

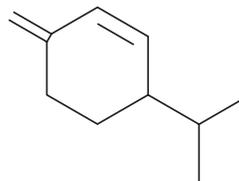
PRODUCT INFORMATION



β -Phellandrene

Item No. 33905

CAS Registry No.: 555-10-2
Formal Name: 3-methylene-6-(1-methylethyl)-cyclohexene
Synonyms: (\pm)- β -Phellandrene, NSC 53044
MF: C₁₀H₁₆
FW: 136.2
Purity: \geq 95%
UV/Vis.: λ_{max} : 232 nm
Supplied as: A neat oil
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

β -Phellandrene is supplied as a neat oil. A stock solution may be made by dissolving the β -phellandrene in the solvent of choice, which should be purged with an inert gas. β -Phellandrene is soluble in the organic solvent chloroform. β -Phellandrene is slightly soluble in ethanol.

Description

β -Phellandrene is a cyclic monoterpene that has been found in various plants, is secreted from live millipedes (*C. hualienensis* and *O. gracilis*), and has apoptotic activity.^{1,2} It is cytotoxic to RAW 264.7 cells when used at concentrations ranging from 50 to 200 $\mu\text{g}/\text{ml}$ and induces apoptosis in the same cells.²

References

1. Ramírez, J., Gilardoni, G., Ramón, E., *et al.* Phytochemical study of the Ecuadorian species *Lepechinia mutica* (Benth.) Epling and high antifungal activity of carnosol against *Pyricularia oryzae*. *Pharmaceuticals (Basel)* **11(2)**, 33 (2018).
2. Taira, J., Tamashiro, M., Naka, K., *et al.* Initial defensive secretory compounds emitted from the live millipede and the induction of apoptotic cell death. *Sci. Rep.* **11(1)**, 8222 (2021).

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, 11/04/2022

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