

## Optimized Settings to Validate the PerkinElmer EnVision®/EnVision® Xcite Microplate Readers with the Transcreener® Fluorescence Intensity Assays.

Meera Kumar<sup>1</sup> and Karen Kleman-Leyer<sup>1</sup>  
<sup>1</sup>BellBrook Labs, Madison, WI, USA

*This Application Note describes the optimal instrument parameters used to validate the PerkinElmer EnVision®/EnVision® Xcite plate readers with the following assays:*

*Transcreener® ADP<sup>2</sup> FI (3013)*  
*Transcreener® GDP FI (3014)*

### Introduction

Transcreener® 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® FI Assays are a single step, competitive immunoassay for direct detection of nucleotides with a red fluorescence intensity (FI) readout. The reagents for all of the assays are a red Tracer bound to a highly-specific monoclonal antibody-quencher conjugate. Nucleotide diphosphate or monophosphate produced by the target enzyme displaces the tracer from the antibody-quencher conjugate, resulting in an increase in fluorescence intensity (Figure 1). The use of a red tracer minimizes interference from fluorescent compounds and light scattering. The Transcreener® FI Assays are designed specifically for HTS with a single addition, mix-and-read format.

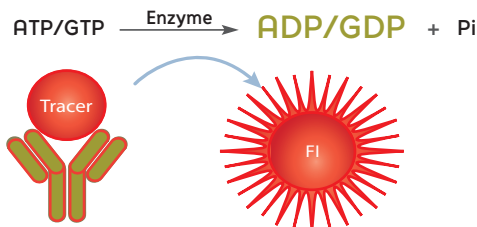


Figure 1. Transcreener® FI Assay Principle

### Validation Criteria

A critical factor in realizing the advantages of the Transcreener® 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® 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® FI Assays, a  $Z' > 0.7$  at 10% conversion of 10  $\mu\text{M}$  ATP was required.*

### EnVision®/EnVision® Xcite Information

- Reads Fluorescence Intensity, Fluorescence Polarization and TR-FRET versions of the Transcreener® Assays.
- Capable of reading 96-, 384- and 1536-well plates.
- Simultaneous dual detection capabilities.
- Interchangeable filters and dichroic modules.



### Instrument Settings

#### Instrument Wavelength Settings (PerkinElmer Catalog #)

Excitation Filter	547/7 nm (2100-5070)
Emission Filters	635/15 nm (2100-5590)
Mirror	Texas Red FP D595 (2100-4190) Single Mirror

#### Optimized EnVision® Settings

Detector Gain 1	750
Detector Gain 2	0
Measurement Height	11.8
Excitation Light (%)	100
Flash Number	Variable

Table 1. Recommended PerkinElmer EnVision® Instrument Settings

## Instrument Setup

The Transcreeper-specific FI mirrors and filters were installed prior to instrument evaluation. Once those components have been installed, proceed with the following steps to optimize the detector gains and Z-height focus:

1. Create a label by replicating an existing label.
2. Associate the installed filters and mirror with the new label in the "General" tab.
3. Create a new protocol by replicating an existing protocol and associate the new label and a plate with the protocol.
4. Add 20  $\mu\text{L}$  of Free Tracer (2 nM) to the four corners of the plate (in the buffer conditions of your enzyme reaction).
5. Run the *Label Optimization Wizard* to optimize the label. Select the appropriate protocol and then select "Plate Dimension", "Measurement Height", and "Detector Gains" to proceed with the *Label Optimization Wizard*.
6. Record the "Measurement Height" value from the "Optimization" tab. Next, delete the optimization. Then, input the *Measurement Height* value into the "General" tab for the label.
7. Run the *Label Optimization Wizard* again. This time, select only "Plate Dimension" for optimization. This allows the instrument to use the correct plate dimensions, while allowing the detector gains to be increased above the reader's recommended settings.

The same measurement settings can be used for subsequent plates as long as the volumes, tracer and concentrations remain the same. A snapshot of the *FI Transcreeper Assay 595 Mirror* screen is shown in Figure 2.

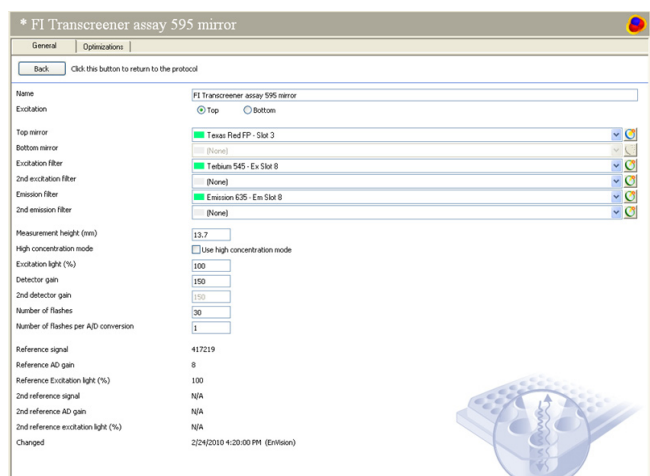


Figure 2. Screen Snapshot of FI Transcreeper Assay 595 Mirror Dialog Box

## Sample FI Standard Curve

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

### Materials

**ATP/ADP Mixture** - 4 mM  $\text{MgCl}_2$ , 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  $\mu\text{M}$ ).

**ADP Detection Mixture** - 1X Stop & Detect Buffer B, 8 nM ADP Alexa594 Tracer, and 10  $\mu\text{g}/\text{mL}$  ADP<sup>2</sup> Antibody-IRDye®QC-1.

**Free Tracer** - 1X Stop & Detect Buffer B and 8 nM ADP Alexa594 Tracer.

**Buffer Blank** - 1X Stop & Detect Buffer B and 10  $\mu\text{g}/\text{mL}$  ADP<sup>2</sup> Antibody-IRDye®QC-1.

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  $\mu\text{L}$  of each ATP/ADP combination across an entire row of a 384-well plate.
2. Add 10  $\mu\text{L}$  of ADP Detection Mix to those rows.
3. Dispense 10  $\mu\text{L}$  of the 10  $\mu\text{M}$  ATP/0  $\mu\text{M}$  ADP combination into row P.
4. Dispense 10  $\mu\text{L}$  of Free Tracer into wells P1-P12.
5. Dispense 10  $\mu\text{L}$  of Buffer Blank into wells P13-P24.

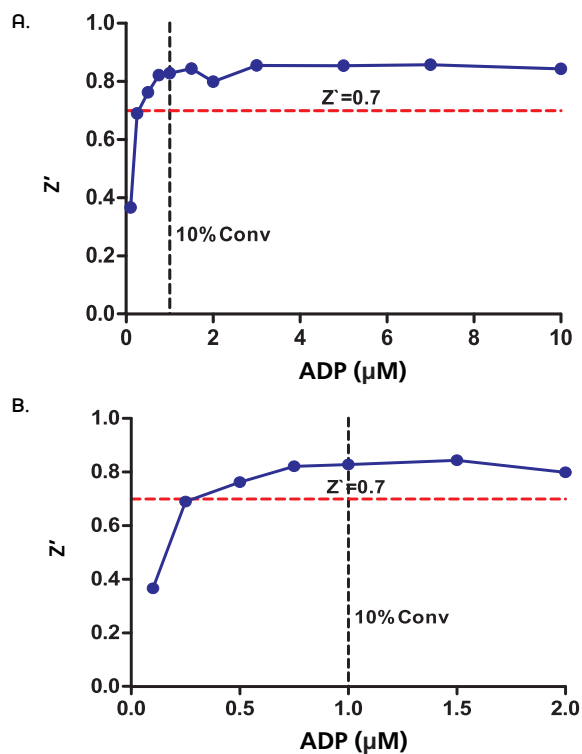


Figure 3. A). Z' values observed in a standard curve mimic conversion of 10  $\mu\text{M}$  ATP to ADP. B). Zoomed view of the 0-2  $\mu\text{M}$  ADP section of the standard curve shows the Z' validation minimal qualification data (red dotted line) and 10% ATP conversion validation point (black dotted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

Assay Performance, 10% Conversion 10 $\mu\text{M}$ ATP						
Flashes	1	5	10	20	50	100
Read Time (minutes)	4:01	4:08	4:11	4:17	4:85	5:69
% CV at 10% ATP Conversion	6.25	3.22	2.40	1.97	1.88	1.65
Z'-Factor at 10% ATP Conversion	0.69	0.84	0.88	0.91	0.92	0.92

Table 2. Assay Performance with Various Instrument Settings

## Conclusions

This application note demonstrates the validation of the PerkinElmer EnVision® and EnVision® Xcite microplate readers for use with the Transcreeper® FI Assays. By utilizing the optimized instrument settings suggested within this Application Note, Z' values > 0.7 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.

EnVision® is a registered trademark of PerkinElmer.

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

IRDye®QC-1 is a registered trademark of LI-COR Biosciences.

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.

