

Methyltransferase, acetyltransferase, kinases, and GTPases can all be measured with Transcreener® assays

Meera Kumar¹, Tom Zielinski¹, Andrew Koop¹, Franka Ganske², and E. J. Dell²
¹BellBrook Labs; ²BMG LABTECH

- The universal transcreener® assays measure the byproducts of common biological enzymatic reactions
- For epigenetic studies, AMP/GMP can be monitored for methyltransferase or acetyltransferase activity
- ADP monitored for kinases, ATPases, or helicases; GDP for G-proteins; and UDP for glycosyltransferases

Introduction

Transcreener® assays from BellBrook Labs offer generic, universal HTS assays for nucleotide detection like ADP, UDP, GDP and AMP. These assays are available in three detection modes - Fluorescent Polarization (FP); Fluorescent Intensity (FI) and Time Resolved Fluorescence Resonance Energy Transfer (TR-FRET).

The PHERAstar series of microplate readers are perfect for all three detection formats. For FP and TR-FRET it uses a unique Simultaneous Dual Emission detection system that measures both emission wavelengths in one read of the microplate. Simultaneous Dual Emission detection not only reduces plate read times by half, it corrects for any signal variations due to differences in well volumes, concentrations, or fluctuations in excitation energy.

In this application note we show the performance of PHERAstar FS for detection of different nucleotides in the Transcreener® FP assay. The instrument can achieve a Z' >0.8 and a Δ mP greater than 100 at 10% conversion for all the Transcreener® assays tested with 10 μ M respective standard curve.

Assay Principle

Transcreener® Fluorescent Polarization Assay

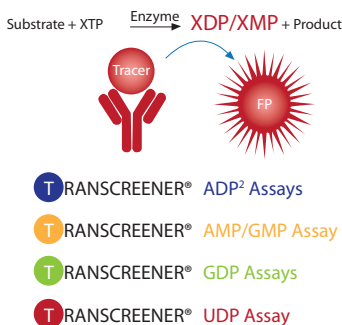


Fig. 1: Transcreener® Fluorescent polarization assays are competitive immunoassays that use a far red tracer bound to a selective antibody for ADP, UDP, GDP and AMP. The tracer is displaced by the above mentioned nucleotide products generated during enzyme reactions. The displaced tracer freely rotates leading to a decrease in fluorescence polarization, relative to the bound tracer.

Materials & Methods

- Transcreener XMP Detection Reagents
- Black 384-well plates from Corning
- PHERAstar® FS microplate reader (BMG LABTECH)

Representative for all XDP Transcreener assays the ADP² assay is explained in more detail. As the ratio of ADP:ATP increases, the proportion of bound tracer versus free tracer decreases, resulting in an overall decrease in mP values. A 15-point standard curve was prepared as described below:

1. Dispense 10 μ L of each ATP/ADP combination across an entire row of the black 384-well plate.
2. Add 10 μ L of ADP Detection Mix to those rows.
3. Dispense 10 μ L of the 10 μ M ATP/0 μ M ADP combination into row P.
4. Dispense 10 μ L of free tracer into wells P1 to P12.
5. Dispense 10 μ L of buffer blank into well P13-P24.

For the FP measurement the Transcreener specific FP modul should be installed. The excitation wavelength is at 590 nm and both emission wavelengths are at 675 nm. A focus and Gain Adjustment should be done, while setting the target mP value to 20 mP using a well that contains only the free tracer. A detailed description of the procedure is given on BellBrookLabs application note: http://www.bellbrooklabs.com/PDFs/ApNote_PHERAstar_FP.pdf

Results & Discussion

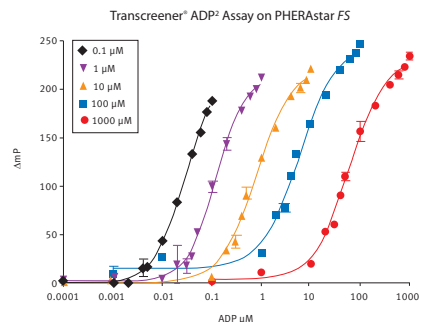


Fig. 2: Graph shows standard curves run on PHERAstar FS using 1000 μ M, 100 μ M, 10 μ M, 1 μ M and 0.1 μ M initial concentrations of ATP using the Transcreener® ADP² assay.



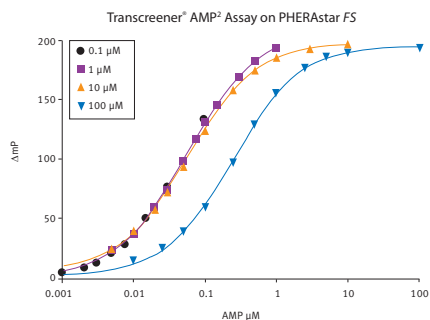


Fig. 3: Graph shows standard curves run on PHERAstar FS using 100 μM, 10 μM, 1 μM and 0.1 μM initial concentrations of ATP using the Transcreener® AMP² assay.

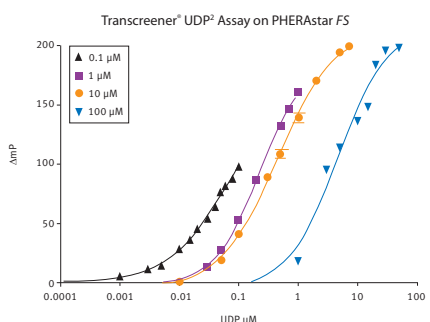


Fig. 4: Graph shows standard curves run on PHERAstar FS using 100 μM, 10 μM, 1 μM and 0.1 μM initial concentrations of UDP Glucuronic Acid using the Transcreener® UDP² assay.

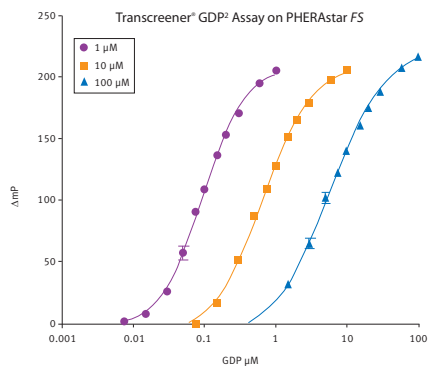


Fig. 5: Graph shows standard curves run on PHERAstar FS using 100 μM, 10 μM and 1 μM initial concentrations of GTP using the Transcreener® GDP assay.

Table 1: Table shows Z' values for 10 μM ATP/ADP; 10 μM ATP/AMP; 10 μM UDP Glucuronic Acid/UDP and 10 μM GTP/GDP generated in PHERAstar FS. The highlighted values show the Z' at 10% conversion demonstrating a superior assay performance in BMG LABTECH's PHERAstar FS.

% Conv	ADP ² Assay	AMP ² Assay	UDP ² Assay	GDP Assay
20	0.88	0.92	0.86	0.86
10	0.85	0.93	0.91	0.86
5	0.82	0.93	0.85	0.81
4	0.77			
3	0.74	0.86	0.83	
2	0.58	0.55	0.57	0.67

Table 2: Read Times for the PHERAstar series of microplate readers in Fluorescence Polarization Mode (10 flashes, different microplate formats).

Fluorescent Polarization Read Times (10 flashes)		
	PHERAstar FS	Reader with No Simultaneous Dual Emission Detection
384-well	0:38	≥ 1:16 (0:38 x 2)
1536-well	1:52	≥ 3:44 (1:52 x 2)
3456-well	9:47	≥ 19:34 (9:47 x 2)

Conclusion

Bellbrook Labs offers four Transcreener assays for direct detection of ADP, AMP/GMP, UDP and GDP.

Excellent Z' values and large signals are shown for low percent conversion with all the four assays using BMG LABTECH's PHERAstar FS.

This application note demonstrates the validation of the BMG LABTECH PHERAstar FS instrument for use with the Transcreener FP Assays. By utilizing the optimized instrument settings suggested within this Application Note, Z' values > 0.7 and ΔmP values > 120 at 10% conversion are achievable.



PHERAstar® FSX

*The PHERAstar FSX is the newest PHERAstar reader.



CLARIOstar®



Omega Series
only in 96-w-format