



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: PolyPIPosomes™

Catalog #	PIP _n	Catalog #	PIP _n
Y-0000	No PIP _n	Y-P034	PI(3,4)P ₂
Y-P000	PI	Y-P035	PI(3,5)P ₂
Y-P003	PI(3)P	Y-P045	PI(4,5)P ₂
Y-P004	PI(4)P	Y-P039	PI(3,4,5)P ₃
Y-P005	PI(5)P		

Composition: PolyPIPosomes™ are prepared in deionized water containing 0.1% sodium azide.

	Mole %	Concentration (mM) in H ₂ O
Polymerizable-PC	65	0.65
Polymerizable-PE	29	0.29
Polymerizable-PIP _n	5	0.05
Polymerizable-biotin-PE	1	0.009
		1 mM total lipid

Specifications and Storage Recommendations: PolyPIPosomes™ are relatively stable at room temperature for short periods of time. For long term storage of up to 6 months, PolyPIPosomes™ can be stored at 4°C. DO NOT FREEZE. Storage in basic buffers (pH > 9.0) will result in slow hydrolysis of the ester chains, and may cause phosphate or acyl migration to occur. Storage in acidic buffers (pH < 4.0) may cause decomposition or phosphate migration.

Field of Interest: PtdInsP_ns compounds are employed as substrates for kinases, phosphatases and binding proteins as described in many publications. PolyPIPosomes have been described in the following publications:

- 1) C. G. Ferguson, *et al*, "Phosphoinositide-Containing Polymerized Liposomes: Stable Membrane-Mimetic Vesicles for Protein-Lipid Binding Analysis" *Bioconjugate Chem* **16**, 2005, 1475.
- 2) S. K Elkin, *et al*, "A PHD Finger Motif in the C Terminus of RAG2 Modulates Recombination Activity" *J. Biol. Chem.*, **280**, 2005, 28701.

Hazardous Properties and Cautions: No hazardous or toxic properties are known for this substance. PolyPIPosomes™ are 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. PolyPIPosomes™ are not listed on the Chemical Inventory of the Toxic Substance Control Act, and are 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. This product contains sodium azide at a concentration of 0.1%. Sodium azide is not classified as a hazardous chemical at this concentration; however, toxicity information regarding sodium azide at this concentration has not been thoroughly investigated—handle with caution. A Material Safety Data Sheet (MSDS) for sodium azide in pure form is available upon request.

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.

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.