

PKCβII Kinase Assay

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

Scientific Background:

PKCβ II is a member of the protein kinase C (PKC) family of serine- and threonine-specific protein kinases that can be activated by calcium and second messenger diacylglycerol and phosphorylate a wide variety of protein targets known to be involved in diverse cellular signaling pathways (1). PKCβ II has been reported to be involved in many different cellular functions such as B cell activation, apoptosis induction, endothelial cell proliferation, and intestinal sugar absorption (2).

1. Greenham, J. et al: Elucidation of the exon-intron structure and size of the human protein kinase C beta gene (PRKCB). *Hum. Genet.* 103: 483-487, 1998.
2. Su, T T. et al: PKC-beta controls I-kappa-B kinase lipid raft recruitment and activation in response to BCR signaling. *Nature Immun.* 3: 780-786, 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.

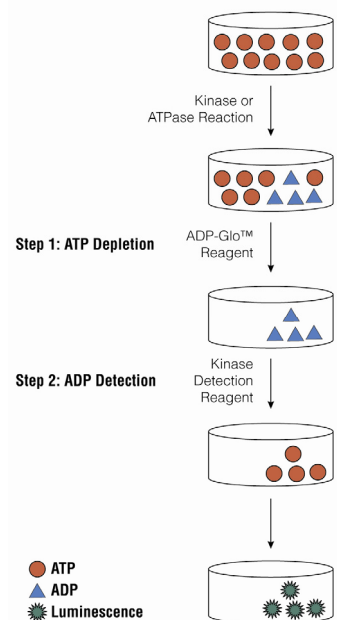


Figure 1. Principle of the ADP-Glo™ Kinase Assay. The ATP remaining after completion of the kinase reaction is depleted prior to an ADP to ATP conversion step and quantitation of the newly synthesized ATP using luciferase/luciferin reaction.

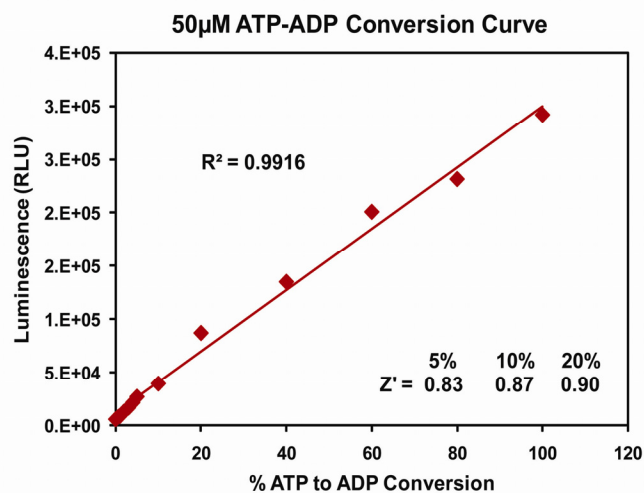


Figure 2. Linearity of the ADP-Glo Kinase Assay. ATP-to-ADP conversion curve was prepared at 50μM ATP+ADP concentration range. This standard curve is used to calculate the amount of ADP formed in the kinase reaction. Z' factors were determined using 200 replicates of each of the % conversions shown.



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

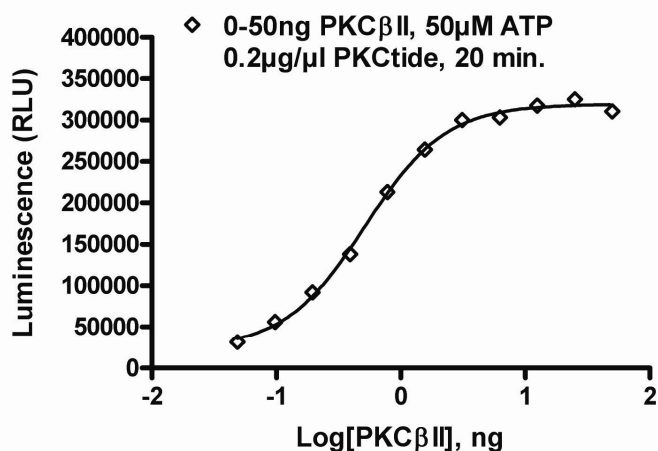
Protocol

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

Table 1. PKC β II 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 β II, ng	12.5	6.3	3.1	1.6	0.8	0.39	0.20	0.10	0.05	0
RLU	303403	284057	300303	264642	213156	138330	92016	56039	31368	5495
S/B	55.2	51.7	54.7	48.2	38.8	25.2	16.7	10.2	5.7	1
% Conversion	100	94.6	100	87.9	70.0	43.9	27.8	15.3	6.7	0

Titration of PKC β II enzyme



Staurosporine Titration

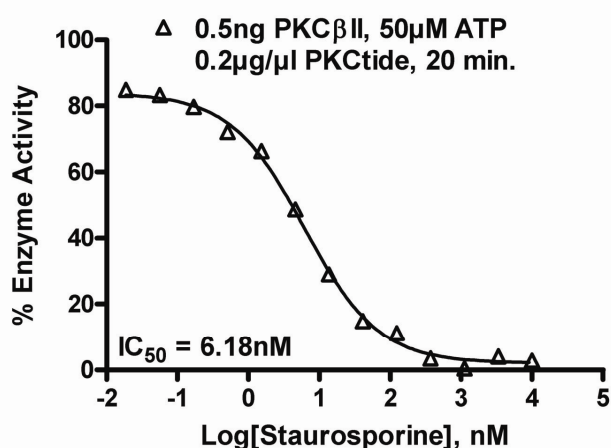


Figure 3. PKC β II Kinase Assay Development: (A) PKC β II enzyme was titrated using 50 μ M ATP and the luminescence signal generated from each of the amounts of the enzyme is shown. (B) Staurosporine dose response was created using 0.5ng of PKC β II to determine the potency of the inhibitor (IC₅₀).

Assay Components and Ordering Information:



Products	Company	Cat.#
ADP-Glo™ Kinase Assay	Promega	V9101
PKC β II Kinase Enzyme System	Promega	V3741
ADP-Glo + PKC β II Kinase Enzyme System	Promega	V9701

PKC β II Kinase Buffer: 40mM Tris, 7.5; 20mM MgCl₂; 0.1mg/ml BSA; 50 μ M DTT; 1 x PKC Lipid activator mix.