

Tracy Worzella<sup>1</sup>, Brad Larson<sup>1</sup>, Karen Kleman-Leyer<sup>1</sup>, Tony A. Klink<sup>1</sup>, Andrew Kopp<sup>1</sup>, Ingo Montenbruck<sup>2</sup>

<sup>1</sup>BellBrook Labs, Madison, WI, USA; <sup>2</sup>Tecan Austria GmbH, Grödig, Austria

## Overview

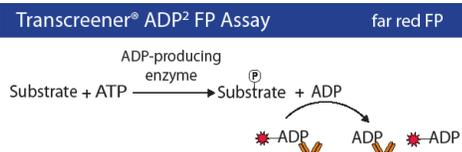
We present instrument validation data and enzyme application data in 384-well format using BellBrook Labs' Transcreener® ADP<sup>2</sup> FP Assay and the Tecan Safire<sup>2</sup>™ microplate reader. The Transcreener ADP<sup>2</sup> Assay is a universal, second generation far red fluorescence polarization assay that detects the ADP nucleotide produced in most ATP-utilizing enzyme reactions. The increased sensitivity of the new ADP<sup>2</sup> antibody allows for detection of ADP at less than 10% ATP consumption, over broader initial ATP concentrations ranging from 0.1 μM to 1,000 μM. Data generated with a standard curve on the Safire<sup>2</sup> met BellBrook Labs' Instrument Validation Program criteria, which include  $Z' > 0.7$  with  $\Delta mP$  shift  $\geq 95$  at 10% conversion of 10 μM ATP to ADP. Optimized instrument settings are provided, allowing the researcher to quickly utilize the assay with their Tecan reader. In addition, we use the Protein Kinase A (PKA) enzyme as a model system to show utility of the assay at ATP concentrations ranging from 0.1 μM to 1,000 μM ATP using one set of reagents.

## Introduction

BellBrook Labs has developed a universal assay technology called Transcreener® HTS Assay Platform that enables seamless incorporation of hundreds of new drug targets into HTS. The Transcreener HTS platform relies on a proprietary fluorescence detection method for group transfer enzymes that enables an entire family of enzymes to be screened using the same detection reagents. The Transcreener ADP<sup>2</sup> FP Assay is a universal, second generation far red fluorescence polarization assay that detects the ADP nucleotide by-product from any ATP-utilizing enzyme reaction. The increased sensitivity of the new ADP<sup>2</sup> antibody allows for detection of ADP at less than 10% ATP consumption, or at initial velocity enzyme reaction rates, and at broader ATP concentrations ranging from 0.1 μM to 1,000 μM. The improvement in sensitivity of the ADP<sup>2</sup> antibody over the previous Transcreener ADP antibody allows the researcher to expand their high throughput screening capabilities to low activity enzymes or those possessing low ATP  $K_m$ .

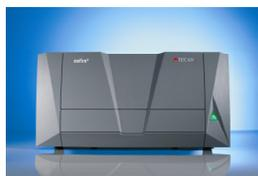
A critical factor in achieving success with an assay signal is to ensure that the microplate reader is compatible with the assay. In this study, we evaluated the Tecan Safire<sup>2</sup> instrument's excitation and emission wavelengths, number of flashes per well, and the delay time between well read using a 10 μM ADP/ATP standard curve. From this curve we determined Z'-factor and mP shift at 10% ATP conversion. We show that data generated with a standard curve in 384-well format on the Safire<sup>2</sup> meet BellBrook Labs' Instrument Validation Program criteria, which include  $Z' > 0.7$  with  $\Delta mP$  shift  $\geq 95$  at 10% conversion of 10 μM ATP. The attributes of the new Transcreener Assay were demonstrated by performing PKA titrations over the range of initial ATP concentration (0.1 μM to 1,000 μM), with the convenience of using one set of Transcreener ADP<sup>2</sup> reagents.

## The Transcreener® ADP<sup>2</sup> FP Assay



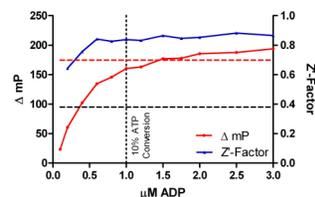
**Figure 1. The Transcreener ADP<sup>2</sup> Assay.** The Transcreener ADP<sup>2</sup> Assay was developed to follow the progress of any enzyme that produces ADP. The Transcreener ADP Detection Mixture comprises an ADP Alexa633 Tracer bound to an ADP<sup>2</sup> antibody. The tracer is displaced by ADP, the invariant product generated during an enzyme reaction. The displaced tracer freely rotates in solution leading to a decrease in fluorescence polarization. Therefore, ADP production is proportional to a decrease in polarization. The far red tracer minimizes interference from fluorescent compounds and light scattering.

## The Tecan Safire<sup>2</sup>™ Microplate Reader



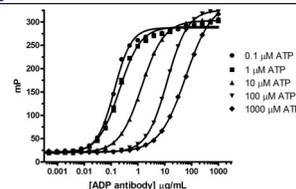
**Figure 2. The Safire<sup>2</sup>™ Microplate Reader.** Safire<sup>2</sup>™ can handle microplate formats with up to 1536-wells with high speed in every detection mode. This reader is a fully modular microplate detection system based on the quad-4 monochromator-quad-4 technology that offers a range of high speed detection techniques. Safire<sup>2</sup>™ eliminates the need for cumbersome filter changes. The autofocus feature automatically adjusts the optics to the assay volume, microplate height or well shape. The modular concept of Safire<sup>2</sup>™ allows upgrades to new detection modes at any time. Detection modules are available for top and bottom fluorescence intensity measurements, fluorescence polarization studies and multi-channel absorbance and luminescence measurements. Quasi-simultaneous readouts in FP allows high speed measurements. For FP, we used the following optimized wavelengths: excitation 635 nm, emission 670/20 nm bandwidth.

## Assay Validation at 10% ATP Conversion



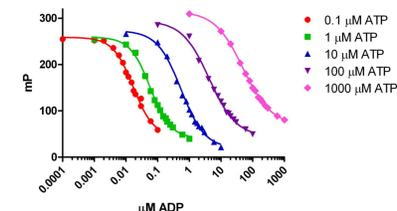
**Figure 3. The Safire<sup>2</sup>™ meets BellBrook Labs' Instrument Validation Criteria.** The validation criteria include:  $Z' \geq 0.7$  at 10% ATP conversion of 10 μM ATP, a  $\Delta mP$  shift  $\geq 95$  mP at 10% conversion and read time  $\leq 5$  minutes. The assay passed validation under the following conditions: 1 flash per well, 50 ms delay between move and read, resulting in a  $Z' = 0.84$ ,  $\Delta mP$  of 161 in 57 seconds.

## ADP<sup>2</sup> Antibody Titrations at Various ATP Concentrations



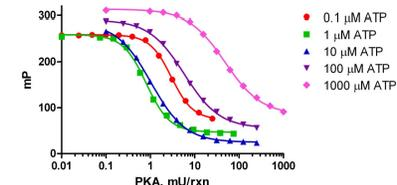
**Figure 4. Antibody titration in the presence of varying amounts of ATP.** The ADP<sup>2</sup> antibody has a finite selectivity for the nucleotide diphosphate (ADP) vs. the nucleotide triphosphate (ATP). The concentration of antibody determines the total assay window and the ADP detection range, and the amount needed is dependent on the ATP concentration in the assay. As a result, antibody must be titrated in the reaction system and ATP concentration desired in the assay. ADP<sup>2</sup> antibody was titrated as shown. The final 20 μL volume consisted of 2 nM ADP Alexa633 tracer, 0.5X Stop & Detect Buffer B, 0.5X enzyme reaction mixture (50 mM HEPES pH 7.5, 2 mM MgCl<sub>2</sub>, 1 mM EGTA, 0.5% DMSO, 0.01% Brij-35, 0.5X ATP), and ADP<sup>2</sup> antibody. The EC<sub>50</sub> antibody concentration is determined, and that concentration used in an ADP/ATP standard curve.  $EC_{50} = ((85/(100-85))^{1/\text{HillSlope}}) * EC_{50}$ .

## ADP/ATP Standard Curves



**Figure 5. ADP/ATP standard curves mimic an enzyme reaction.** In an enzyme reaction, ATP concentration decreases as ADP concentration increases; adenine concentration remains constant. Standard curves were run using the EC<sub>50</sub> antibody concentration. The final concentration in the 20 μL reaction consisted of 2 nM ADP Alexa633 tracer, 0.5X Stop & Detect Buffer B, and 0.5X enzyme reaction mixture (25 mM HEPES pH 7.5, 2 mM MgCl<sub>2</sub>, 1 mM EGTA, 0.5% DMSO, 0.005% Brij-35, 0.5X ATP). The EC<sub>50</sub> antibody concentrations for each ATP concentration are as follows: 0.4 μg/mL ADP<sup>2</sup> antibody (0.1 μM ATP); 1 μg/mL ADP<sup>2</sup> antibody (1 μM ATP); 9 μg/mL ADP<sup>2</sup> antibody (10 μM ATP); 58 μg/mL ADP<sup>2</sup> antibody (100 μM ATP); 535 μg/mL ADP<sup>2</sup> antibody (1,000 μM ATP).

## PKA Titrations



**Figure 6. PKA titrations at varying concentrations of ATP.** PKA serial titrations (1:2) were performed in a 10 μL reaction with the ATP concentrations indicated above and kemptide substrate at 5X ATP concentration. Reactions were incubated for one hour at room temperature. An equal volume of detection reagent was then added and reactions equilibrated for one hour before reading. The concentration of components in the final 20 μL reaction were as noted previously. Enzyme concentrations that resulted in 10% ATP conversion (interpolated from an ADP/ATP standard curve at each ATP concentration) were as follows: 3 mU/rxn (0.1 μM ATP); 1 mU/rxn (1 μM ATP); 2 mU/rxn (10 μM ATP); 16 mU/rxn (100 μM ATP); 125 mU/rxn (1,000 μM ATP).

## Conclusions

- For optimal performance of the Transcreener ADP<sup>2</sup> Assay on the Tecan Safire<sup>2</sup> reader, we recommend the following: excitation wavelength 635 nm, emission wavelength 670 nm with a 20 nm bandwidth.
- The Safire<sup>2</sup> passes BellBrook Labs' Validation Program under the following conditions: 1 flash, 50 ms delay between move and read. These settings yielded a  $Z' > 0.7$  in 57 seconds.
- The Transcreener ADP<sup>2</sup> FP Assay can be used with a wide range of initial ATP concentrations, simplifying assay optimization and reducing reagent cost.
- The Transcreener Assay can detect ADP production from an enzyme reaction within initial velocity rates ( $\leq 10\%$  ATP conversion) using ATP concentrations from 0.1 μM to 1,000 μM.

For more information on Transcreener validation on Tecan readers, or for a copy of this poster, please visit Tecan at booth # 205.

©2008 BellBrook Labs. All rights reserved.  
Transcreener is a registered trademark of BellBrook Labs.  
Transcreener is a patented technology of BellBrook Labs.  
AlexaFluor® is a registered trademark of Molecular Probes, Inc. (Invitrogen).  
BellBrook Labs 5500 Nobel Drive, Suite 250, Madison, WI 53711  
866.313.7881 or 608.443.2400