



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: PI(3,4,5)P₃ Grip™

Product Number: G-3901

Quantity: 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.

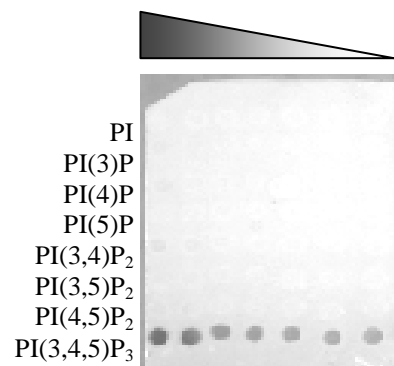
Source: Recombinant Grp1 PH domain GST-tagged protein

M.W. of Protein: 45 kDa including GST-tag

Purity: > 80% by SDS-PAGE

Specificity: Recognizes PI(3,4,5)P₃

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



PI(3,4,5)P₃ Grip
On PIP Array (P-6100)

For PIP Strips™, PIP MicroStrips™, PIP Array™:

Recommended concentration = 0.25 – 0.5 µg/mL for PIP Strips™ and PIP Arrays™, and 0.1 – 0.2 µg/mL for MicroStrips™. Reconstitute protein in blocking buffer (PBS-T 0.1% (v/v) Tween-20 with 3% BSA for best results). Add solution directly to previously blocked PIP Strips™, PIP Arrays™, or PIP MicroStrips™. 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 = 0.4 - 1 µg/mL. Add 100 µL/well of protein solution to PIP Plate. Detect with anti-GST, enzyme conjugated secondary antibody.

Reference: 1. 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 Ipl1 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
MultiPIP Grip™	LL5α	G-9901	All phosphoinositides
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