



Salt Lake City, Ut 84158-0537
 Telephone 866-588-0455
 Fax 801-588-0497
 Email: echelon@echelon-inc.com
 http://www.echelon-inc.com

Product Name: PIP Strips™

Technical Data Sheet

Product Number: P-6001

For research use only
 Not intended or approved for
 diagnostic or therapeutic use

General Description: PIP Strips™ are 2 x 6 cm hydrophobic membranes that have been spotted with 15 different biologically active lipids. These membranes can be used to determine lipid-protein interactions, through a simple lipid-protein overlay experiment. This allows researchers a convenient way to determine if their protein of interest interacts with one or more of the bound lipids.

Storage: Store at 2-8 °C. Product is moisture and light sensitive.

Format: The membrane has a diagonal cut on its top left corner and Ponceau S* staining to assist in orientation of the strip. See template below for location of lipids.

Membrane Template:

Lysophosphatidic Acid (LPA)	○	○	Sphingosine-1-phosphate (S1P)
Lysophosphocholine (LPC)	○	○	PtdIns(3,4)P ₂
PtdIns	○	○	PtdIns(3,5)P ₂
PtdIns(3)P	○	○	PtdIns(4,5)P ₂
PtdIns(4)P	○	○	PtdIns(3,4,5)P ₃
PtdIns(5)P	○	○	Phosphatidic Acid (PA)
Phosphatidylethanolamine (PE)	○	○	Phosphatidylserine (PS)
Phosphatidylcholine (PC)	○	○	Blank

Suggested Usage: See support protocol “basic Protocol for PIP strips™, PIP Array™, PIP Microstrips™, and SphingoStrips™” on our website www.echelon-inc.com

References: Dowler, S., Currie, R.A., Downes, P.C., and Alessi, D.R. DAPP1: a dual adaptor for phosphotyrosine and 3-phosphoinositides *Biochemical Society.J.* **342**, 7-12 (1999).
 Dowler, S., Kular, G., and Alessi, R.D., Protein lipid overlay assay, *Sci STKE*, 2002, L6. (2002).

*Final concentration of 0.1% (v/v) Ponceau S was added for accuracy during spotting.

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.

Product Name: PIP Array™

Technical Data Sheet

Product Number: P-6100

For research use only

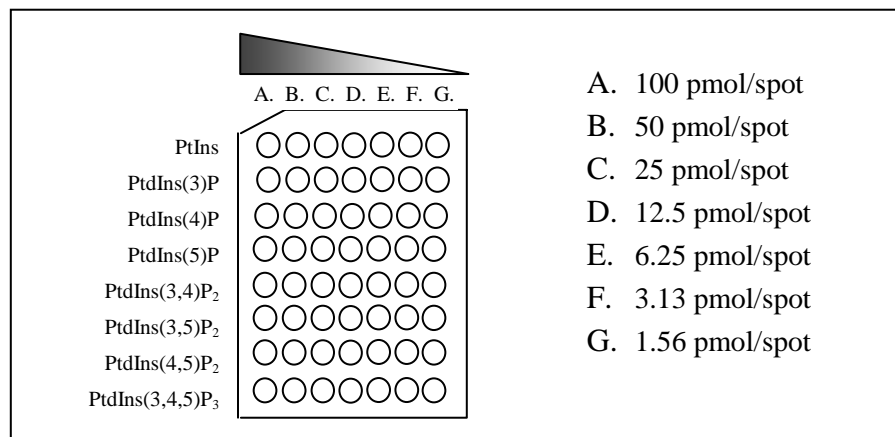
Not intended or approved for diagnostic or therapeutic use

General Description: PIP Array™ are 4 x 5 cm hydrophobic membranes that have been spotted with a concentration gradient of all eight phosphoinositides. These membranes can be used to determine lipid-protein interactions through a simple protein-lipid-overlay experiment. This allows researchers a convenient way to determine if their protein of interest interacts with one or more of the bound lipids and the degree of binding.

Storage: Store at 2-8 °C. Product is moisture and light sensitive.

Format: The membrane has a diagonal cut on its top left corner and Ponceau S* staining to assist in orientation of the strip. See template below for location of lipids.

Membrane Template:



Suggested Usage: See support protocol “basic Protocol for PIP strips™, PIP Array™, PIP Microstrips™, and SphingoStrips™” on our website www.echelon-inc.com

References: Dowler, S., Currie, R.A., Downes, P.C., and Alessi, D.R. DAPP1: a dual adaptor for phosphotyrosine and 3-phosphoinositides *Biochemical Society.J.* **342**, 7-12 (1999).
Dowler, S., Kular, G., and Alessi, R.D., Protein lipid overlay assay, *Sci STKE*, 2002, L6. (2002).

*Final concentration of 0.1% (v/v) Ponceau S was added for accuracy during spotting.



P.O. Box 58537
Salt Lake City, Ut 84158-0537
Telephone 866-588-0455
Fax 801-588-0497
Email: echelon@echelon-inc.com
<http://www.echelon-inc.com>

Product Name **LL5- α (MultiPIP Grip™)**

Document Type

Product Number: **G-1100**

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Quantity: 2.5 μ g

Storage: Store pellets at -20°C before reconstitution.
After reconstitution: Short term storage: Protein is stable at $+4^{\circ}\text{C}$
for 24 hours.
Long term storage: Aliquot and store at -20°C .
For best results, we recommend reconstituting the protein
immediately before use.

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

Protein Size: 49.95 kD (24.95 kD LL5- α + 25 kD GST-tag)

Purity: > 80% (Purity determined via SDS-PAGE)

Specificity: Recognizes all phosphoinositides, interacts more strongly with bis-
and tris- phosphates.

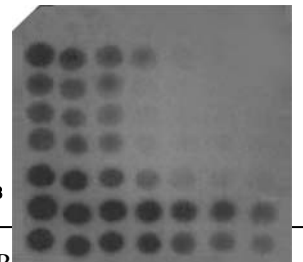
Suggested Usage: As a control reagent for PIP Strips™, PIP MicroStrips™, PIP
Array™, or PIP Plates™.

For PIP Strips™, PIP MicroStrips™,
PIP Array™:
Recommended concentration = 0.5 μ g/mL

Add solution directly to
previously blocked PIP Strips™,
PIP MicroStrips™, or PIP Array™.

LL5 α (MultiPIP Grip™)
PIP array at 0.5 μ g/mL

PtdIns
PtdIns(3)P
PtdIns(4)P
PtdIns(5)P
PtdIns(3,5)P₂
PtdIns(4,5)P₂
PtdIns(3,4)P₂
PtdIns(3,4,5)P₃



Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications, are covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our products at busdev@echelon-inc.com.

TDS G-1100 Rev. 5 (4/20/05)

For PIP Plate™.

1.0 µg/mL is the suggested
2.0 concentration of the stock solution.

Then add 50-100 µL/well of
stock 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 Ip1 and Tih1 The Journal of Biological Chemistry (2002) 277 (51) 49935-49944

Related Products:

- | | |
|---|---------------------------|
| 1. Grp1 (PIP₃ Grip™)
Specificity for PI(3,4,5)P ₃ | (Catalog # G-1200) |
| 2. PLCδ (PIP₂ Grip™)
Specificity for PI(4,5)P ₂ | (Catalog # G-1000) |