



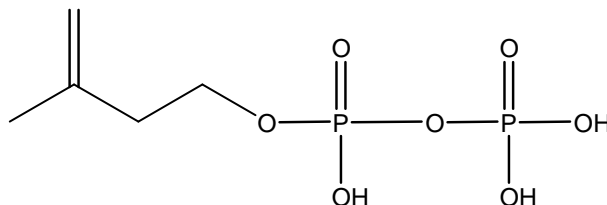
Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

*For research use only*

**Product Name:**  
**ISOPENTENYL DIPHOSPHATE (IPP)**

**Product Number:** I-0050



Provided as	Chemical Formula	MW (g/mol)	Solubility	CAS Number
Tris-ammonium salt	C <sub>5</sub> H <sub>21</sub> N <sub>3</sub> O <sub>7</sub> P <sub>2</sub>	297.18	>1 mg/ml H <sub>2</sub> O or MeOH	116057-53-3

**PACKAGING:** Isopentenyl diphosphate (IPP) is provided as the lyophilized tris-ammonium salt.

**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. The samples should be kept at 0°C or kept 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. Isopentenyl diphosphate 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 MSDS for Organic Phosphates in Solution available on request.

Isopentenyl diphosphate 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.

Rohdich, F., Hecht, S., Gartner, K., Asam, P., Krieger, C., Amslinger, S., Arigoni, D., Bacher, A., & Eisenrich, W. Studies on the nonmevalonate terpene biosynthetic pathway; metabolic role of IspH (LytB) protein. *Proc. Natl. Acad. Sci. USA* **99**, 1158-63. (2002).

Burke, C. & Croteau, R. Interaction with the small subunit of geranyl diphosphate synthase modifies the chain length specificity of geranylgeranyl diphosphate synthase to produce geranyl diphosphate. *J Biol Chem.* **277**, 3141-9. (2002).