



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:	Biotinylated Anti-PI(3,4)P₂ Antibody Mouse Monoclonal IgG
Product Number:	Z-B034
Quantity:	25 µg
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:	0.5 mg/mL
How Supplied:	Clear solution, diluted to the indicated concentration in PBS, pH 7.4, 0.02% Azide.
Antibody Subclass:	IgG3
Suggested Use:	<p>Z-B034 is from the same lot of purified antibody (Z-P034) which recognizes PtdIns(3,4)P₂, biotinylated PtdIns(3,4)P₂, or fluorescent PtdIns(3,4)P₂ at the following concentrations:</p> <p><u>0.5 µg/mL</u> for ELISA (PIP Plate™) and Lipid Overlay (PIP Strip™) assays, <u>0.25 to 1 µg/mL</u> for AlphaScreen™ and, <u>250 µg/mL</u> for Fluorescent Polarization assays.</p> <p>This biotinylated product has been tested to confirm the presence of biotin covalently linked to mouse IgG in cellular immunofluorescence experiments. Other <i>in vitro</i> and cellular applications are possible using this antibody, but have not been verified by Echelon Biosciences.</p> <p><i>Note: 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.</i></p>
Specificity:	Z-P034 reacts primarily with the head group of the indicated phosphoinositide, 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.