



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)P Grip™

Product Number: G-0302

Quantity: 2.5 µg, 10 µg, or 50 µg as 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 2.5 µg size, reconstitute the protein in assay buffer immediately before use. For multiple uses, reconstitute 10 µg and 50 µg protein with 50 µL and 250 µL ddH₂O respectively, aliquot the obtained 200 µg/mL protein stock solution and store at -80 °C for up to 1 month.

Source: Recombinant p40 PX domain GST-tagged protein

M.W. of Protein: 43.6 kDa including GST-tag

Purity: > 90% by SDS-PAGE

Specificity: Phosphatidylinositol 3-phosphate, PI(3)P

Applications: As a PI(3)P detector or as a control reagent for PIP Strips™, PIP Array™, PIP Beads™, or PIP Plates™.

For PIP Strips™, and PIP Array™:

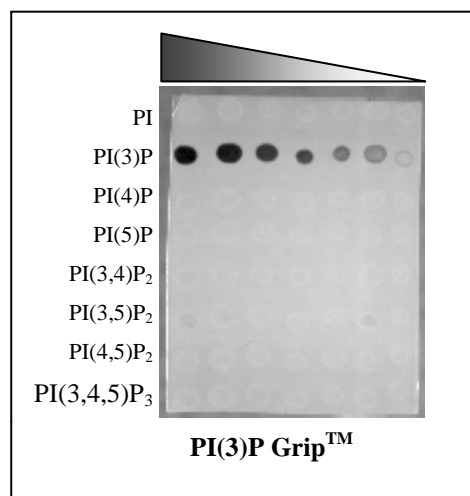
Recommended concentration = 1 µg/mL

Reconstitute 2.5 µg protein in blocking buffer (PBS-0.1% (v/v) Tween-20 with 3% fatty acid free BSA for best results). Add solution directly to previously blocked PIP Strips™ or PIP Array™. Detect with anti-GST, enzyme-conjugated secondary antibody. See protocol for PIP Strip products for more information.

For PIP / Cova Plate™:

Recommended concentration = 1 - 2 µg/mL

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



References:

1. Chris D. Ellison, Simon Andrews, Len R. Stephens, and Phill T. Hawkins. The PX domain: a new phosphoinositide-binding module. *Journal of Cell Science* (2002) 115: 1099-1105.
2. Robert V. Stahelin, Aura Burian, Karol S. Bruzik, Diana Murray, and Wonhwa Cho. Membrane Binding Mechanisms of the PX Domains of NADPH Oxidase p40phox and p47phox. *Journal of Biological Chemistry* (2003) 278 (16) 14469-14479.

Related Products:

1. Anti-PtdIns(3)P Antibody (cat# **Z-P003**)
2. Other PIP Grips™: PI(4)P Grip™ (cat# **G-0402**), PI(4,5)P₂ Grip™ (cat# **G-4501**), PIP₃ Grip™ (cat# **G-3901**), and MultiPIP Grip™ (cat# **G-9901**)
3. PI(3)P Mass ELISA kit (cat# **K-3300**)
4. Class III PI3-K Activity ELISA kit (cat# **K-3000**)
5. PI(3)P (cat# **P-3004**, **P-3008**, **P-3008a**, and **P-3016**)
6. Biotin labeled PI(3)P (cat# **C-03B6**) and GloPIPs™ Biotin labeled PI(3)P (cat# **C-03B6a**)
7. BODIPY FL-PI(3)P (cat# **C-03F6**) and GloPIPs™ BODIPY FL-PI(3)P (cat # **C-03F6a**)
8. BODIPY® TMR-PI(3)P (cat# **C-03M6**) and GloPIPs™ BODIPY® TMR-PI(3)P (cat# **C-03M6a**)
9. PIP Beads™ PI(3)P (Catalog # **P-B003a**) and PI(3)P PolyPIPosome (cat# **Y-P003**)
10. PIP Strips™ (cat# **P-6001**) and PIP Array™ (cat# **P-6100**)
11. PIP Plate (cat# **H-6203**, **H-6300**, **H-6403**, and **H-6500**) and PIP Microarray plate (cat# **M-1000**)

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.