



Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

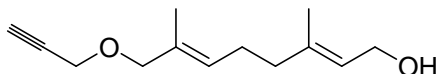
Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

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

Product Number: S-0110



Molecular Formula: C₁₃H₂₀O₂ **MW:** 208.3 **CAS:** 109334-38-5

Solubility: Chloroform, dichloromethane, ethyl acetate,

Storage: -20°C or below

Field of Interest: 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" *Chem. Biol. Drug Des.* **2010**, 76, 460-471 2) B.P. Duckworth, *et al.* "Selective Labeling of Proteins by Using Protein Farnesyltransferase" *ChemBioChem*, **2007**, 8, 98-105.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. Propargyl-geraniol 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.