

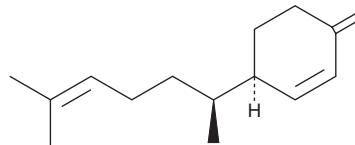
PRODUCT INFORMATION



(-)-β-Sesquiphellandrene

Item No. 35749

CAS Registry No.: 20307-83-9
Formal Name: 3R-[(1S)-1,5-dimethyl-4-hexen-1-yl]-6-methylene-cyclohexene
MF: C₁₅H₂₄
FW: 204.4
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

(-)-β-Sesquiphellandrene is supplied as an oil. A stock solution may be made by dissolving the (-)-β-sesquiphellandrene in the solvent of choice, which should be purged with an inert gas. (-)-β-Sesquiphellandrene is slightly soluble in chloroform.

Description

(-)-β-Sesquiphellandrene is a sesquiterpene that has been found in *Z. officinale* and has antiviral and anticancer activities.^{1,2} It reduces rhinovirus IB virus replication in a plaque reduction assay (IC₅₀ = 0.44 μM).¹ (-)-β-Sesquiphellandrene is cytotoxic to HCT116 cells when used at a concentration of 10 μM.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. Tyagi, A.K., Prasad, S., Yuan, W., et al. Identification of a novel compound (β-sesquiphellandrene) from turmeric (*Curcuma longa*) with anticancer potential: Comparison with curcumin. *Invest. New Drugs* **33**(6), 1175-1186 (2015).

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, 12/13/2022

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