

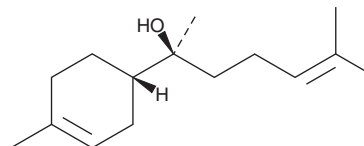
PRODUCT INFORMATION



(-)- α -Bisabolol

Item No. 9002495

CAS Registry No.: 23089-26-1
Formal Name: (α S,1S)- α ,4-dimethyl- α -(4-methyl-3-penten-1-yl)-3-cyclohexene-1-methanol
MF: $C_{15}H_{26}O$
FW: 222.4
Purity: $\geq 90\%$
Supplied as: A neat oil
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

(-)- α -Bisabolol is supplied as a neat oil. A stock solution may be made by dissolving the (-)- α -bisabolol in the solvent of choice, which should be purged with an inert gas. (-)- α -Bisabolol is soluble in the organic solvent ethanol.

Description

(-)- α -Bisabolol is a sesquiterpene alcohol that has been found in the essential oils of several aromatic plants, including *C. sativa*, *C. indica*, and *C. sativa/C. indica* hybrid strains, with diverse biological activities.¹⁻⁵ It inhibits the growth of *L. amazonensis* promastigotes and amastigotes ($\text{IC}_{50}\text{s} = 8.07$ and $4.29 \mu\text{g/ml}$, respectively).² It also reduces the growth of *L. infantum* and *L. donovani* amastigotes ($\text{IC}_{50}\text{s} = 56.9$ and $39.4 \mu\text{M}$, respectively) with a cytotoxic concentration (CC_{50}) value of greater than $1,000 \mu\text{M}$ in L929 cells.³ *In vivo*, (-)- α -bisabolol (200 mg/kg) reduces parasite load in the liver and spleen in a mouse model of visceral leishmaniasis. (-)- α -Bisabolol (200 mg/kg) reduces the size of ethanol-induced lesions, decreases malondialdehyde (MDA) production, and increases superoxide dismutase (SOD) activity in gastric mucosa in mice.⁴ It also decreases serotonin-, carrageenan-, and dextran-, but not histamine-, induced paw edema as well as acetic-acid induced writhing and paw-licking time in the formalin test in mice, indicating anti-inflammatory and antinociceptive activity.⁵

References

1. Elzinga, S., Fischedick, J., Podkolinski, R., *et al.* Cannabinoids and terpenes as chemotaxonomic markers in cannabis. *Nat. Prod. Chem. Res.* **3**(4), (2015).
2. Rottini, M.M., Amaral, A.C., Ferreira, J.L., *et al.* *In vitro* evaluation of (-)- α -bisabolol as a promising agent against *Leishmania amazonensis*. *Exp. Parasitol.* **148**, 66-72 (2015).
3. Corpas-López, V., Morillas-Márquez, F., Navarro-Moll, M.C., *et al.* (-)- α -Bisabolol, a promising oral compound for the treatment of visceral leishmaniasis. *J. Nat. Prod.* **78**(6), 1202-1207 (2015).
4. Rocha, N.F., Oliverira, G.V., Araújo, F.Y., *et al.* (-)- α -Bisabolol-induced gastroprotection is associated with reduction in lipid peroxidation, superoxide dismutase activity and neutrophil migration. *Eur. J. Pharm. Sci.* **44**(4), 455-461 (2011).
5. Rocha, N.F., Rios, E.R., Carvalho, A.M., *et al.* Anti-nociceptive and anti-inflammatory activities of (-)- α -bisabolol in rodents. *Naunyn Schmiedeberg's Arch. Pharmacol.* **384**(6), 525-533 (2011).

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, 01/10/2023

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