



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: PIP Beads™

One ml of agarose beads containing 10 nanomoles of bound PIP_n. Beads are in PBS buffer. Each order of PIP Beads™ includes a small amount of control beads.

Product Numbers

P-B001- DiC₆ PtdIns

P-B003- DiC₆ PI(3)P

P-B004 - DiC₆ PI(4)P

P-B005- DiC₆ PI(5)P

P-B034- DiC₆ PI(3,4)P₂

P-B035- DiC₆ PI(3,5)P₂

P-B045- DiC₆ PI(4,5)P₂

P-B345- DiC₆ PI(3,4,5)P₃

Product Storage: Store at 2-8°C. Product is moisture and light sensitive. Do not freeze.

Phospholipid source: Phosphoinositides and inositol phosphates are from Echelon Biosciences, Inc.

Suggested Usage: lipid-protein pull-down

Reference: Rao, V. et al. Expression Cloning of Protein Targets for 3-phosphorylated phosphoinositides. *J. Biol. Chem.* **274**(53): 37893-37900 (1999).

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.