



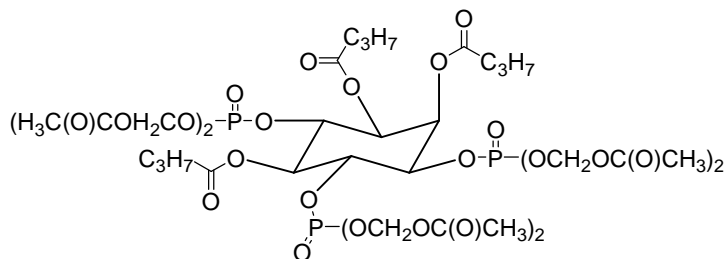
Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Product Name: 2,3,5-Tri-O-butyryl-Ins(1,4,6)P₃/AM, Membrane Permeant Ins(1,4,6)P₃

Product Number: Q-M0146



Chemical Formula	MW (g/mol)	Solubility
C ₃₆ H ₅₇ O ₃₀ P ₃	1062.75	< 5mg/ml, DMSO, toluene

Storage: Membrane permeant inositol polyphosphates are stable for at least one year when stored as a syrup, protected from moisture, at -20 °C. After reconstitution, solutions should be stored at -20 °C between uses.

Field Of Interest: Membrane permeant inositol polyphosphates are employed in cell signaling experiments to introduce exogenous inositol polyphosphates across cell membranes. Some examples of applications are listed below. See the following page for an example protocol.

- 1) M. Vajanaphanich *et al.* "Long-term uncoupling of chloride secretion from intracellular calcium levels by Ins(3,4,5,6)P₄" *Nature*, **1994**, 371, 711-714.
- 2) M.T. Rudolf *et al.* "A membrane-permeant, bioactivatable derivative of Ins(1,3,4)P₃ and its effect on Cl(-)-secretion from T84 cells" *Bioorg. Med. Chem.Lett.*, **1998**, 8, 1857-1860.
- 3) X. Yang *et al.* "Inositol 1,3,4-trisphosphate acts in vivo as a specific regulator of cellular signaling by inositol 3,4,5,6-tetrakisphosphate." *J. Biol. Chem.*, **1999**, 274, 18973-18980.

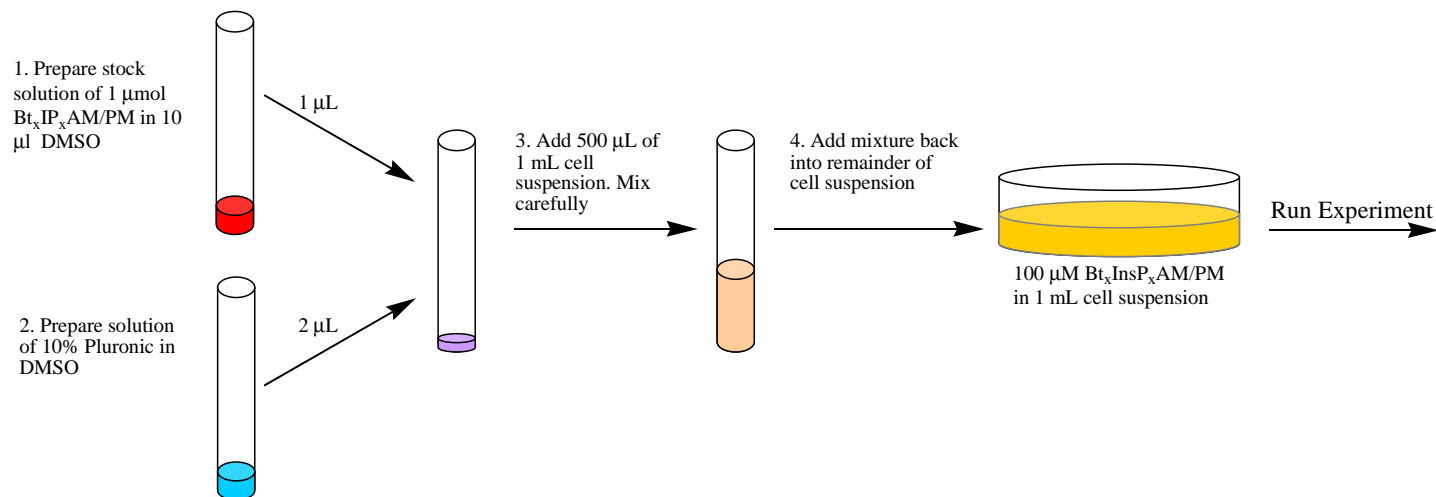
Hazardous Properties and Cautions: No hazardous or toxic properties are known for this substance. 2,3,5-Tri-O-butyryl-Ins(1,4,6)P₃/AM is not listed as a hazardous waste or as a Toxic Chemical subject to release reporting under the Emergency Planning and Community Right-to-Know Act. For solutions containing methanol or other solvents, see MSDS for Phosphoinositides in Solution available on request. 2,3,5-Tri-O-butyryl-Ins(1,4,6)P₃/AM 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.

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.

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.

Example protocol for preparing final concentration of 100 μ M membrane permeant inositol phosphate (Bt_xInsP_xAM/PM) in 1 mL of cell suspension.



- 1) Add 10 μ L of DMSO to 1 μ mol Bt_xIP_xAM/PM to prepare 100 mM stock solution
- 2) Prepare a solution of 10% Pluronic[®] in DMSO
- 3) Add 1 μ L of Bt_xIP_xAM/PM and 2 μ L of Pluronic[®] solution to a vial.
- 4) Add 500 μ L of your cell suspension to the vial and mix carefully
- 5) Add this mixture back into your remaining cell suspension and mix carefully. Proceed with planned experiment.

Notes

- 1) The cell suspension should not contain more than 1% DMSO
- 2) The stock solution of Bt_xIP_xAM/PM in DMSO can be frozen for later use but **do not** freeze the solution containing Bt_xIP_xAM/PM **and** Pluronic[®].