

PRODUCT INFORMATION



Fangchinoline

Item No. 29243

CAS Registry No.: 436-77-1
Formal Name: (4a*S*,16a*S*)-3,4,4a,5,16a,17,18,19-octahydro-12,21,26-trimethoxy-4,17-dimethyl-16*H*-1,24:6,9-dietheno-11,15-metheno-2*H*-pyrido[2',3':17,18][1,11]dioxacycloeicosino[2,3,4-*ij*]

Synonyms: isoquinolin-22-ol
7-O-Demethyltetrandrine,
Hanfangchin B, NSC 77036

MF: C₃₇H₄₀N₂O₆

FW: 608.7

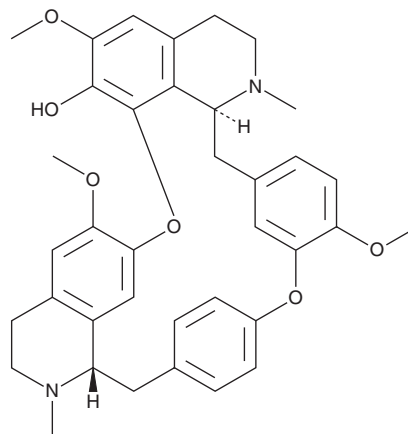
Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Stephaniae tetrandrae radix*



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

Laboratory Procedures

Fangchinoline is supplied as a crystalline solid. A stock solution may be made by dissolving the fangchinoline in the solvent of choice, which should be purged with an inert gas. Fangchinoline is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of fangchinoline in these solvents is approximately 30 mg/ml.

Fangchinoline is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fangchinoline should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fangchinoline 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

Fangchinoline is an alkaloid that has been found in *S. tetrandrae* and has diverse biological activities, including anti-inflammatory, neuroprotective, and anticancer properties.¹⁻³ It inhibits the activity of sheep COX and decreases IL-6-induced proliferation of MH60 mouse hybridoma cells with IC₅₀ values of 129 and 3.7 μM, respectively.¹ Fangchinoline (1 and 10 μM) reduces cyanide-induced increases in secreted glutamate levels and cell death in primary neonatal rat cerebellar granule neurons.² It decreases proliferation of PC3 human prostate cancer cells by 63 and 86% when used at concentrations of 20 and 30 μM, respectively.³ Fangchinoline reduces tumor growth in a PC3 mouse xenograft model when administered at a dose of 5 mg/kg per day for 12 days.

References

1. Choi, H.S., Kim, H.S., Min, K.R., *et al.* Anti-inflammatory effects of fangchinoline and tetrandrine. *J. Ethnopharmacol.* **69**(2), 173-179 (2000).
2. Cho, S.O. and Seong, Y.H. Protective effect of fangchinoline on cyanide-induced neurotoxicity in cultured rat cerebellar granule cells. *Arch. Pharm. Res.* **25**(3), 349-356 (2002).
3. Wang, C.D., Huang, J.G., Gao, X., *et al.* Fangchinoline induced G1/S arrest by modulating expression of p27, PCNA, and cyclin D in human prostate carcinoma cancer PC3 cells and tumor xenograft. *Biosci. Biotechnol. Biochem.* **74**(3), 488-493 (2010).

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, 10/12/2022

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