

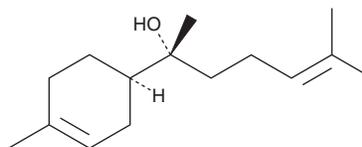
PRODUCT INFORMATION



(±)- α -Bisabolol

Item No. 31397

CAS Registry No.: 515-69-5
Formal Name: (α R,1R)-*rel*- α ,4-dimethyl- α -(4-methyl-3-penten-1-yl)-3-cyclohexene-1-methanol
Synonym: DL- α -Bisabolol
MF: C₁₅H₂₆O
FW: 222.4
Purity: \geq 95% (mixture of isomers)
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

(±)- 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 organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (±)- α -bisabolol in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

(±)- α -Bisabolol is a sesquiterpene alcohol that has been found in *M. chamomilla* and has anticancer activity.^{1,2} It induces apoptosis in, and is cytotoxic to, T67 and U87 glioma cells when used at concentrations ranging from 3.5 to 10 μ M.¹ (±)- α -Bisabolol is also cytotoxic to primary acute lymphoblastic leukemia (ALL) and acute myeloid leukemia (AML) cells (IC₅₀s = 14-65 μ M).²

References

1. Cavalieri, E., Mariotto, S., Fabrizi, C., *et al.* α -Bisabolol, a nontoxic natural compound, strongly induces apoptosis in glioma cells. *Biochem. Biophys. Res. Commun.* **315**(3), 589-594 (2004).
2. Cavalieri, E., Rigo, A., Bonifacio, M., *et al.* Pro-apoptotic activity of α -bisabolol in preclinical models of primary human acute leukemia cells. *J. Transl. Med.* **9**, 45 (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, 11/28/2022

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