

# PKC $\epsilon$ Kinase Assay

By Juliano Alves, Ph.D., Said A. Goueli, Ph.D., and Hicham Zegzouti, Ph.D., Promega Corporation

## Scientific Background:

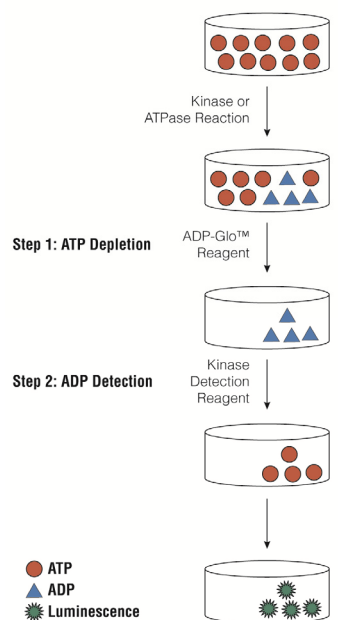
PKC $\epsilon$  is a member of the protein kinase C (PKC) family of serine- and threonine-specific protein kinases that can phosphorylate a wide variety of protein targets known to be involved in diverse cellular signaling pathways. PKC $\epsilon$  is involved in many different cellular functions, such as neuron channel activation, cardioprotection from ischemia (1), heat shock response, as well as insulin exocytosis. Knockout studies in mice suggest that this kinase is important for lipopolysaccharide (LPS)-mediated signaling in activated macrophages and may control anxiety-like behavior (2).

1. Chen, C H. et al: Cardioprotection from ischemia by a brief exposure to physiological levels of ethanol: role of epsilon protein kinase C. *Proc. Nat. Acad. Sci.* 96: 12784-12789, 1999.
2. Hodge, C W. et al: Decreased anxiety-like behavior, reduced stress hormones, and neurosteroid supersensitivity in mice lacking protein kinase C-epsilon. *J. Clin. Invest.* 110: 1003-1010, 2002.

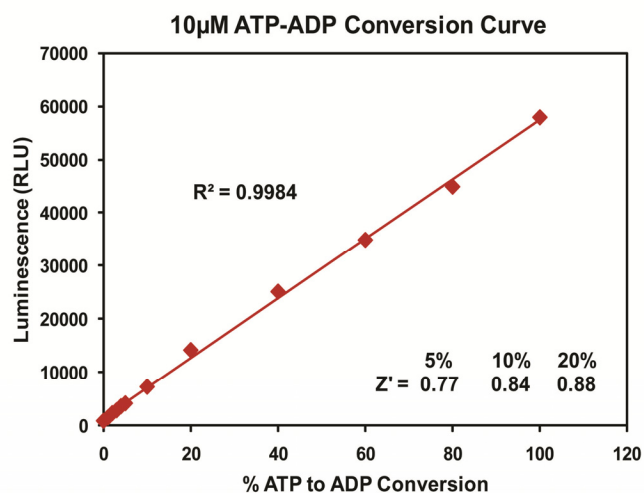
## ADP-Glo™ Kinase Assay

### Description

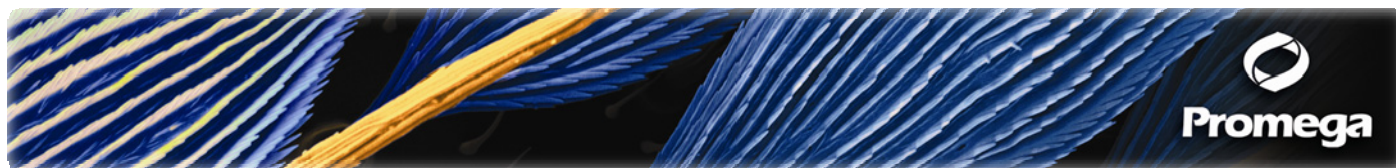
ADP-Glo™ Kinase Assay is a luminescent kinase assay that measures ADP formed from a kinase reaction; ADP is converted into ATP, which is converted into light by Ultra-Glo™ Luciferase (Fig. 1). The luminescent signal positively correlates with ADP amount (Fig. 2) and kinase activity (Fig. 3A). The assay 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 10µM 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.



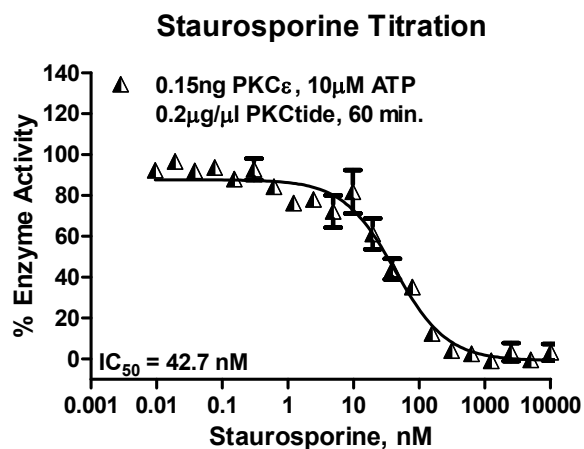
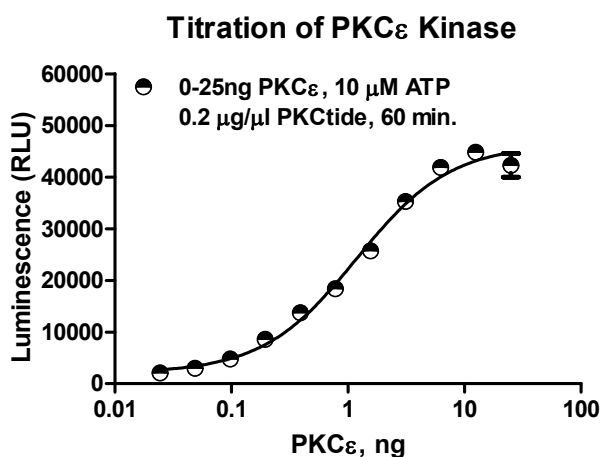
For detailed protocols on conversion curves, kinase assays and inhibitor screening, see *The ADP-Glo™ Kinase Assay Technical Manual #TM313*, available at [www.promega.com/tbs/tm313/tm313.html](http://www.promega.com/tbs/tm313/tm313.html)

## 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 60 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 $\epsilon$  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 $\epsilon$ , ng	25	12.5	6.3	3.1	1.6	0.8	0.4	0.2	0.1	0.05	0.02	0
RLU	44572	44500	41871	35260	25712	18363	13717	8591	4781	2975	2062	293
S/B	152	152	143	121	88	63	47	29	16	10	7	1
% Conversion	98	98	92	77	55	39	28	17	8	4	2	0



**Figure 3. PKC $\epsilon$  Kinase Assay Development.** (A) PKC $\epsilon$  enzyme was titrated using 10 $\mu$ M ATP and the luminescence signal generated from each of the amounts of the enzyme is shown. (B) Staurosporine dose response was created using 0.15ng of PKC $\epsilon$  to determine the potency of the inhibitor (IC<sub>50</sub>).

Assay Components and Ordering Information:		Promega	SignalChem <small>Specialists in Signaling Proteins</small>
Products	Company	Cat.#	
ADP-Glo™ Kinase Assay	Promega	V9101	
PKC $\epsilon$ Kinase Enzyme System	Promega	V4036	
ADP-Glo™ + PKC $\epsilon$ Kinase Enzyme System	Promega	V4037	

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