

Class III PI3-Kinase Kit

K-3000 (96 tests)

Support: echelon@echelon-inc.com

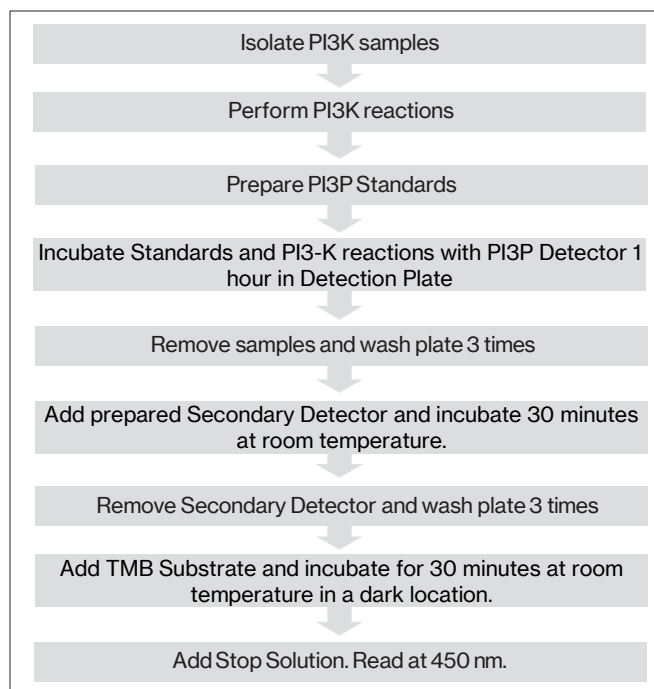
Description: 96-well ELISA Assay for Detection of PI(3)P

Storage: Store kit part 1 at 4°C. Store kit part 2 at -20°C.

Materials Provided

Catalog #	Description	Quantity
K-3001	PI(3)P Detection Plate, 12 x 8-strip well	1 plate
K-3002	PI Substrate, diC8, MW = 608.6 g/mol	2 vials
K-3003	PI(3)P Standard, diC8, MW = 732.5 g/mol	1 vial
K-3004	5x PI(3)P Detection Buffer (5x DB)	1 bottle
K-EDTA	100 mM EDTA, pH 8	1 vial
K-TBST	10x TBS-T Buffer	1 bottle
K-3305	PI(3)P Detector	2 vials
K-DIL3	5x Diluent	1 vial
K-SEC2	Secondary Detector	1 vial
K-TMB1	TMB Solution	1 bottle
K-STOPt	1 N H2SO4 Stop Solution	1 bottle
---	Colored 96-well U-bottom plate	1 plate
---	Clear acetate sheet, 1 side adhesive	3 sheets

Quick Protocol



Additional Materials Provided by User

- Source of purified PI3-Kinase
- 2X Kinase reaction buffer (see Assay Note 1)
- Absorbance plate reader capable of reading at 450 nm
- Plate shaker

This kit and all non-radioactive, competitive assays for determining phosphoinositide-3-kinase (PI3-K) activity are protected by Echelon Biosciences Inc. U.S. Patent 7,067,269. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. For inquiries, email echelon@echelon-inc.com.

Background

PI3-Kinases (PI3-K) are grouped into three classes according to their structural homology and in vivo lipid substrate preference. Class III PI3-K uses phosphatidylinositol (PI) as substrate to generate PI(3)P product. While Class I PI3-Ks are well-recognized for their role in cell growth and division, the physiological roles of the Class II and III enzymes are still emerging. Typically, experiments to measure PI3-K activity have involved phosphorylation of a phosphoinositide substrate using 32P, then extraction of radioactive products, and separation using thin-layer chromatography or HPLC. The assay plate method developed by Echelon Biosciences allows the user to determine PI3-K activity, using either recombinant or immunoprecipitated enzyme, in a standard ELISA format, eliminating the need for radioactivity, and thin layer chromatography or HPLC.

Assay Design

Echelon's Class III PI3K Activity Assay is a competitive ELISA in which the signal is inversely proportional to the amount of PI(3)P produced. After the PI3-K reactions are complete and quenched, reaction products are diluted and added to the PI(3)P-coated microplate, for competitive binding to a PI(3)P detector protein. The amount of PI(3)P detector bound to the plate is determined through colorimetric detection. This ELISA is specific to PI(3)P detection. As low as 1 pmol PI(3)P can be detected in a 100 μ L detection mixture, corresponding to a sensitivity of 10 nM.

Assay Notes

1. The composition of the 2x kinase reaction buffer should be determined by the user. Two sample recipes for 2x kinase reaction buffer are listed below for human Vps34 (PIK3C3). For the best result, prepare the 2x kinase reaction buffer freshly on day of assay. To offset potential interference, make sure to dilute PI(3)P standards with Detection Buffer Blend (DBB) as shown in assay protocol.
 - a. Recipe #1: 100 mM HEPES pH 7.5, 300 mM NaCl, and 2 mM CHAPS, 10 mM MnCl₂, 2 mM DTT and 100 μ M ATP.
 - b. Recipe #2: 20 mM Tris pH 8, 200 mM NaCl, 2 mM EDTA, 20 mM MnCl₂, and 100 μ M ATP.
2. The user should optimize the amount of PI3-Kinase by doing an enzyme titration. For purified enzyme, we suggest 5 nM PI3-K enzyme as a starting point for optimization.
3. Increase the PI Substrate if low or no PI(3)P product is detected. We suggest including a reaction without PI3-K enzyme as a "No Enzyme" control; and a reaction without PI as a "No Substrate" control.
4. It's important to dilute PI3-K reaction 8-fold into 2x PI(3)P detection buffer. Greater concentrations of reaction buffer will interfere with the PI(3)P detection.
5. Unused strip wells should be removed from the plate, sealed in a plastic bag, and stored at 4°C for later use.
6. Never let the detection plate dry out after the assay has started. Always prepare the next solution before discarding the current one from the detection plate.

Reagent Preparation

Place a vial of PI(3)P Detector (K-3305) and the vial of Secondary Detector (K-SEC2) on ice. Bring PI Substrate (K-3002), PI(3)P Standard (K-3003), Detection Plate (K-3001) and other reagents to room temperature prior to use.

1. Prepare 2 mL Detection Buffer Blend (DBB) by adding 800 μ L of 5x PI(3)P Detection Buffer (K-3004), 50 μ L of 100 mM EDTA (K-EDTA), and 125 μ L of 2x Kinase Reaction Buffer to 1,025 μ L ddH₂O.
2. Combine entire bottle of 10x TBS-T buffer (K-TBST) with 180 mL ddH₂O for 200 mL TBS-T buffer.
3. Prepare 1x Diluent for entire plate, dilute 1.4 mL of 5x Diluent (K-DIL3) with 5.6 mL ddH₂O. Leave 1x Diluent at room temperature prior to use. Prepare only the volume needed for the day of assay.

Assay Protocol

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

Kinase Reaction

1. Isolate or prepare PI3-Kinase according to user's standard protocol. For human cells and tissue samples, it's recommended to use the anti-hVps34 antibody (cat# Z-R015) to immunoprecipitate human Vps34 enzyme. Please refer to Z-R015 Technical Data Sheet (TDS) for the IP protocol.
2. Prepare a 500 μ M PI stock by adding 99 μ L of ddH₂O to a vial of 30 μ g PI Substrate (K-3002). Vortex the vial at maximum speed for 30 seconds. Spin down vial. Place vial at room temperature. After use, store PI stock solution at -20°C for up to 3 months.
3. We suggest the following for setting up PI3-K reactions in the colored incubation plate:
 - a. For each PI3-K reaction, add the following to a single well of the 96-well incubation plate:
 - i. 4 μ L of 500 μ M PI substrate (2 nmol, see assay note 3)
 - ii. 12.5 μ L of 2x kinase reaction buffer (see assay note 1).
 - iii. Enough ddH₂O to bring the total volume (including PI3-Kinase) to 25 μ L
 - iv. Volume of PI3-Kinase to start the reaction. Seal the plate with a plate sealer (see assay note 2).

Note: If using immunoprecipitated enzyme bound to beads (in a centrifuge tube), add 12.5 μ L of 2x kinase reaction buffer to the beads, 8.5 μ L ddH₂O, then add 4 μ L of 500 μ M PI substrate to start the reaction. Scale up the reaction volume to cover all the beads if needed.

4. Let the kinase reaction proceed. Reaction conditions are to be determined by the user. Typically, reactions are 0.5-3 hours at desired temperature: room temperature, 30°C or 37°C. Do not shake vigorously.
5. Quench the kinase reactions by adding 5 μ L of 100 mM EDTA (K-EDTA) to each well of 25 μ L reaction. Then dilute each quenched reaction with 90 μ L

ddH₂O. If the enzyme is bound to beads or otherwise immobilized, the quenched enzyme reaction (120 µL) can be separated from the beads with centrifugation before being transferred to the colored incubation plate for ELISA. These reactions can be stored at -20°C.

6. Add 80 µL of 5x PI(3)P Detection Buffer (K-3004) to each 120 µL reaction mixture to bring the final volume to 200 µL ([PI] + [PI(3)P] = 10 µM total lipid in Detection Buffer Blend). This will provide enough sample for triplicate assay points.

ELISA Assay

1. Prepare PI(3)P Standard solutions.
 - a. Add 410 µL of DBB to the vial of PI(3)P Standard (K-3003) for a concentration of 10 µM. Vortex at maximum speed for 30 seconds. Spin down. Place the vial at room temperature. Store unused portion of 10 µM PI(3)P at -20°C for up to 2 weeks.
 - b. Further dilute 10 µM PI(3)P 4-fold serially in DBB for concentrations of 2.5 µM through 9.8 nM, respectively. Refer to Table 1.
2. Reconstitute PI(3)P Detector (K-3305) by adding 200 µL 1x Diluent directly to a vial of PI(3)P Detector. Place the vial of reconstituted detector on ice for a few minutes before use in step 4. Do not vortex PI(3)P Detector. Proceed immediately to the next step. Reconstituted PI(3)P Detector is not stable. Do not retain stock for second use.

Table 1: PI(3)P Standards

[PI(3)P]	Dilution Factor	PI(3)P, Volume and Source	DBB	PI(3)P per 50 µL
10 µM	-	10 µM	-	500 pmol
2.5 µM	4x	60 µL of 10 µM	180 µL	125 pmol
625 nM	4x	60 µL of 2.5 µM	180 µL	31.25 pmol
156 nM	4x	60 µL of 625 nM	180 µL	7.81 pmol
39.1 nM	4x	60 µL of 156 nM	180 µL	1.95 pmol
9.8 nM	4x	60 µL of 39.1 nM	180 µL	0.49 pmol

Table 2: Suggested Detection Plate Layout

Row	Standards and Controls			PI3K Reaction Samples (from step 3 of Kinase Reaction Section)								
	1	2	3	4	5	6	7	8	9	10	11	12
A	500 pmol PI(3)P			Reaction #1			Reaction #9			Reaction #17		
B	125 pmol PI(3)P			Reaction #2			Reaction #10			Reaction #18		
C	31.25 pmol PI(3)P			Reaction #3			Reaction #11			Reaction #19		
D	7.81 pmol PI(3)P			Reaction #4			Reaction #12			Reaction #20		
E	1.95 pmol PI(3)P			Reaction #5			Reaction #13			Reaction #21		
F	0.49 pmol PI(3)P			Reaction #6			Reaction #14			Reaction #22		
G	NL (No Lipid)			Reaction #7			Reaction #15			Reaction #23		
H	Blank			Reaction #8			Reaction #16			Reaction #24		

3. Add PI(3)P standards and PI3-K samples to Detection Plate (K-3001). See table 2, suggested plate layout.
 - a. Add 50 µL DBB to No Lipid (NL) wells, G1-G3 and to Blank wells H1-H3.
 - b. Add 50 µL PI(3)P standard dilutions (step 1) to wells 1-3 of rows A through F.
 - c. Add 50 µL of each PI3-K reaction mixture from incubation plate (or centrifuge tube) to triplicate wells. All wells in use should have 50 µL.
4. Add PI(3)P Detector to Detection Plate (K-3001)
 - a. Invert the vial of reconstituted PI(3)P detector (from step 2) 10 times to mix. Spin down vial. Pipet up and down multiple times then dilute. For 96-wells use 190 µL of the detector stock solution with 5.7 mL of 1x Diluent. Only prepare what is needed for the day of use. Mix by inverting the tube. Do not vortex.
 - b. Add 50 µL of diluted PI(3)P detector to all wells except blank wells, H1-H3. All wells in use should have 100 µL.
 - c. Add 50 µL 1x Diluent to Blank wells, H1-H3.
 - d. Seal plate and incubate on a plate shaker (250-300 rpm) at room temperature for exactly 1 hour. Caution: Prolonged incubation time will reduce the assay signal strength.
5. Discard solution from Detection Plate and wash plate 3 times with TBS-T 200 µL/well.
6. Dilute Secondary Detector (K-SEC2) 1:100 with TBS-T. Prepare enough for current assay only. For entire plate, dilute 120 µL Secondary Detector with 12 mL TBS-T and mix gently by inverting the tube 10 times. Discard the last TBS-T wash from plate and add 100 µL diluted Secondary Detector to each well. Seal plate and incubate for 30 minutes at room temperature on a plate shaker.

