

Flexible and targeted inhibitor profiling using the luminescent ADP-Glo™ kinase assay platform

Zegzouti Hicham, Alves Juliano, Hsiao Kevin, Vidugiriene Jolanta and Goueli Said

Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711, USA

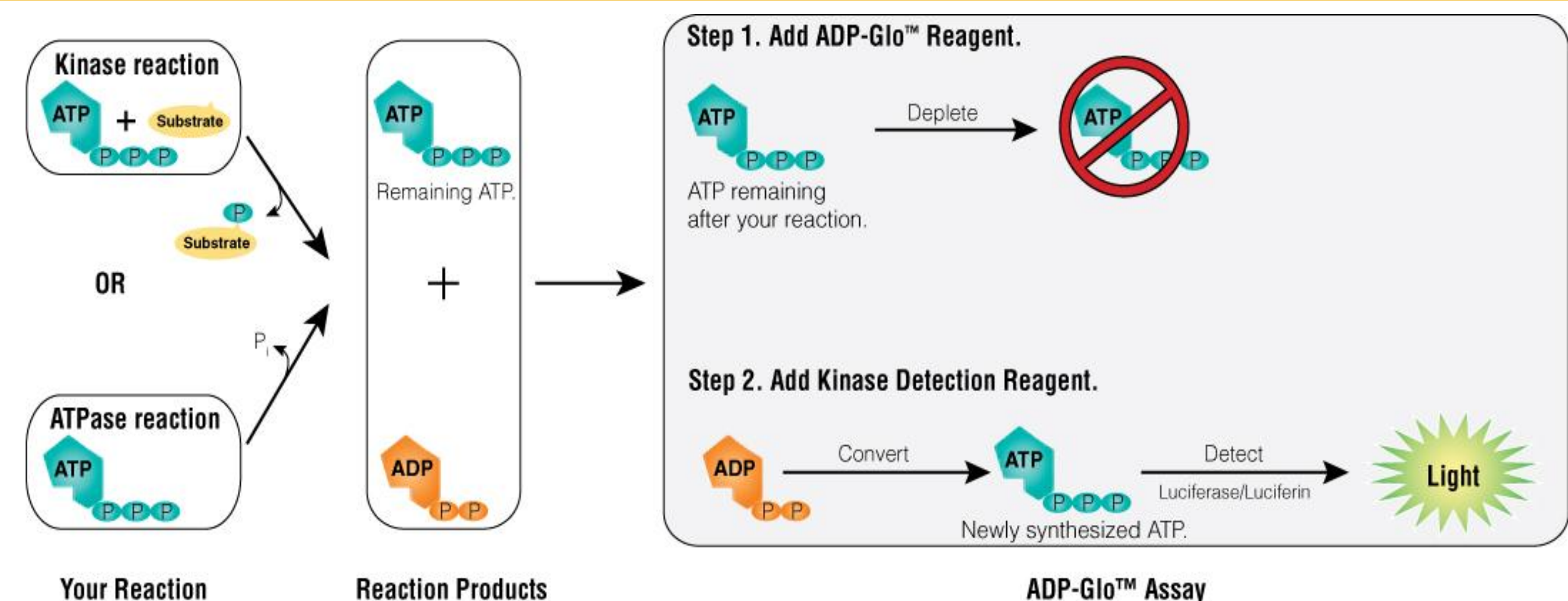


2012

1. Abstract

Profiling kinase inhibitors is a critical step during drug development. Knowing the inhibitory profile of a compound across a broad collection of kinases can be important for better understanding its biological activity, obviating any off-target activities, and in some cases identifying new targets may lead to novel therapeutic applications. Unlike an HTS phase, where a target-specific kinase assay is used to identify hit compounds, during profiling a robust universal assay is needed to assess the selectivity and potency of the inhibitor on multiple kinases from different classes. Often, these kinases use substrates with different chemical structures and target-specific screening assays are not suitable or can be cost prohibitive. We previously introduced the luminescent ADP-Glo™ kinase assay that measures kinase activity by quantifying the amount of ADP produced during the enzymatic reaction. This technology addressed all the needs of kinase screening, mode of action (MOA) studies and profiling using one assay format during the drug discovery process. Here, we show that we can achieve the sensitivity and robustness required for profiling the different kinase families covering the human kinome with one platform. The ADP-Glo™ assay was previously validated with more than hundred kinases and is now optimized for use with a large panel of complete Kinase Enzyme Systems (KES) that span different families of the human kinome. We also demonstrate the profiling of different kinase inhibitors using newly designed kinase strips. Each multi-well strip contains a group of kinases from the same family. The kinase stock volumes were standardized in a way that, when all the kinases were diluted into the kinase reaction, the kinases generated 5–10% ATP to ADP conversion. The substrate stocks were standardized in similar fashion and are located in a second strip at corresponding positions. We show that using the new kinase profiling strips and an optimized protocol, we could easily generate selectivity profiles using small or large kinase panels, as well as detecting compound promiscuity towards members of a single kinase subfamily or different subfamilies of the kinome. The fact that ADP-Glo™ platform offers so many positive attributes makes it an ideal assay not only for primary and secondary screening but also for profiling compounds in a cost-effective manner using one single platform.

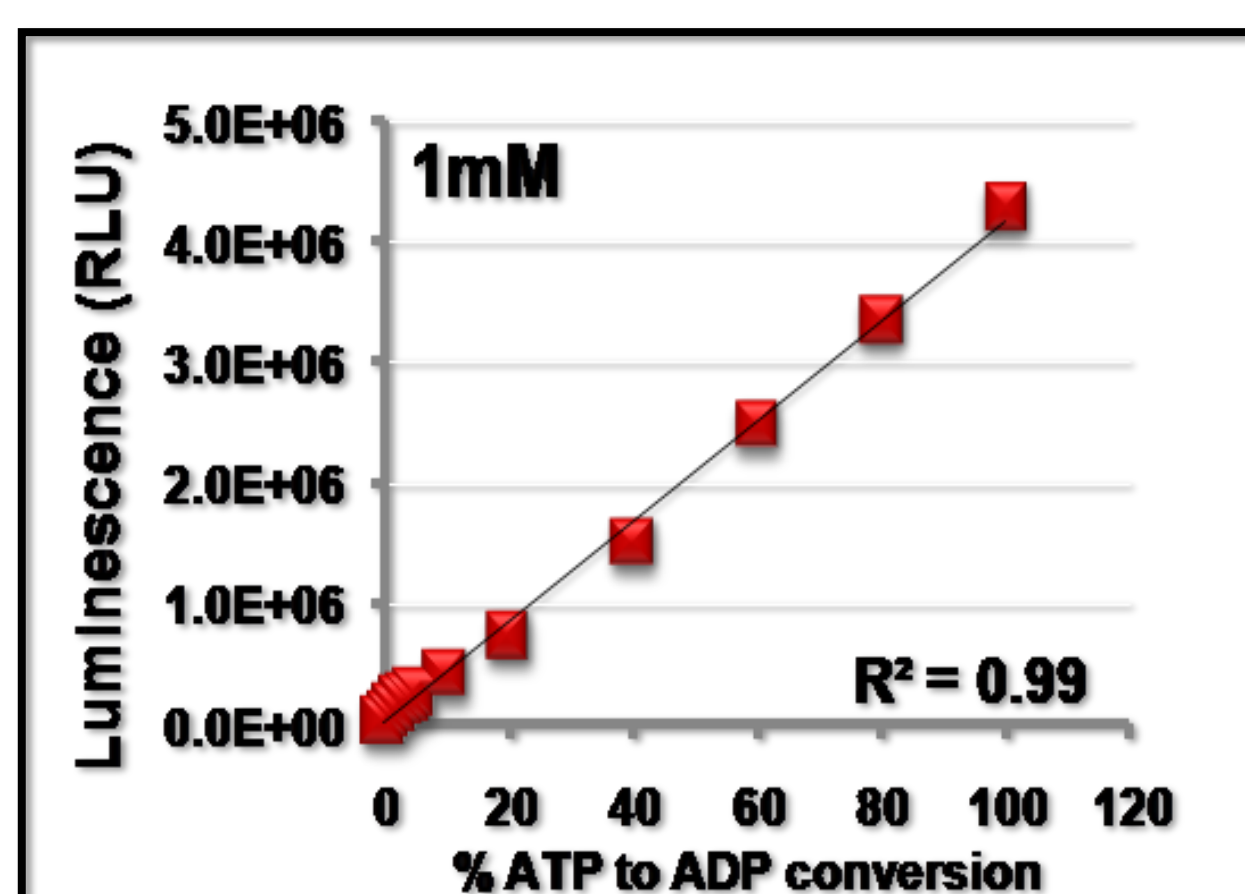
2. ADP-Glo™ is a positive detection assay for product formation



- Step 1: Depletion of unconsumed ATP after the kinase reaction.
- Step 2: ADP is converted into ATP that is detected via a luciferase/luciferin reaction.
- Luminescent signal is proportional to ADP produced and the kinase activity.

3. ADP-Glo™ can be used with a broad range of ATP concentrations

ADP detection up to 1mM ATP in the reaction



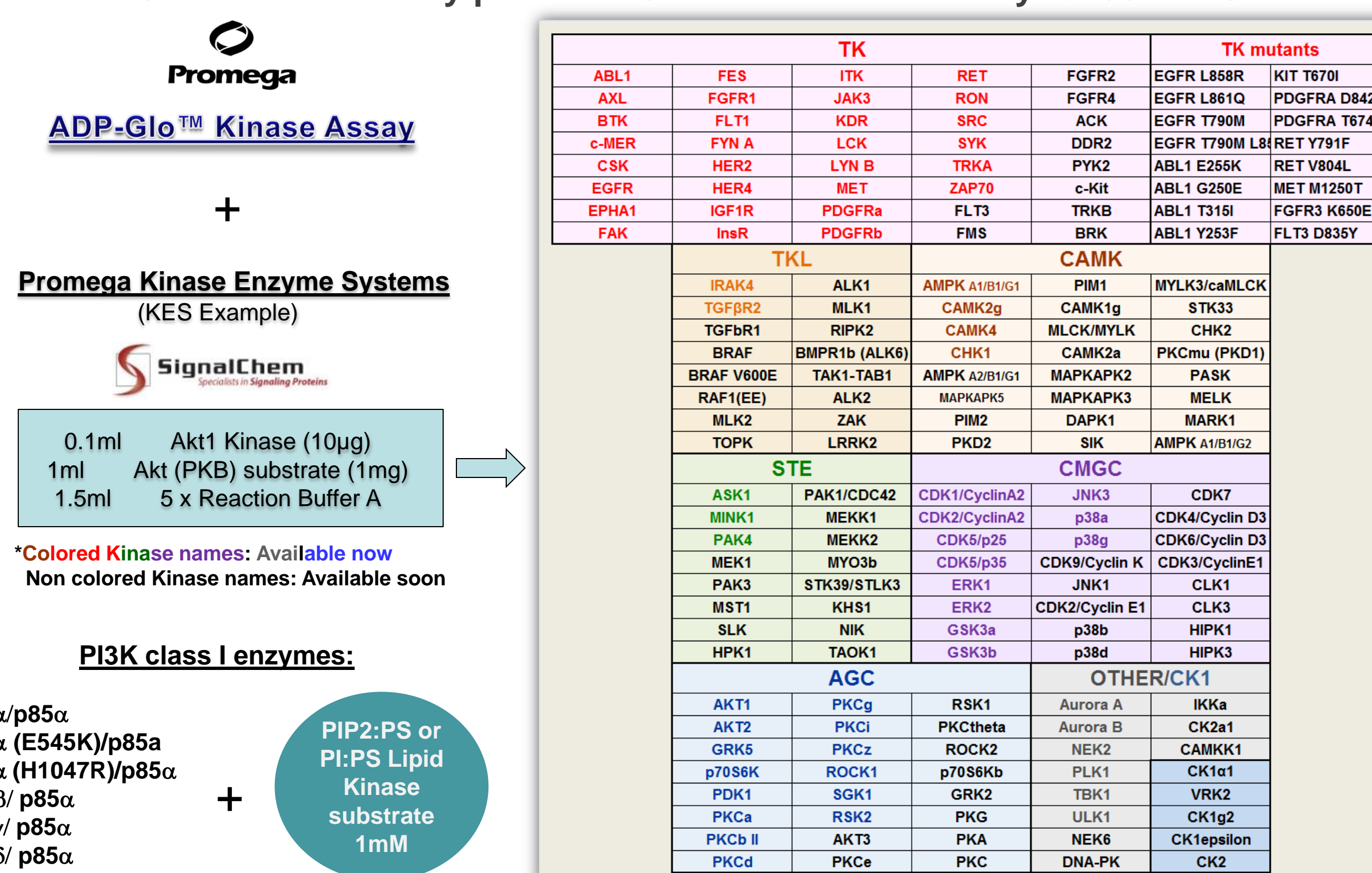
More sensitivity with Promega new ultrapure ATP compared to other ATP sources

	ATP-ADP ranges								
	1mM			100µM			10µM		
% ADP in ATP + ADP mixture	20%	10%	5%	20%	10%	5%	20%	10%	5%
Promega	45	23	13	54	32	18	37	20	10
Competitor S	26	13	8	35	21	11	31	16	8
Competitor T	13	7	4	15	10	5	16	9	5
Competitor G	12	6	4	16	10	6	18	10	6

- Sensitivity: ADP-Glo™ generates wide dynamic range with Z' > 0.7
- Linearity: ADP-Glo™ Kinase assay detects up to 1mM ADP in a linear fashion
- ADP-Glo™ sensitivity is highly improved with the new Ultra Pure ATP

4. ADP-Glo™ is a universal assay validated with large kinase panel spanning the kinome

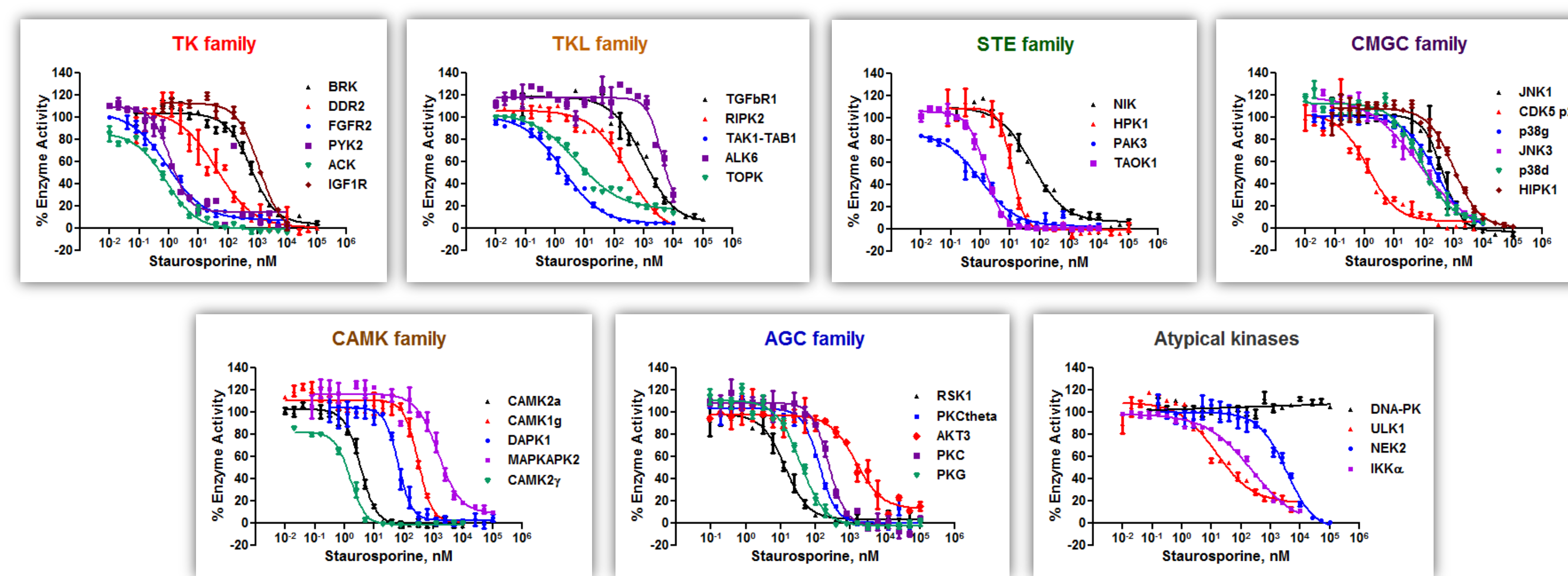
ADP-Glo™ Kinase Assay platform: Universal Kinase Assay & 150+ KES*



Larger kinase panel for easy profiling and selectivity assays

5. Profiling kinase inhibitors with ADP-Glo™ kinase assay

Staurosporine selectivity profiles of small kinase family panels



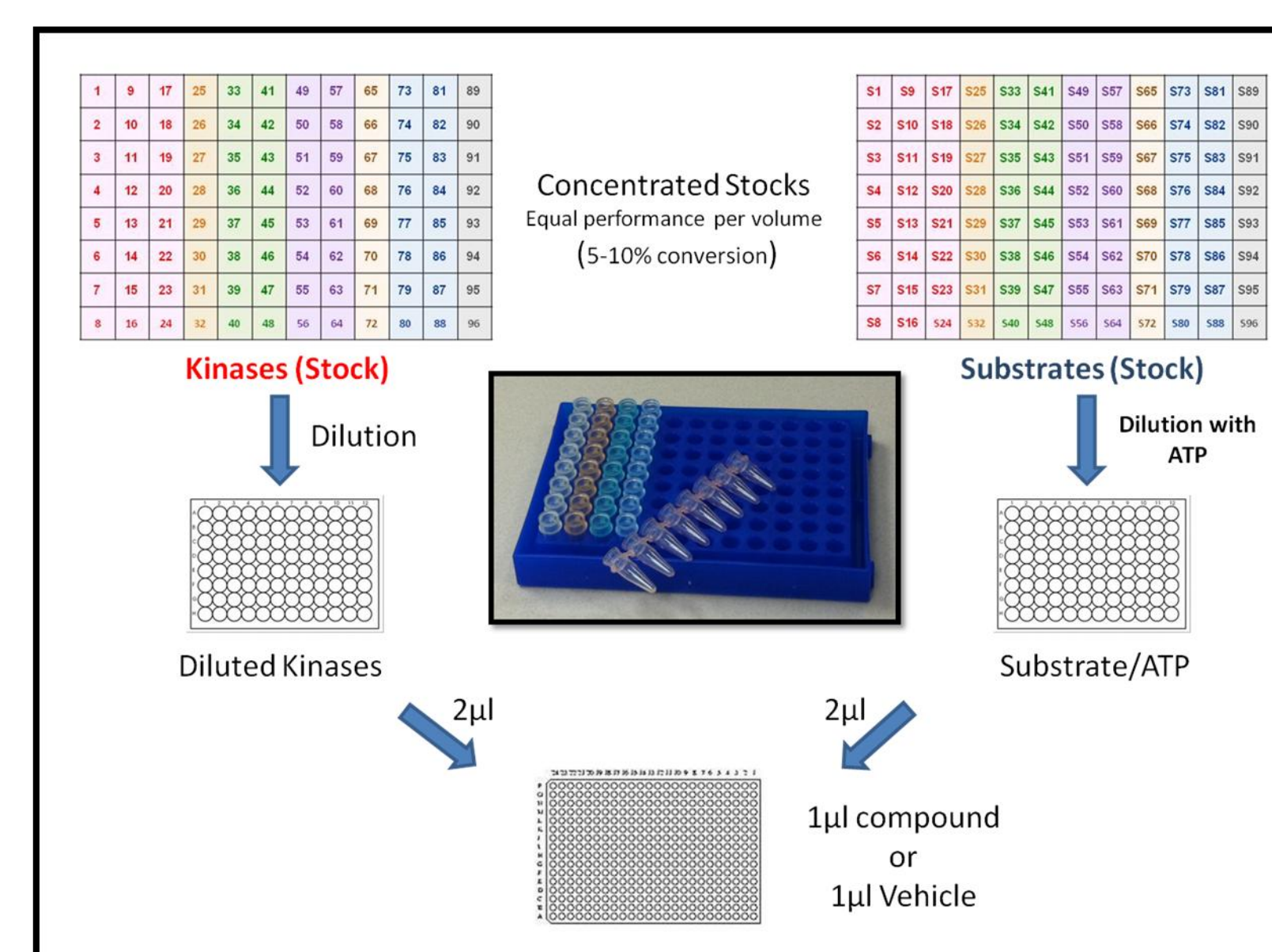
"One Assay for all", ADP-Glo™ is ideal for profiling inhibitors against large or small kinase panels

6. Kinase strips for flexible and targeted inhibitor profiling

Important kinase targets organized in strips by kinase families

Strip	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	
TK	ABL1	FES	RET	FGFR2	EGFR L858R	IGT T870I	TKL	ALK1	AMPK α1/β1	PKA	PKCα	PKCβ	PKCγ	PKCδ	PKCε	PKCζ	PKCη	PKCθ	PKCι	PKCκ	PKCλ
TKL	ABL1	FES	RET	FGFR2	EGFR L858R	IGT T870I	TKL	ALK1	AMPK α1/β1	PKA	PKCα	PKCβ	PKCγ	PKCδ	PKCε	PKCζ	PKCη	PKCθ	PKCι	PKCκ	PKCλ
STE	ASK1	PAK1	CDK1	JNK1	CDK2	CDK3	CDK5	CDK6	CDK7	CDK8	CDK9	CDK10	CDK11	CDK12	CDK13	CDK14	CDK15	CDK16	CDK17	CDK18	CDK19
AGC	AKT1	PKCα	PKCβ	PKCγ	PKCδ	PKCε	PKCζ	PKCη	PKCθ	PKCι	PKCκ	PKCλ	PKCμ	PKCν	PKCξ	PKCο	PKCπ	PKCρ	PKCσ	PKCτ	PKCυ
Other	AMPK α1/β1	PKA	PKCα	PKCβ	PKCγ	PKCδ	PKCε	PKCζ	PKCη	PKCθ	PKCι	PKCκ	PKCλ	PKCμ	PKCν	PKCξ	PKCο	PKCπ	PKCρ	PKCσ	PKCτ

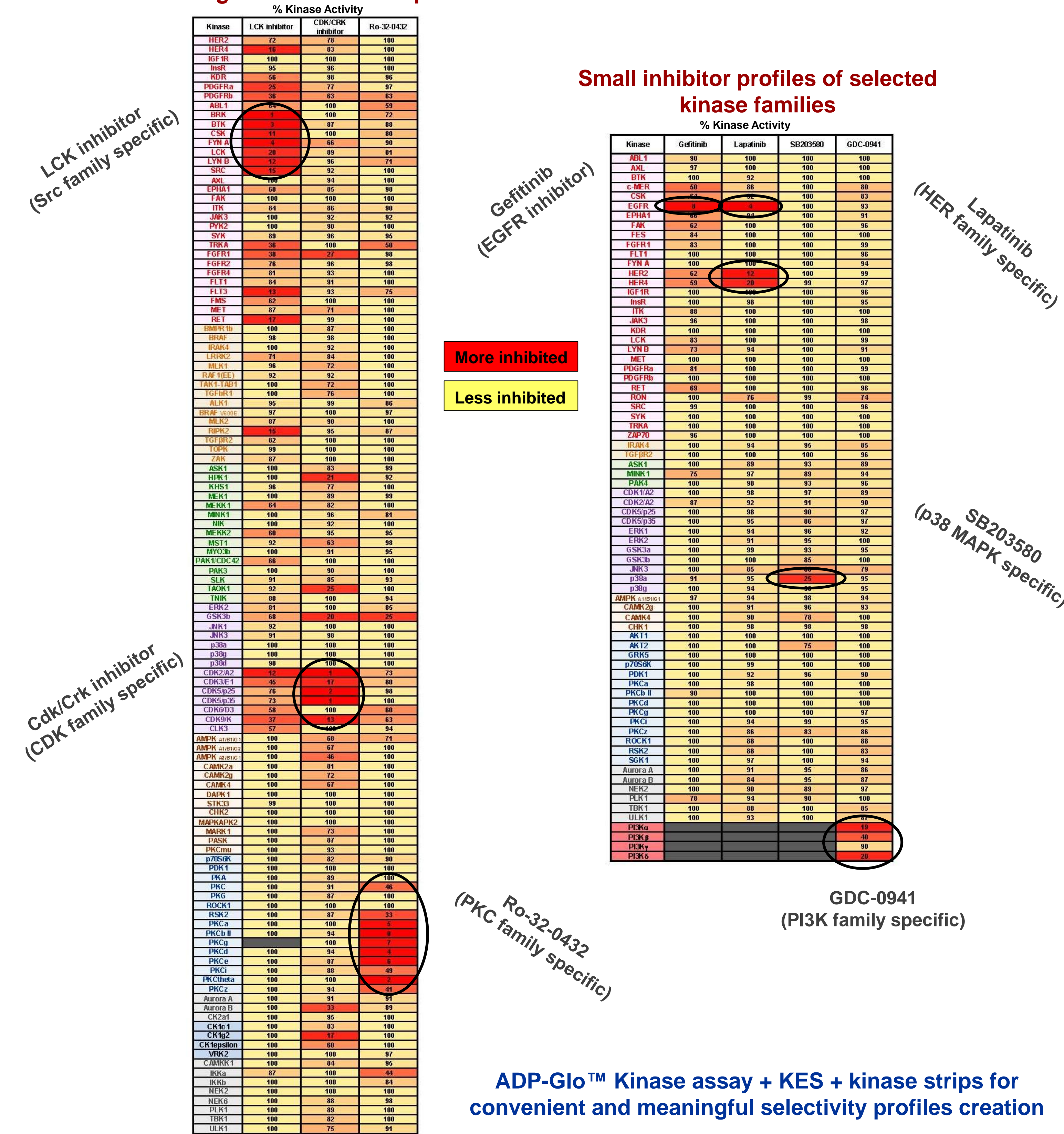
Streamlined profiling protocol



www.promega.com/kinase

7. Determination of inhibitor profiles of different kinase families with ADP-Glo™ platform

Inhibitor selectivity profiles created using all the kinase strips



8. General features of ADP-Glo™ platform

- High sensitivity assay: less than 1% ATP conversion detected with more than 2.5 signal to background ratio with new Ultra Pure ATP
- High dynamic range: High signal to background at low % ATP to ADP conversion allows use of lower amount of enzyme during HTS or profiling
- Homogenous, non radioactive and antibody Free
- Robust assay (Z' higher than 0.7)
- Universal assay: Any ADP-producing enzyme (Kinase or ATPase). Ideal for inhibitor profiling with easy protocol.
- Larger kinase panel: Complete Kinase Enzyme Systems for easy profiling and selectivity assays
- Convenient profiling concept: Strips can be used in a mix and match fashion to profile representatives of a single kinase family or multiple families with the same assay platform

More info: hicham.zegzouti@promega.com