

## Optimized Settings to Validate the BMG LABTECH PHERAstar Plus/FS HTS Microplate Readers with the Transcreener® Fluorescence Intensity Assays.

Meera Kumar<sup>1</sup>, Brad Larson<sup>1</sup>, Franka Ganske<sup>2</sup> and EJ Dell<sup>2</sup>  
<sup>1</sup>BellBrook Labs, Madison, WI, USA; <sup>2</sup>BMG Labtech, Durham, NC, USA.

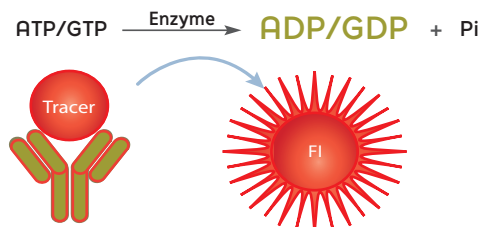
*This Application Note describes the optimal instrument parameters used to validate the BMG LABTECH PHERAstar Plus/FS plate readers with the following assays:*

*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.



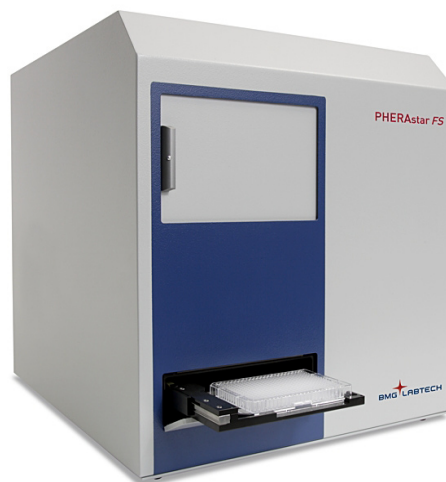
**Figure 1. Transcreener® FI Assay Principle**

### 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 µM ATP/ADP standard curve (24 replicates), as standard curves of this type mimic enzyme reactions. Starting with 10 µM ATP, ADP was added in increasing amounts and ATP is decreased proportionately, maintaining a total adenine nucleotide concentration of 10 µM. 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 µM ATP was required.*

### PHERAstar Plus/FS Information

- Eight detection modes, separate measurement electronics, 5 PMTs, and assay specific optic modules.
- Equipped to read 6- to 1536- well plates.
- SDE decreases read times and increases accuracy as measured by %CV and Z' values.



### Instrument Settings

| Instrument Wavelength Settings |                                                                 |
|--------------------------------|-----------------------------------------------------------------|
| Optical Module                 | Transcreener FI Application Specific Module (EXC: 580/EMS: 620) |
| Optimized PHERAstar Settings   |                                                                 |
| Detector Gain 1                | Should be Adjusted                                              |
| Measuring Height               | 11.3 mm                                                         |
| Positioning Delay              | 0.1 sec                                                         |
| Flash Number                   | 30-100                                                          |

**Table 1. Recommended PHERAstar Instrument Settings**

## Instrument Setup

To install the Transcreeper-specific FI Optical Module, refer to <http://www.bmg-labtech.com/technology/optic-modules.cfm> for a detailed protocol. Once the Transcreeper-specific Optical Module has been installed, proceed with the following steps to optimize the detector gains and Z-height focus:

1. Select the "Focus and Adjustment" tab from the *Measurement* screen.
2. For optimization purposes, select a well containing *only Free Tracer* from the plate layout.
3. Select "Focus Adjustment" and "Gain Adjustment".
4. Select "Start Adjustment" to begin the optimization process.

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 *Fluorescence Intensity* screen is shown in Figure 2.

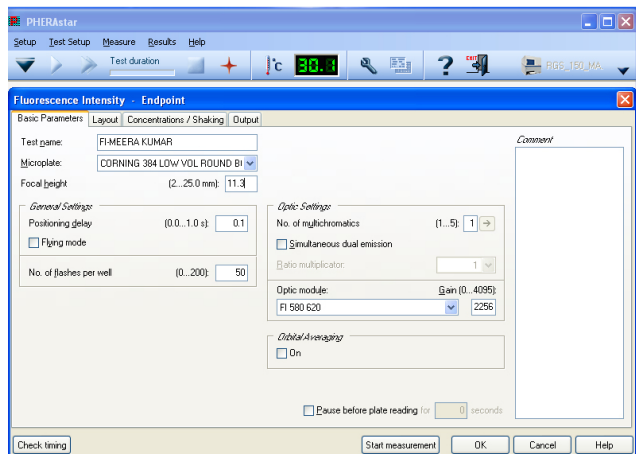


Figure 2. Screen Snapshot of Start Measurement Dialog Box

## 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 PHERAstar Plus HTS 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, 8 nM ADP Alexa594 Tracer, and 10 μg/mL ADP<sup>2</sup> Antibody-IRDye<sup>®</sup>QC-1.

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

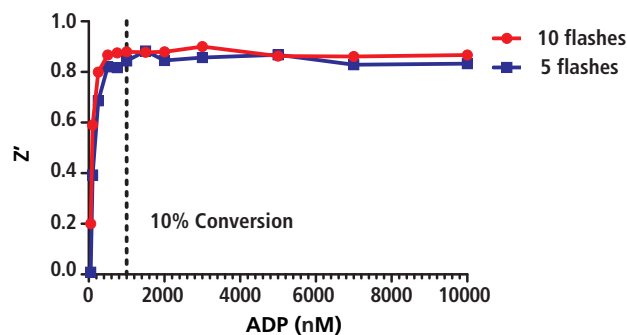
**Buffer Blank** - 1X Stop & Detect Buffer B and 10 μg/mL ADP<sup>2</sup> Antibody-IRDye<sup>®</sup>QC-1.

For a detailed procedure on how to prepare a standard curve, please refer to the appropriate Transcreeper<sup>®</sup> 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.

A.



B.

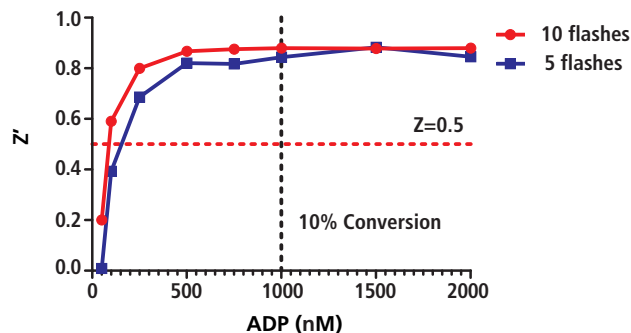


Figure 3. A). Z' values observed in a standard curve mimic conversion of 10 μM ATP to ADP. B). Zoomed view of the 0-2 μM ADP section of the standard curve shows the Z' validation minimal qualification data (red dotted line) and 10% ATP conversion validation point (black dotted line). Plate reader set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

Assay Performance at 10% Conversion 10 μM ATP

| Flashes                         | 1    | 5    | 10   | 20   | 50   |
|---------------------------------|------|------|------|------|------|
| Read Time (minutes)             | 1:10 | 1:18 | 1:25 | 1:39 | 2:19 |
| % CV at 10% ATP Conversion      | 7.01 | 3.34 | 2.73 | 1.92 | 2.23 |
| Z'-Factor at 10% ATP Conversion | 0.69 | 0.85 | 0.88 | 0.91 | 0.91 |

Table 2. Assay Performance with Various Instrument Settings

## Conclusions

This application note demonstrates the validation of the BMG LABTECH PHERAstar Plus and PHERAstar FS instruments for use with the Transcreeper<sup>®</sup> FI Assays. By utilizing the optimized instrument settings suggested within this Application Note, Z' values > 0.7 at 10% conversion 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.

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.

