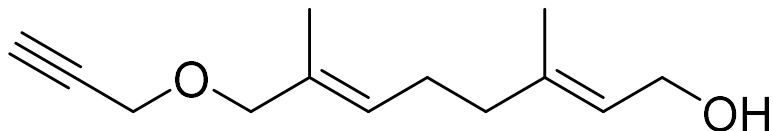


Propargyl-geraniol

Catalog number: S-0110



Molecular Formula: C₁₃H₂₀O₂

MW: 208.3

CAS#: 109334-38-5

Alternate Name: (E,E)-3,7-Dimethyl-8-(2-propargyloxy)-2,6-octadien-1-ol

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

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

Background: The alkyne-modified isoprenoids were developed by the Distefano group at the University of Minnesota as chemical reporters of protein prenylation. The isoprenoids readily penetrate mammalian cells and are used as substrates by prenyltransferases. The pendant alkyne can be derivatized with an azide-containing tag, such as a fluorophore or biotin, via a Cu(I) catalyzed “Click” reaction allowing the prenylated proteins to be identified and quantified.

References: 1) A.J. DeGraw, et al. “Evaluation of alkyne-modified isoprenoids as chemical reporters of protein prenylation” (2010) Chem. Biol. Drug Des. 76, 460-47, doi: 10.1111/j.1747-0285.2010.01037.x
2) B.P. Duckworth, et al “Selective Labeling of Proteins by Using Protein Farnesyltransferase” (2007) ChemBioChem 8, 98-105. DOI: 10.1002/cbic.200600340

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 2, 5/13/24– **For research use only.** Not intended for diagnostic or therapeutic use.

