

Application Note

Introduction

Transcreener® is a universal, high throughput biochemical assay based on detection of nucleotides, which are formed by thousands of cellular enzymes — many of which catalyze the covalent regulatory reactions that are central to cell signaling and are high value targets in drug discovery. The advantages of the Transcreener® HTS Assay Platform over existing assay methods include the following:

Universality

The detection of invariant nucleotide reaction product means that a single set of detection reagents can be used for all of the enzymes in a family and all acceptor substrates.

Far Red Fluorescence Intensity Detection

Use of far red shifted dyes with a simple relative fluorescence output greatly reduces interference and particulate-based light scattering from fluorescent compounds.

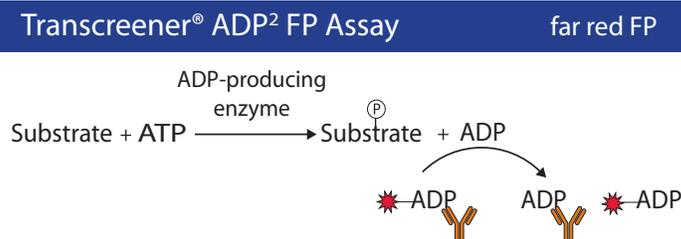
Sensitivity

High affinity antibodies enable robust detection of low levels of substrate conversion (<10%) with less enzyme than other methods.

Assay Principle

The Transcreener ADP² assay is a simple one-step homogenous detection assay. Transcreener ADP Detection Mixture, comprised of an ADP Alexa633 Tracer bound to an ADP² Antibody, is added to an equal volume of enzyme reaction mix. Enzymatically generated ADP displaces the tracer resulting in a decrease in fluorescence polarization. Standard curves are generated using varying concentrations of ATP and ADP to mimic the conversion of ATP to ADP during the course of an enzyme reaction.

Transcreener® ADP² FP Assay Principle



Materials

Instrument: Microplate Reader

Microplates: Corning® 384 Well Low Volume Black Round Bottom PS NBS™ Microplate (Product #3676)

Reagent	Kit/Component Catalog #
Transcreener® ADP ² FP Assay	3010-1K
ADP Alexa 633 Tracer	2009
Stop & Detect Buffer B, 10X	2027
ADP ² Antibody	2051
5 mM ADP	2052
5 mM ATP	2053
Buffer Components	
500 mM EGTA	Not Provided
1000 mM HEPES	
500 mM MgCl ₂	
1% Brij-35	
100% DMSO	

Protocol

Enzyme Reaction:

The enzyme was titrated in 7.5 μ L reaction volume in the presence of substrate in 4 mM MgCl₂, 2 mM EGTA, 50 mM HEPES, pH 7.5, 1% DMSO, and 0.01% Brij-35. 2.5 μ L of ATP was added to the enzyme / substrate mix to start the kinase reaction. After incubating at room temperature for a specific period of time, the kinase reaction was stopped by addition of 10 μ L of detection mix that consisted of 4nM tracer, 1X Stop and Detect buffer B and ADP² antibody. The concentration of the antibody depended on the ATP concentration and was determined from the linear equation. For 100 μ M ATP, the antibody concentration was 109 μ g/mL.

Standard Curve Reaction:

Standard curve was set up by adding 10 μ L of ATP/ADP mixes in 4 mM MgCl₂, 2 mM EGTA, 50 mM HEPES, pH 7.5, 1% DMSO, and 0.01% Brij-35. The total adenine concentration was maintained at 100 μ M. After incubating at room temperature for one hour, 10 μ L of detection mix was added which consisted of 4 nM tracer, 1X Stop and Detect buffer B and ADP² antibody. The concentration of the antibody depended on the ATP concentration and was determined from the linear equation. For 100 μ M ATP, the antibody concentration was 109 μ g/mL.

Controls

Free Tracer:

The Free Tracer consists of 20 μL of 0.5X Stop & Detect Buffer B, and 2 nM ADP Alexa633 Tracer.

Buffer Blank:

The Buffer Blank consists of 10 μL of the Kinase buffer plus 10 μL of 1X Stop & Detect Buffer B, and 109 $\mu\text{g}/\text{mL}$ ADP2 Antibody.

Final concentration of components in 20 μL reaction:

2 mM MgCl_2 , 1 mM EGTA, 25 mM HEPES (pH 7.5), 0.5% DMSO, 0.005% Brij-35, 0.5X Stop & Detect Buffer B (25 mM HEPES, pH 7.5, 200 mM NaCl, 10 mM EDTA, and 0.01% Brij-35), 2 nM ADP Alexa633 Tracer, and 109 $\mu\text{g}/\text{mL}$

Initial Reaction Kinetics

An appropriate working concentration of PKA was determined by running an enzyme titration in the presence of saturating concentrations of Kemptide and ATP. This experiment was run to determine the initial rates of the enzyme. PKA was titrated in the presence of 100 μM Kemptide and 100 μM ATP in a total reaction volume of 10 μL . The reactions were stopped at different time intervals, for a period of one hour by adding 1X Stop and Detect Buffer B with 4 nM Alexa 633 tracer and antibody concentration at 109 $\mu\text{g}/\text{mL}$ as determined from the linear equation in the technical manual. A standard curve of 100 μM ATP/ADP was run in the same plate. The raw mP values were converted to ADP based on the standard curve. The data indicated that 6 mU/Rxn of PKA was sufficient to produce a ΔmP 145 and produced 6 μM ADP at the end of one hour, which corresponds to less than 10% conversion of ADP.

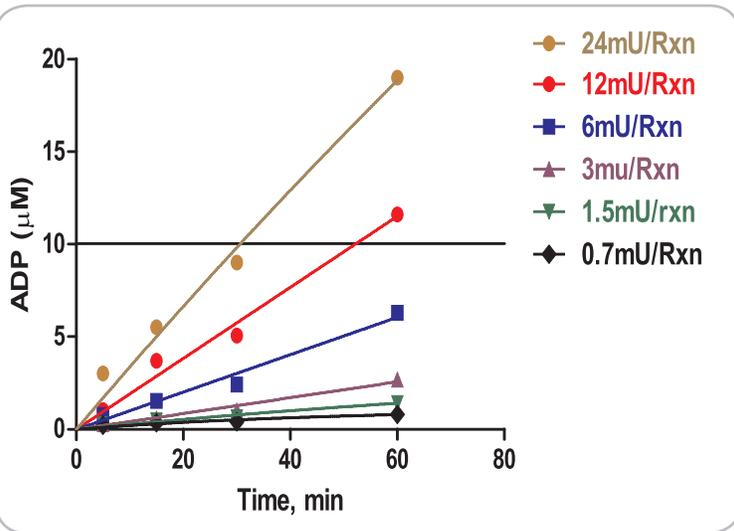


Figure 1. Initial rates of PKA enzyme in triplicates of 20 μL reactions containing varying amounts of PKA, 100 μM ATP and 100 μM Kemptide showing the amount of product formed with time.

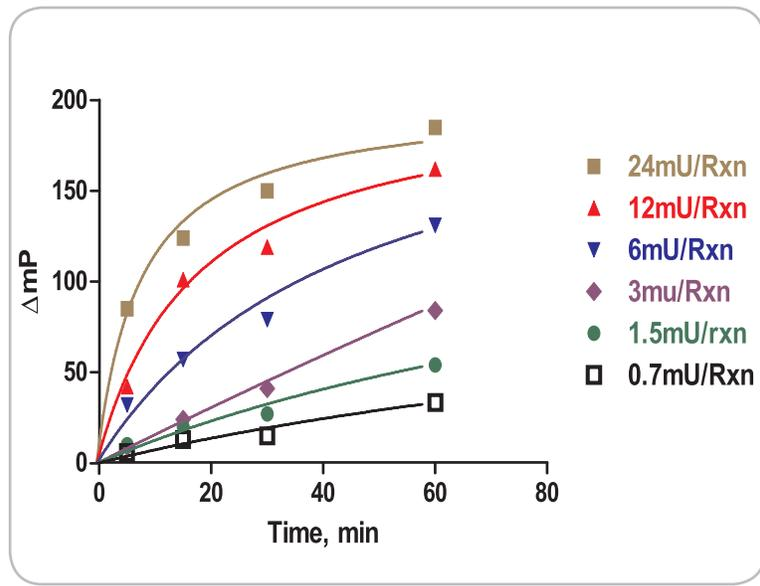


Figure 2. Initial rates of PKA enzyme in triplicates of 20 μL reactions containing varying amounts of PKA, 100 μM ATP and 100 μM Kemptide showing the change in polarization with time.

K_m Determinations

K_m values were determined for Kemptide and ATP using the Transcreener ADP² FP assay. The K_m of Kemptide was determined using saturating concentrations of ATP and varying the acceptor substrate, Kemptide concentration. 10 μL enzyme reactions containing increasing concentrations of kemptide, 6 mU of PKA/ reaction and 100 μM ATP was run. The reactions were stopped at time=0, 5, 15, 30, 45 and 60 minutes by adding 10 μL of 1X stop and Detect buffer B with 4nM tracer and 109 $\mu\text{g}/\text{mL}$ of ADP² antibody as determined from the linear equation. A 100 μM ATP/ADP standard curve was run on the same plate. The raw mP values were converted to ADP based on the standard curve. The initial reaction velocities (V) were calculated for each substrate concentration by finding the slopes of product formed vs. time during the first 60 minutes of the reaction. K_m was determined by plotting the (V) vs. substrate concentration and applying non linear regression analysis using the Michaelis –Menton equation. The K_m of Kemptide was determined to be 2 μM .

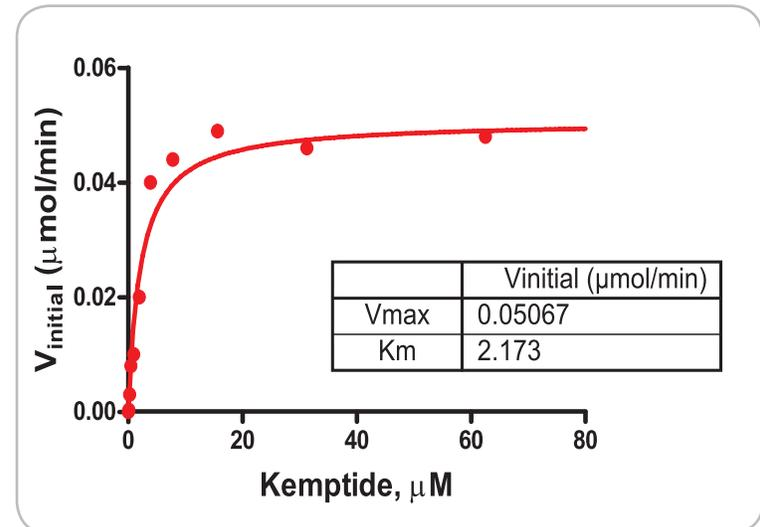


Figure 3. The plot shows Michaelis-Menten non linear regression analysis of initial rates vs peptide concentration. Reactions were triplicates of 20 μL reactions with 6 mU PKA/Rxn, 100 μM ATP with increasing concentrations of Kemptide.

The K_m of ATP was determined using saturating concentrations of Kemptide and varying the ATP concentration. 10 μL enzyme reactions containing different ATP amounts, 6 mU of PKA/reaction and 100 μM Kemptide was run. The reactions were stopped at time- 0, 5, 15, 30, 45 and 60 minutes by adding 10 μL of 1X stop and Detect buffer B with 4 nM tracer and ADP2 antibody, whose concentration depended on the amount of ATP used for each point and was determined from the linear equation. Standard curves for each of the ATP concentration was run on the same plate to convert the raw mP values to determine the amount of ADP formed. The initial reaction velocities (V) were calculated for each substrate concentration by finding the slopes of product formed vs. time during the first 60 minutes of the reaction. K_m was determined by plotting the (V) vs. substrate concentration and applying non linear regression analysis using the Michaelis –Menton equation. The K_m of ATP was determined to be 23 μM .

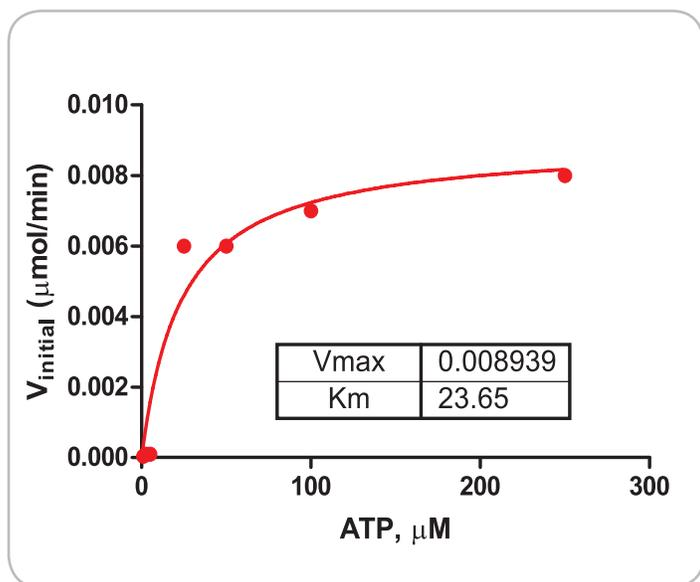


Figure 4. The plot shows Michaelis-Menten non linear regression analysis of initial rates vs ATP concentration. Reactions were triplicates of 20 μL reactions with 6 mU PKA/Rxn, 100 μM Kemptide with increasing concentrations of ATP.

Summary

K_m of Kemptide was determined to be 2 μM .

K_m of ATP was determined to be 23 μM .

For inhibitor screening we recommend using the K_m of ATP and saturating concentrations of acceptor substrate.

Additional Information

Related Products

Transcreener [®] ADP ² FP Assay.....	3010-1K
Transcreener [®] ADP ² TR-FRET Red Assay.....	3011-1K
Transcreener [®] ADP ² FI Assay.....	3013-1K
Transcreener [®] AMP/GMP Assay.....	3006-1K
Transcreener [®] UDP Assay.....	3007-1K
Transcreener [®] GDP FP Assay.....	3009-1K
Transcreener [®] GDP FI Assay.....	3014-1K
Transcreener [®] AMP ² /GMP ² Assay.....	3015-1K
Transcreener [®] UDP ² Assay.....	3018-1K
Transcreener [®] EPIGEN Methyltransferase Assay.....	3017-1K

Technical Information

For technical information, please contact
Meera Kumar, Applications Scientist

608 · 443 · 2400

866 · 313 · 7881

meera.kumar@bellbrooklabs.com

Ordering Information

Please contact BellBrook Labs for product pricing.
Custom quotes are available for orders of 10,000 wells or more.

Phone orders:
608 · 443 · 2400
866 · 313 · 7881

Fax orders:
608 · 441 · 2967

Email orders:
info@bellbrooklabs.com

Transcreener[®] HTS Assay Platform is a patented technology of BellBrook Labs. Transcreener[®] is a registered trademark of BellBrook Labs. AlexaFluor[®] is a registered trademark of Molecular Probes, Inc (Invitrogen).

©2010 BellBrook Labs. All rights reserved.
Rev. 062510