

## 1.0 INTRODUCTION

PDK1 (3-Phosphoinositide-Dependant Protein Kinase-1) activates the conventional PKCs and PKC $\zeta$  (zeta) through phosphorylation at Threonine 500 in the 'activation loop'. PDK1 also phosphorylates Protein Kinase B (PKB/Akt) at Threonine 308 in the presence of phosphatidylinositol-3,4,5-trisphosphate. Active PKB/Akt inactivates Glycogen Synthase Kinase-3 (GSK3), eventually leading to the dephosphorylation and activation of glycogen synthase and the stimulation of glycogen synthesis (1). Because of the role PDK1 plays in insulin-induced glycogen synthesis and PKC activation, it is potentially an important target for metabolic drug research. Conventional threonine kinase assays are tedious, utilize radioactive reagents, and are not easily automated or converted to a high-throughput format for drug screening. PanVera's PDK1 assay kit is a major advance because it is simple, sensitive, non-radioactive, homogeneous and formatted for high-throughput screening. This assay kit contains all the materials necessary to analyze PDK1-like kinase activity via PDK1 phosphorylation using fluorescence polarization (FP) as the detection method. The PDK1 kits contain reagent volumes sufficient for one hundred (PanVera Part No. P2884) and one thousand (PanVera Part No. P2885) 100  $\mu$ L assays, respectively. These kits are formatted for use with either a multi-well plate instrument capable of measuring FP in 96-well plates or a single-well FP instrument, such as the Beacon® 2000 System (PanVera Corporation). These kits will also work in 384- and 1536-well plates by using reduced reagent volumes.

## 2.0 THEORY

### 2.1 What is Fluorescence Polarization?

Fluorescence polarization is a technique for monitoring molecular interaction in a homogenous environment at equilibrium. The binding of a small, fluorescent molecule to another, larger molecule can be quantified by a change in the rate of rotation of the fluorescent molecule. Plane-polarized light is used to excite fluorescent molecules in solution; if the excited molecule remains stationary during the period of excitation (4 nanoseconds for fluorescein) the emitted light will remain highly polarized. However, if the excited molecule can tumble or rotate during this period, the emitted light will be depolarized (2, 3).

Since FP is a measure of this tumbling rate, it is directly related to the molecular volume of the fluorescent molecule. An increase in molecular volume, due to biological processes such as receptor-ligand binding, antibody-antigen binding, DNA hybridization or DNA-protein binding, or a decrease in molecular volume due to enzymatic degradation or dissociation, can be measured directly by FP.

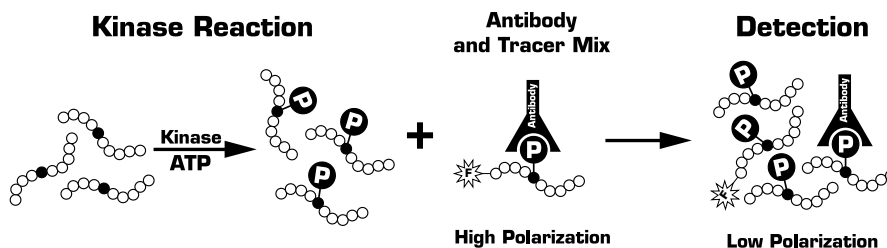
The measured FP value is the weighted average of the FP values of the individual bound and free fluorescent molecules and is therefore a direct measure of the fraction bound. These data can be handled exactly the same way as traditional radioligand-binding assay data. Fluorescence polarization values are plotted against the logarithm of the receptor concentration, resulting in the familiar saturation-binding isotherm.

For more information, see our on-line Fluorescence Polarization Applications Guide at:

<http://www.panvera.com/tech/appguide/index.html>

### 2.2 Assay Theory

Fluorescence polarization can be exploited in a PDK1 assay because fluorescein-labeled phosphopeptides (F-phosphopeptides, also referred to as the "tracer") will compete with phosphopeptides or phosphoproteins generated during a kinase reaction for binding to anti-phosphothreonine peptide-specific antibodies. When there are no kinase reaction products present, a significant portion of the F-phosphopeptide tracer will be bound by the anti-phosphothreonine peptide-specific antibody, resulting in a high polarization value. However, after a PDK1 reaction has occurred, phosphorylation can be detected because the anti-phosphothreonine peptide-specific antibodies will bind to both the F-phosphopeptide tracer and the non-fluorescent reaction products, decreasing the amount of bound tracer and thus decreasing the fluorescence polarization value of the sample. This is illustrated below. If enough competitor phosphopeptide is generated during the reaction, the fluorescent tracer can be completely displaced from the anti-phosphothreonine peptide-specific antibodies and the emitted light will be totally depolarized. Thus, the change in FP is directly related to PDK1 kinase activity.





# Threonine Kinase Assay Kit, PDK1

## Protocol

Part # P2884, P2885

Lit. # L0520 Rev. 07/01

Page 2 of 6

### 3.0 SAFETY PRECAUTIONS

Normal precautions exercised in handling laboratory reagents should be followed. All reagents in this kit are considered non-hazardous according to 29 CFR 1910.1200. The chemical, physical, and toxicological properties of these products may not have been thoroughly investigated. We recommend the use of gloves, lab coats, eye protection and a fume hood when working with any chemical reagents.

### 4.0 KIT COMPONENTS

#### 4.1 Materials Provided

Description	Composition	P2884		P2885	
		Size	Part #	Size	Part #
Anti-PDK1 Antibody, 4X	Anti-phosphothreonine peptide-specific antibody in 20 mM Tris (pH 7.5), 0.02% NaN <sub>3</sub>	2.5 mL	P2914	25 mL	P2915
PDK1 Tracer, 10X	Fluorescein-labeled phosphopeptide in 20 mM Tris (pH 7.5), 0.02% NaN <sub>3</sub>	1 mL	P2916	10 mL	P2917
PDK1 Competitor	5 $\mu$ M phosphopeptide in 20 mM Tris (pH 7.5), 0.02% NaN <sub>3</sub>	500 $\mu$ L	P2918	500 $\mu$ L	P2918
Kinase Quench Buffer	500 mM EDTA (pH 8.0)	1 mL	P2825	10 mL	P2832
FP Dilution Buffer	20 mM Tris (pH 7.5), 0.02% NaN <sub>3</sub>	5 mL	P2838	50 mL	P2839

#### 4.2 Materials Available Separately

- PDK1 Substrate Peptide, Non-phosphorylated, 100  $\mu$ M (PanVera Part No. P2925, 1 mL)

#### 4.3 Materials Required but not Supplied

- Kinase
- ATP
- Kinase Reaction Buffer
- Multi-well plates suitable for fluorescence polarization or 6mm borosilicate tubes
- Multi-well fluorescence polarization instrument or a Beacon® 2000 Fluorescence Polarization instrument with suitable 485 nm excitation and 530 nm emission interference filters
- Pipeting devices
- Reagent reservoir(s)
- Laboratory timer
- Beacon® FP One-Step Standardization Kit (PanVera Part No. P2581). This kit is recommended to validate instrument performance and may be required as a polarization standard to calibrate certain instruments.

### 5.0 STORAGE AND STABILITY

The 4X PDK1 Anti-pThr Antibody should be stored at  $-20^{\circ}\text{C}$ . Once thawed, it should be stored at  $+4^{\circ}\text{C}$ . The 10X PDK1 Tracer and the PDK1 Competitor should be stored at  $-20^{\circ}\text{C}$ , thawed when needed, and then refrozen. Avoid subjecting these reagents to repeated freeze/thaw cycles. The FP Dilution Buffer and Kinase Quench Buffer can be thawed upon receipt and stored at  $20-30^{\circ}\text{C}$ . All reagents are stable for 6 months from the date of receipt.

## 6.0 DETERMINING THE IC<sub>50</sub> FOR THE PDK1 COMPETITOR

The following instructions for testing and validating the kit are optional, but highly recommended for a first-time user. Generating a standard curve demonstrates the type of data that can be expected during the enzymatic reaction. It is also recommended that you generate this type of standard curve using the threonine-phosphorylated form of the peptide substrate you intend to use for your PDK1 reaction. Once the IC<sub>50</sub> for the threonine-phosphorylated form of the substrate is known, one can use these data to determine (based on the turnover rate of your enzyme) how much enzyme you will need, how much substrate your PDK1 reaction will require, and how long the reaction should be run.

### 6.1 96-well Plate Format

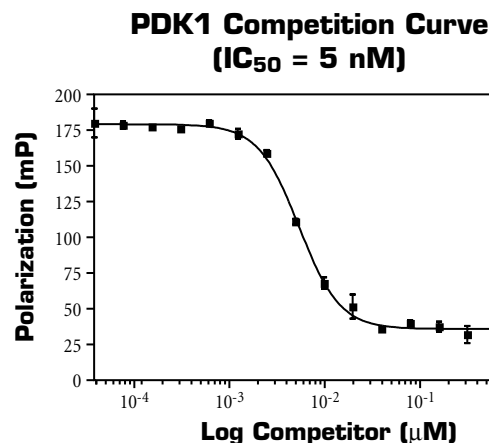
1. In a 96-well plate suitable for fluorescence polarization, pipette 100 µL of 5 µM PDK1 Competitor (**VORTEXED**) into well A1.
2. Dispense 50 µL FP Dilution Buffer into wells A2 to A12 and wells B1 to B6 (17 wells total).
3. Perform a 2-fold serial dilution of the PDK1 Competitor by transferring and mixing 50 µL from well A1 into well A2.
4. Repeat this process (from well A2 to A3, then from well A3 to A4 etc.) until well B6 is reached (18 wells total).
5. Discard 50 µL from well B6. There should be 50 µL in every well from A1 to B6.
6. **VORTEX** 10X PDK1 Tracer.
7. Add 25 µL of the 4X PDK1 Antibody and 10 µL of the 10X PDK1 Tracer to every well from A1 to B6 (18 wells total). Add 15 µL of FP Dilution Buffer to each well. (**Do not pre-mix the tracer and antibody** or displacement by the phosphopeptide competitor will not occur.) This step will reduce the concentration of all of the detection components to 1X and reduce the concentration of the PDK1 Competitor in each well by 2-fold (*i.e.*, well A1 now has a competitor concentration of 2.5 µM, while B6 has a competitor concentration of approximately 0.02 nM).
8. Mix, cover the multi-well plate to reduce evaporation and protect the reagents from light, and incubate for 1 hour at room temperature.
9. Measure the polarization values of your samples, following the procedures required by your FP instrument. We recommend that you use 100 µL FP Dilution Buffer as your blank and the Beacon® FP One-Step Standardization Kit (Panvera Part No. P2581) for your low and high polarization standards, if necessary. We recommend that you may also run controls for the antibody/tracer mixture (without competitor) and for the tracer alone to determine the low and high polarization values for the assay itself. See Section 6.3.
10. Perform non-linear regression on a semi-log plot of the data (mP vs. log [Competitor]). The IC<sub>50</sub> of the competitor control should be less than 20 nM. Sample data (n = 4) generated on a TECAN Polarion (Research Triangle Park, NC) appear in Section 6.3.

### 6.2 Single-tube Format

1. Label eighteen borosilicate tubes #1 through #18. Pipette 100 µL of 5 µM PDK1 Competitor (**VORTEX**) into tube #1.
2. Dispense 50 µL FP Dilution Buffer into tube #2 through #18 (17 tubes total).
3. Perform a 2-fold serial dilution of the PDK1 Competitor by transferring and mixing 50 µL from tube #1 into tube #2.
4. Repeat this process (from tube #2 to #3, then from tube #3 to #4 etc.) until tube #18 is reached.
5. Discard 50 µL from tube #18. There should now be 50 µL in all 18 tubes.
6. **VORTEX** 10X PDK1 Tracer.
7. Add 25 µL of the 4X PDK1 Antibody and 10 µL of the 10X PDK1 Tracer to every tube. Add 15 µL of FP Dilution Buffer to each tube. (**Do not pre-mix the tracer and antibody** or displacement by the phosphopeptide will not occur.) This step will reduce the concentration of all of the detection components to 1X and reduce the concentration of the PDK1 Competitor in each tube by 2-fold (*i.e.*, tube #1 now has a competitor concentration of 2.5 µM, while tube #18 will have a competitor concentration of approximately 0.02 nM).
8. Mix, cover the tubes to reduce evaporation and protect the reagents from light, and incubate for 1 hour at room temperature.
9. Measure the polarization values of your samples, following the procedures required by your FP instrument. It is recommended that you use 100 µL FP Dilution Buffer as your blank and the Beacon® FP One-Step Standardization Kit (Panvera Part No. P2581) for your low and high polarization standards, if necessary. We recommend that you also run controls for the antibody/tracer mixture (without competitor) and for the tracer alone to determine the low and high polarization values for the assay itself. See Section 6.3.
10. Perform non-linear regression on a semi-log plot of the data (mP vs. log [Competitor]). The IC<sub>50</sub> of the competitor control should be less than 20 nM. Sample data (n = 4) generated on a TECAN Polarion (Research Triangle Park, NC) appear in Section 6.3.

### 6.3 Standard Curve Controls

To demonstrate the minimum polarization value of this detection system under the standard curve conditions, add 10  $\mu$ L of the 10X PDK1 Tracer to 90  $\mu$ L of the FP Dilution Buffer in a single well of a multi-well plate or borosilicate tube. The minimum polarization value represents conditions in which the tracer has been completely displaced from the antibody (*i.e.*, total phosphorylation of the substrate by a kinase). To demonstrate the maximum polarization value of this detection system under the standard curve conditions, add 10  $\mu$ L of the 10X PDK1 Tracer and 25  $\mu$ L of the 4X Anti-PDK1 Antibody in 65  $\mu$ L of the FP Dilution Buffer. The high polarization value represents conditions in which the tracer is completely bound by the antibody (*i.e.*, no phosphorylated substrate has been generated by a kinase that can compete for binding sites and displace the tracer from the antibody).



## 7.0 PDK1 THREONINE KINASE ASSAY

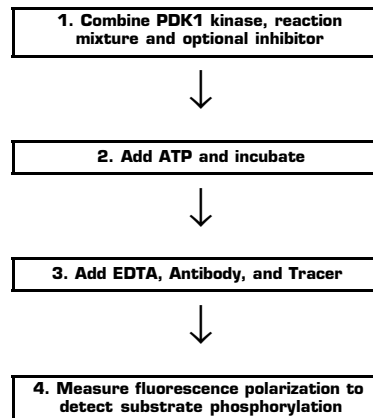
### 7.1 PDK1 Threonine Kinase Assay Outline

Each complete PDK1 Threonine Kinase Assay Outline reaction must contain an appropriate buffer system, enzyme, ATP, substrate, PDK1 Tracer, and Anti-PDK1 Antibody. EDTA is then used to quench kinase activity.

There are two modes of analysis for this assay: kinetic and end-point. The kinetic mode is useful for rapidly optimizing enzyme concentration, reaction time, substrate concentration, and buffer conditions. The end-point assay mode is useful when screening kinase inhibitors or activators.

In the end-point mode (Section 7.4), four general steps are required. First, the reaction mixture, including the substrate (PanVera offers PDK1 Substrate Peptide, Part No. P2925) and optional inhibitor, is combined. ATP is then added to start the reaction. Following incubation, stop the kinase reaction with EDTA (Kinase Quench Buffer), then add anti-phosphothreonine antibodies and tracer **separately** to initiate the competition for antibody binding. Finally, polarization values are measured to determine the extent of PDK1 phosphorylation. If a kinetic experiment is performed (Section 7.3), the components of Steps 1 and 3 (without EDTA) are combined first and then Steps 2 and 4 are executed simultaneously.

**Note:** The following protocol is configured for use with 96-well plates and a multi-well FP instrument. However, the protocol can also be performed in 384- or 1536-well plates, or borosilicate tubes if you are using a single-tube FP instrument such as the Beacon® 2000 Instrument.



### 7.2 Experimental Controls

To demonstrate the minimum polarization value of this detection system under the conditions of your reaction, **DO NOT** add the Anti-PDK1 Antibody to one reaction as described in Step 2 of Section 7.3 or Step 1 of Section 7.4.2. Substitute this volume with FP Dilution Buffer. This is referred to as the “No Antibody” control throughout the remainder of this protocol. The “No Antibody” control will indicate the minimum polarization value of the reaction system, which simulates a kinase reaction that has completely displaced the fluorescent phosphopeptide tracer from the anti-phosphothreonine peptide-specific antibody. If a time-course experiment is performed, the polarization value should decrease with time.

To demonstrate the maximum polarization value of this detection system under the conditions of your reaction, add in at least one tube or well, all of your kinase reaction and detection components, **except the ATP**, which is added in Step 3 of Section 7.3 or Step 2 of Section 7.4.1. Substitute this volume with FP Dilution Buffer. Perform the rest of the experiment as described. This sample is referred to as the “No ATP” control throughout the remainder of this protocol. The “No ATP” control should have a high polarization value since no competitor phosphopeptide was produced during the PDK1 reaction to displace the phosphopeptide tracer from the antibody.

If kinase inhibitors are to be tested or screened, we recommend that you initially serially dilute the inhibitors in 100% DMSO or another solvent. Perform a 10-fold dilution of the inhibitors once from 100X (in 100% DMSO) to 10X (in 10% DMSO) using kinase reaction buffer

as the diluent, followed by another 10-fold dilution into the PDK1 reaction(s), producing a final concentration of 1X (in 1% DMSO). Include a “zero” inhibitor sample (“vehicle” or buffer control) containing the appropriate amount of DMSO or another solvent to measure its effect on the PDK1 reaction. If kinase inhibitors were titrated into the reaction, the polarization value should remain high as the kinase inhibitor concentration is increased, while low or zero concentrations of PDK1 inhibitor will have low polarization values, indicating that the PDK1 reaction was able to proceed largely (or totally) uninhibited.

### 7.3 Kinetic Assay

In the kinetic assay, all of the reaction and detection reagents (FP Dilution Buffer, ATP, substrate, Anti-PDK1 Antibody, and the PDK1 phosphopeptide tracer) are present during the reaction phase of the assay. Phosphorylation can be monitored in real-time as the reaction proceeds; the polarization value of the sample will be reduced as phosphothreonine-containing peptides are generated and compete with the phosphopeptide tracer for binding to the antibody.

1. In a volume of less than 50  $\mu$ L, set-up a PDK1 reaction without ATP. The final reaction volume (including ATP) will be 50  $\mu$ L prior to the addition of the detection reagents. A typical reaction will require a protein threonine kinase, an appropriate buffer system, peptide or protein substrate, and ATP (which is added last to initiate the reaction).
2. Once you have dispensed the reactants into a multi-well plate or borosilicate tube, add 25  $\mu$ L of the 4X Anti-PDK1 Antibody and 10  $\mu$ L of the 10X PDK1 Tracer (**VORTEXED**), to each sample. **DO NOT** pre-mix the tracer and antibody or displacement by the phosphopeptides generated during your kinase reaction *will not* occur. Add 15  $\mu$ L of FP Dilution Buffer to each well. Preincubate each sample at your desired reaction temperature.
3. Add an ATP solution of the appropriate concentration to each sample to start the reaction. The final volume should be brought to 100  $\mu$ L with water. A “No ATP” control (substitute water or buffer for ATP) should be included in every experiment.
4. Quickly transfer your multi-well plate or borosilicate tube to your FP instrument (operating at your preferred reaction temperature) and begin reading the fluorescence polarization values of your sample(s). Continue to measure the polarization values of your samples as long as necessary.

### 7.4 End-point Assay

The end-point PDK1 assay is divided into two phases: the reaction phase and the detection phase. The reaction phase is performed similar to a traditional radioactive kinase assay, except that no radioactive label is required. The concentrations of substrate, ATP and enzyme required should be determined experimentally.

For the end-point assay, a kinase reaction mix is prepared in a volume of less than 50  $\mu$ L, so that the final reaction volume (including ATP and Kinase Quench Buffer) will equal 50  $\mu$ L. The reaction is started by the addition of ATP, incubated for the desired length of time, then quenched with EDTA. The final volume of the mixture should be 50  $\mu$ L. Then, 25  $\mu$ L of the 4X Anti-PDK1 Antibody, 10  $\mu$ L of the 10X PDK1 Tracer (**VORTEXED**), and 15  $\mu$ L of FP Dilution Buffer are added to each well. This addition reduces the concentration of each component of the kit to its final 1X concentration. The polarization values are then read and the data analyzed.

It should be noted that the reaction can be run in any volume, keeping in mind that the quenched kinase reaction will be diluted with the addition of antibody and tracer. The final volume must be less than the maximum volume of the multi-well plate you are using.

#### 7.4.1 Reaction Phase

1. In a volume of less than 50  $\mu$ L, set-up a PDK1 reaction without ATP. A typical reaction will require a protein kinase, an appropriate buffer system and peptide substrate.
2. Once you have dispensed the reaction(s) into a multi-well plate or borosilicate tube, add ATP (at a concentration appropriate for your reaction) to each well to start the reaction. The total volume should now be 50  $\mu$ L. Please note that a “No ATP” control (substitute water or buffer for ATP) should be included in every experiment.
3. Incubate at your preferred reaction temperature. Incubation time will vary depending on the amount of kinase used, as well as the turnover rate of your kinase. We recommend running an initial time-point experiment to determine the optimal incubation length.
4. Quench the kinase reaction(s) at the end of the incubation. We recommend that you add EDTA to a final concentration of 10 mM and bring the final volume of each quenched PDK1 reaction to 50  $\mu$ L. After the addition of EDTA, mix and incubate for 5 minutes at room temperature. You may use more or less EDTA (or another quenching reagent) for your specific kinase. EDTA at a concentration greater than 20 mM is known to interfere with the anti-phosphothreonine antibody/phosphopeptide binding.
5. Proceed to the Detection Phase of the experiment in Section 7.4.2.



# Threonine Kinase Assay Kit, PDK1

Protocol

Part # P2884, P2885

Lit. # L0520 Rev. 07/01

Page 6 of 6

## 7.4.2 Detection Phase

1. After quenching the PDK1 reaction, add 25  $\mu$ L of the 4X Anti-PDK1 Antibody, 10  $\mu$ L of the 10X PDK1 Tracer (**VORTEXED**), and 15  $\mu$ L of FP Dilution Buffer to each well. **DO NOT** pre-mix the tracer and antibody or displacement by the phosphopeptides generated during your kinase reaction *will not* occur.
2. Cover the wells or tubes to reduce evaporation and protect them from light, then incubate for 1 hour at room temperature.
3. Measure the fluorescence polarization value of each reaction following the procedures required by your FP instrument. It is recommended that you use 100  $\mu$ L FP Dilution Buffer as your blank and the Beacon® FP One-Step Standardization Kit (PanVera Part No. P2581) for your low and high polarization standards, if necessary.

## 8.0 REFERENCES

1. Vanhaesebroeck, B. and Alessi, D.R. (2000) *Biochem. J.* **346**:561-76.
2. Checovich, W.C., Bolger, R.E. and Burke, T. (1995) *Nature* **375**:254-6.
3. Jameson, D.M. and Sawyer, W.H. (1995) *Meth. Enzymol.* **246**:283-300.

\*United States and International Patents Pending

The performance of this product is guaranteed for six months from the date of purchase if stored and handled properly.

"The Protein Company" is a trademark of PanVera Corporation.

Beacon and CoreHTS are registered trademarks of PanVera Corporation.

© 2001, PanVera Corporation. All rights reserved.

PURCHASER SHALL USE THE PRODUCTS PURCHASED HEREUNDER (the "Products") SOLELY FOR THE PURPOSE OF CONDUCTING INTERNAL RESEARCH AT ITS ORGANIZATION ("RESEARCH"). Purchaser will not sell, transfer, disclose or otherwise provide access to the Products to any third party. Purchaser agrees that Purchaser shall not have the right to authorize any third party to use or sell any Products or derivatives thereof. Unless a license has been executed between Purchaser and PanVera explicitly providing otherwise, Purchaser will not: (i) reformulate or create derivatives of the Products; or (ii) use the Products for providing services to a third party (e.g. screening or profiling); use the Products in a diagnostic process; or (iii) or use the Products in a quality control or quality assurance process for the manufacture of a product for sale.