

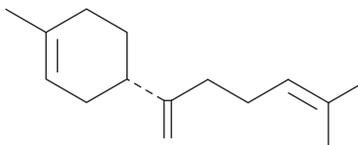
PRODUCT INFORMATION



β-Bisabolene

Item No. 43567

CAS Registry No.: 495-61-4
Formal Name: (4S)-1-methyl-4-(5-methyl-1-methylene-4-hexen-1-yl)-cyclohexene
Synonyms: (-)-β-Bisabolene, (S)-β-Bisabolene
MF: C₁₅H₂₄
FW: 204.4
Purity: ≥85%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/Mandarin oil



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

Laboratory Procedures

β-Bisabolene is supplied as a neat oil. A stock solution may be made by dissolving the β-bisabolene in the solvent of choice, which should be purged with an inert gas. β-Bisabolene is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

β-Bisabolene is a sesquiterpene that has been found in *Z. officinale* and has antiviral and neuroprotective activities.^{1,2} It reduces rhinovirus IB virus replication in a plaque reduction assay (IC₅₀ = 14.3 μg/plate).¹ β-Bisabolene (31.25-125 μg/ml) protects against decreases in viability, reduces production of reactive oxygen species (ROS), and decreases levels of COX-2, nitric oxide (NO), TNF-α, and IL-1β in HMC3 human microglial cells in an *in vitro* model of ischemia induced by glucose-oxygen deprivation.²

References

1. Denyer, C.V., Jackson, P., and Loakes, D.M. Isolation of antirhinoviral sesquiterpenes from ginger (*Zingiber officinale*). *J. Nat. Prod.* **57**(5), 658-662 (1994).
2. Zhang, X., Gao, L., Wang, J., Zhang, W., *et al.* Study of the ameliorative effect of β-Bisabolene on ischemic stroke via COX-2 with the Keap1/Nrf2 and MAPK pathways. *Eur. J. Pharmacol.* **1001**, 177773 (2025).

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, 10/29/2025

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