

# Echelon Biosciences Inc.

## PI(4,5)P2 Grip™ (PLC-δ1-PH)

G-4501

Support: echelon@echelon-inc.com

### Description:

N-terminal GST-tagged recombinant PLC-δ1 PH domain expressed in E. coli.

### Properties:

**Size** – 2.5 µg, 10 µg, & 50 µg

**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 ddH<sub>2</sub>O respectively, aliquot the obtained 200 µg/mL protein stock solution and store at -80°C for up to 1 month.

**M.W. of Protein** – 42.3 kDa including GST tag

**Purity** – >85% by SDS-PAGE

**Specificity** – recognizes PI(4,5)P2

### Applications:

As a PI(4,5)P2 detector or as a control reagent for PIP Strips™, Membrane Lipid Strips™, PIP Array™, or PIP Plates™.

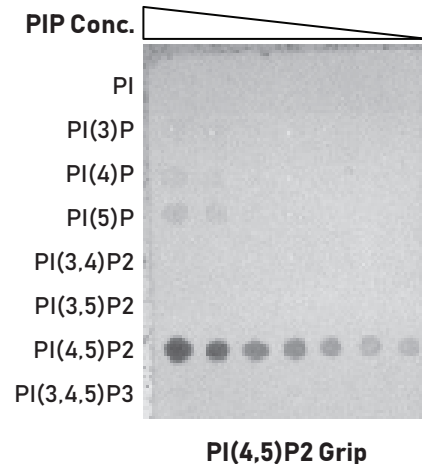
For PIP Strips™, Membrane Lipid Strips™, PIP Array™: recommended concentration = 0.5 - 1.0 µg/mL.

Reconstitute the 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™, PIP Array™, or Membrane Lipid Strips™. Detect with anti-GST, enzyme-conjugated secondary antibody. See protocol for PIP Strip products for more information.

For PIP / Cova Plate™: recommended concentration = 1 µg/mL.

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

### Data: PIP Array (Cat. # P-6100)



### References:

1. Barnett KC, Coronas-Serna JM, Zhou W, Ernandes MJ, Cao A, Kranzusch PJ, et al. (2019) Phosphoinositide Interactions Position cGAS at the Plasma Membrane to Ensure Efficient Distinction between Self- and Viral DNA. *Cell*. 2019;176(6):1432-46.
2. Xu J, Kozlov G, McPherson PS, Gehring K. A (2018) PH-like domain of the Rab12 guanine nucleotide exchange factor DENND3 binds actin and is required for autophagy. *The Journal of Biological Chemistry*. 293(12):4566-74.

### Related Products:

Products	Catalog Number
<b>Reagents</b>	
PIP Beads™ PI(4,5)P2	P-B045a
Flow PIPs: Multiplex Beads	P-FB01
PolyPIPosome	Y-P045
<b>Lipids</b>	
PI(4,5)P2	P-4504, P-4508, P-4516
BODIPYFL-PI(4,5)P2	C-45B6
BODIPYTMR-PI(4,5)P2	C-45M6
<b>Antibody &amp; Enzyme</b>	
Purified Anti-PI(4,5)P2	Z-P045
Acites Anti-PI(4,5)P2	Z-A045
Biotin Anti-PI(4,5)P2	Z-B045
Fluorescein Anti-PI(4,5)P2	Z-G045
<b>Assays</b>	
PIP Strips	P-6001
Membrane Lipid Strips	P-6002
PIP Array	P-6100
PI(4,5)P2 Mass ELISA	K-4500

Technical Data Sheet Rev. 3, 08-01-19 - For research use only. Not intended or approved for diagnostic or therapeutic use.



Echelon Biosciences Inc.

Ph: 866-588-0455

Fax: 801-588-0497

Echelon-inc.com