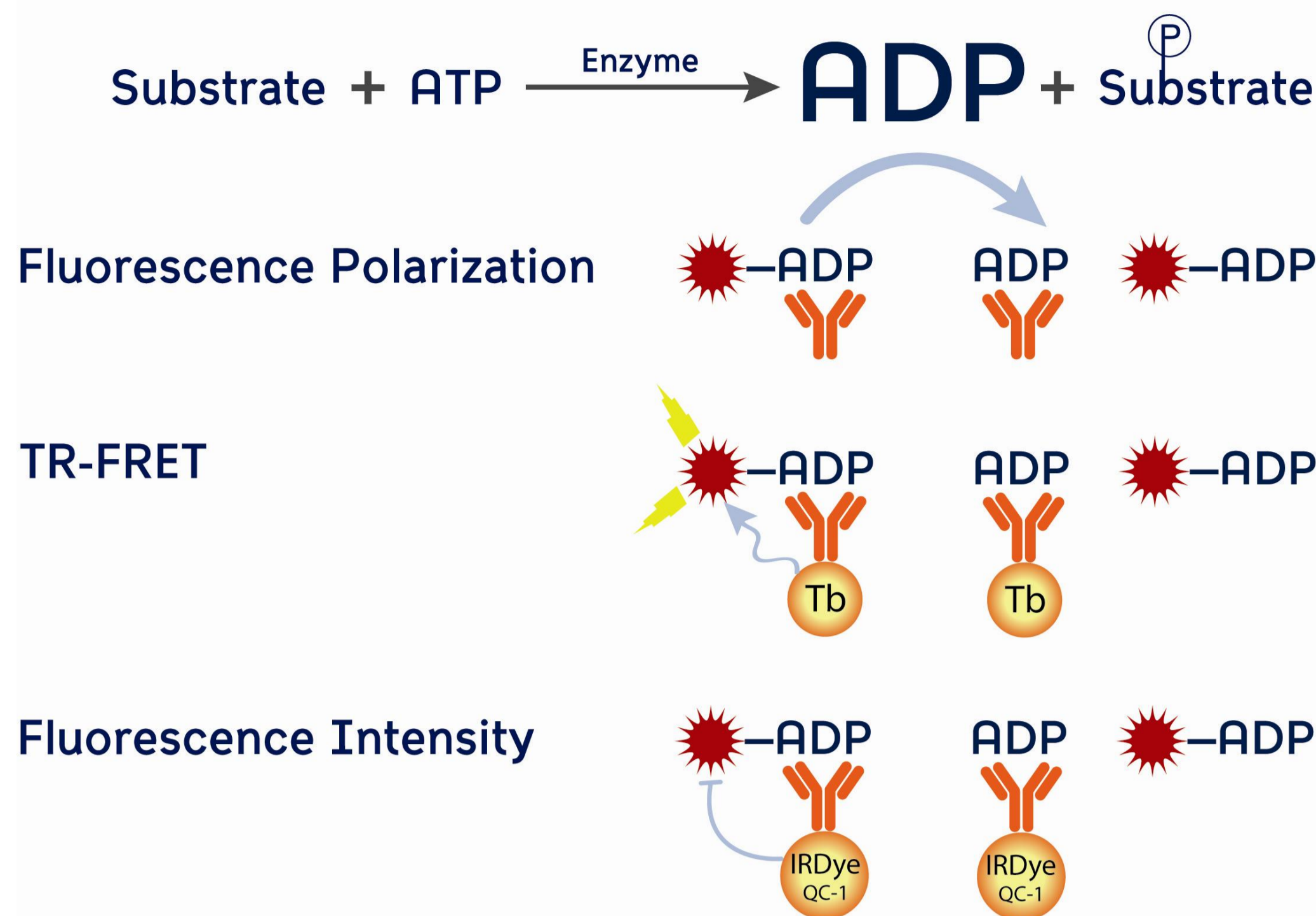


Abstract

The Transreener ADP² Assay is a single addition, homogenous assay method that directly measures ADP formation in kinase or ATPase reactions. The first generation Transreener ADP assay, introduced in 2006 as a far red fluorescence polarization (FP) assay, was recently improved by incorporating a new antibody (ADP²) that enables a 20-fold increase in sensitivity. Together these Transreener ADP assays have been validated for high throughput screening with diverse targets including protein, lipid and carbohydrate kinases, carboxylases, and chaperonin ATPases. To extend the instrument compatibility for the ADP² assay, we have now formatted it for two additional red fluorescence detection modes, time resolved Forster resonance energy transfer (TR-FRET) and fluorescence intensity (FI). Here we showcase the sensitivity and robustness of the three Transreener assays and compare them with other nucleotide detection assay methods. With each of the Transreener assay formats, Z' values greater than 0.7 were observed at less than 10% ATP conversion over the full range of initial ATP concentrations (0.1 μM to 1 mM ATP). Similar EC₅₀ values and inhibitor profiles for PKA were also observed with the three Transreener detection modes demonstrating the equivalence of the assays for inhibitor profiling.

A TRANSCREENER[®] ADP² Assays



B

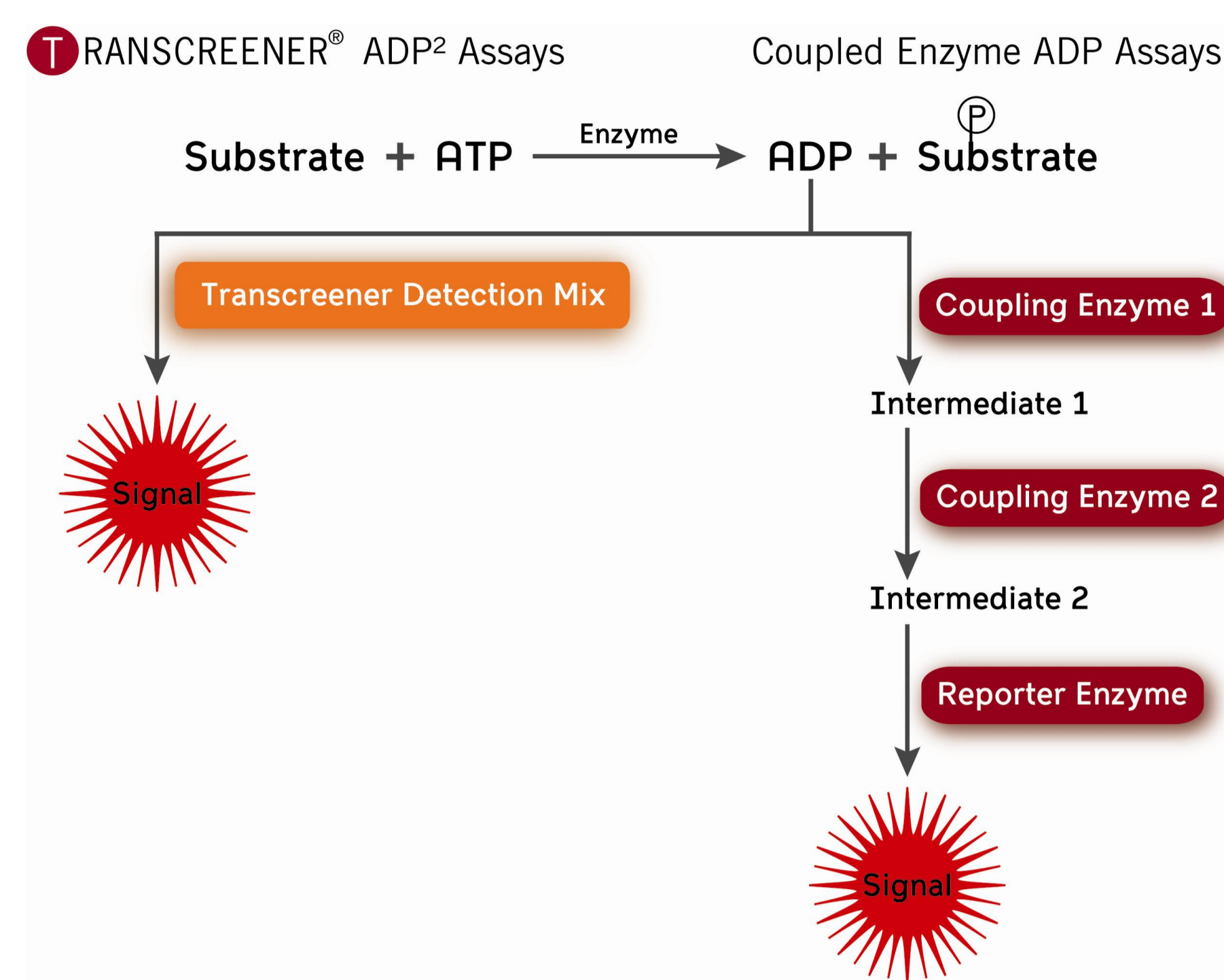


Figure 1. The Transreener ADP² Assays are competitive immunoassays that directly measure enzymatically generated ADP. A) Transreener is the simplest, most direct ADP detection method available: binding of ADP to antibody displaces a tracer, causing a change in its fluorescence properties. The assay has been formatted for three detection modes: Fluorescence polarization, TR-FRET and Fluorescence intensity. B) Other ADP or ATP detection methods use coupling enzymes to convert ADP to a product that can be detected as well as reporter enzymes, to generate a signal. Each coupling and reporter enzyme is a potential target for the compounds being screened, which increases the risk of false positives or of missing a hit. Some coupled methods require extra liquid addition and incubation steps, which complicates assay automation.

Sensitive Initial Velocity Detection over a Broad Range of ATP

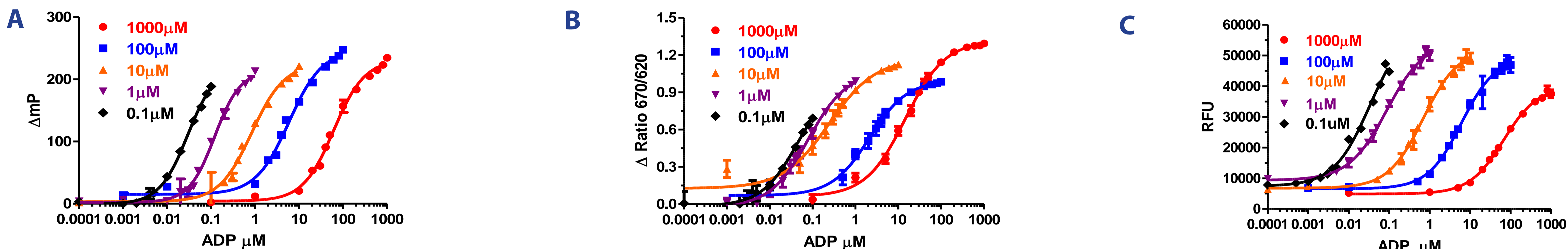


Figure 2. Standard curves mimicking the conversion of ATP to ADP. Figure 2A, 2B and 2C show ADP/ATP standard curves for a broad range of initial ATP concentrations (0.1 μM to 1000 μM ATP) using the FP, TR-FRET and FI assay respectively. ADP was increased as total ATP + ADP was held constant. Antibody was adjusted based on a linear relationship between EC₈₅ and initial ATP. All assays were performed in 384-well plates (n=24) and read on the Tecan Safire2TM (FP), BMG Labtech PHERAstar Plus (TR-FRET), or the Perkin Elmer EnVision (FI) multimode plate readers.

Greater Sensitivity than Luciferase Based ATP or ADP Detection

	1 μM ATP/ADP standard curve		10 μM ATP/ADP standard curve		100 μM ATP/ADP standard curve	
	Z' at 10% ATP Conv	*LLD (μM)	Z' at 10% ATP Conv	*LLD (μM)	Z' at 10% ATP Conv	*LLD (μM)
Transreener ADP ² FP	0.86	0.020 ± 0.07	0.85	0.01 ± 0.12	0.89	1 ± 0.3
Transreener ADP ² TR-FRET	0.71	0.03 ± 0.06	0.72	0.1 ± 0.09	0.72	1 ± 0.25
Transreener ADP ² FI	0.92	0.030 ± 0.01	0.88	0.05 ± 0.04	0.92	0.5 ± 0.37
Luc- ADP Detection Assay	ND	0.4 ± 0.87	0.3	0.5 ± 0.32	0.62	5 ± 0.67
Luc-ATP Depletion Assay	ND	0.25 ± 0.4	ND	1.5 ± 0.3	0.52	7 ± 0.56

Table 1. Z' values and Lower Limits of Detection for ADP in the Presence of Excess ATP

Each of the Transreener ADP² Assays show robust Z' at low ADP concentrations (10% ATP conversion) with a 20-fold better lower limit of detection to luciferase ADP detection methods. The Transreener ADP² Assays consistently show Z' > 0.7 at 10% ATP conversion for 1 μM, 10 μM and 100 μM ATP/ADP standard curves (shown highlighted). A Z' > 0.5 is considered a robust assay. *LLD=Lower Limit of Detection, defined as the concentration of ADP that generates Z' > 0.

More Sensitive ADP Detection Minimizes Enzyme Usage

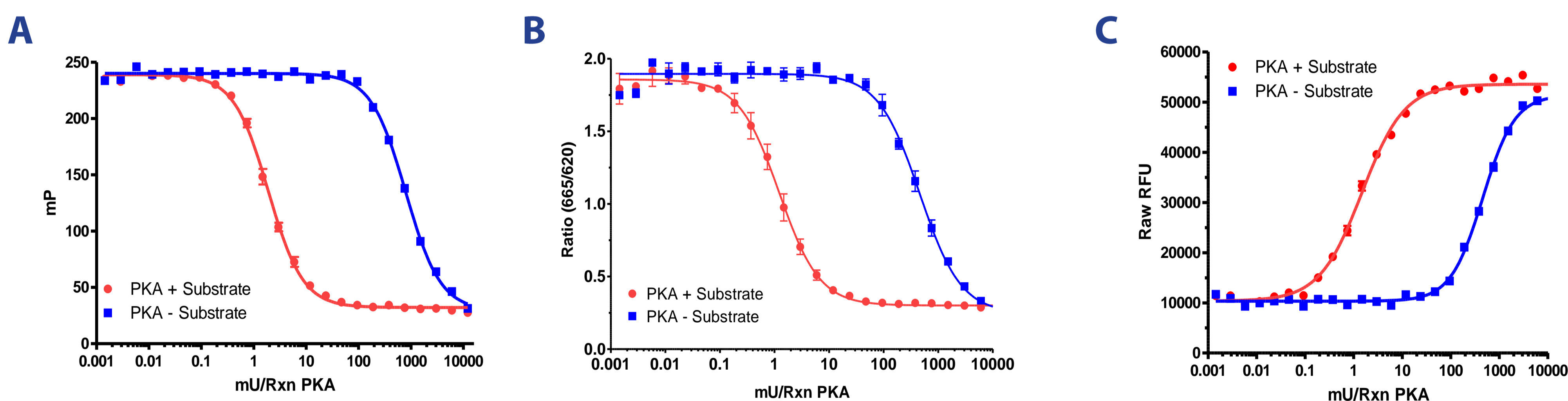


Figure 3. PKA Enzyme Titrations. Excellent signal is demonstrated at low [enzyme] for each of the assay readouts. PKA generates 400-times more ADP in the presence of substrate (productive phosphorylation) than in its absence (intrinsic ATP hydrolysis). A) FP EC₅₀ values = 1.9 and 793 mU/RXN with and without substrate. B) TR-FRET EC₅₀ values = 1.2 and 458 mU/RXN with and without substrate. C) FI EC₅₀ values = 1.5 and 454 mU/RXN with and without substrate. Detection Mixture (10 μL) was added and reactions equilibrated for one hour before reading the final 20 μL assay volume.

Accurate Inhibitor Pharmacology with All Three Readouts

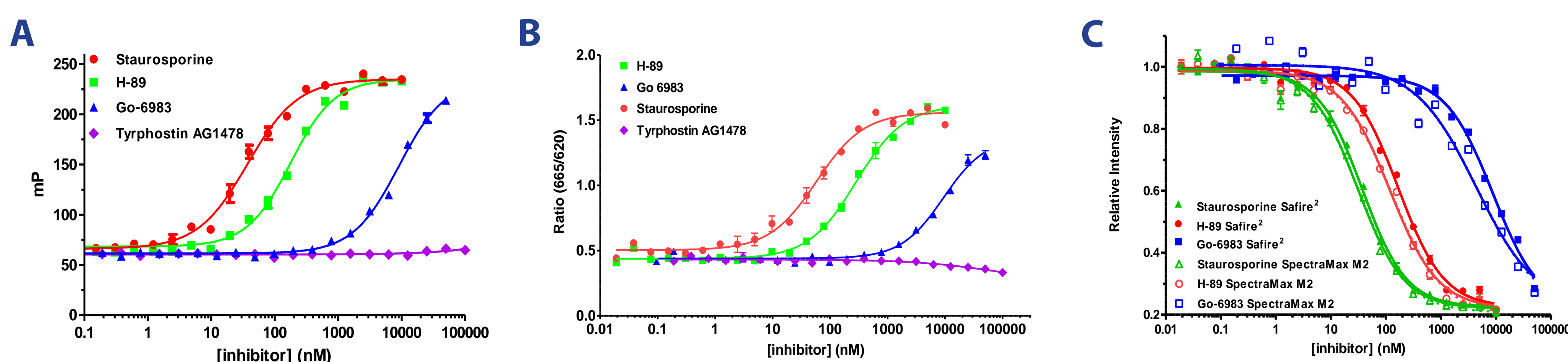


Figure 4. Inhibition curves with known inhibitors and PKA. Dose dependency curves were generated for Stauroporine, H-89, and Go-6983 (known PKA inhibitors) by serially titrating into the PKA reaction conditions. A) For FP, IC₅₀ values for stauroporine, H-89 and Go-6983 were 37 nM, 179 nM and 9.0 μM respectively. B) For TR-FRET, IC₅₀ values for stauroporine, H-89 and Go-6983 were 58 nM, 283 nM and 9.3 μM respectively. C) Similar IC₅₀ values are shown using both instruments (Safire2/SpectraMax M2) for stauroporine, H-89 and Go-6983 were 37/32 nM, 165/124 nM and 9.0/4.7 μM respectively.

Conclusions

- 1) All three ADP² Assays are capable of robust, initial velocity ADP detection (≥ 9:1 ATP:ADP ratio) with Z' values greater than 0.7 using total (ATP + ADP) concentrations ranging from 0.1 μM to 1,000 μM.
- 2) The lower limits of detection for ADP in the presence of 1 μM total (ATP + ADP) are 20 nM, 30 nM and 30 nM for the FP, TR-FRET and FI assays, respectively.
- 3) Compared to luciferase-based ATP and ADP detection, the Transreener ADP² Assays exhibited significantly higher Z' values for initial velocity detection and greater sensitivity (LLD) at all (ATP + ADP) levels tested.
- 4) Sensitivity of detection of PKA enzyme activity was similar with the three Transreener ADP² Assays: EC₅₀ values of 1.9, 1.2, and 1.5 mU per 10 μL enzyme reaction were observed for the FP, TR-FRET and FI assays respectively.
- 5) Similar potencies were observed for known inhibitors of PKA using all three ADP² assays.

Acknowledgements: Funded in part by NIH SBIR grants R44CA110535 and R44GM073290