

PI3-Kinase Fluorescence Polarization Activity Assay

K-1100 (384 tests)

Support: echelon@echelon-inc.com

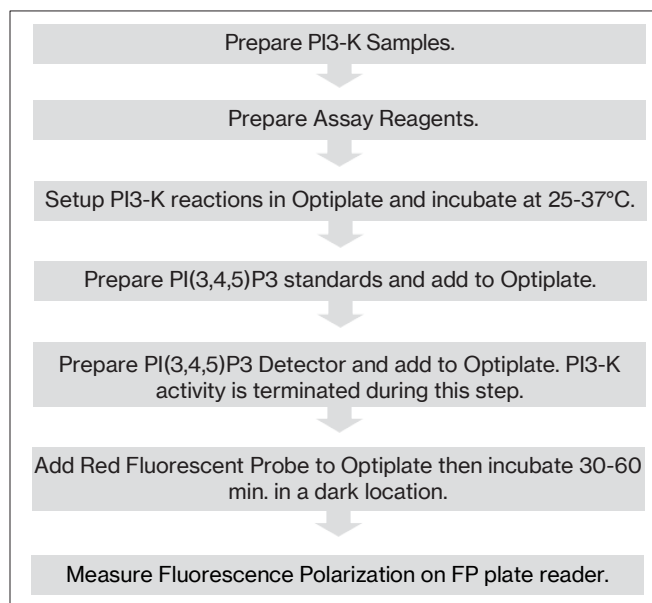
Description: The PI3-Kinase Activity Fluorescence Polarization Assay determines PI3-K activity through quantification of its product, PI(3,4,5)P3 (PIP3).

Storage: Upon receipt, store the assay kit at -20°C . Some components are light sensitive. Under proper storage conditions, this product is stable for at least 6 months from date of receipt. Opened and reconstituted solutions are less stable. All components and solutions should be protected from excessive light and heat.

Materials Provided

Catalog #	Description	Quantity
K-1101L	PI(4,5)P2 Substrate	69 μg
K-1003s	PI(3,4,5)P3 Standard	0.88 μg
K-1103	PI(3,4,5)P3 Detector	1 pellet
K-1104	Red Fluorescent Probe	50 μL , 2.5 μM
K-1105	Detector Diluent	10 mL
K-KBZ	5X KBZ buffer	4 mL
K-ATP1	10 mM ATP	50 μL
K-DTT1	50 μmol DTT (DL-dithiothreitol)	1 pellet
---	OptiPlate-384 F	1 plate
---	Plate Sealers	1 seal

Quick Protocol



Additional Materials Provided by User:

- Source of immunoprecipitated or purified PI 3-Kinase. The use of crude lysates is not recommended.
- Fluorescence plate reader equipped for Fluorescence Polarization using red fluorophores with appropriate filters (550 nm excitation/580 nm polarizing emission filters)

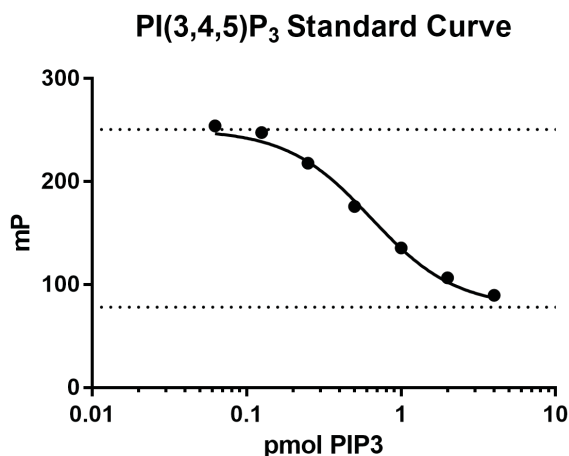
Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. This kit and all non-radioactive, competitive assays for determining phosphoinositide-3-kinase (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. We welcome inquiries about licensing the use of our trademarks and technologies at echelon@echelon-inc.co

Background

The production of PI(3,4,5)P3 from PI(4,5)P2 by PI3-Kinases (PI3-K) is important in multiple cell signaling pathways. Typically, experiments to measure PI3-K activity have involved phosphorylation of a phosphoinositide substrate using 32P, extraction of radioactive products, and separation using thin-layer chromatography. The fluorescence polarization PI3-K activity assay developed by Echelon Biosciences, Inc. allows the user to determine PI3-K activity using a homogenous mix and read format, eliminating the need for radioactivity, organic solvents, and thin layer chromatography. The assay can be adapted for HTS applications.

Assay Design

The assay is a competitive assay in which the degree of polarization (mP) of the fluorescent PI(3,4,5)P3 probe is inversely proportional to the amount of PI(3,4,5)P3 produced by PI3-K activity. After the PI3-K reactions are complete, reaction products are mixed with a PI(3,4,5)P3 detector protein and the fluorescent PI(3,4,5)P3 probe. Polarization (mP) values decrease as probe binding to the PI(3,4,5)P3 detector is displaced by PI(3,4,5)P3 produced by enzymatic activity and the amount of unbound fluorescent probe in the mixture increases. The graph (below) shows a PI(3,4,5)P3 standard curve using sigmoidal dose-response (variable slope) curve fit. The Top and Bottom of the standard curve have been constrained to Enzyme Only control and Probe Alone control signals, respectively.



Assay Notes

1. We suggest running duplicate or triplicate assay points for each enzyme reaction.
2. The 5x KBZ buffer has been tested with PI3-K isoforms (α , β , γ) and has shown to improve PI3-K activity. If you have your own PI3-K reaction buffer, it can be substituted in the assay where the protocol says "Reaction Buffer".
3. The amount of enzyme to use per 100 pmol of substrate will vary according to your individual experiment. Whether you are using purified PI3-K or enzyme immunoprecipitated from cell extracts, you

will need to try reactions using different amounts of enzyme to determine the optimum amount. In our testing, we found 10 ng of purified recombinant PI3-K α per point, or enzyme immunoprecipitated from cell lysates containing 5 ug cellular protein (approximately 50,000 cells), to be sufficient. A support protocol for immunoprecipitation of PI3-K from cell lysates is provided below.

4. The assay was developed using a Fusion Alpha Universal Microplate reader equipped for Fluorescence Polarization. The sensitivity of the assay and the amount of substrate, detector, and fluorescent probe required for each assay point may vary depending on the specific fluorescence polarization detection system you are using.
5. The provided OptiPlate-384 F assay plate works well with PI3K reactions. If you choose to use a different plate, please test PI3K reactions in the assay to make sure the well surface will not inhibit PI3-K reactions. Corning plates with NBS surface are not recommended.
6. The concentrations of DTT, ATP, and PI(4,5)P2 substrate suggested are based on our experience using recombinant PI3K α . The assay conditions used in your enzyme reaction can affect your enzyme activity and the activity of a potential inhibitor. You may want to titrate DTT, ATP, and/or PI(4,5)P2 substrate to determine the optimum conditions for your experiments.
7. Suggested controls include:
 - Buffer Only: 5 μ L of 2X Reaction Buffer, 5 μ L of enzyme diluent, 15 μ L of Detector Diluent. This control may be required as FP Blank for plate reading in FP mode.
 - Probe Alone: 5 μ L of 2X Reaction Buffer, 5 μ L of enzyme diluent, 10 μ L of Detector Diluent, and 5 μ L of 50 nM Probe.
 - No enzyme: 5 μ L of 2X Reaction Buffer, 5 μ L of enzyme diluent, 10 μ L of PI(3,4,5)P3 Detector, and 5 μ L of 50 nM Probe.

Reagent Preparation

Place PI(3,4,5)P3 Detector (K-1103), Red Fluorescent Probe (K-1104), ATP (K-ATP1), and DTT (K-DTT1) on ice. Equilibrate lipids (K-1101L and K-1003s) to room temperature before use. Leave buffers (K-KBZ, K-1105) and assay plate at room temperature. Unless indicated otherwise, all steps are performed at room temperature. Centrifuge each vial briefly prior to opening its cap.

- **PI(3,4,5)P3 Standard (K-1003s)** Add 22.4 μ L of dH2O to the vial of PI(3,4,5)P3 for a 40 μ M solution; vortex to fully reconstitute; spin down and leave vial at room temperature prior to use. Enough standard is provided for approximately 10 separate dilution series. Store frozen at -20°C after use for up to 6 months. Multiple freeze-thaw cycles do not affect stability.
- **DTT (K-DTT1)** Add 50 μ L of dH2O to the vial of DTT for a 1 M solution and vortex to fully

reconstitute; spin down and keep on ice. Once reconstituted DTT is less stable, store frozen at –20°C after use for up to 1 month.

- **ATP (K-ATP1)** One vial with 50 µL of 10 mM ATT. Store frozen at –20°C after use.
- **PI(4,5)P2 Substrate (K-1101L)** Add 100 µL of dH2O to the vial of PI(4,5)P2 for a 800 µM solution; vortex to fully reconstitute; spin down and leave vial at room temperature prior to use. There is substrate for up to 800 assay points using 100 pmol per assay point (Final concentration in reaction: 10 µM). Store frozen at –20°C after use for up to 6 months.
- **2X Reaction Buffer** Dilute 5x KBZ buffer (K-KBZ) 2.5-fold in dH2O and supplement with 4 mM DTT, 20 µM PI(4,5)P2 substrate, and 50 µM ATP. Only prepare the amount needed. Make it fresh before each use. Store remaining 5x KBZ at 4°C. For 2 mL of 2X Reaction Buffer add:
 - 800 µL 5x KBZ Buffer
 - 8 µL 1M DTT
 - 10 µL 10 mM ATP
 - 1,132 µL dH2O
 - 50 µL 800 µM PI(4,5)P2 solution
- **1X Reaction Buffer** Prepare a 1X reaction buffer for preparation of the PI(3,4,5)P3 standards by adding 250 µL of the 2X Reaction Buffer to 250 µL of dH2O or your enzyme diluent.
- **PI(3,4,5)P3 Detector (K-1103)** Each pellet contains enough PI(3,4,5)P3 detector for 400 assay points. Reconstitute with 4.25 mL of Detector Diluent (K-1105) for a 250 nM stock. Do not vortex. Mix gently by pipetting and keep solution on ice. Aliquot and store frozen at –80°C. The reagent is stable for several weeks through multiple freeze-thaw cycles.
- **Red Fluorescent Probe (K-1104)** Prior to use, dilute 2.5 µM Red Fluorescent Probe (K-1104) 50-fold in Detector Diluent (K-1105) for a 50 nM working solution. Store remaining Red Fluorescent Probe frozen at –20°C. Multiple freeze-thaw cycles do not affect stability. IMPORTANT: Minimize exposure of this reagent to light.
- **PI3-Kinase Reaction**
 - Conditions for enzyme activity will depend on the characteristics and source of the enzyme used in each specific application. The following protocol has been used at Echelon to detect the activity of a recombinant 6xHis-tagged PI3-K α and is

given as a guideline. Reaction Buffers and working solutions of Substrate, Enzyme, and Inhibitors should be freshly made prior to each experiment. Substrate concentration in the 1X reaction buffer is 10 µM.

- Set up PI3-kinase reactions in 10 µL volume per assay point. For best result, set up reactions directly in the wells of the provided OptiPlate.
- For each 10 µL reaction add the following:
 - 5 µL PI3-kinase enzyme. If inhibitors are to be added to the reaction they should be added with the enzyme and a 5 min pre-incubation of the enzyme with the inhibitor is suggested. Note: We suggest a “no enzyme” control.
 - 5 µL 2X Reaction Buffer.
- Incubate at 25 to 37°C for appropriate time, depending on the activity of your enzyme. The exact amount of enzyme and conditions of incubation will vary with different enzyme preparations and will need to be optimized for each specific application.
- In our hands, each assay point uses approximately 10 ng of 6xHis-tagged recombinant PI3-K α per 100 pmol of PI(4,5)P2 substrate, with 1 hour incubation at 37°C.

Protocol for Fluorescence Polarization Assay

Please read this entire section and the Assay Notes section before beginning the assay.

1. Prepare six 2-fold serial dilutions of PI(3,4,5)P₃ from the 4 µM stock using 1X Reaction Buffer, as outlined in the table below. Each standard will use 10 µL/well of each dilution (0.625, 1.25, 2.5, 5, 10, 20, and 40 pmol PI(3,4,5)P₃). Also set up a standard containing 10 µL 1X Reaction Buffer alone with no PI(3,4,5)P₃ competitor. It is suggested that the standards are run in duplicate or triplicate. Prepare only what is needed for each individual experiment.
2. Add reagents in the order below for 25 µL/well total volume in the provided Optiplate.
 - a. 10 µL of enzyme reactions, controls, or PI(3,4,5)P₃ standards in each well.

Table 1, PIP3 Standards

Concentration of PIP3	PIP3 / 10µL	100 µM stock or previous dilution	1X Reaction Buffer
4.0 µM	40 pmol	8 µL (40 µM PI(3,4,5)P3 Stock Soln.)	72 µL
2.0µM	20 pmol	40 µL (4 µM Solution)	40 µL
1.0 µM	10 pmol	40 µL (2 µM Solution)	40 µL
0.5 µM	5 pmol	40 µL (1 µM Solution)	40 µL
0.25 µM	2.5 pmol	40 µL (0.5 µM Solution)	40 µL
0.125 µM	1.25 pmol	40 µL (0.25 µM Solution)	40 µL
0.0625 µM	0.625 pmol	40 µL (0.125 µM Solution)	40 µL
0 µM	0 pmol	-	40 µL

Note: we suggest running a Probe Alone and Buffer Only control with each experiment. See assay notes for more information.

- b. 10 µL of 250 nM stock of PI(3,4,5)P₃ Detector.

Note: Addition of the PI(3,4,5)P₃ Detector solution quenches the kinase reaction. For time-course experiments, add the Detector solution at desired time points to terminate enzyme activity. After all reactions are quenched, add the Red Fluorescent Probe solution to proceed with detection.

- c. 5 µL of 50 nM Red Fluorescent Probe working solution.
3. Tap plate to mix gently. Seal plate and incubate in a dark location for 30 - 60 minutes. Incubations may be as long as six hours with minimal effect on final measurements. Measure fluorescence polarization using an appropriate instrument and filter set compatible with TAMRA dyes. (550 nm excitation/580 nm polarizing emission filters). Values obtained for enzyme reactions can be compared to the standard curve to determine conversion of substrate to PI(3,4,5)P₃.

Support Protocol: Immunoprecipitation of PI3-Kinase

1. Grow cells to 80% confluence in 10 cm dishes.
2. Induce quiescence by incubating overnight in serum-free medium containing 0.5% insulin-free BSA.
3. Remove medium and stimulate cells with 100 nM insulin for 10 minutes at 37°C.
4. Remove solution and place cells on ice. Add 10 mL per dish of ice-cold Buffer A (137 mM NaCl, 20 mM Tris-HCl, pH 7.4, 1 mM CaCl₂, 1 mM MgCl₂, and 0.1 mM sodium orthovanadate). Rinse three times with this solution.
5. Remove Buffer A and add 1 mL of ice-cold Lysis Buffer (Buffer A plus 1% NP-40 and 1 mM PMSF). Keep plates on ice for 20 minutes.
6. Scrape cells from dish, transfer to 1.5 mL microfuge tubes. Centrifuge for 10 minutes at high speed to sediment insoluble material.
7. Transfer supernatant to new tubes, add 5 µL of anti-PI3 kinase antibody (Millipore, catalog # 06-195) to each tube. Incubate for one hour at 4°C with gentle rotation.
8. Add 60 mL of a 50% slurry of Protein A-agarose beads in PBS to each tube. Incubate with mixing for one hour at 4°C.
9. Collect immunoprecipitated enzyme by centrifuging 5 seconds, and wash with freshly prepared buffers as follows:
 - Three times with Buffer A plus 1% NP-40
 - Three times with 0.1 M Tris-HCl, pH 7.4; 5 mM LiCl, and 0.1 mM sodium orthovanadate.
 - Twice with TNE (10 mM Tris-HCl, pH 7.4, 150 mM NaCl, 5 mM EDTA) containing 0.1 mM sodium orthovanadate.

10. Remove last wash as completely as possible. Wash twice with 1X KBZ and proceed immediately with kinase reactions as described in the Basic Protocol.

References

1. Miller MS, Maheshwari S, McRobb FM, Kinzler KW, Mario Amzel L, Vogelstein B, et al. Identification of allosteric binding sites for PI3K1± oncogenic mutant specific inhibitor design. *Bioorganic & Medicinal Chemistry*. 2017.
2. Heffron TP, Heald RA, Ndubaku CO, Wei B, Augustin M, Do S, et al. The Rational Design of Selective Benzoxazepin Inhibitors of the α-Isoform of Phosphoinositide 3-Kinase Culminating in the Identification of (S)-2-((2-(1-isopropyl-1H-1,2,4-triazol-5-yl)-5,6-dihydrobenzo[f]imidazo [1,2-d][1,4]oxazepin-9-yl)oxy)propanamide (GDC-0326). *Journal of medicinal chemistry*. 2016.
3. Safina BS, Baker S, Baumgardner M, Blaney PM, Chan BK, Chen Y-H, et al. Discovery of Novel PI3-Kinase d Specific Inhibitors for the Treatment of Rheumatoid Arthritis: Taming CYP3A4 Time-Dependent Inhibition. *Journal of medicinal chemistry*. 2012;55(12):5887-900.
4. Dehnhardt CM, Mansour Tarek S. Lead Optimization of N-3-Substituted 7-Morpholinotriazolopyrimidines as Dual Phosphoinositide 3-Kinase/Mammalian Target of Rapamycin Inhibitors: Discovery of PKI-402. *J Med Chem*. 2009.
5. Ma H, Deacon S, Horiuchi K. The Challenge of Selecting Protein Kinase Assays for Lead Discovery Optimization. *Expert Opinion on Drug Discovery*. 2008;3(6):607-21.
6. Carson JD, Van Aller G, Lehr R, Sinnamon RH, Kirkpatrick RB, Auger KR, et al. Effects of oncogenic p110α subunit mutations on the lipid kinase activity of phosphoinositide 3-kinase. *Biochem J*. 2008;409(2):519-24.
7. Nagoshi T, Matsui T, Aoyama T, Leri A, Anversa P, Li L, et al. PI3K rescues the detrimental effects of chronic Akt activation in the heart during ischemia/reperfusion injury. *J Clin Invest*. 2005;115(8):2128-38.
8. Drees BE, Weipert A, Hudson H, Ferguson CG, Chakravarty L, Prestwich GD. Competitive fluorescence polarization assays for the detection of phosphoinositide kinase and phosphatase activity. *Comb Chem High Throughput Screen*. 2003;6(4):321-30.

Related Product

Products	Catalog Number
Assays	
PI3-Kinase Activity ELISA: Pico	K-1000s
Class III PI3K Elisa Kit	K-3000
PIP3 Binding Protein	
Anti-PtdIns(3,4,5)P ₃ IgM	Z-P345
PI(3,4,5)P ₃ Grip (GRP1-PH)	G-3901
Lipids and inhibitors	
PIP3	P-3908, P-3916
Wortmannin	B-0222
LY294002	B-0294

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