



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.

PIP₃ Grip™ (Grp1)

Product Number: G-1200

Quantity: 10 µg (2 pellets x 5 µg)

Storage: Store pellets at -20 °C before reconstitution.
After reconstitution: Short term storage: Protein is stable at +4 °C for 12-24 hours.
Long term storage: Aliquot and store at -20 °C.
For best results, we recommend reconstituting the protein immediately before use.

Source: Recombinant Grp1 PH domain GST-tagged protein

Protein Size: 46.8 kD (21.8 kD Grp1 + 25 kD GST-tag)

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

Specificity: Recognizes PI(3,4,5)P₃

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

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

0.005 µg/mL binds PI(3,4,5)P₃

0.5 µg/mL binds all PIPs

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

For PIP Plate™:

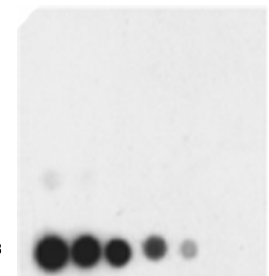
Recommended concentration = 1.0 µg/mL

Add 50-100 µL/well of protein solution to PIP plate.
Detect with anti-GST, enzyme-conjugated secondary antibody.

PIP₃ Grip™ (Grp1)

PIP Array at 0.005 µg/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



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. **MultiPIP Grip™ (LL5α)** (Catalog # G-1100)
Recognizes all phosphoinositides, interacts more strongly with bis- and tris- phosphates.
2. **PIP₂ Grip™ (PLCδ)** (Catalog # G-1000)
Specificity for PI(4,5)P₂

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-1200 Rev: 3 (12/6/05)