

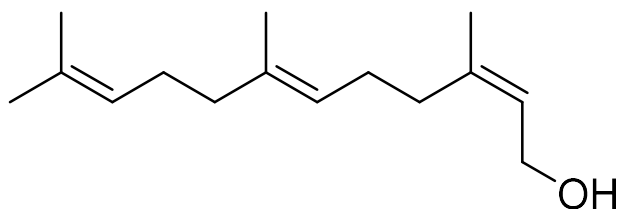
2Z,6E-Farnesol

Catalog number: S-0180

Molecular Formula: C₁₅H₂₆O

MW: 222.3

CAS#: 3790-71-4



Alternate Name: (2Z,6E)-3,7,11-trimethyl-2,6,10-dodecatriene-1-ol, *cis,trans*-farnesol

Solubility: Chloroform, ethyl acetate, hexane, and most organic solvents

Storage and Handling: Store dry at -20 °C. Stock solutions should be stored frozen (-20 °C or below).

Background: Isoprenoids and isoprenols are key metabolic intermediates for biosynthesis of steroids, including cholesterol and hormones that regulate human reproduction or fertility, and posttranslational modifications of proteins in signal transduction pathways for cell division and regulation of vesicular trafficking. Farnesol is an acyclic sesquiterpene alcohol and the 2E,6E-isomer being the most common in nature. 2Z,6E-Farnesol is a metabolite in the sesquiterpenoid and triterpenoid biosynthesis, formed by farnesol 2-isomerase from 2E,6E-farnesol.

References: 1) Panayiotis Anastasis, Isabel Freer, Christopher Gilmore, Hugh Mackie, Karl Overton, Douglas Picken, and Stephen Swanson. (1984) "Biosynthesis of γ -bisabolene in tissue cultures of *Andrographispaniculata*." *Canadian Journal of Chemistry*. **62**(10): 2079-2088. doi:10.1139/v84-355
2) Thulasiram HV, Poulter CD. (2006) "Farnesyl diphosphate synthase: the art of compromise between substrate selectivity and stereoselectivity. *J Am Chem Soc*. 128(49):15819-23. doi: 10.1021/ja065573b.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

Technical Data Sheet, Rev 1, 5/13/24– **For research use only.** Not intended for diagnostic or therapeutic use.

