

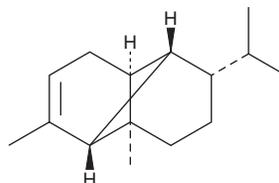
PRODUCT INFORMATION



α -Copaene

Item No. 25770

CAS Registry No.: 3856-25-5
Formal Name: (1R,2S,6S,7S,8S)-1,3-dimethyl-8-(1-methylethyl)-tricyclo[4.4.0.0^{2,7}]dec-3-ene
Synonym: (-)-Copaene
MF: C₁₅H₂₄
FW: 204.4
Purity: \geq 95%
Supplied as: An oil
Storage: -20°C
Stability: \geq 2 years
Item Origin: Plant/Angelica oil



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

Laboratory Procedures

α -Copaene is supplied as an oil. A stock solution may be made by dissolving the α -copaene in the solvent of choice, which should be purged with an inert gas. α -Copaene is slightly soluble in chloroform and methanol.

Description

α -Copaene is a sesquiterpene found in a variety of plants, including *Cannabis*, that has antiproliferative and antioxidant properties.^{1,2} It inhibits proliferation of primary rat neurons and N2a rat neuroblastoma cells in a concentration-dependent manner and cultured human lymphocytes when used at concentrations greater than 100 mg/L.^{2,3} α -Copaene (25 mg/L) increases total antioxidant capacity in primary rat neurons and N2a neuroblastoma cells, as well as cultured human lymphocytes, when used at concentrations of 50 and 100 mg/L.

References

1. Marchini, M., Charvoz, C., Dujourdy, L., *et al.* Multidimensional analysis of cannabis volatile constituents: Identification of 5,5-dimethyl-1-vinylbicyclo[2.1.1]hexane as a volatile marker of hashish, the resin of *Cannabis sativa* L. *J. Chromatogr. A* **1370**, 200-215 (2014).
2. Türkez, H., Çelik, K., and Toğar, B. Effects of copaene, a tricyclic sesquiterpene, on human lymphocytes cells *in vitro*. *Cytotechnology* **66(4)**, 597-603 (2014).
3. Turkez, H., Togar, B., Tatar, A., *et al.* Cytotoxic and cytogenetic effects of α -copaene on rat neuron and N2a neuroblastoma cell lines. *Biologia* **69(7)**, 936-942 (2014).

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, 07/08/2021

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