



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: MultiPIP Grip™ (LL5α-PH)

Product Number: G-9901

Quantity: 2.5 µg lyophilized protein

Storage: Store lyophilized protein at -80 °C or -20 °C. Protein will be stable for at least one year from date of shipment. For best results, reconstitute the protein in assay buffer immediately before use.

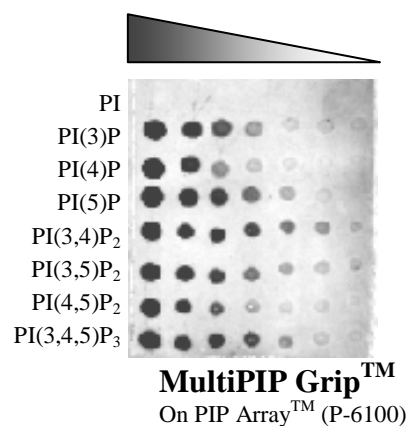
Source: Recombinant LL5-α PH domain GST-tagged protein

Protein Size: 50 kD including GST-tag

Purity: > 80% by SDS-PAGE

Specificity: Recognizes all phosphoinositides

Suggested Use: As a control reagent for PIP Strips™, MicroStrips™, PIP Array™, PIP Beads™, or PIP Plates™.



For PIP Strips™, PIP Microstrips™, PIP Arrays™:

Recommended concentration = 0.5 – 1.0 µg/mL for PIP Strips™ and PIP Arrays, and 0.1 µg/mL – 0.25 µg/mL for MicroStrips™. Reconstitute protein in blocking buffer (TBS-T 0.1% (v/v) Tween-20 with 3% fatty acid free BSA for best results). Add solution directly to previously blocked PIP Strips™, PIP MicroStrips™, or PIP Arrays™. Detect with anti-GST or enzyme conjugated secondary antibody. See protocol for PIP Strip™ products for more information at www.echelon-inc.com

For PIP / Cova Plate™:

Recommended concentration = 1 - 3 µg/mL. Add 100 µL/well of protein solution to PIP plate. Detect with anti-GST, enzyme conjugated secondary antibody.

Reference: Anjana Saxena, Pavel Morozov, Dale Frank, Raymond Musalo, Mark A. Lemmon, Edward Y. Skolnik, and Benjamin Tycko. Phosphoinositide Binding by the Pleckstrin Homology Domains of Ip1 and Tih1 The Journal of Biological Chemistry (2002) 277 (51) 49935-49944

Related Products: Lipid-Protein Interaction tools:
GloPIPs™, PIP Beads™, PolyPIPosomes™, PIP Strips™, PIP Arrays™, PIP Plates™
Anti-lipid antibodies (see website)

Other PIP Grips™:

PIP Grip™	Name	Catalog	Specificity
PI(3)P Grip™	p40PX	G-0302	Recognizes PI(3)P
PI(4)P Grip™	SidC_3C	G-0402	Recognizes PI(4)P
PI(4,5)P ₂ Grip™	PLCδ	G-4501	Recognizes PI(4,5)P ₂
PI(3,4,5)P ₃ Grip™	Grp1	G-3901	Recognizes PI(3,4,5)P ₃

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 G-9901 Rev: 1 (05/19/08)