



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

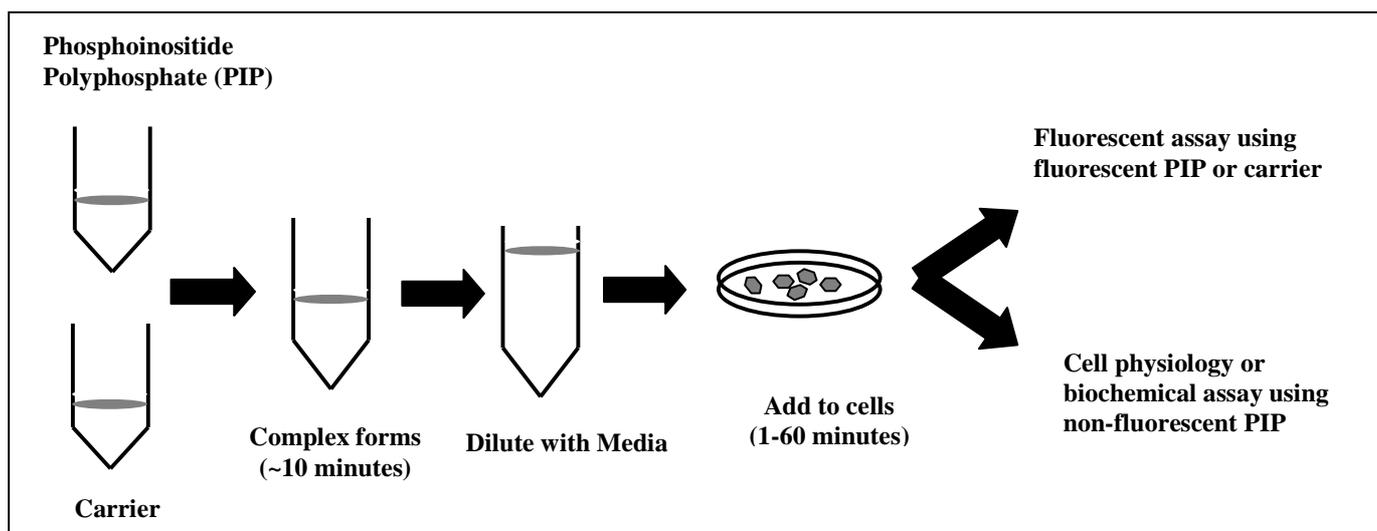
Technical Data Sheet

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Shuttle PIP™ Intracellular delivery of phosphoinositides

General Protocol

Add carriers at a 1 to 1 molar ratio with phosphoinositides (1:1) for 5-15 minutes in a test tube at room temperature. Brief vortex mixing and/or bath sonication may help complexes dissolve completely. Once formed, the complex is diluted to the desired final concentration in media or PBS containing cells (either adherent or suspended). *Note, we recommend final concentrations between 0.1 and 50 μ M, but the optimal concentration for a given experimental system should be determined by each investigator.* Depending on the cell type, the phosphoinositide, and the system of study; significant uptake of fluorescent phosphoinositide or change in cell physiology is seen within 1 to 60 minutes. Slower-developing effects may require repeated application of carriers and phosphoinositides.



Product Test

Kits were tested using fluorescent phosphoinositides and carriers on NIH/3T3 cells imaged with a Nikon TS100 microscope. A successful test shows intracellular staining. Briefly, carriers and fluorescent phosphoinositides, 13 μ L combined, were incubated in a test tube at a 1 to 1 molar ratio (~100 μ M final concentration each) for 10 minutes at room temperature. The complex was diluted with 12 μ L of medium and then added to 100 μ L media covering adherent cells grown in an 8-well chamber slide on glass cover-slips. After 30-60 minutes, the dye-containing media was replaced with PBS and images were collected. The final shuttle and PIP concentrations on cells were both 10 μ M. See Ozaki et.al. (2000) Intracellular delivery of phosphoinositides and inositol phosphates using polyamine carriers *Proc Natl Acad Sci U S A* **97** (21)11286-11291. Please visit our web site or contact our technical service department (1-866-588-0455) for additional questions.

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 P-9199 Rev: 7 (1/16/04)