

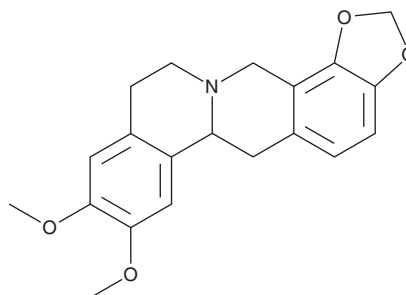
PRODUCT INFORMATION



Tetrahydroepiberberine

Item No. 36177

CAS Registry No.: 38853-67-7
Formal Name: 6,6a,11,14-tetrahydro-8,9-dimethoxy-12H-benzo[a]-1,3-benzodioxolo[4,5-g]quinolizine
Synonym: Sinactine
MF: C₂₀H₂₁NO₄
FW: 339.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Sinomenium acutum*



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

Laboratory Procedures

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

Description

Tetrahydroepiberberine is an alkaloid originally isolated from *P. rhoeas* that has diverse biological activities.^{1,2} It is active against the fungus *C. albicans* (MIC = 8 µg/ml), as well as the bacteria *E. coli*, *P. aeruginosa*, *P. mirabilis*, *K. pneumoniae*, *A. baumannii*, *S. aureus*, and *B. subtilis* (MICs = 32-64 µg/ml, respectively).² Tetrahydroepiberberine (16 µg/ml) inhibits the cytopathogenic effect of parainfluenza-3 (PI-3) virus in Vero cells. It decreases the viability of HaCaT keratinocytes (IC₅₀ = 150 µM).¹

References

1. Hijazi, M.A., Aboul-Ela, M., Bouhadir, K., *et al.* Cytotoxic activity of alkaloids from *Papaver rhoeas* growing in Lebanon. *Rec. Nat. Prod.* **11(2)**, 211-216 (2017).
2. Orhana, I., Ozcelik, B., Karaoğlu, T., *et al.* Antiviral and antimicrobial profiles of selected isoquinoline alkaloids from *Fumaria* and *Corydalis* species. *Z. Naturforsch. C. J. Biosci.* **62(1-2)**, 19-26 (2007).

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, 12/19/2022

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