



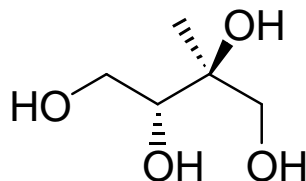
Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

*For research use only*

**Product Name:** 2-C-Methyl-erythritol (ME)

**Product Number:** I-M051A



Provided as	Chemical Formula	MW (g/mol)	Solubility
Oil	C <sub>5</sub> H <sub>12</sub> O <sub>4</sub>	136.15	>1 mg/ml MeOH

**PACKAGING:** 2-C-Methyl-erythritol is available in sealed vials as an oil.

### SPECIFICATIONS AND STORAGE RECOMMENDATIONS:

Provided as a racemic mixture. Purity not less than 95% (by weight) when shipped. For optimum shelf life, store below -20°C.

### FIELD OF INTEREST:

2-C-Methyl-erythritol (ME) is a substrate for the first pathway specific enzymes in the biosynthesis of thiamine diphosphate (vitamin B1), pyridoxol phosphate (vitamin B6), and the nonmevalonate family of isoprenoid compounds recently discovered in bacteria and plant chloroplasts. ME was prepared via chemical synthesis from racemic precursors.

### HAZARDOUS PROPERTIES AND CAUTIONS:

No hazardous or toxic properties are known for this substance. 2-C-Methyl-erythritol (ME) 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.

2-C-Methyl-erythritol (ME) 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

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.

TDS I-M051A Rev: 1 (1/14/04)

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.