

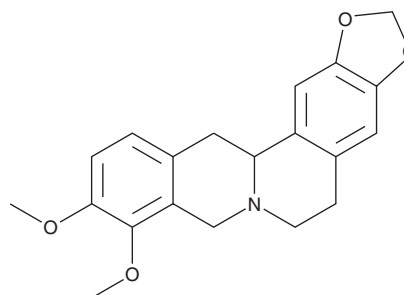
PRODUCT INFORMATION



Tetrahydroberberine

Item No. 33157

CAS Registry No.: 522-97-4
Formal Name: 5,8,13,13a-tetrahydro-9,10-dimethoxy-6H-benzo[g]-1,3-benzodioxolo[5,6-a]quinolizine
Synonyms: Canadine, NSC 36351, NSC 94918, (±)-Tetrahydroberberine, Xanthopuccine
MF: C₂₀H₂₁NO₄
FW: 339.4
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Phellodendron chinense*



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

Laboratory Procedures

Tetrahydroberberine is supplied as a crystalline solid. A stock solution may be made by dissolving the tetrahydroberberine in the solvent of choice, which should be purged with an inert gas. Tetrahydroberberine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of tetrahydroberberine in these solvents is approximately 1 mg/ml.

Tetrahydroberberine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tetrahydroberberine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Tetrahydroberberine has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Tetrahydroberberine is an isoquinoline alkaloid that has been found in *C. turttschaninovii* and has diverse biological activities.¹⁻⁴ It inhibits iron- and ascorbate-induced lipid peroxidation in rat liver microsomes (IC₅₀ = 13.7 μM).¹ Tetrahydroberberine binds to dopamine D₂ receptors and the serotonin (5-HT) receptor subtype 5-HT_{1A} (K_is = 0.83 and 4.2 μM, respectively).² It inhibits potassium currents induced by acetylcholine (Item No. 23829) in primary rat hippocampal CA1 pyramidal neurons (IC₅₀ = 13 μM).³ Tetrahydroberberine stimulates mouse C12C12 myoblast differentiation in a concentration-dependent manner.⁴

References

1. Ubeda, A., Montesinos, C., Payá, M., *et al.* Antioxidant action of benzyloisoquinoline alkaloids. *Free Radic. Res. Commun.* **18**(3), 167-175 (1993).
2. Lee, T.H., Kim, K.H., Lee, S.O., *et al.* Tetrahydroberberine, an isoquinoline alkaloid isolated from corydalis tuber, enhances gastrointestinal motor function. *J. Pharmacol. Exp. Ther.* **338**(3), 917-924 (2011).
3. Wu, J. and Jin, G.Z. Tetrahydroberberine inhibits acetylcholine-induced K⁺ current in acutely dissociated rat hippocampal CA1 pyramidal neurons. *Neurosci. Lett.* **222**(2), 115-118 (1997).
4. Lee, H., Lee, S.-J., Bae, G.-U., *et al.* Canadine from *Corydalis turttschaninovii* stimulates myoblast differentiation and protects against myotube atrophy. *Int. J. Mol. Sci.* **18**(12), 2748 (2017).

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