

PI(4)P 5-Kinase (PIP5KI) Activity Assay

K-5700 (96 wells)

Support: echelon@echelon-inc.com

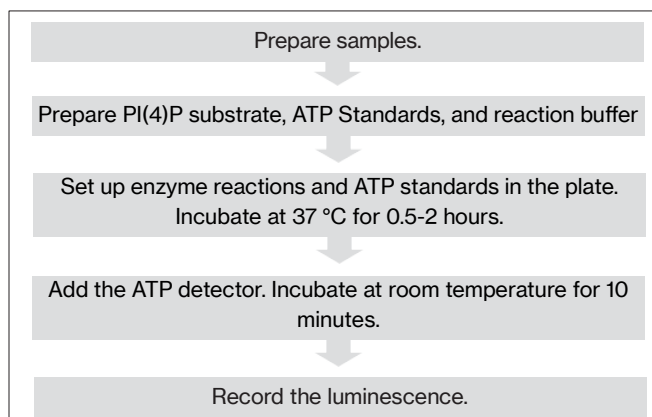
Description: The PI(4)P 5-Kinase (PIP5K1) Activity Assay is an ATP depletion assay which quantifies the remaining ATP levels in solution following the kinase reaction. The assay can be used to quantify PIP5K1 activity in vitro or in cell lysate.

Storage: Upon receipt, store at -20°C. Under proper storage conditions, the kit components should remain stable for at least 6 months from date of receipt. Allow the reagents to warm to room temperature before opening vials.

Materials Provided

Catalog #	Description	Quantity
K-5701	PI(4)P Substrate (522 µmol)	1 vial
K-KBZ	5x KBZ Buffer (4 mL)	1 bottle
K-5403	Diluent (150 µL)	1 vial
K-CF01	Kinase Cofactor (100 µL)	1 vial
K-LUMa	ATP Detector (5 mL)	1 bottle
K-ATP1	10 mM ATP (50 µL)	1 vial
K-DTT1	50 µmol dried dithiothreitol (DTT) powder	1 vial
---	96-well round bottom white plate	1 each
---	Microtiter Plate Seal	1 each

Quick Protocol



Additional Materials Provided by User

- Source of purified PI(4)P 5-Kinase as positive control (optional)
- Plate reader capable of reading luminescence in a 96-well microtiter plate
- Microcentrifuge tubes (0.5-1.5 mL)
- Multichannel pipettes

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 exposures 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

Type I PI(4)P 5-Kinase (PIP5KI) catalyzes the phosphorylation of phosphatidylinositol-4-phosphate [PI(4)P] to phosphatidylinositol-4,5-bisphosphate [PI(4,5)P₂]. PIP5KI has been shown to play a role in several cellular functions including gene transcription, actin cytoskeleton reorganization, secretion and endocytosis. It is thought that PIP5KI activity is responsible for the majority of the PI(4,5)P₂ found in the cell which is an important second messenger for many cellular processes.

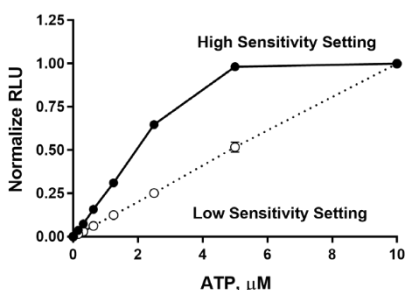
Assay Design

Echelon's PI(4)P 5-Kinase Activity Assay is an ATP depletion assay which quantifies the remaining ATP levels in solution following the kinase reaction. The assay reaction is prepared by the addition of PIP5KI (enzyme), PI(4)P (substrate) in reaction buffer and initiated by the addition of ATP. The reaction is then stopped by adding the ATP detector after the chosen reaction time. The luminescent signal is inversely correlated with the kinase activity. Echelon's PI(4)P 5-Kinase Activity Assay is tested using SignalChem's PIP5KIA (P16-10AG) kinase.

Assay Kit Notes

1. The ATP detector's (K-LUMa) linear range of detection may vary between instruments. Depending on the instrument capacity and sensitivity setting, the optimal range and sensitivity for the ATP detector (K-LUMa) is up to 100 μ M (Figure 1).

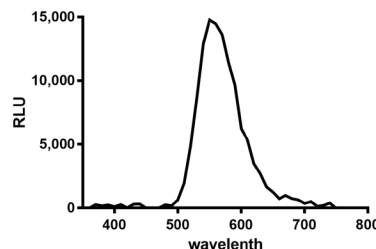
Figure 1 | Detection System Linearity



2. The luminescent signal, or light, collected by your plate reader is not restricted to a particular wavelength. When setting up your instrument, select the option on your reader that allows the instrument to collect all wavelengths. For most plate readers this setting is labeled "all wavelengths". For BMG Labtech plate readers the "lens" setting will read all wavelengths. Please refer to your plate reader's manual and/or the manufacturer of your reader if you have questions on how to set up your specific instrument. If your instrument does not have this setting, or is limited to a certain filter, choose a broad filter centered at 550 nm

(Figure 2). When the luminescent signal is read at one wavelength, a lower signal may be observed.

Figure 2 | Wavelength Emission Spectra of ATP detector; K-LUMa



3. The buffer of your sample may interfere with the luminescent output. Sample buffers should always be included in the ATP standard curve and as a "buffer only" background control.
4. Two commonly used lysis buffers (figure 3) and common lysis buffer components (figure 4) were tested in the assay. Since some buffers and lysis buffer components affect the assay, we suggest using RIPA lysis buffer. Alternatively, cells can be lysed with sonication and freeze thaw cycles in the complete reaction buffer (step 3 of assay protocol). If you plan to use your own lysis buffer, we suggest testing it in the assay for interference before running samples. If your lysis buffer has a strong effect on the assay, substitute lysis buffer or dilute sample with complete reaction buffer.

Figure 3 | Lysis Buffer Effects

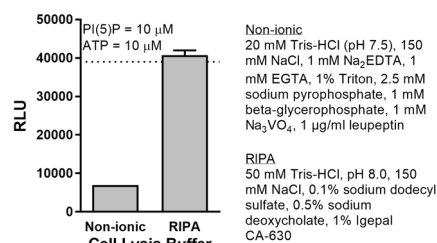
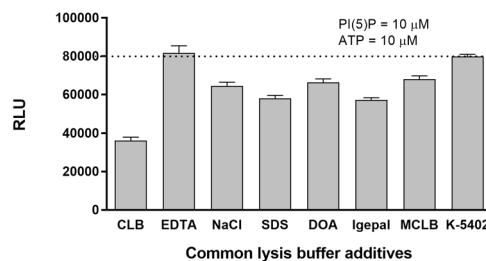


Figure 4 | Lysis Buffer Effects



5. Optimization of the kinase reactions by enzyme concentration, reaction time and temperature may be required.
6. Kinase reactions can be run at either 37°C or at room temperature. Reactions run at 37°C should be allowed

- to equilibrate to room temperature before the ATP detector is added to avoid a temperature gradient.
- The provided 5X KBZ Buffer (K-KBZ) and supplements (DTT, Kinase Cofactor) must be used for enzyme reactions.
 - The amount of kinase to use per reaction will vary according to your individual experiment. Whether you are using purified enzyme or cell lysate, you will need to test reactions using different amounts of enzyme to determine the optimum conditions. When using purified PIP5KI from SignalChem (Cat# P16-10AG), enzyme concentrations of 0.5-2 ng/ μ L are suggested as a starting point. In testing, we found that 5-10 μ g of cell lysate was sufficient to detect significant kinase activity.
 - Sample Reaction: 10 μ L samples (PIP5KI) + 10 μ L PI(4)P + 20 μ L ATP = 40 μ L reaction.
ATP Standards: 10 μ L complete reaction buffer or sample buffer of your choice + 10 μ L PI(4)P + 20 μ L ATP standard dilution = 40 μ L solution.

Assay Protocol

Please read this entire section and the assay notes section before beginning the assay.

- Bring all reagents to room temperature before use. Place enzyme samples and ATP on ice until use. All steps are performed at room temperature unless noted.
- Add 50 μ L deionized water to 50 μ mol dried DTT (K-DTT1). Vortex to mix. This is the 1 M DTT stock. Keep on ice. The 1 M DTT stock is stable for at least 1 month when stored at -20 $^{\circ}$ C.
- Prepare the complete reaction buffer by diluting the 5X KBZ Buffer (K-KBZ) to 1X and supplement with 1:1000 dilution of DTT stock and 1:500 dilution of Kinase Cofactor (K-CF01). Dilute only what is needed. The 1X KBZ buffer, DTT and Kinase Cofactor are not stable as complete reaction buffer.
- Prepare a 5 mM stock of PI(4)P Substrate by adding 105 μ L of the provided diluent (K-5403) to the vial of PI(4)P Substrate (K-5701). Vortex solution for a full two minutes followed by a 5 minute incubation at 37 $^{\circ}$ C to fully resuspend the lipid. From the 5 mM stock prepare a 4X PI(4)P working solution by diluting the 5 mM PI(4)P stock in the complete reaction buffer (step 3). Suggested PI(4)P concentration is 100 μ M (4X at 400 μ M).
- Prepare a 2X ATP solution by diluting the 10 mM ATP stock (K-ATP1) in the complete reaction buffer (step 3). Suggested ATP concentration is 10 μ M (2X at 20 μ M).

- Prepare a 2-fold dilution series of ATP standards by diluting 10 mM ATP stock (K-ATP1) in 0.5 mL centrifuge tubes. A sample of a 10 μ M ATP 2X dilution standard curve is shown below (Table 1).

Table 1

Final ATP Conc. in reaction (1X)	2 X ATP Solutions	ATP Stock (K-ATP1) or previous dilution	Complete Reaction Buffer (step 3)
10 μ M	20 μ M	1 μ L of K-ATP1	499 μ L
5 μ M	10 μ M	100 μ L of 20 μ M soln.	100 μ L
2.5 μ M	5 μ M	100 μ L of 10 μ M soln.	100 μ L
1.25 μ M	2.5 μ M	100 μ L of 5 μ M soln.	100 μ L
0.625 μ M	1.25 μ M	100 μ L of 2.5 μ M soln.	100 μ L
0.3125 μ M	0.625 μ M	100 μ L of 1.25 μ M soln.	100 μ L
0.15625 μ M	0.3125 μ M	100 μ L of 0.625 μ M soln.	100 μ L

- Using the provided 96-well plate, add 10 μ L of the 4X PI(4)P solution per well. Then, add 10 μ L of the kinase samples per well. For the ATP standard wells, add 10 μ L of the complete reaction buffer or sample buffer. Suggested plate layout in duplicates is shown below (Table 2).
- Start the reaction by adding 20 μ L of the 2X ATP solution to the kinase sample wells. Add 20 μ L of the 2-fold ATP dilution series (prepared in step 6) to the standard wells. Cover with plate seal. Mix briefly by tapping the plate or by plate shaker. Incubate the reaction at 37 $^{\circ}$ C. Incubation times may vary depending on the PIP5KI activity within samples. Suggested incubation time – 30 minutes to 2 hours. Note – The last 10 minutes of the incubation should be done at room temperature to allow the plate temperature to equilibrate.
- After incubation, add 40 μ L of ATP Detector (K-LUMa) per well. Mix briefly by tapping the plate or by plate shaker. Incubate at room temperature for at least 10 minutes to stabilize the luminescent signal. Protect from light.
Note: if K-LUMa shows precipitate centrifuge at 1,000 RPM for 15 minutes.

Data Analysis

Kinase activity can be determined by quantifying the remaining ATP levels in the reaction buffer. An example of an ATP standard curve and PIP5KI titration are shown below.

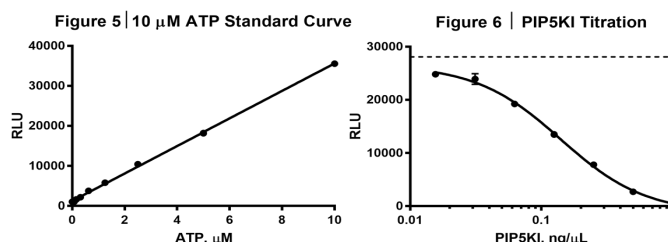


Table 2

	1	2	3	4	5	6	7	8	9	10	11	12
A	10 μ M ATP	10 μ M ATP	Sample 1	Sample 1	Sample 9	Sample 9	Sample 17	Sample 17	Sample 25	Sample 25	Sample 33	Sample 33
B	5 μ M ATP	5 μ M ATP	Sample 2	Sample 2	Sample 10	Sample 10	Sample 18	Sample 18	Sample 26	Sample 26	Sample 34	Sample 34
C	2.5 μ M ATP	2.5 μ M ATP	Sample 3	Sample 3	Sample 11	Sample 11	Sample 19	Sample 19	Sample 27	Sample 27	Sample 35	Sample 35
D	1.25 μ M ATP	1.25 μ M ATP	Sample 4	Sample 4	Sample 12	Sample 12	Sample 20	Sample 20	Sample 28	Sample 28	Sample 36	Sample 36
E	0.63 μ M ATP	0.63 μ M ATP	Sample 5	Sample 5	Sample 13	Sample 13	Sample 21	Sample 21	Sample 29	Sample 29	Sample 37	Sample 37
F	0.31 μ M ATP	0.31 μ M ATP	Sample 6	Sample 6	Sample 14	Sample 14	Sample 22	Sample 22	Sample 30	Sample 30	Sample 38	Sample 38
G	0.16 μ M ATP	0.16 μ M ATP	Sample 7	Sample 7	Sample 15	Sample 15	Sample 23	Sample 23	Sample 31	Sample 31	Sample 39	Sample 39
H	0 μ M ATP	0 μ M ATP	Sample 8	Sample 8	Sample 16	Sample 16	Sample 24	Sample 24	Sample 32	Sample 32	Sample 40	Sample 40

Support Protocol

Cell Lysate Preparation

The following support protocol for cell lysate preparation has been validated for use in the PIP5KI Activity Assay. Further optimization of this protocol and/or different protocols can be used depending on the needs and experience of the user.

Materials Needed	Company	Catalog Number
1.5 mL Centrifuge Tubes	N/A	N/A
RIPA Lysis Buffer (50 mM Tris-HCl pH 8.0, 150 mM NaCl, 0.1% SDS, 1% Igepal CA-630, 0.5% Na deoxycholate)	Sigma	R0278
Protease Inhibitor Cocktail	Sigma	P8340
PBS	N/A	N/A

This protocol is written for a 100 mm dish of HEK 293 cells (90% confluent) for about 10-15 mg/mL of cellular protein. To maximize kinase activity; keep all solutions ice cold, carry out all reactions on ice or at 4 °C, and use a centrifuge that is equilibrated at 4°C.

Reagent Preparation

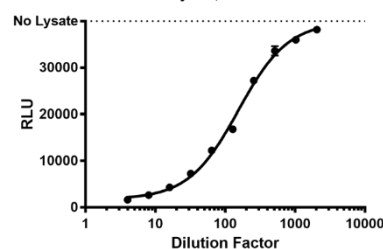
Lysis Buffer: Use a commercially available RIPA Lysis Buffer or prepare according to the recipe listed in the materials. Add the protease inhibitor cocktail fresh at a 1:100 dilution. Place buffer on ice and chill until ice cold.

Cell Lysis

- Place cell culture dish on ice and wash cells twice with ice cold PBS.
- Add 0.5 mL of ice cold Lysis Buffer to cells. Scrape cells and transfer mixture into a cooled 1.5 mL centrifuge tube.

- Incubate cells for 15 minutes with constant agitation at 4°C.
- Centrifuge cells for 10 minutes at 14,000 x g to pellet cells.
- Transfer supernatant to a fresh, cooled 1.5 mL centrifuge tube and freeze immediately at -80°C or proceed with the kinase reactions.

Figure 7 | Cell Lysate Titration
HEK 293 Lysate, 2 Hour Incubation



Note: ATP, DTT, Kinase Cofactor and PI(4)P Substrate levels should be kept consistent when setting up kinase reactions containing cell lysate. Depending on how much lysate is added to the reaction it may be necessary to supplement the lysate with these components.

Related Products

Products	Catalog Number
Substrates	
diC16 PI(4)P	P-4016
diC8 PI(4)P	P-4008
Assays	
PI(5)P 4-Kinase Activity Assay	K-5400
PI3-Kinase Activity ELISA: Pico	K-1000s
PTEN Activity ELISA	K-4700