

PI(4)P Mass ELISA Assay Kit

K-4000E (96 tests)

Support: echelon@echelon-inc.com

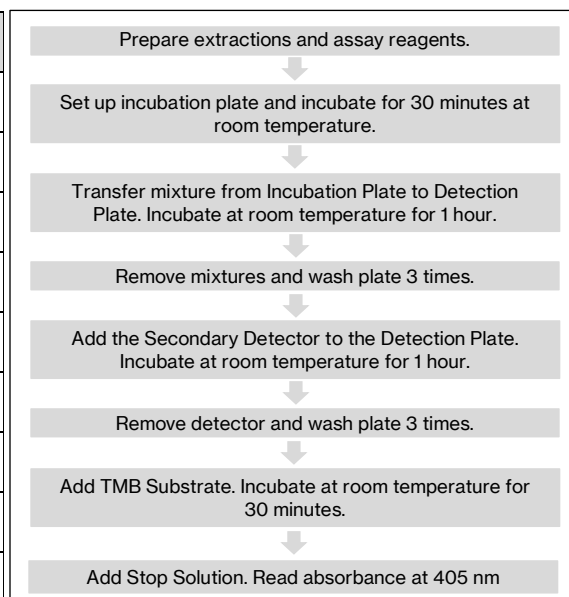
Description: 96-well ELISA assay for the detection and quantification of PI(4)P from cells.

Storage: The kit comes in two parts with different storage requirements. Upon receipt store Kit Part 1 at 4°C and Kit Part 2 at -80°C. Store prepared reagents as indicated in the protocol.

Materials Provided

Catalog #	Description	Quantity
K-4005	PI(4)P Detector	1 vial
K-4001E	PI(4)P Detection Plate	1 plate
K-4006	PI(4)P Standard (diC16)	1 vial
K-DIL2	Diluent (lyophilized pellet)	1 bottle
K-PBST2	10X PBS-T Buffer	1 bottle, 20 mL
K-SEC2	Secondary Detector	1 vial, 300 µL
K-TMB1	TMB Substrate	1 bottle, 12 mL
K-STOPt	1 N H2SO4 Stop Solution	1 bottle, 8 mL
Incubation Plate	Natural 96-well Polypropylene Plate	1 plate
---	Plate Seals Clear acetate sheet, 1 side Adhesive	3 seals

Quick Protocol



Additional Materials Provided by User

- Extracted PI(4)P samples (See Support Protocol PI(4)P extraction at the end of this document)
- Buffers and solvents for PI(4)P extractions: Trichloroacetic Acid, EDTA, Methanol, Chloroform, and 12 N HCl
- Pipettes (capable of delivering between 5 and 1,000 µL with appropriate tips)
- Absorbance microplate reader capable of reading at 405 nm
- Vacuum dryer

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposure or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

Background

Phosphatidylinositol 4-phosphate (PtdIns(4)P, PI(4)P) is the most abundant monophosphorylated phosphoinositide found in mammalian cells¹. It is produced by PtdIns 4-Kinases (PI4K) which phosphorylate the D-4 position of the inositol ring of PtdIns. PI(4)P is then converted to PtdIns(4,5)P₂ by PtdIns(4)P 5-Kinases.

Assay Design

Echelon's PI(4)P Mass ELISA kit is designed to detect and quantify PI(4)P by means of a competitive ELISA format, eliminating the need for radioactivity and thin layer chromatography. The PI(4)P Mass ELISA directly detects PI(4)P over all other phosphoinositides at relevant biological levels.

The PI(4)P Mass assay is a competitive ELISA in which the signal is inversely proportional to the amount of PI(4)P measured. The PI(4)P extractions are added to the Incubation Plate with PI(4)P Detector. This mixture is then transferred to the PI(4)P Detection Plate for competitive binding. A peroxidase-linked secondary detector and colorimetric detection is used to detect the amount of PI(4)P Detector binding to the plate. The assay is sensitive to 0.7 pmol PI(4)P.

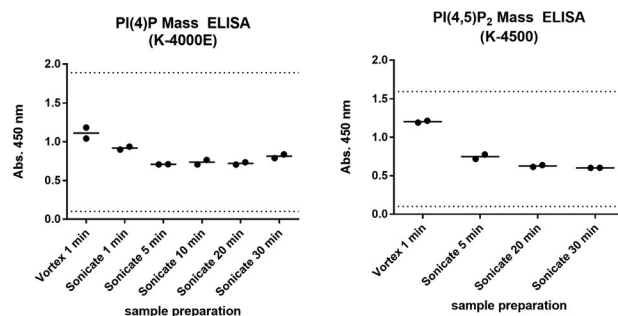
Disclaimer

The PIP Mass Assays are used to quantify the total amount of the specific lipid extracted from cells. However, since the lipids are substrates for enzymes in multiple pathways the data obtained may not correlate with what has been observed with isolated enzyme reactions or visualized with immunohistochemistry.

Assay Notes

1. The incubation steps for this assay require a plate shaker. If a plate shaker is unavailable, the incubation steps can be run for two hours without shaking. A reduction in signal and some loss in sensitivity may be observed.
2. The PI(4)P Detection Plate is composed of 12 8-well strips. Unused strips should be removed from the plate frame and stored in a clean sealable plastic bag at 4°C. Save the plate frame after assay for future use of the remaining strip wells.
3. Never let the detection plate dry out after the ELISA assay has started. Always prepare the next solution needed before discarding the current one from wells in use.
4. For lipid extraction questions please see "PIP Mass FAQ: PIP Mass Customer Information Sheet" located on the webpage for this product. This document can also be requested at echelon@echelon-inc.com.
5. It can be difficult to reproduce conditions of sonication; due to variation in the number of vials between

batches, temperature of the water bath, and sonicator tuning. The suggested sonication time of 5-10 min was developed using a water bath sonicator. The results observed with your water bath sonicator may be different. It is suggested that your sonicator is tested with PI(4)P extraction samples for day-to-day variation and time dependent consistency. If you are running other lipid mass assays you may want to consider dissolving your lipids in the same manner. How you dissolve the lipid will affect how it goes into solution and can cause inconsistencies in your data if it is not held constant.



Protocol for the detection of PI(4)P

Please read the entire Assay Procedure, Lipid Extraction Protocol, and Assay Notes before beginning assay. Bring PI(4)P Standard (K-4006), PI(4)P extraction samples, Buffers (K-PBST2, K-DIL2), TMB Solution (K-TMB1), 1 N H₂SO₄ Stop Solution (K-STOPT), and Detection Plate (K-4001E) to room temperature before use. Place PI(4)P Detector (K-4005) and Secondary Detector (K-SEC2) on ice until use.

Buffer Preparation

1. PBS-T Buffer: Prepare by adding the entire bottle of 10X PBS-T Buffer (K-PBST2) to 180 mL DI water. Mix. Keep at room temperature.
2. Working Diluent: Prepare by adding 12 mL DI water to the bottle of Diluent (K-DIL2). Vortex. Let sit for 5 minutes. Vortex. Keep at room temperature until use. For long term storage keep Working Diluent at -20°C.

PI(4)P Mass ELISA General Protocol

1. Prepare PI(4)P extraction samples and PI(4)P Standards
 - a. Make sure PI(4)P Standard (K-4006), PI(4)P extraction samples, and PBS-T buffer are at room temperature before proceeding to next step.
 - b. Add 558 µL of PBS-T Buffer to the vial of PI(4)P standard (K-4006). Vortex for 1 min at moderate speed to resuspend lipid. Follow with centrifugation. This is the 450 pmol/well standard. Further dilute 450 pmol/well standard

in 0.5 mL tubes, using the dilution table below. The 450 pmol/well standard can be stored at -20°C for 1 month and can withstand one freeze/defrost cycle.

PI(4)P Standard curve dilution table			
Row	pmol/well	Vol. stock solution or previous dilution	Vol. PBS-T
B	450.00	200 µL 450 pmol/well Standard	0 µL
C	150.00	100 µL 450 pmol/well Standard	200 µL
D	50.00	100 µL previous dilution	200 µL
E	16.66	100 µL previous dilution	200 µL
F	5.55	100 µL previous dilution	200 µL
G	1.85	100 µL previous dilution	200 µL

- c. To the PI(4)P extraction samples add 200 µL PBS-T Buffer then sonicate 5-10 min in a room temperature water bath. Do not add ice to water bath. See assay notes section for more information on sample preparation.
2. Prepare working PI(4)P Detector (K-4005) in Working Diluent
 - a. Add 50 µL of dH₂O to the vial of PI(4)P Detector

Incubation Plate/PI(4)P Detection Plate layout

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	Blank	ES #1	ES #1	ES #9	ES #9	ES #17	ES #17	ES #25	ES #25	ES #33	ES #33
B	450 pmol	450 pmol	ES #2	ES #2	ES #10	ES #10	ES #18	ES #18	ES #26	ES #26	ES #34	ES #34
C	150 pmol	150 pmol	ES #3	ES #3	ES #11	ES #11	ES #19	ES #19	ES #27	ES #27	ES #35	ES #35
D	50 pmol	50 pmol	ES #4	ES #4	ES #12	ES #12	ES #20	ES #20	ES #28	ES #28	ES #36	ES #36
E	16.6 pmol	16.6 pmol	ES #5	ES #5	ES #13	ES #13	ES #21	ES #21	ES #29	ES #29	ES #37	ES #37
F	5.55 pmol	5.55 pmol	ES #6	ES #6	ES #14	ES #14	ES #22	ES #22	ES #30	ES #30	ES #38	ES #38
G	1.85 pmol	1.85 pmol	ES #7	ES #7	ES #15	ES #15	ES #23	ES #23	ES #31	ES #31	ES #39	ES #39
H	0 pmol	0 pmol	ES #8	ES #8	ES #16	ES #16	ES #24	ES #24	ES #32	ES #32	ES #40	ES #40

ES = PI(4)P extraction sample

4. Seal plate with plate sealer and incubate on plate shaker for 30 minutes at room temperature.
5. Transfer from Incubation Plate to PI(4)P Detection Plate.
 - a. From each well of the Incubation Plate transfer 100 µL to the corresponding well of the PI(4)P Detection Plate (K-4001E). For best results, mix each well 3-6 times with pipette before transferring.
 - b. Seal plate with plate sealer and incubate on plate shaker for 2 hours at room temperature.
6. Discard protein/lipid solution and wash PI(4)P Detection Plate 3 times with 200 µL/well PBS-T Buffer. Leave last wash in plate and proceed to next step.
7. Prepare and add Secondary Detector to PI(4)P

(K-4005) for a 10 µg/mL stock. Keep on ice for 1 minute to fully dissolve protein. Mix vial by flicking with finger. Do not vortex. The 10 µg/mL PI(4)P Detector stock solution can be stored for up to 3 months at -20°C.

- b. Further dilute the PI(4)P Detector by adding 40 µL of the 10 µg/mL stock (prepared in step 2a) to 7 mL Working Diluent. Keep at room temperature and proceed immediately to next step.
3. Set up the Incubation Plate using plate layout as guide
 - a. Add 60 µL/well prepared PI(4)P Standards (step 1b) to standard wells.
 - b. Add 60 µL/well prepared PI(4)P extraction samples (step 1c) to extraction sample wells.
 - a. Add 60 µL/well PBS-T Buffer to Blank and 0 pmol wells.
 - c. Add 60 µL/well prepared 0.05 µg/mL PI(4)P Detector (step 2b) to all wells except Blank control.
 - d. Add 60 µL/well working Diluent to Blank control.

Detection Plate.

- a. Add 200 µL of Secondary Detector (K-SEC2) and 3 mL Diluent to 9 mL PBS-T Buffer. Mix well.
 - c. Discard PBS-T Buffer from PI(4)P Detection Plate and add 100 µL/well prepared Secondary Detector (step 7a) to all wells of PI(4)P Detection Plate.
 - d. Seal plate with plate sealer and incubate on plate shaker for 1 hour at room temperature.
8. Discard Secondary Detector and wash PI(4)P Detection Plate 3 times with 200 µL/well PBS-T Buffer. Remove last wash from plate and proceed to next step.
 9. Develop plate
 - a. Add 100 µL/well TMB (K-TMB1) to PI(4)P Detection Plate and develop for 30 minutes at

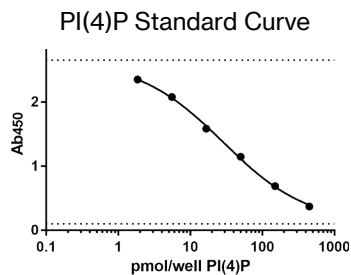
room temperature in a dark location.

- b. Stop reaction by adding 50 μL /well 1 N H_2SO_4 Stop Solution (K-STOPt) and read at 450 nm with an absorbance plate reader.

Data Analysis

Generate a best fit curve for the PI(4)P standards and interpolate relative sample values (See figure 1. for example). The final PI(4)P concentration, for each sample, should include any dilutions made to the sample before it is added to the Incubation Plate.

Figure 1
PI(4)P standard curve was generated using non-linear regression analysis with GraphPad software. A log(agonist) vs. response -- Variable slope (four parameters) analysis was utilized. For best results, constrain standard top and bottom using the 0 pmol PI(4)P and Blank control.



Support Protocol: Lipid Extraction

The PI(4)P extraction protocol was verified with 2×10^6 HL-60 cells. The number of cells necessary for PI(4)P quantification needs to be determined for each cell type. Larger or smaller amounts of cells require proportional adjustments of volumes. If you have never run lipid extractions or have little experience with the reagents listed below, please read the "PIP Mass FAQ: PIP Mass Customer Information Sheet" before running extractions. The FAQ can be found on the webpage of this product, or it can be requested at echelon@echelon-inc.com.

Solutions for Extraction

1. 0.5 M TCA: For 50 mL, dissolve 4.08 g TCA (Trichloroacetic Acid) in dH_2O and bring volume to 50 mL.
2. 5% TCA with 1 mM EDTA: For 50 mL, dissolve 2.5 g TCA in dH_2O , add 100 μL 0.5 M EDTA, and bring volume to 50 mL with dH_2O .
3. MeOH:CHCl₃ (2:1): For 60 mL, add 40 mL MeOH to 20 mL CHCl₃
 - a. Measure CHCl₃ with a glass pipette. Pure CHCl₃ may dissolve plasticware.
 - b. MeOH: CHCl₃ (2:1) should be prepared in an amber glass bottle. This solution is not stable long term and should be used within a month of preparation. It's safe to use plasticware to transfer this solution.
4. MeOH:CHCl₃:HCl (80:40:1): For 60 mL, combine 40 mL MeOH, 20 mL CHCl₃, and 0.5 mL 12 N HCl
 - a. Measure CHCl₃ with a glass pipette. Pure CHCl₃

may dissolve plasticware.

- b. MeOH:CHCl₃:HCl (80:40:1) should be prepared in an amber glass bottle. This solution is not stable long-term and should be used within a month of preparation. It's safe to use plasticware to transfer this solution.
 - c. Use 12 N concentrated 36% - 38% HCl. Do not use diluted acid.
5. 0.1 N HCl: For 50 mL, add 0.42 mL 12 N HCl to 50 mL dH_2O .

Extraction of PI(4)P from cells

1. Collect Cells
 - a. For adherent cells in a 75 cm² flask, remove medium by gentle aspiration and immediately add 5 mL ice cold 0.5 M TCA. Incubate cells on ice for 5 minutes. Scrape the cells from the flask with additional 0.5 M TCA if needed and transfer to a 15 mL centrifuge tube. Centrifuge at 3,000 RPM (approximately 900-1,000 RCF) for 7 minutes at 4°C. Discard the supernatant. The remaining steps are performed at room temperature.
 - b. For non-adherent cells in a 75 cm² flask, collect cells into 15 mL centrifuge tube, spin the cells down, decant media, add 5 mL ice cold 0.5 M TCA and vortex. Incubate cells on ice for 5 minutes. Centrifuge at 3000 RPM (approximately 900-1,000 RCF) for 7 minutes at 4°C. Discard the supernatant. The remaining steps are performed at room temperature.
2. Wash Pellet
Add 3 mL 5% TCA/ 1 mM EDTA to the pellet. Vortex for 30 seconds. Centrifuge at 3000 RPM for 5 minutes. Discard the supernatant. Repeat wash one more time.
3. Extract Neutral Lipids
Add 3 mL MeOH : CHCl₃ (2:1) and vortex for 10 minutes at room temperature. Centrifuge at 3000 RPM for 5 minutes, discard the supernatant. Repeat neutral lipids extraction one more time. A small white pellet should be visible after this step.
4. Extract Acidic Lipids
Add 750 μL MeOH:CHCl₃:1HCl (80:40:1) and vortex for 25 minutes at room temperature. Centrifuge at 3000 RPM for 5 minutes. Transfer supernatant to a new 2 mL vial. Discard pellet.
5. Phase Split
To supernatant from step 4, add 250 μL of CHCl₃ and 450 μL of 0.1 N HCl. Vortex for 30 seconds. Centrifuge at 3000 RPM for 5 minutes to separate organic and aqueous phases. Disregard any excess cellular debris that may appear between the two layers. Collect 0.5 mL of the organic (lower) phase with a positive displacement pipette, into a new 1.5 mL vial and dry in a

vacuum dryer (45-60 minutes). Dried lipid can be stored at -20°C for up to 12 months. The dried lipid should not be visible. If there is a visible substance at the end of this step, it is most likely cell debris that was not eliminated in the extraction. Do not attempt to dissolve the cell debris in the PBS-T buffer (step 1c). This black to yellow substance should also be avoided when pipetting the lipid extraction samples into the PI(4)P Mass ELISA wells.

References

PI(4)P Background References

1. Lemmon, M. A. 2008. Membrane recognition by phospholipid-binding domains. *Nat. Rev. Mol. Cell Biol.* 9,99-111.

Lipid Extraction Protocol References

2. Grey, H. Olsson, I. H. Batty, L. Priganica, and C. P. Downes, Nonradioactive methods for the assay of phosphoinositid 3-kinases and phosphoinositide phosphatases and selective detection of signaling lipids in cell and tissue extracts, *Analytical Biochemistry* 313(2003) 234-245.
3. Costa et al., Measurement of PIP3 Levels Reveals an Unexpected Role for p110β in Early Adaptive Responses to p110α-Specific Inhibitors in Luminal Breast Cancer, *Cancer Cell* (2015).

K-4000E References

4. Zhang L, Li J, Zhang P, Gao Z, Zhao Y, Qiao X, et al. PI4KIIα regulates insulin secretion and glucose homeostasis via a PKD-dependent pathway. *Biophysics Reports*. 2018:1-14.

5. Arellanes-Robledo J, Reyes-Gordillo K, Ibrahim J, Leckey L, Shah R, Raj Lakshman M. Ethanol Targets Nucleoredoxin/Dishevelled Interactions and Stimulates Phosphatidylinositol 4-phosphate Production In vivo and In vitro. *Biochemical pharmacology*. 2018.
6. Choi S, Hedman AC, Sayedyahosseini S, Thapa N, Sacks DB, Anderson RA. Agonist-stimulated phosphatidylinosi-tol-3,4,5-trisphosphate generation by scaffolded phosphoinositide kinases. *Nat Cell Biol.* 2016;16(12):1324-35.
7. Choi JY, Park S. Role of protein kinase A and class II phosphatidylinositol 3-kinase C2I² in the downregulation of KCa3.1 channel synthesis and membrane surface expression by lyso-globotriaosylceramide. *Biochemical and biophysical research communications*. 2016.
8. Esteban-Pretel G, MarÃn MP, Romero AM, Timoneda J, Ponsoda X, BallestÃn RI, et al. Polyphosphoinositide Metabolism and Golgi Complex Morphology in Hippocampal Neurons in Primary Culture is Altered by Chronic Ethanol Exposure. *Alcohol and Alcoholism*. 2012.
9. Reiss S, Rebhan I, Backes P, Romero-Brey I, Erfle H, Matula P, et al. Recruitment and activation of a lipid kinase by hepatitis C virus NS5A is essential for integrity of the membranous replication compartment. *Cell Host Microbe*. 2011;9(1):32-45.
10. Berger KL, Kelly SM, Jordan TX, Tartell MA, Randall G. Hepatitis C virus stimulates the phosphatidylinositol 4-kinase III alpha-de-pendent phosphatidylinositol 4-phosphate production that is essential for its replication. *Journal of virology*. 2011.

Related Products

Products	Catalog Number
Other PIP Mass Assays	
PIP ₃ Mass ELISA	K-2500s
PI(3)P Mass ELISA	K-3300
PI(3,4)P ₂ Mass ELISA	K-3800
PI(4,5)P ₂ Mass ELISA	K-4500
Activity Assays	
PI3-Kinase Activity ELISA	K-1000s
PTEN Activity ELISA	K-4700
PI(4)P 5-Kinase Activity Assay	K-5700
PI(4)P Binding Proteins and Antibodies	
PI(4)P Grip	G-0402
Purified Anti-PtdIns(4)P IgM	Z-P004

Please visit our website at www.echelon-inc.com for more phospholipid products.