 (1992).

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, 12/12/2022

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