



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.

## PIP<sub>2</sub> Grip™ (PLC-δ)

- Product Number:** G-1000
- Quantity:** One pellet of 10.5 µg or 2.5 µg of PLC-δ1
- Storage:** Store pellets at -20 °C before reconstitution.  
For best results, we recommend reconstituting the protein immediately before use.
- Source:** Recombinant PLC-δ1 PH domain GST-tagged protein
- Protein Size:** 42.4 kD (17.4 kD PLC-δ1 + 25 kD GST-tag)
- Purity:** > 60% (Determined by SDS-PAGE)
- Specificity:** Recognizes PI(4,5)P<sub>2</sub>
- Suggested Usage:** As a control reagent for PIP Strips™, PIP MicroStrips™, PIP Array™, or PIP Plates™.

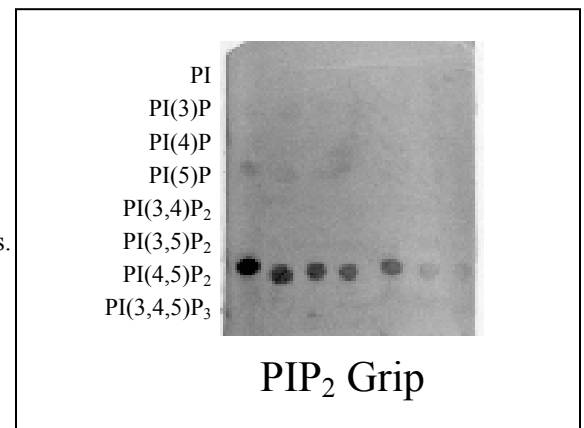
For PIP Strips™, PIP MicroStrips™, PIP Array™:

Reconstitute 1 pellet in 10 mL blocking buffer (PBS-1% Milk for best results).  
1.05 µg/mL for the 10.5 µg pellet and 0.25 µg/mL for the 2.5 µg pellet produce similar results.  
Add solution directly to previously blocked PIP Strips™, PIP MicroStrips™, PIP Array™, or Membrane Lipid Strips.

For PIP / Cova Plate™:

Recommended concentration = 1.0 µg/mL

Add 50-100 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 Ip1 and Tih1 The Journal of Biological Chemistry (2002) 277 (51) 49935-49944

- Related Products:**
- MultiPIP Grip™ (LL5α)** (Catalog # G-1100)  
Recognizes all phosphoinositides, interacts more strongly with bis- and tris- phosphates.
  - PIP<sub>3</sub> Grip™ (Grp1)** (Catalog # G-1200)  
Specificity for PI(3,4,5)P<sub>3</sub>

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-1000 Rev: 7 (12/29/05)