

# PKC $\delta$ Kinase Assay

By Hicham Zegzouti, Ph.D., Jolanta Vidugiriene, Ph.D., and Said A. Goueli, Ph.D., Promega Corporation

## Scientific Background:

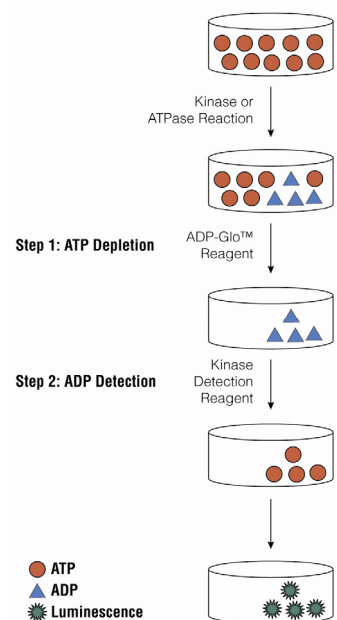
Protein kinase C delta (PKC $\delta$ ) is a member of the protein kinase C (PKC) family of serine-threonine kinases. It is a 79 kD protein kinase that shows strict dependence on the presence of phospholipids, but shows no activation by Ca<sup>2+</sup> (1). Phosphatidylinositol is the most potent activator of PKC delta. Northern blot analysis indicated that PKCdelta is widely distributed in almost all the tissues and is a major isoform of PKC expressed in hemopoietic cells (2). PKCdelta is involved in fundamental cellular functions regulated by diacylglycerols and mimicked by phorbol esters.

1. Leibersperger, H. et al: Immunological demonstration of a calcium-unresponsive protein kinase C of the delta-type in different species and murine tissues. Predominance in epidermis. *J Biol Chem.* 1991 Aug 5;266(22):14778-84.
2. Mischak, H. et al: Mouse protein kinase C-delta, the major isoform expressed in mouse hemopoietic cells: sequence of the cDNA, expression patterns, and characterization of the protein. *Biochemistry.* 1991 Aug 13;30(32):7925-31.

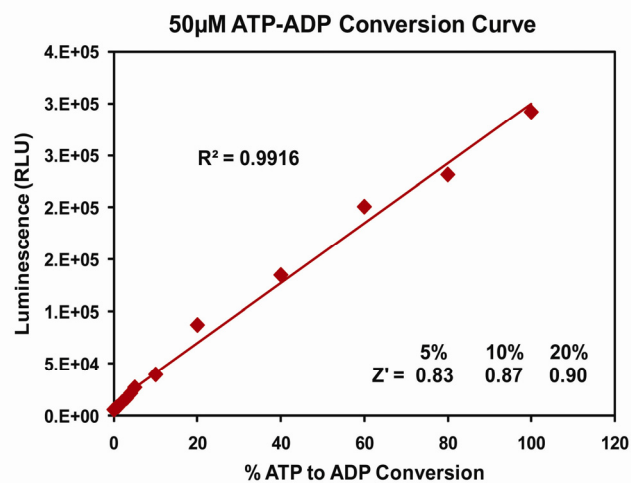
## ADP-Glo™ Kinase Assay

### Description

ADP-Glo™ Kinase Assay is a luminescent kinase assay that measures ADP formed from a kinase reaction; ADP is converted into ATP, which is converted into light by Ultra-Glo™ Luciferase (Fig. 1). The luminescent signal positively correlates with ADP amount (Fig. 2) and kinase activity (Fig. 3A). The assay is well suited for measuring the effects chemical compounds have on the activity of a broad range of purified kinases—making it ideal for both primary screening as well as kinase selectivity profiling (Fig. 3B). The ADP-Glo™ Kinase Assay can be used to monitor the activity of virtually any ADP-generating enzyme (e.g., kinase or ATPase) using up to 1mM ATP.



**Figure 1. Principle of the ADP-Glo™ Kinase Assay.** The ATP remaining after completion of the kinase reaction is depleted prior to an ADP to ATP conversion step and quantitation of the newly synthesized ATP using luciferase/luciferin reaction.



**Figure 2. Linearity of the ADP-Glo Kinase Assay.** ATP-to-ADP conversion curve was prepared at 50µM ATP+ADP concentration range. This standard curve is used to calculate the amount of ADP formed in the kinase reaction. Z' factors were determined using 200 replicates of each of the % conversions shown.



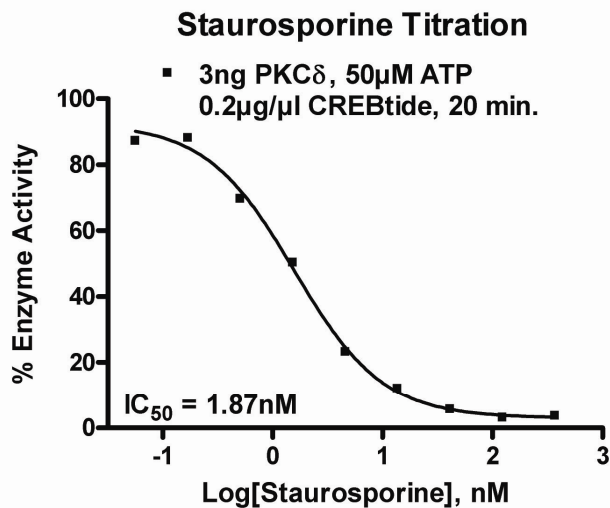
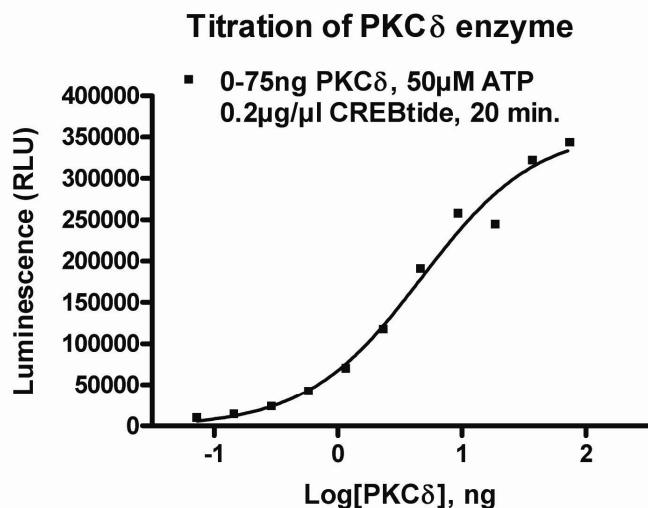
For detailed protocols on conversion curves, kinase assays and inhibitor screening, see *The ADP-Glo™ Kinase Assay Technical Manual #TM313*, available at [www.promega.com/tbs/tm313/tm313.html](http://www.promega.com/tbs/tm313/tm313.html)

## Protocol

- Dilute enzyme, substrate, ATP and inhibitors in Kinase Buffer.
- Add to the wells of 384 low volume plate:
  - 1  $\mu$ l of inhibitor or (5% DMSO)
  - 2  $\mu$ l of enzyme (defined from table 1)
  - 2  $\mu$ l of substrate/ATP mix
- Incubate at room temperature for 20 minutes.
- Add 5  $\mu$ l of ADP-Glo™ Reagent
- Incubate at room temperature for 40 minutes.
- Add 10  $\mu$ l of Kinase Detection Reagent
- Incubate at room temperature for 30 minutes.
- Record luminescence (Integration time 0.5-1second).

**Table 1. PKC $\delta$  Enzyme Titration.** Data are shown as relative light units (RLU) that directly correlate to the amount of ADP produced. The correlation between the % of ATP converted to ADP and corresponding signal to background ratio is indicated for each kinase amount.

PKC $\delta$ , ng	18.75	9.38	4.69	2.34	1.17	0.59	0.29	0
RLU	244149	257044	190643	117157	69432	41688	23097	4365
S/B	56	59	44	27	16	10	5	1
% Conversion	90	95	69	40	22	11	3	0



**Figure 3. PKC $\delta$  Kinase Assay Development:** (A) PKC $\delta$  enzyme was titrated using 50 $\mu$ M ATP and the luminescence signal generated from each of the amounts of the enzyme is shown. (B) Staurosporine dose response was created using 3ng of PKC $\delta$  to determine the potency of the inhibitor (IC<sub>50</sub>).

### Assay Components and Ordering Information:



#### Products

ADP-Glo™ Kinase Assay  
PKC $\delta$  Kinase Enzyme System  
ADP-Glo + PKC $\delta$  Kinase Enzyme System

#### Company

Promega  
Promega  
Promega

#### Cat.#

V9101  
V3401  
V9721

PKC $\delta$  Kinase Buffer: 40mM Tris, 7.5; 20mM MgCl<sub>2</sub>; 0.1mg/ml BSA; 50 $\mu$ M DTT; 1 x PKC Lipid activator mix.