

## Optimized Settings to Validate the Tecan Infinite<sup>®</sup> M1000/Safire<sup>2™</sup>\* HTS Microplate Readers with the Transcreener<sup>®</sup> Fluorescence Polarization Assays.

Meera Kumar<sup>1</sup>, Karen Kleman-Leyer<sup>1</sup> and Christian Oberdanner<sup>2</sup>  
<sup>1</sup>BellBrook Labs, Madison, WI, USA; <sup>2</sup>Tecan Austria GmbH, Grödig, Austria

*This Application Note describes the optimal instrument parameters used to validate the Tecan Infinite<sup>®</sup> M1000/Safire<sup>2™</sup> plate readers with the following assays:*

*Transcreener<sup>®</sup> ADP<sup>2</sup> FP (3010) Transcreener<sup>®</sup> AMP/GMP (3006)  
 Transcreener<sup>®</sup> UDP FP (3007) Transcreener<sup>®</sup> GDP FP (3009)*

### Introduction

Transcreener<sup>®</sup> HTS is a universal, high throughput biochemical assay platform based on the 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 of great value as targets in drug discovery.

The Transcreener<sup>®</sup> FP Assays are a single step, competitive immunoassay for direct detection of nucleotides with a far red fluorescence polarization (FP) readout. The reagents for all of the assays are a far red Tracer bound to a highly-specific monoclonal/polyclonal antibody. Nucleotide diphosphate or monophosphate produced by the target enzyme displaces the tracer from the antibody, leading to increased rotational freedom and results in a decrease in polarization (Figure 1). The use of a far red tracer minimizes interference from fluorescent compounds and light scattering. The Transcreener<sup>®</sup> FP Assays are designed specifically for HTS with a single addition, mix-and-read format.

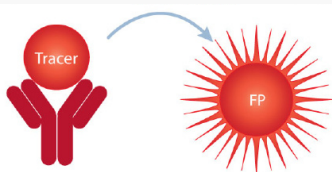


Figure 1. Transcreener<sup>®</sup> FP Assay Principle

### Validation Criteria

A critical factor in realizing the advantages of the Transcreener<sup>®</sup> HTS assays is the correct setup of the microplate reader used for data readout. Proper selection of filters, dichroics, gain and flashes can impact the instrument's sensitivity for any given assay. The key instrument parameters for Transcreener<sup>®</sup> HTS assay performance were identified by running a 10  $\mu\text{M}$  ATP/ADP standard curve (24 replicates), as standard curves of this type mimic enzyme reactions. Starting with 10  $\mu\text{M}$  ATP, ADP was added in increasing amounts and ATP is decreased proportionately, maintaining a total adenine nucleotide concentration of 10  $\mu\text{M}$ . The flash numbers were varied to determine the requirements for a  $Z' > 0.5$ . *In order to validate an instrument for use with the Transcreener<sup>®</sup> FP Assays, a  $Z' > 0.7$  and a  $\Delta mP > 120$  at 10% conversion of 10  $\mu\text{M}$  ATP were required.*

### Infinite<sup>®</sup> M1000/Safire<sup>2™</sup> Information

- Quad-4 Monochromators<sup>™</sup> technology.
- Z-Focusing in Top Reading Modes - Fluorescence Intensity, Fluorescence Polarization, and Luminescence.
- User-Selected Frequency Modes (patented) - Choose between "High Sensitivity" and "High Speed" settings to adjust for individual assay needs.
- Fast Ratiometric Measurements - Immediate changes in wavelength in (TR)-FRET based assays and instant change of polarization direction for FP measurements.
- Free Definition of Plate Formats - Up to 1,536 wells.
- Top & Bottom Reading - Variety of microplate formats in the fluorescence mode.
- Modular Design - Allows upgrades to new detection modes at any time if further applications are required.
- Optional Stacker Module - Batch processing of up to 50 microplates.
- i-Control<sup>™</sup> Software - Convenient workflow-oriented reader control.



### Instrument Settings

Instrument Wavelength Settings	
EXC Wavelength and Bandwidth	635/5 nm
EMS Wavelength and Bandwidth	680/20 nm
Optimized Infinite <sup>®</sup> M1000 Settings	
Gain	Optimal
Z-Position	Calculate from Well containing Fluor
Settle Time	50 msec
Flash Number	1-20

Table 1. Recommended Tecan Infinite<sup>®</sup> M1000 Instrument Settings

\* Safire<sup>2™</sup> Instrument No Longer Commercially Available.

## Instrument Setup

1. Open the Tecan i-control™ software, connect to the Infinite® M1000 reader, and open a new method file.
2. Select the appropriate labware/plate from the "Plate Definition" dropdown menu, and select the portion of the labware/plate to be read.
3. Select "Fluorescence Polarization" from the "Measurements" menu.
4. Select "635 (5 nm)" from the "Excitation" dropdown menu and "670" for "Emission (20.0 nm for Bandwidth)".
5. To allow the instrument to calculate the proper gain based on each well, select "Optimal" for the Gain. If the plate is being read more than once, select "Manual" for the Gain and input the value that the instrument previously calculated.
6. To set the measurement blank, select "Change" from the "Measurement" section and select the wells containing the buffer blank.

The same measurement settings can be used for subsequent plates as long as the volumes, tracer and concentrations remain the same.

## Sample FP Standard Curve

As the ratio of ADP:ATP increases, the proportion of bound tracer vs. free tracer decreases, resulting in an overall decrease in mP values. Assay plates containing the 15-point standard curve were read on the Tecan Infinite® M1000 Microplate Reader.

### Materials

**ATP/ADP Mixture** - 4 mM MgCl<sub>2</sub>, 2 mM EGTA, 50 mM HEPES, pH 7.5, 1% DMSO, 0.01% Brij-35, and ATP/ADP (combined to a constant adenine concentration of 10 μM).

**ADP Detection Mixture** - 1X Stop & Detect Buffer B, 4 nM ADP Alexa633 Tracer, and 14.8 μg/mL ADP<sup>2</sup> Antibody.

**Free Tracer** - 1X Stop & Detect Buffer B and 4 nM ADP Alexa633 Tracer.

**Buffer Blank** - 1X Stop & Detect Buffer B and 14.8 μg/mL ADP<sup>2</sup> Antibody.

For a detailed procedure on how to prepare a standard curve, please refer to the appropriate Transcreeper® Technical Manual ([http://www.bellbrooklabs.com/transcreeper\\_hts\\_assays.html](http://www.bellbrooklabs.com/transcreeper_hts_assays.html)).

### Method

1. Dispense 10 μL of each ATP/ADP combination across an entire row of a 384-well plate.
2. Add 10 μL of ADP Detection Mix to those rows.
3. Dispense 10 μL of the 10 μM ATP/0 μM ADP combination into row P.
4. Dispense 10 μL of Free Tracer into wells P1-P12.
5. Dispense 10 μL of Buffer Blank into wells P13-P24.

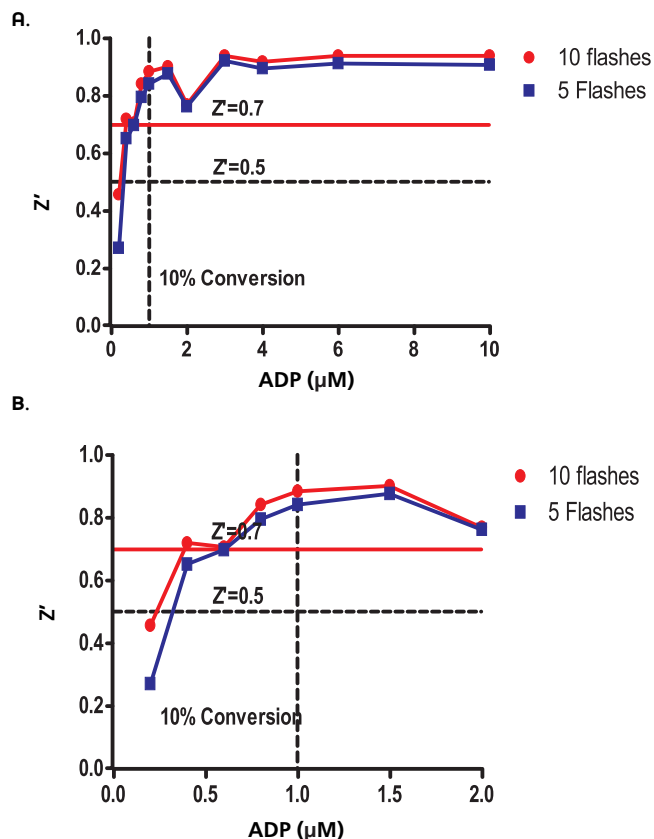


Figure 3. A). Z' values observed in a standard curve mimic conversion of 10 μM ATP to ADP. B). Zoomed view of the 0-2 μM ADP section of the standard curve shows the Z' validation minimal qualification data (red dotted line) and 10% ATP conversion validation point (black dotted line). Plate reader set at 10 or 5 flashes.

Assay Performance, 10% Conversion 10 μM ATP						
Flashes	1	3	5	10	15	20
Read Time (minutes)	0:19	0:25	0:32	0:47	1:03	1:18
Δ mP at 10% ATP Conversion	117	121	122	124	125	125
Z'-Factor at 10% ATP Conversion	0.61	0.74	0.71	0.89	0.89	0.93

Table 2. Assay Performance with Various Instrument Settings

## Conclusions

This application note demonstrates the validation of the Tecan Infinite® M1000 and Safire<sup>2</sup>™ microplate readers for use with the Transcreeper® FP Assays. By utilizing the optimized instrument settings suggested within this Application Note, Z' values > 0.7 and ΔmP > 120 are achievable.

## Additional Information

### Ordering Information

Please visit [www.bellbrooklabs.com](http://www.bellbrooklabs.com) or contact BellBrook Labs for pricing for the Transcreener® HTS Assays. Custom quotes are available for bulk orders.

#### Phone Orders:

608.443.2400

866.3137881

#### Fax Orders:

608.441.2967

#### Email Orders:

[info@bellbrooklabs.com](mailto:info@bellbrooklabs.com)

### Related Products

Transcreener® ADP <sup>2</sup> FP Assay.....	3010-1K
Transcreener® ADP <sup>2</sup> FI Assay.....	3013-1K
Transcreener® ADP <sup>2</sup> TR-FRET Red Assay.....	3011-1K
Transcreener® AMP/GMP FP Assay.....	3006-1K
Transcreener® UDP FP Assay.....	3007-1K
Transcreener® GDP FP Assay.....	3009-1K
Transcreener® GDP FI Assay.....	3014-1K

### Technical Information

For technical information, please contact:

#### **Meera Kumar, Applications Scientist**

Tel: 608.443.2400

Toll-Free: 866.313.7881

Email: [meera.kumar@bellbrooklabs.com](mailto:meera.kumar@bellbrooklabs.com)

## References & Notes

Transcreener® HTS Assay Platform is a patented technology of BellBrook Labs.

Transcreener® is a registered trademark of BellBrook Labs.

Infinite® is a registered trademark and Safire™ is a trademark of Tecan.

AlexaFluor® is a registered trademark of Molecular Probes, Inc (Invitrogen).

The Transcreener® product line is the subject of U.S. Patent No. 7,332,278, 7,355,010 and 7,378,505 issued. U.S. Patent Application Nos. 11/353,500, 11/958,515 and 11/958,965, U.S. Divisional Application 12/029,932, and foreign equivalents licensed to BellBrook Labs.

