

Tracy Worzella<sup>1</sup>, Brad Larson<sup>1</sup>, Karen Kleman-Leyer<sup>1</sup>, Tony A. Klink<sup>1</sup>, Xavier Amouretti<sup>2</sup>, Peter Banks<sup>2</sup>  
<sup>1</sup>BellBrook Labs, Madison, WI, USA; <sup>2</sup>BioTek Instruments, Winooski, VT, USA

## Overview

We present data showing the validation of BellBrook Labs' Transcreener® ADP<sup>2</sup> FP Assay on the BioTek Synergy™ 2 and Synergy™ 4 plate readers. The Transcreener ADP<sup>2</sup> 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 assay is amenable to volume scaling, as shown by data generated in 96-well and 384-well microplate formats. Data generated with a standard curve in both microplate formats resulted in Z' values  $\geq 0.6$ . Further, the 96-well assay using the standard PMT and the 384-well assay using the red-shifted PMT met BellBrook Labs' Instrument Validation Program criteria, which include Z' > 0.7 with  $\Delta mP$  shift  $\geq 95$  at 10% conversion of 10  $\mu M$  ATP to ADP. Further enzyme studies demonstrate that 96- and 384-well plate formats yield excellent Z'-factor results with Protein Kinase A (PKA) at low % ATP conversion. Known PKA inhibitors show comparable potency in both 96- and 384-well plate formats, and between different Synergy plate readers. Although the standard PMT Synergy does not meet validation criteria in 384-well plate format, comparable inhibitor potency results are obtained between the standard and red-shifted PMT instruments.

## Introduction

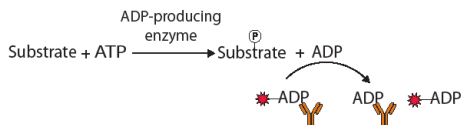
BellBrook Labs has developed a universal assay technology called Transcreener® HTS Assay Platform that enables seamless incorporation of hundreds of new drug targets into HTS. The Transcreener HTS platform relies on a proprietary fluorescence detection method for group transfer enzymes that enables an 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 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, or at initial velocity enzyme reaction rates, and at broader ATP concentrations ranging from 0.1  $\mu M$  to 1,000  $\mu M$ .

A critical factor in achieving success with an assay is to ensure that the detection instrumentation is compatible with the reagents being used. Assay reagents must also demonstrate scalability from larger volume to smaller volume wells to accommodate higher throughput processing for the transition from assay development to screening. We present data showing the validation of BellBrook Labs' Transcreener ADP<sup>2</sup> FP Assay on the BioTek Synergy™ 2 and Synergy™ 4 plate readers. We evaluated different filter sets and number of flashes per well using a 10  $\mu M$  ADP/ATP standard curve in both 96- and 384-well format. From the curve we determined Z'-factor and mP shift at 10% ATP conversion at different settings.

Attributes of the new Transcreener Assay were demonstrated in both well formats by performing PKA titrations, Z'-factor determinations, and inhibitor potency studies. Results indicate that researchers can have confidence in both reagent and reader performance regardless of which well format or detection instrumentation they are using.

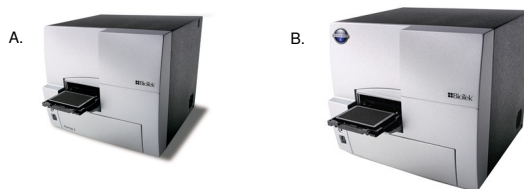
## The Transcreener® ADP<sup>2</sup> FP Assay

### Transcreener® ADP<sup>2</sup> FP Assay far red FP



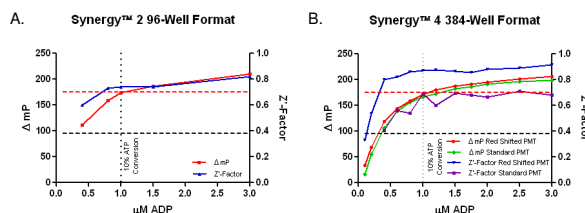
**Figure 1. The Transcreener ADP<sup>2</sup> Assay.** The Transcreener ADP<sup>2</sup> Assay was developed to follow the progress of any enzyme that produces ADP. The Transcreener ADP Detection Mixture comprises an ADP Alexa633 Tracer bound to an ADP<sup>2</sup> antibody. The tracer is displaced by ADP, 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 BioTek™ Synergy 2 and 4 Multi-Mode Microplate Readers



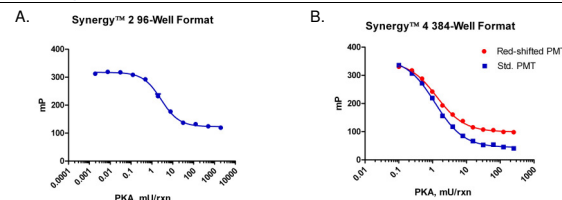
**Figure 2. The Synergy 2 and Synergy 4 Microplate Readers.** The Synergy 2 and 4 plate readers can be used in well formats from 6- to 1536-well plates. Both readers have 2 broad spectrum light sources for fluorescence. Gen5 Data Analysis Software comes standard with both readers. The Synergy 2 (A) can be customized with combinations of monochromator, filters and dichroic mirrors for optimal performance in all detection modes. The Synergy 4 (B) contains Hybrid Technology comprising a high-performance filter system with flexible monochromator detection. For optimal FP performance, an excitation filter at 620/40 nm was used, combined with an emission filter at 680/30 nm and 660 nm cutoff dichroic mirror. The instrument comes with a standard PMT, or optional red-shifted PMT detector which we tested in this study.

## Assay Validation at 10% ATP Conversion



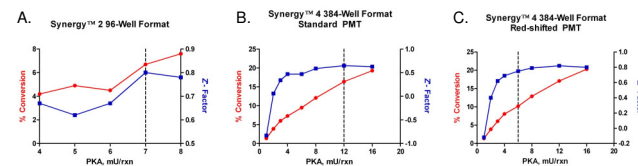
**Figure 3. The Synergy 2 and 4 meet BellBrook Labs' Instrument Validation Criteria.** The validation criteria include: Z'  $\geq 0.7$  at 10% conversion of 10  $\mu M$  ATP, a  $\Delta mP$  shift  $\geq 95$  mP at 10% conversion and read time  $\leq 5$  minutes. The 96-well format assay passed validation under the following conditions: 15 flashes per well, Z' = 0.74,  $\Delta mP$  of 174 in 1 minute, 39 seconds. The 384-well format assay passed validation with the red-shifted PMT under the following conditions: 3 flashes per well, Z' = 0.81,  $\Delta mP$  of 168 in 1 minute, 52 seconds. The 384-well format assay using the standard PMT did not meet the validation criteria of the Program, but results yielded acceptable data by many standards: 12 flashes per well, Z'  $\geq 0.6$ ,  $\Delta mP$  of 166 in 4 minutes, 11 seconds.

## PKA Enzyme Titrations



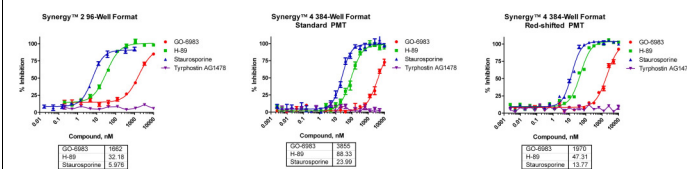
**Figure 4. PKA titrations in 96- and 384-well plate formats show comparable EC<sub>50</sub>.** PKA was titrated (1:2) and reactions were performed in a 10  $\mu L$  assay volume with 10  $\mu M$  ATP and 50  $\mu M$  kemptide substrate, and incubated for one hour. Detection Mixture was added and reactions equilibrated for one hour before reading. The final 100  $\mu L$  (96-well) and 20  $\mu L$  (384-well) assay volume with the Detection Mixture consisted of 2 nM ADP Alexa633 tracer, 0.5X Stop & Detect Buffer B, 8.5  $\mu g/mL$  ADP<sup>2</sup> antibody, 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, 5  $\mu M$  ATP and 25  $\mu M$  kemptide). EC<sub>50</sub> for 96-well assay was 2.8 mU/rxn. EC<sub>50</sub> for 384-well assay was 1.3 mU/rxn (red-shifted PMT), and 1.2 mU/rxn (standard PMT).

## Z'-Factor at Low % ATP Conversion



**Figure 5. Excellent Z' values achieved at low % ATP conversion.** Z'-factor experiments were run to confirm % conversion achieved in the PKA titration. Enzyme concentrations were chosen as indicated on the graphs above, and reactions were performed under the same conditions used for the enzyme titration. Z'-factor was determined by comparing wells with enzyme to wells without enzyme. Enzyme concentrations chosen for subsequent inhibitor titration experiments achieved a Z'  $\geq 0.7$ , highlighted by the dashed lines above. 96-well format (A): 7 mU/rxn achieved a Z' = 0.8 at 7% conversion. 384-well format, standard PMT (B): 12 mU/rxn achieved Z' = 0.7 at 16% conversion. 384-well format, red-shifted PMT (C): 6 mU/rxn achieved Z' = 0.8 at 10% conversion.

## PKA Inhibitor Potency Study



**Figure 6. Potency of PKA inhibitors are comparable between well formats.** GO-6983, H-89, and staurosporine (known PKA inhibitors) and Tyrphostin AG1478 (a no inhibition control for PKA) were serially titrated into the PKA reaction following the assay conditions noted previously. Results show that the potency of compound inhibition (nM) of the known inhibitors are all comparable between the different well formats tested. There is also good agreement between the inhibitor results obtained in 384-well format using the standard and red-shifted PMT.

## Conclusions

- For optimal performance of Transcreener ADP<sup>2</sup> Assay on BioTek microplate readers, we recommend the following: 620/40 nm excitation filter, 680/30 nm emission filter, 660 nm cutoff dichroic mirror.
- The Synergy 2 and 4 pass BellBrook Labs' Validation Program in 96-well format under the following conditions: standard PMT, 15 flashes/well to yield Z' > 0.7 in 1 minute, 39 seconds.
- The Synergy 2 and 4 pass BellBrook Labs' Validation Program in 384-well format under the following conditions: red-shifted PMT, 3 flashes/well to yield Z' > 0.7 in 1 minute, 52 seconds.
- Studies with a PKA model system in 96- and 384-well plate formats show comparable results on the Synergy 2 and 4 microplate readers.

For more information on Transcreener validation on BioTek readers, or for a copy of this poster, please visit BioTek at booth # 443.

©2008 BellBrook Labs. All rights reserved.  
 Transcreener is a registered trademark of BellBrook Labs.  
 Synergy 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