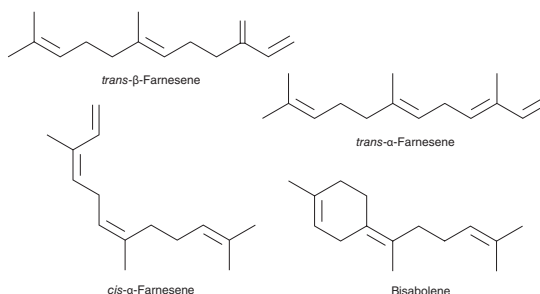


# PRODUCT INFORMATION

## Farnesene (mixed isomers)

Item No. 26891

MF:  $C_{15}H_{24}$   
 FW: 204.4  
 UV/Vis.:  $\lambda_{max}$ : 229, 280 nm  
 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

Farnesene (mixed isomers) is supplied as a neat oil. A stock solution may be made by dissolving the farnesene (mixed isomers) in the solvent of choice, which should be purged with an inert gas. Farnesene (mixed isomers) is miscible in organic solvents such as ethanol, DMSO, and dimethyl formamide.

### Description

Farnesene (mixed isomers) is a mixture of isomers of the sesquiterpene hydrocarbon farnesene, which has been found in *G. hirsutum* essential oils.<sup>1</sup> This product may contain *trans*-β-farnesene, *cis*-α-farnesene, *trans*-α-farnesene, and bisabolene, among others.

### Reference

1. Minyard, J.P., Tumlinson, J.H., Thompson, A.C., *et al.* Constituents of the cotton bud. Sesquiterpene hydrocarbons. *J. Agric. Food Chem.* **14**(4), 332-336 (1966).

#### 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, 03/04/2024

#### 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