

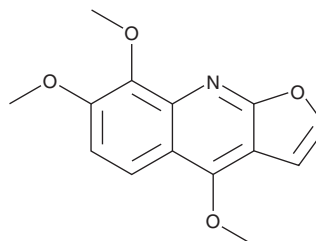
PRODUCT INFORMATION



Skimmianine

Item No. 34643

CAS Registry No.: 83-95-4
Formal Name: 4,7,8-trimethoxy-furo[2,3-b]quinoline
Synonyms: β -Fagarine, NSC 94654, NSC 217986
MF: C₁₄H₁₃NO₄
FW: 259.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 250 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/Skimmia japonica



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

Laboratory Procedures

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

Description

Skimmianine is a furoquinoline alkaloid that has been found in *Z. nitidum* and has diverse biological activities.¹⁻⁴ It inhibits acetylcholinesterase (AChE; IC₅₀ = 8.6 μ g/ml for the human enzyme).¹ Skimmianine is active against *P. falciparum* (IC₅₀ = 48.2 μ g/ml), as well as *L. gongylophorus*, a symbiotic fungus of the agricultural pest *A. sexdens* (leaf-cutting ant).^{2,3} It reduces ear edema induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in mice (ED₄₅ = 0.75 mg/ear).⁴

References

1. Yang, Z.-d., Zhang, D.-b., Ren, J., *et al.* Skimmianine, a furoquinoline alkaloid from *Zanthoxylum nitidum* as a potential acetylcholinesterase inhibitor. *Med. Chem. Res.* **21**, 722-725 (2011).
2. Khalid, S.A., Farouk, A., Geary, T.G., *et al.* Potential antimalarial candidates from African plants: An in vitro approach using *Plasmodium falciparum*. *J. Ethnopharmacol.* **15(2)**, 201-209 (1986).
3. Biavatti, M.W., Vieira, P.C., da Silva, M.F.d.G.F., *et al.* Biological activity of quinoline alkaloids from *Raulinoa echinata* and X-ray structure of flindersiamine. *J. Braz. Chem. Soc.* **13(1)**, 66-70 (2002).
4. García-Argáez, A.N., Ramírez Apan, T.O., Parra Delgado, H., *et al.* Anti-inflammatory activity of coumarins from *Decatropis bicolor* on TPA ear mice model. *Planta Med.* **66(3)**, 279-281 (2000).

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, 10/06/2022

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