



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(4)P Grip (SidC_3C)

Product Number: G-0402

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: N-terminal GST-tagged *Legionella pneumophila* SidC protein, C-terminal fragment.

Protein Size: 62 kDa (including GST-tag)

Purity: > 80% by SDS-PAGE

Specificity: Phosphatidylinositol 4-phosphate with long carbon chain, PI(4)P

Suggested Use: As a PI(4)P detector or as a control reagent for PIP Strips, PIP Array.

For PIP Strips or PIP Array:

Recommended concentration = 0.5 µg/mL

Reconstitute protein in blocking buffer (PBS-0.1% (v/v)

Tween-20 with 3% fatty acid free BSA for best results).

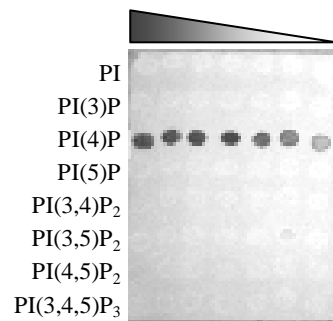
Add solution directly to previously blocked PIP Strips or

PIP Array. Detect with anti-GST, enzyme conjugated

secondary Antibody. See protocol for PIP Strip

membrane type products on our website at [www.echelon-](http://www.echelon-inc.com)

[inc.com](http://www.echelon-inc.com).



PI(4)P Grip

References:

1. Stefan S. Weber, Curdin Ragaz, Katrin Reus, Yves Nyfeler, and Hubert Hilbi. *Legionella pneumophila* Exploits PI(4)P to Anchor Secreted Effector Proteins to the Replicative Vacuole. *PLoS Pathogens* (2006) 2 (5) e46.
2. Curdin Ragaz, Holger Pietsch, Simon Urwyler, André Tiaden, Stefan S. Weber and Hubert Hilbi. *The Legionella pneumophila phosphatidylinositol-4 phosphate-binding type IV substrate SidC recruits endoplasmic reticulum vesicles to a replication-permissive vacuole*. *Cellular Microbiology* (2008) 10(12):2416-33

Related Products:

1. Anti-PtdIns(4)P Antibody (cat# **Z-P004**)
2. Other PIP Grips: PI(3)P Grip (cat # **G-0302**), PI(4,5)P₂ Grip (cat# **G-4501**), PIP₃ Grip (cat# **G-3901**), and MultiPIP Grip (cat# **G-9901**)
3. PI(4)P Mass Strip Kit (cat# **K-4000**)
4. PI(4)P (cat# **P-4004**, **P-4008**, **P-4008a**, **P-4016**) and GroPI(4)P (cat# **P-4000**)
5. Biotin labeled PI(4)P (cat# **C-04B6**) and GloPIPs Biotin labeled PI(4)P (cat# **C-04B6a**)
6. BODIPY FL-PI(4)P (cat# **C-04F6**) and GloPIPs BODIPY FL-PI(4)P (cat# **C-04F6a**)
7. BODIPY® TMR-PI(4)P (cat# **C-04M6**) and GloPIPs BODIPY® TMR-PI(4)P (cat# **C-04M6a**)
8. PIP Beads PI(4)P (cat# **P-B004a**) and PI(4)P PolyPIPosome (cat# **Y-P004**)
9. PIP Strips (cat# **P-6001**), PIP Array (cat# **P-6100**), and Membrane Lipid Strips (cat# **P-6002**)

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.