



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:** Purified Anti-PI(3,4)P<sub>2</sub> Antibody  
Mouse Monoclonal IgG

**Product Number:** Z-P034

**Quantity:** 10µg      25µg      50µg      125µg

**Storage:** Store at 4 °C for up to 30 days. Aliquot and store at -20 °C if longer storage is necessary. Avoid repeated freeze/thaw cycles.

**Concentration:** 1 mg/mL

**How Supplied:** Clear solution, diluted to the indicated concentration in PBS, pH 7.4.

**Antibody Subclass:** IgG3

**Suggested Use:** Z-P034 recognizes PtdIns(3,4)P<sub>2</sub>, biotinylated PtdIns(3,4)P<sub>2</sub>, or fluorescent PtdIns(3,4)P<sub>2</sub> at the following concentrations:  
0.5 µg/mL for ELISA and Protein Lipid Overlay (PIP Strip™) assays,  
0.1 to 1 µg/mL for AlphaScreen™ and,  
250 µg/mL for Fluorescent Polarization assays.  
Other *in vitro* and cellular applications are possible using this antibody, but have not been verified by Echelon Biosciences. *These suggested uses are intended as a starting point for optimization, and the optimal antibody concentration and other experimental conditions for each application should be determined by the user.*

**Specificity:** Z-P034 reacts primarily with the head group of the indicated phosphoinositide (of synthetic or natural origin), and demonstrates low cross-reactivity with other phosphoinositides or phospholipids depending on the assay format. Not all antibodies are validated for use in all types of assays. Please check the updated technical data sheets on our web site or call our customer service department (toll free, 1-866-588-0455) for antibody application information and 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.