

TGFβ2 Kinase Assay

By Kevin Hsiao, M.S., Hicham Zegzouti, Ph.D., Jolanta Vidugiriene, Ph.D., and Said A. Goueli, Ph.D., Promega Corporation

Scientific Background:

TGFβ2 is a member of the TGFβ receptor subfamily and is a ser/thr protein kinase. TGFβ2-induced protein phosphorylation plays a key role in signal transduction that leads to mitogenic responses (1). The TGFβ2 receptor transmits signals from the cell surface to the nucleus and provides instructions for making transforming growth factor (TGF)-beta type II receptor. Mutations in TGFβ2 gene have been associated with Marfan Syndrome, Loeys-Deitz Aortic Aneurysm Syndrome, and the development of various types of tumors (2).

1. Cheng, N. et al. Enhanced hepatocyte growth factor signaling by type II transforming growth factor-beta receptor knockout fibroblasts promotes mammary tumorigenesis." *Cancer Res.* 2007;67(10):4869-77.
2. Sakai, H. et al. Comprehensive genetic analysis of relevant four genes in 49 patients with Marfan syndrome or Marfan-related phenotypes. *Am. J. Med. Genet.* 2006; 140(16):1719-25.

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.

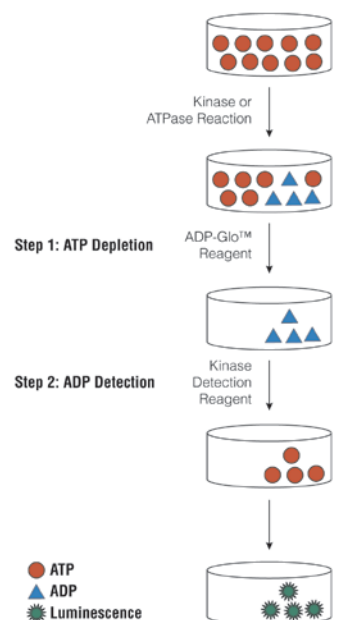


Figure 1. Principle of the ADP-Glo™ Kinase Assay. The ATP remaining after completion of the kinase reaction is depleted prior to an ADP to ATP conversion step and quantitation of the newly synthesized ATP using luciferase/luciferin reaction.

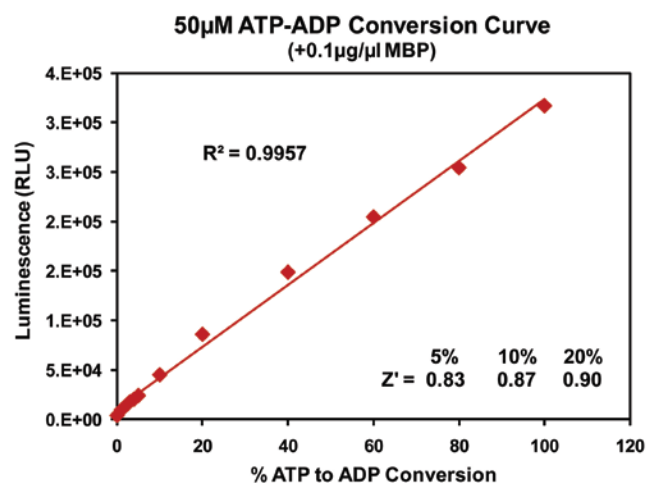


Figure 2. Linearity of the ADP-Glo Kinase Assay. ATP-to-ADP conversion curve was prepared at 50μM ATP+ADP concentration range. This standard curve is used to calculate the amount of ADP formed in the kinase reaction. Z' factors were determined using 200 replicates of each of the % conversions shown.

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 120 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. TGF β R2 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.

TGF β R2 , ng	200	100	50	25	12.5	6.3	0
RLU	360099	183498	88199	35400	21956	11836	4918
S/B	73	37.3	17.9	7.2	4.5	2.4	1
% Conversion	102	50	22	7.0	3.1	0.14	0

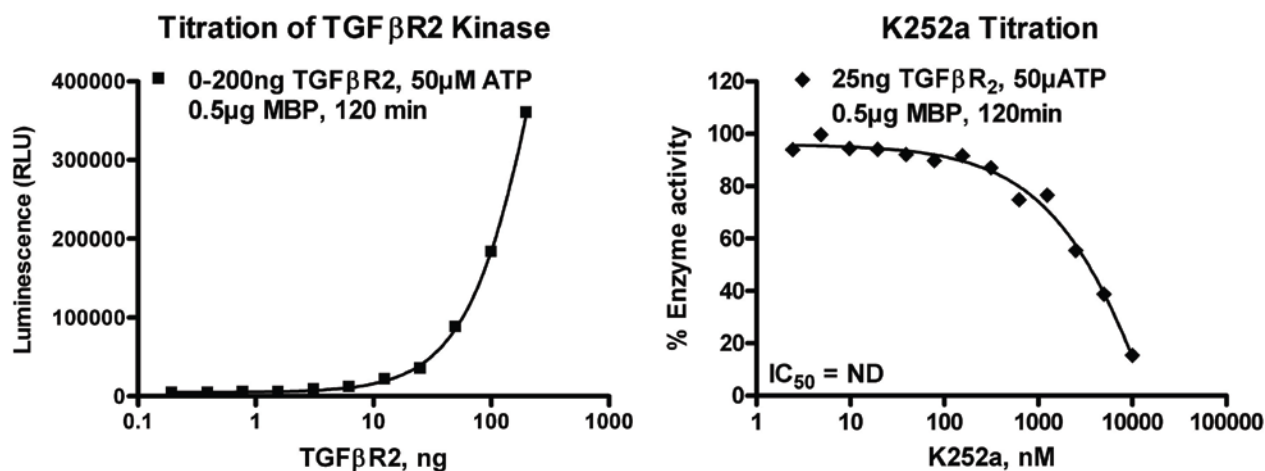


Figure 3. TGF β R2 Kinase Assay Development. (A) TGF β R2 enzyme was titrated using 50 μ M ATP and the luminescence signal generated from each of the amounts of the enzyme is shown. (B) K252a dose response was created using 25ng of TGF β R2 to determine the potency of the inhibitor (IC₅₀).

Assay Components and Ordering Information:



Products

ADP-Glo™ Kinase Assay
 TGF β R2 Kinase Enzyme System
 ADP-Glo™ + TGF β R2 Kinase Enzyme System

Company

Promega
 Promega
 Promega

Cat.#

V9101
 V3931
 V8301

TGF β R2 Kinase Buffer: 40mM Tris, 7.5; 20mM MgCl₂; 0.1mg/ml BSA; 50 μ M DTT.