



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:**  
Shuttle PIP™ Kits  
Intracellular delivery of phosphoinositides

**Product Number:** P-9005

### **Kit Contents:**

#### *Phosphoinositides*

<u>Catalog #</u>	<u>Description</u>	<u>Molecular Weight</u>	<u>Quantity</u>
P-5016	PtdIns(5)P diC <sub>16</sub>	956.96	100 µg
C-05F16	*BODIPY® FL-PtdIns(5)P,C <sub>16</sub>	1,372.88	25 µg

#### *Carriers*

P-9C3	Carrier 3	1,551	3 x 50 nmoles
-------	-----------	-------	---------------

### **Storage and Handling:**

Certain kit components are moisture and light sensitive. Store unopened kit for up to one year frozen at -20°C protected from moisture and light. Reconstitute phosphoinositides and carriers in aqueous buffers or media, and store at 4°C for up to 3 months. Multiple freeze thawing is not recommended. Note: Vortex mixing, brief bath sonication, and addition of small amounts of methanol, ethanol, or DMSO may facilitate complete dissolution of phosphoinositides. *Phosphate buffers are not recommended and may alter complex formation between carriers and phosphoinositides.* We do not recommend storing carriers and PIPs together as complexes.

Carriers are shipped lyophilized and are hygroscopic, so avoid moisture. Reconstitute carriers with water or buffered aqueous solutions, then store at 4°C for up to 3 months. Multiple freezing and thawing is not recommended.

### **References:**

1. Ozaki, S.; DeWald, D. B.; Shope, J. C.; Chen, J.; Prestwich, G. D. (2000) Intracellular delivery of phosphoinositides and inositol phosphates using polyamine carriers. *Proc Natl Acad Sci U S A* **97**(21)11286-11291.
2. Scheid, M.P., M. Huber, et. al. (2002). Phosphatidylinositol (3,4,5)P<sub>3</sub> is essential but not sufficient for protein kinase B (PKB) activation; Phosphatidylinositol (3,4)P<sub>2</sub> is required for PKB phosphorylation at Ser-473: studies using cells from SH2-containing inositol-5-phosphatase knockout mice. *J Biol Chem* **277**(11): 9027-35.
3. Weiner, O.D., Neilsen, P.O., Prestwich, G.D., Kirschner, M.W., Cantley, L.C. and Bourne, H.R. (2002). A PtdInsP<sub>3</sub>- and Rho GTPase-mediated positive feedback loop regulates neutrophil polarity. *Nat Cell Biol* **4** 509-512.

\* **BODIPY® FL** has maximal excitation at 505 nm and maximal emission at 513 nm

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.