

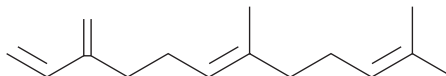
PRODUCT INFORMATION



trans-β-Farnesene

Item No. 31502

CAS Registry No.: 18794-84-8
Formal Name: (6E)-7,11-dimethyl-3-methylene-1,6,10-dodecatriene
Synonyms: (E)-β-Farnesene, β-Farnesene
MF: C₁₅H₂₄
FW: 204.4
Purity: ≥98%
UV/Vis.: λ_{max}: 225 nm
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

trans-β-Farnesene is supplied as a neat oil. A stock solution may be made by dissolving the *trans*-β-farnesene in the solvent of choice, which should be purged with an inert gas. *trans*-β-Farnesene is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of *trans*-β-farnesene in these solvents is approximately 30 mg/ml.

Description

trans-β-Farnesene is a volatile sesquiterpene hydrocarbon that has been found in aphids and has alarm pheromone and kairomonal attractant activities.^{1,2} It is produced by aphids in response to predator or parasite attack.¹ *trans*-β-Farnesene (0.2 ng) induces dispersion of 50% of *M. persicae* aphids from a cluster. *trans*-β-Farnesene (>2 μg) also acts as a kairomone, attracting larvae and adults of *A. bipunctata*, a predator of aphids.²

References

1. Montgomery, M.E. and Nault, L.R. Comparative response of aphids to the alarm pheromone (E)-β-Farnesene. *Entomol. Exp. Appl.* **22(3)**, 236-242 (1977)
2. Francis, F., Lognay, G., and Haubruge, E. Olfactory responses to aphid and host plant volatile releases: (E)-β-farnesene an effective kairomone for the predator *Adalia bipunctata*. *J. Chem. Ecol.* **30(4)**, 741-755 (2004).

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, 11/21/2022

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