

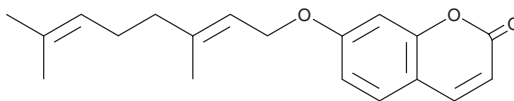
PRODUCT INFORMATION



Auraptene

Item No. 14000

CAS Registry No.: 495-02-3
Formal Name: 7-[[[(E)-3,7-dimethyl-2,6-octadien-1-yl]oxy]-2H-1-benzopyran-2-one
Synonym: 7-Geranyloxy coumarin
MF: C₁₉H₂₂O₃
FW: 298.4
Purity: ≥98%
UV/Vis.: λ_{max}: 323 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Auraptene is supplied as a crystalline solid. A stock solution may be made by dissolving the auraptene in the solvent of choice, which should be purged with an inert gas. Auraptene is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of auraptene in DMF is approximately 14 mg/ml and approximately 12 mg/ml in ethanol and DMSO.

Auraptene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, auraptene should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Auraptene has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Auraptene is a coumarin derived from citrus plants that bears a geranyloxy moiety at its C-7. It has anti-inflammatory, anti-carcinogenic, anti-bacterial, neuroprotective, and hepatoprotective activities.¹⁻⁵ It inhibits leukocyte activation and induces phase II enzymes during the initiation phase of carcinogenesis.¹ When examined for its potential use in Alzheimer's disease treatment, auraptene was shown to inhibit β-secretase (BACE1) activity with an IC₅₀ value of 345.1 μM.⁶

References

1. Murakami, A., Nakamura, Y., Tanaka, T., *et al.* Suppression by citrus auraptene of phorbol ester- and endotoxin-induced inflammatory responses: Role of attenuation of leukocyte activation. *Carcinogenesis* **21(10)**, 1843-1850 (2000).
2. Krishnan, P., Yan, K.J., Windler, D., *et al.* Citrus auraptene suppresses cyclin D1 and significantly delays N-methyl nitrosourea induced mammary carcinogenesis in female Sprague-Dawley rats. *BMC Cancer* **9**, (2009).
3. Takeda, K., Utsunomiya, H., Kakiuchi, S., *et al.* Citrus auraptene reduces *Helicobacter pylori* colonization of glandular stomach lesions in Mongolian gerbils. *J. Oleo Sci.* **56(5)**, 253-260 (2007).
4. Furukawa, Y., Watanabe, S., Okuyama, S., *et al.* Neurotrophic effect of Citrus auraptene: Neuritogenic activity in PC12 cells. *Int. J. Mol. Sci.* **13**, 5338-5347 (2012).
5. Sahebkar, A. Citrus auraptene: A potential multifunctional therapeutic agent for nonalcoholic fatty liver disease. *Ann. Hepatol.* **10(4)**, 575-577 (2011).
6. Marumoto, S. and Miyazawa, M. Structure-activity relationships for naturally occurring coumarins as β-secretase inhibitor. *Bioorg. Med. Chem.* **20(2)**, 784-788 (2012).

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, 10/18/2022

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