



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

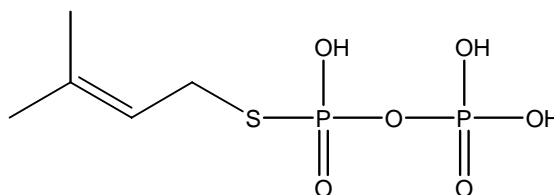
*Not intended or approved for
diagnostic or therapeutic use.*

Product Name:

DIMETHYLALLYL S-THIOLODIPHOSPHATE

Product Number:

I-S051



Provided as	Chemical Formula	MW (g/mol)	Solubility
Tris-ammonium salt	C ₅ H ₂₁ N ₃ O ₆ P ₂ S	313.25	>1mg/ml H ₂ O or MeOH

PACKAGING:

Dimethylallyl S-thiolodiphosphate is available as the lyophilized solid.

SPECIFICATIONS AND STORAGE RECOMMENDATIONS:

Purity not less than 95% (by weight for solids and as percent of stated concentration for solutions) when shipped. Organic phosphates vary in stability and shelf life. Samples should be kept at 0°C or at -70°C for prolonged storage.

FIELD OF INTEREST:

Isoprenoid diphosphates 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.

HAZARDOUS PROPERTIES AND CAUTIONS:

No hazardous or toxic properties are known for this substance. Dimethylallyl S-thiolodiphosphate is not listed as a hazardous waste or as a Toxic Chemical subject to release reporting under the Emergency Planning and Community Right-to-Know Act. For solutions containing methanol or other solvents, see

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.

MSDS for Organic Phosphates in Solution available on request.

Dimethylallyl S-thiolodiphosphate is not listed on the Chemical Inventory of the Toxic Substance Control Act, and 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. Purchase 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.

Phan, R.M. and Poulter, C.D. Synthesis of (S)-Isoprenoid Thiodiphosphates as substrates and inhibitors. *J. Org. Chem.* **66**(20), 6705-10. (2110).

Phan, R.M. and Poulter, C.D. Synthesis of Geranyl S-Thiolodiphosphate. A new alternative substrate/inhibitor for prenyltransferases. *Organic Letters* **2**(15), 2287-89. (2000).