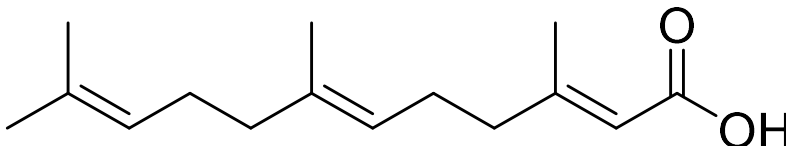


(*E,E*)-Farnesoic Acid

Catalog number: S-0151



Molecular Formula: C₁₅H₂₄O₂

MW: 236.18

CAS#: 462-11-3

Alternate Name: (2*E*,6*E*)-3,7,11-trimethyldodeca-2,6,10-trienoic acid

Solubility: Soluble in most organic solvents, poor water solubility

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

Background: Farnesoic acid is the substrate of farnesoic acid O-methyltransferase which produces the crustacean reproductive hormone, methyl farnesoate (MF). MF is responsible for enhancing reproductive maturation, maintaining juvenile morphology, and influencing male sex determination. Farnesoic acid also inhibits filament formation in *C. albicans*.

References: 1) Chung, S.-C., T.-I. Kim, et al. (2010). "Candida albicans PHO81 is required for the inhibition of hyphal development by farnesoic acid." *FEBS letters* 584(22): 4639.
2) Zhao, N., J. Guan, et al. (2007). "Molecular cloning and biochemical characterization of indole-3-acetic acid methyltransferase from poplar." *Phytochemistry* 68(11): 1537.
3) Li, Y., G. C. Unnithan, et al. (2003). "Stimulation of JH biosynthesis by the corpora allata of adult female *Aedes aegypti* in vitro: effect of farnesoic acid and *Aedes* allatotropin." *J Exp Biol* 206(Pt 11): 1825-32.

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 6, 5/13/24– **For research use only.** Not intended for diagnostic or therapeutic use.

