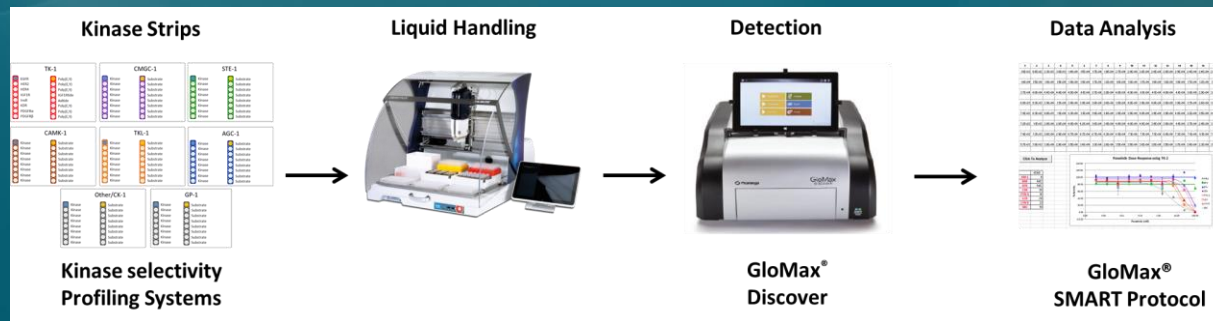
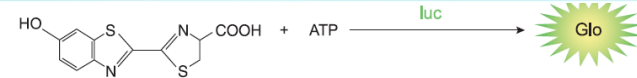


Streamlining Kinase Profiling Workflow by Integrating Bioluminescent Kinase Strips with Multimode Detection System

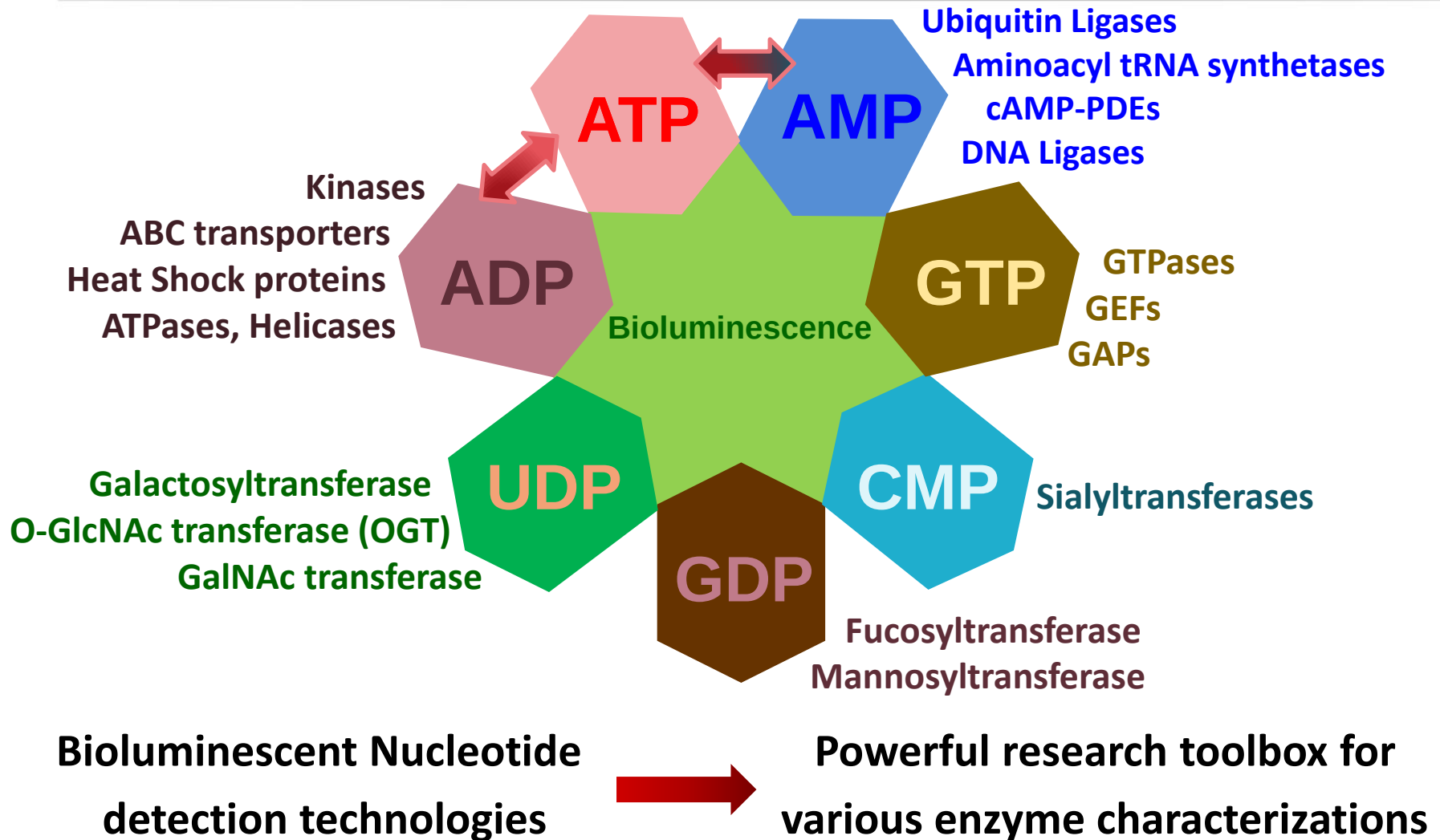
Hicham Zegzouti, Ph.D.

Assay Design Group, Promega, Madison WI

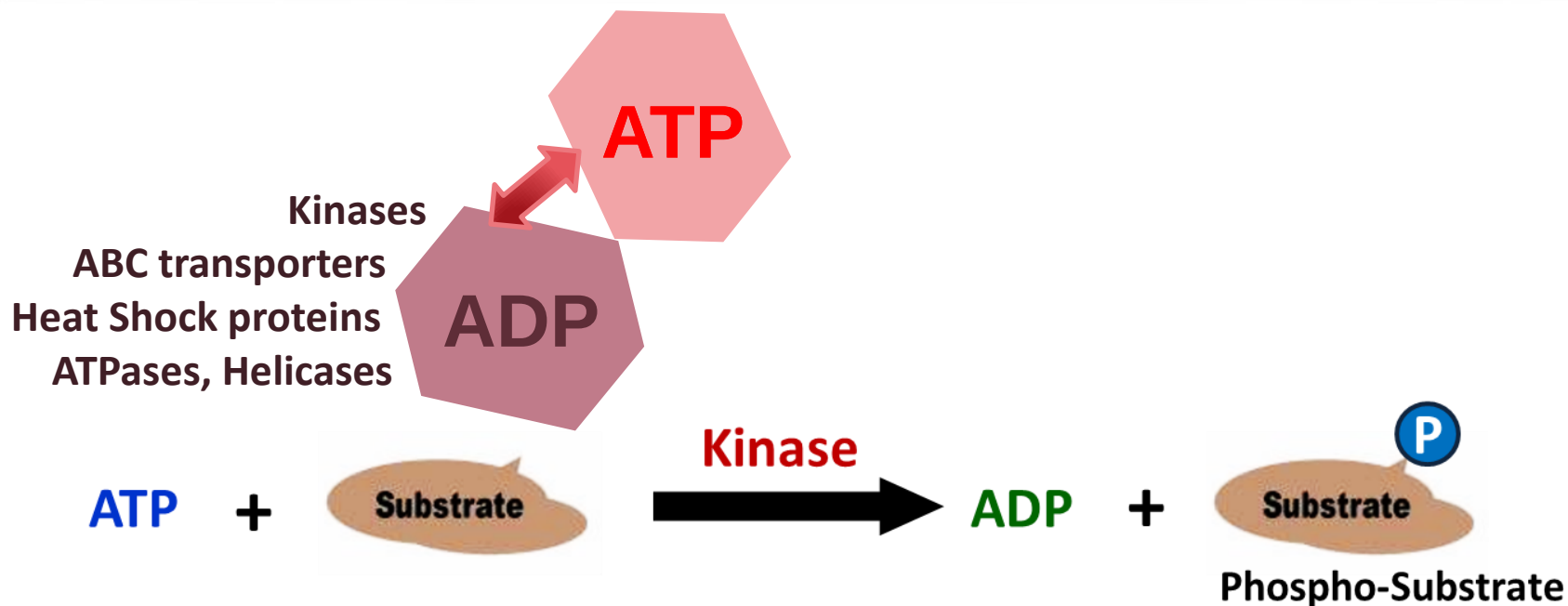




Nucleotides enzyme targets



Nucleotides enzyme targets: **ADP**



Kinases play a critical role in Human biology

Kinases are directly involved in many diseases



Kinases became the largest target in the drug discovery market

Importance of Kinases in Drug Discovery

The Facts:

US Marketed Small Molecule Kinase Inhibitors

Table 1. Targeted Therapeutics in Cancer.^a

Gene	Genetic Alteration	Tumor Type	Therapeutic Agent
Receptor tyrosine kinase			
<i>EGFR</i>	Mutation, amplification	Lung cancer, glioblastoma	Gefitinib, erlotinib
<i>ERBB2</i>	Amplification	Breast cancer	Lapatinib
<i>FGFR1</i>	Translocation	Chronic myeloid leukemia	PKC412, BIBF-1120
<i>FGFR2</i>	Amplification, mutation	Gastric, breast, endometrial cancer	PKC412, BIBF-1120
<i>FGFR3</i>	Translocation, mutation	Multiple myeloma	PKC412, BIBF-1120
<i>PDGFRA</i>	Mutation	Glioblastoma, gastrointestinal stromal tumor	Sunitinib, sorafenib, imatinib
<i>PDGFRB</i>	Translocation	Chronic myelomonocytic leukemia	Sunitinib, sorafenib, imatinib
<i>ALK</i>	Mutation or amplification	Lung cancer, neuroblastoma, anaplastic large-cell lymphoma	Crizotinib
<i>c-MET</i>	Amplification	Gefitinib-resistant non-small-cell lung cancer, gastric cancer	Crizotinib, XL184, SU11274
<i>IGF1R</i>	Activation by insulin-like growth factor II ligand	Colorectal, pancreatic cancer	CP-751,871, AMG479
<i>c-KIT</i>	Mutation	Gastrointestinal stromal tumor	Sunitinib, imatinib
<i>FLT3</i>	Internal tandem duplication	Acute myeloid leukemia	Lestaurtinib, XL999
<i>RET</i>	Mutation, translocation	Thyroid medullary carcinoma	XL184
Non-receptor tyrosine kinase			
<i>ABL</i>	Translocation (BCR-ABL)	Chronic myeloid leukemia	Imatinib
<i>JAK2</i>	Mutation (V617F), translocation	Chronic myeloid leukemia, myeloproliferative disorders	Lestaurtinib, INCB018424
<i>SRC</i>	Overexpression	Non-small-cell lung cancer; ovarian, breast cancer; sarcoma	KX2-391, dasatinib, AZD0530
Serine-threonine-lipid kinase			
<i>BRAF</i>	Mutation (V600E)	Melanoma; colon, thyroid cancer	SB-590885, PLX-4032, RAF265, XL281
Aurora A and B kinases	Overexpression	Breast, colon cancer; leukemia	MK-5108 (VX-689)
Polo-like kinases	Overexpression	Breast, lung, colon cancer; lymphoma	BI2536, GSK461364
<i>MTOR</i>	Increased activation	Renal-cell carcinoma	Temsirolimus (CCI-779), BEZ235
<i>PI3K</i>	PIK3CA mutations	Colorectal, breast, gastric cancer; glioblastoma	BEZ235

Diseases targeted:

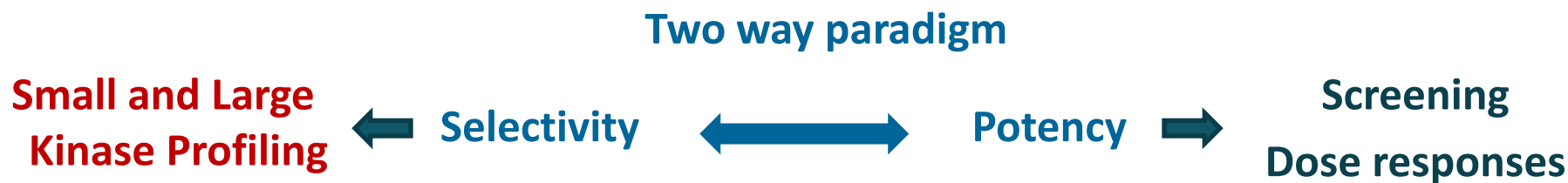
- **Oncology,**
- **inflammation,**
- **Cardiovascular,**
- **Neurological (AD, PD),**
- **Diabetes...**

McDermott, U., Downing, J. R. and Stratton, M. R. (2011) *N Engl J Med* **364**, 340-50.

All Kinase Inhibitors are not equal

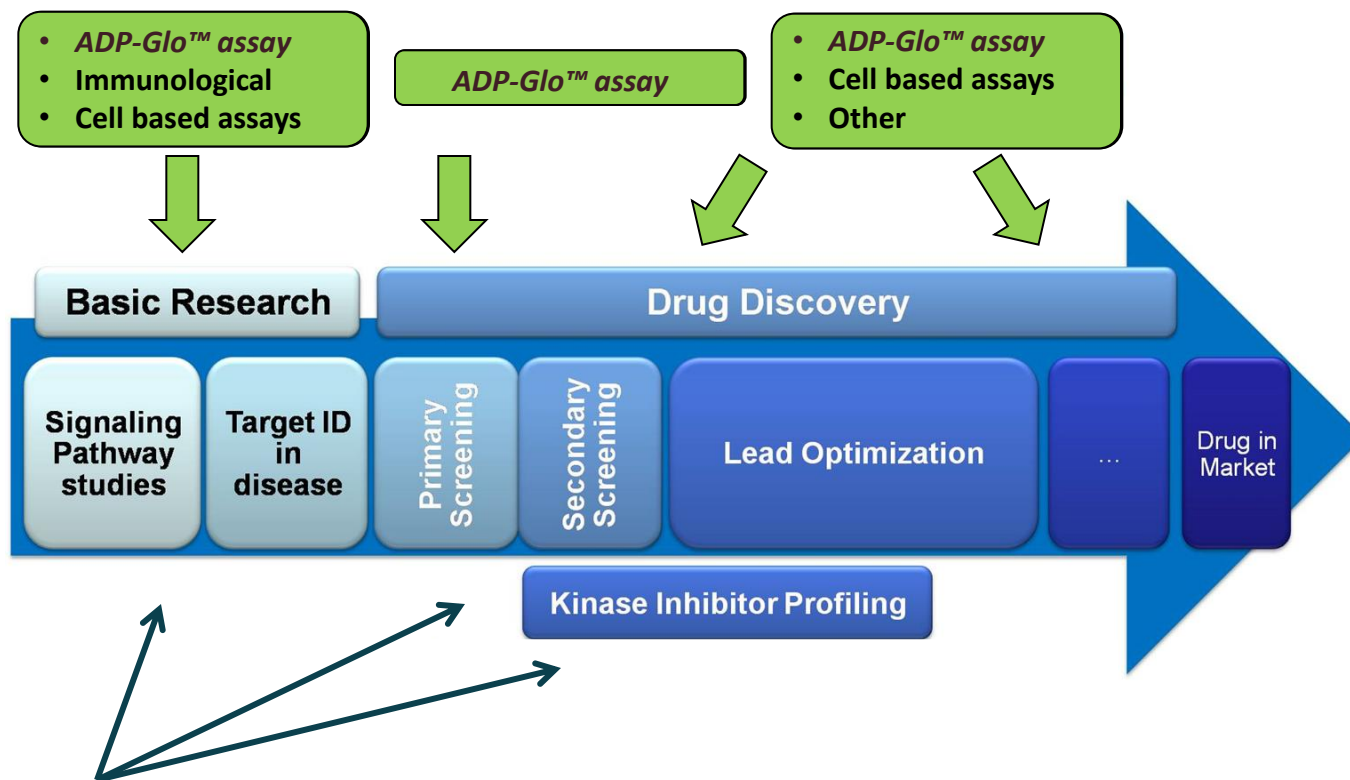
(Different trends of screenings)

- ATP competitive (**low selectivity**)
- ATP non competitive or Allosteric (**more selective**)
- DFG out inhibitors (**inactive form, more selective**)
- DFG in inhibitors (**active form**)
- Type 1 ½ (**Back pocket of ATP site, improve selectivity**)
- DFG in after HTS then optimize to DFG out



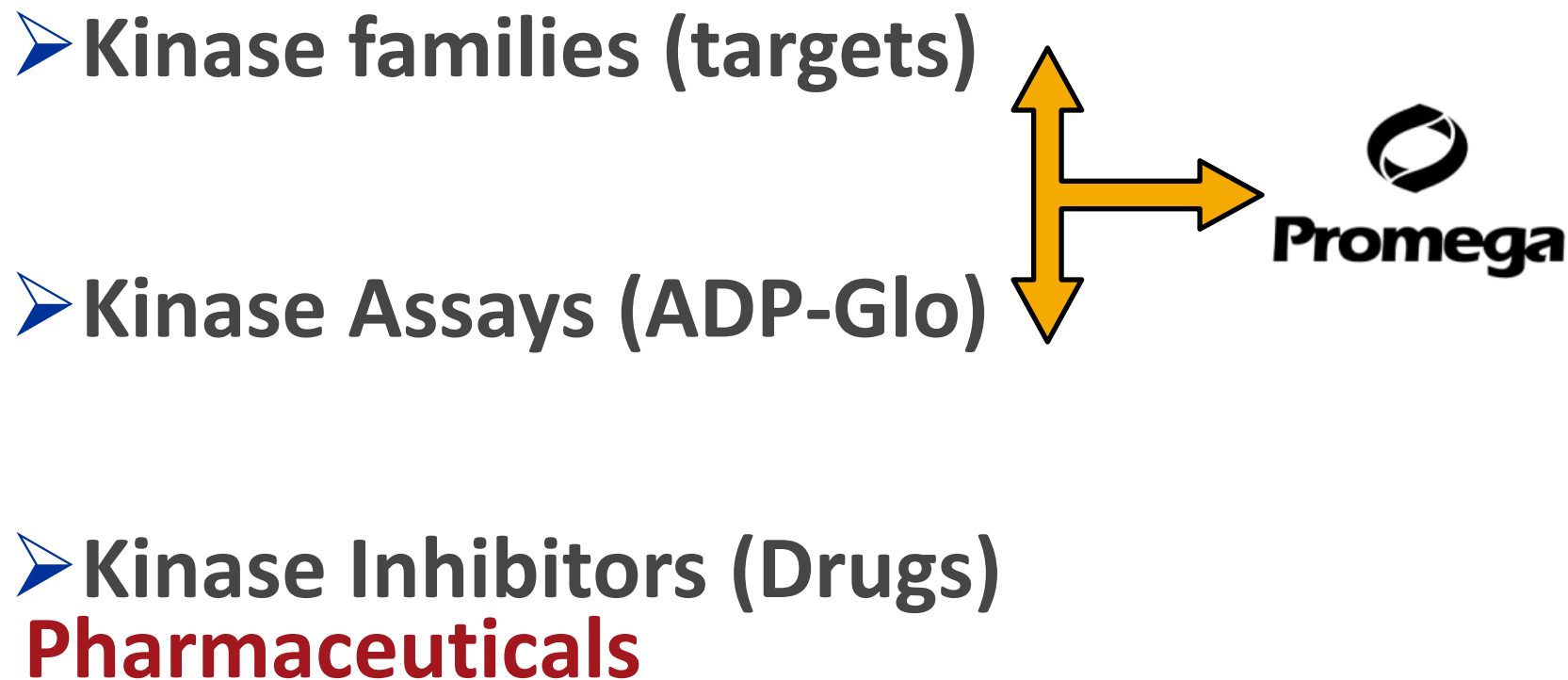
Kinase Studies in Basic Research and Drug Discovery

Kinase Studies and Screening Approaches

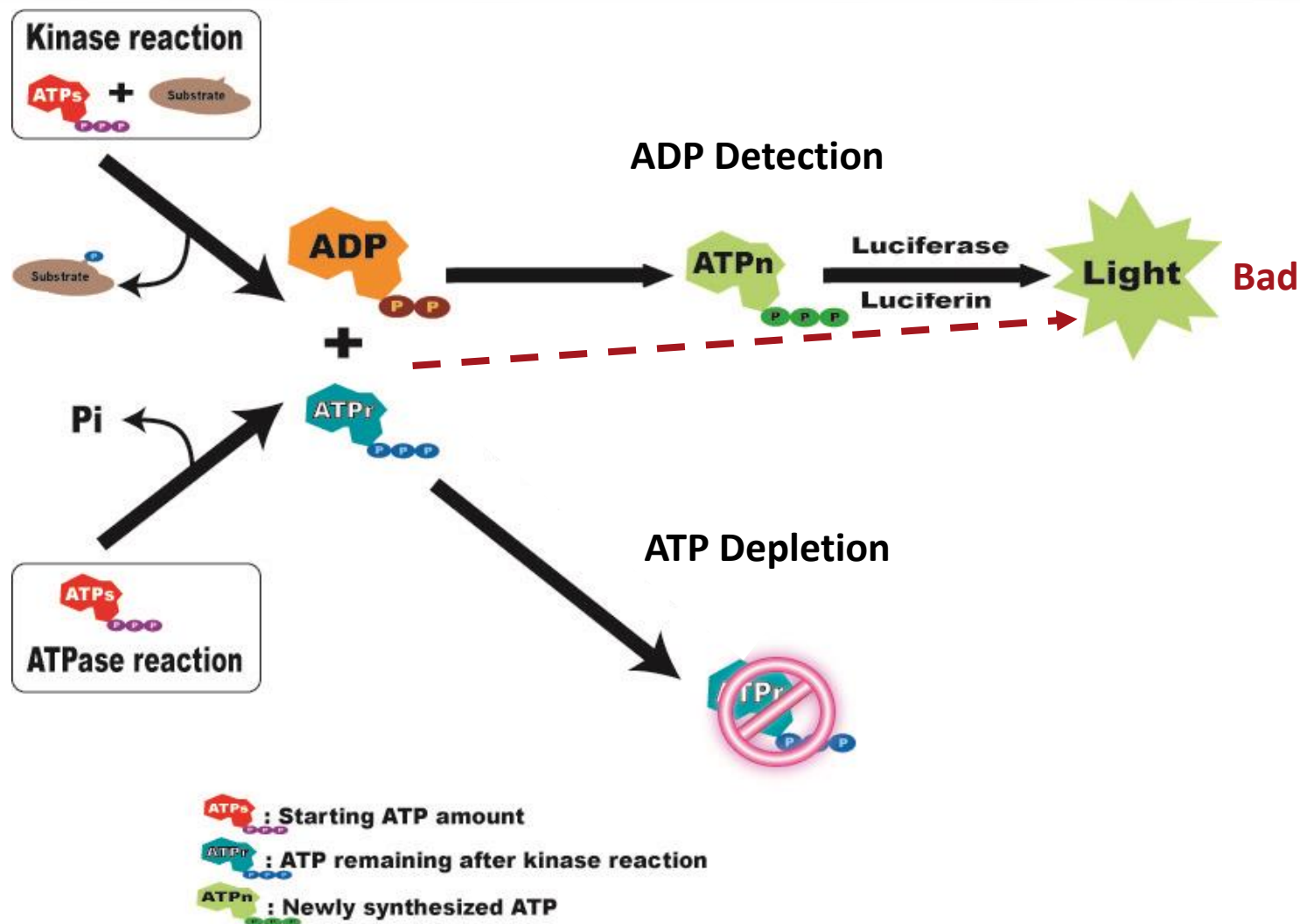


Need for a Universal Kinase Assay that can be applied to all types of Kinase Studies

Kinase Research: parts of the equation



Luminescent ADP detection assay principle



ADP-Glo Assay Format- 1:1:2

384-well plate

5µl kinase reaction

+

5µl ADP-Glo Reagent

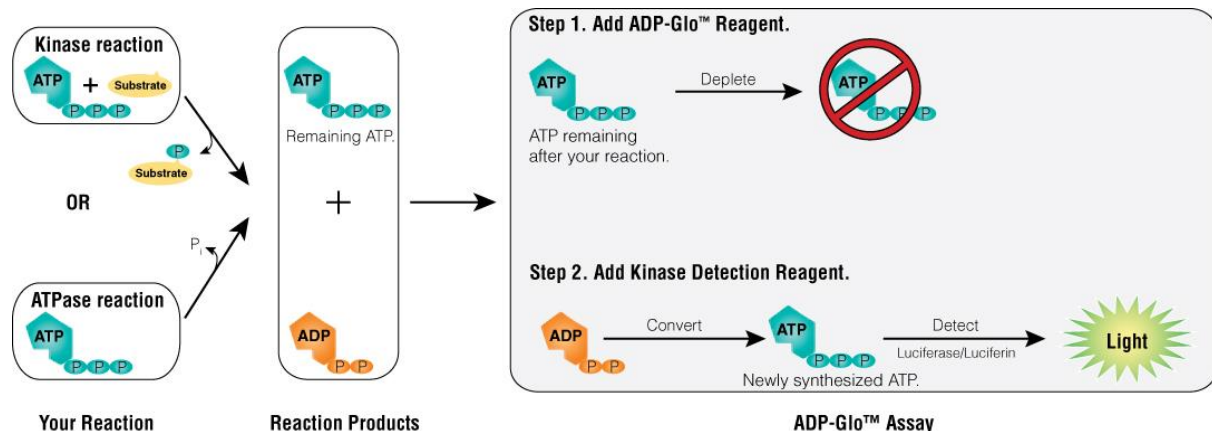
40 min. Incubation

+

10µl Kinase Detection Reagent

30-60 min. Incubation

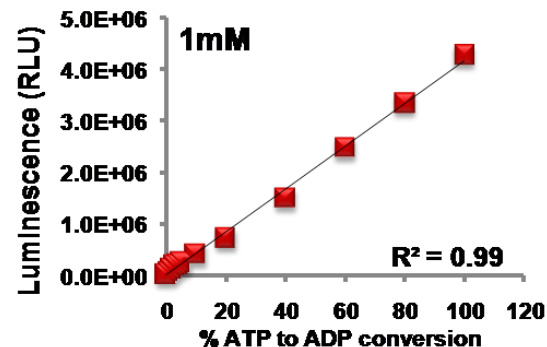
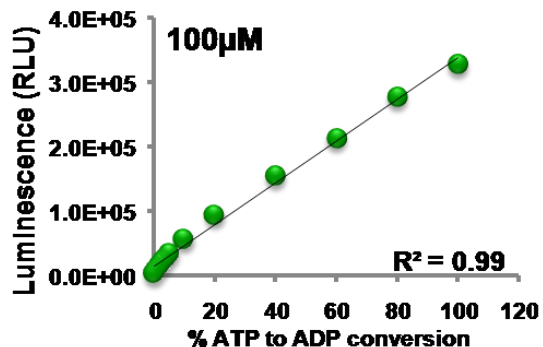
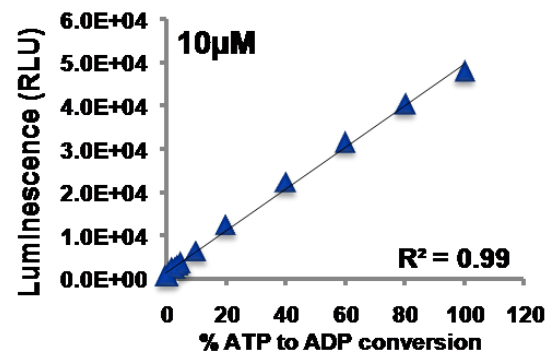
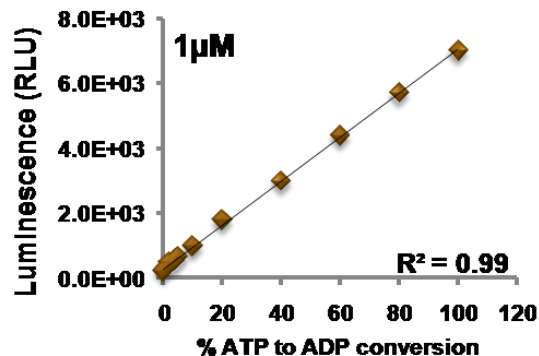
Record Luminescence



White Plates	Format 1:1:2 (µl)	
1,536-well	2.5/2.5/5	
384-well	5/5/10	10/10/20
96-well	25/25/50	50/50/100

Linearity of the ADP-Glo assay

ADP conversion curves at different ADP/ATP concentrations



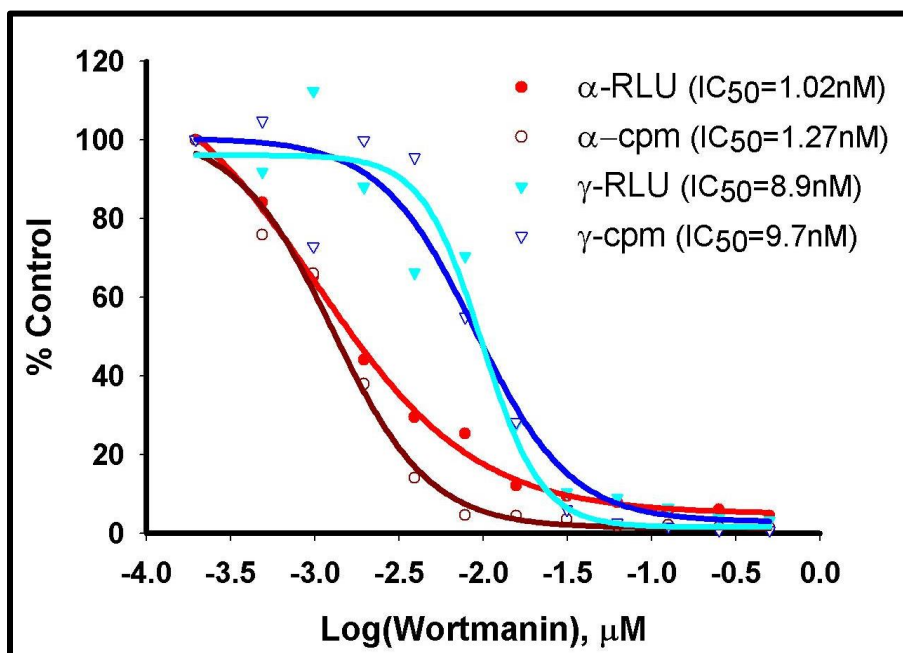
ADP-Glo can be used virtually with any ATP concentration

Sensitivity of the luminescent ADP detection assay

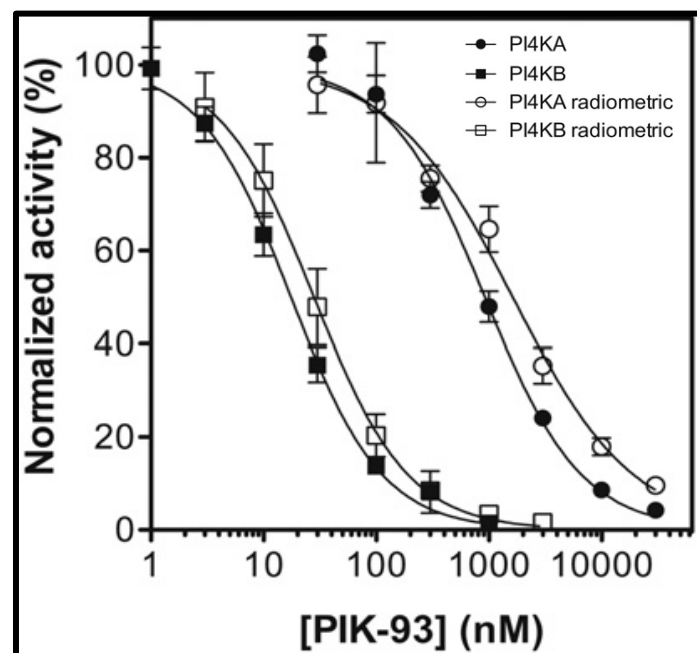
ADP-Glo™ Assay Performance Using Four Different Sources of ATP.									
	ATP-ADP ranges								
	1mM			100μM			10μM		
% ADP in ATP + ADP mixture	20%	10%	5%	20%	10%	5%	20%	10%	5%
Promega	45	23	13	54	32	18	37	20	10
Competitor S	26	13	8	35	21	11	31	16	8
Competitor T	13	7	4	15	10	5	16	9	5
Competitor G	12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

ADP-Glo sensitivity is a proven fact

Comparison between ADP Glo and Radioactivity Assay



[Vidugiriene J, et al \(2009\)](#). Evaluating the utility of a bioluminescent ADP-detecting assay for lipid kinases. [Assay Drug Dev Technol](#). 7(6):585-97.

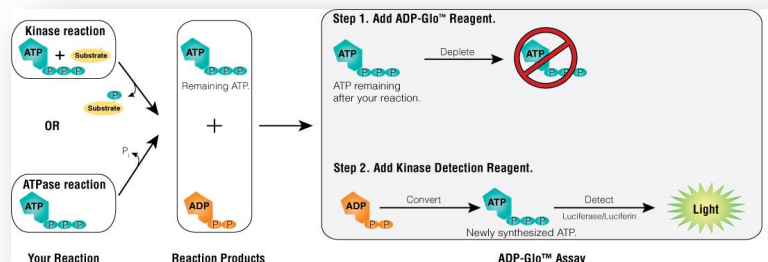


[Tai AW, Bojjireddy N, Balla T. \(2011\)](#) A homogeneous and nonisotopic assay for phosphatidylinositol 4-kinases. [Anal Biochem](#). 417(1):97-102.

Specific activity and response to known inhibitors determined by ADP-Glo correlated well with data from radiometric assay

One assay platform - many applications

ADP-Glo is a Universal *in vitro* Biochemical Assay for all types of Kinase Studies

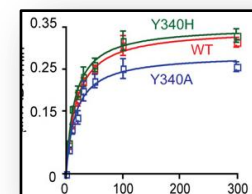


ADP-Glo Assay Platform

High-Throughput Screening



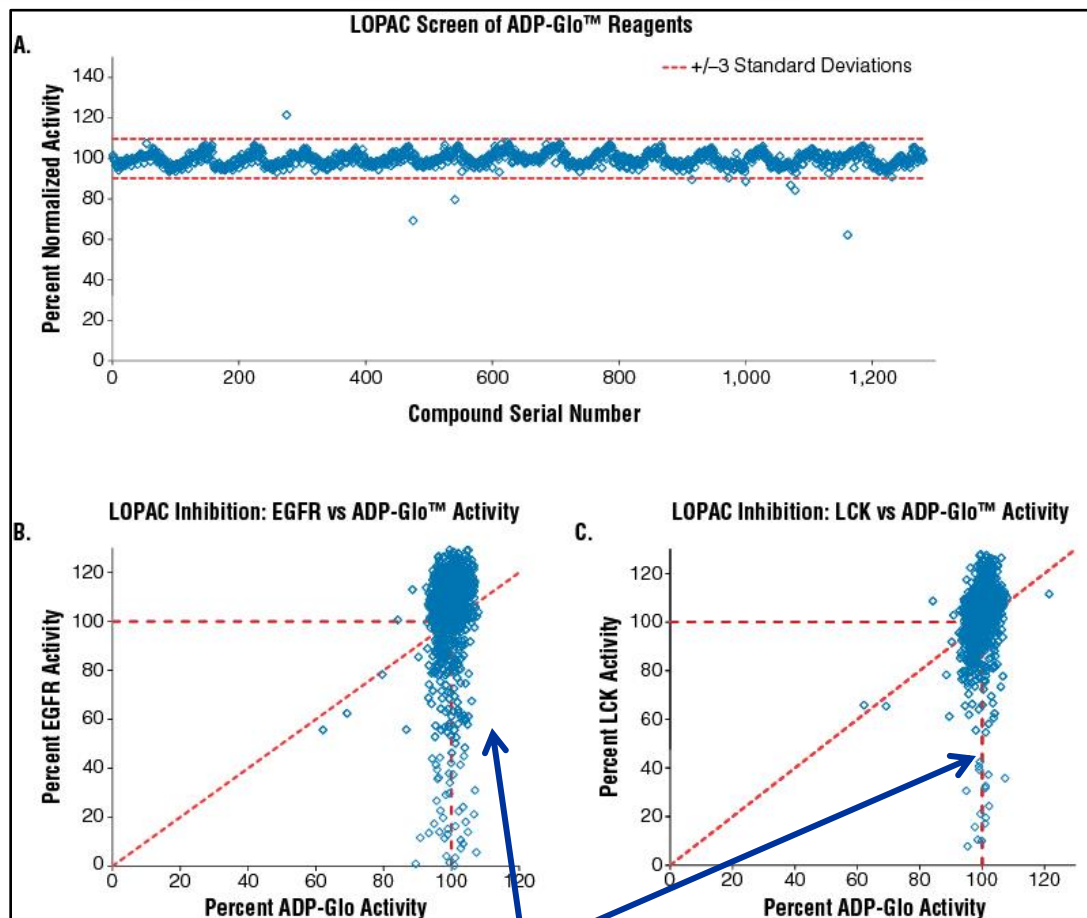
Mode of action studies



Kinase inhibitor profiling

A heatmap showing kinase inhibitor profiling. The table has 10 columns and 20 rows. The columns are labeled with kinase names: PKA, PKB, PKC, PKD, PKF, PKG, PKH, PKI, PKJ, PKL. The rows are labeled with inhibitor names: 1, 2, 3, 4, 5, 6, 7, 8, 9, 10, 11, 12, 13, 14, 15, 16, 17, 18, 19, 20. The cells are colored red or yellow, indicating the level of inhibition.

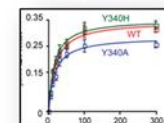
ADP-Glo is resistant to chemical interference



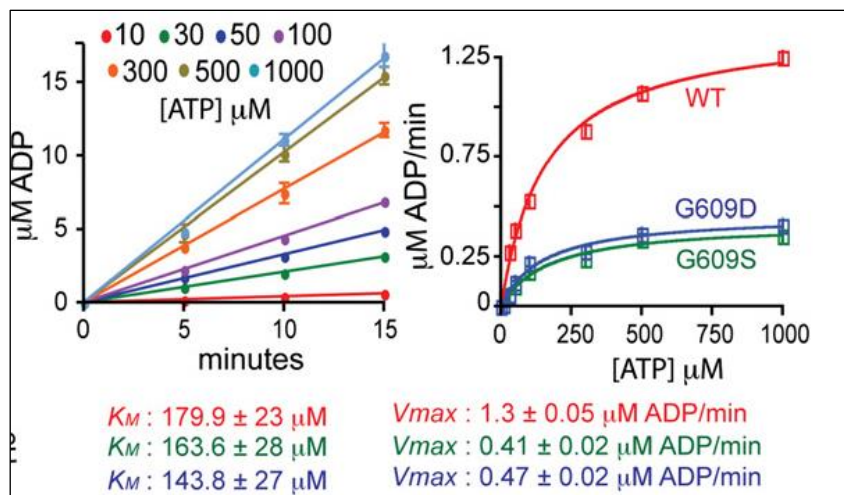
Identification of true **Kinase inhibitors** vs **false positives** using
Luminescence Kinase assay

Determination of Biochemical Values Using ADP-Glo™ Assay

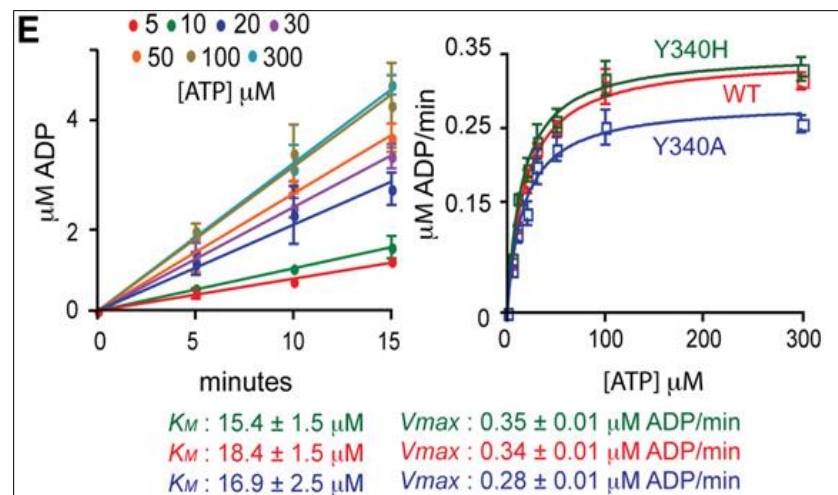
Mode of action studies



C-Src



Haspin

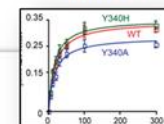


Balzano D et al., (2011). A general framework for inhibitor resistance in protein kinases. *Chem Biol.* 18(8):966-75.

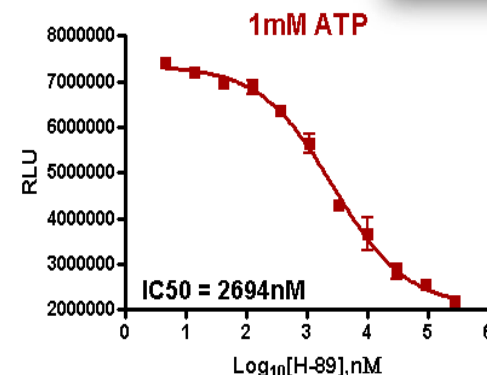
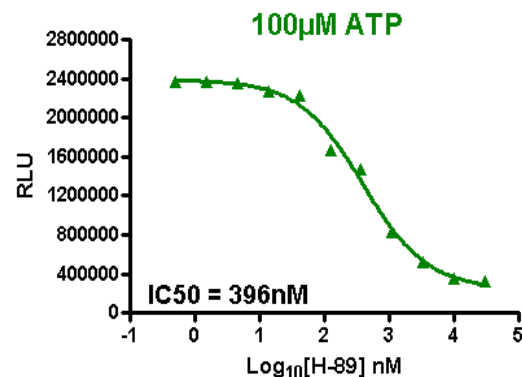
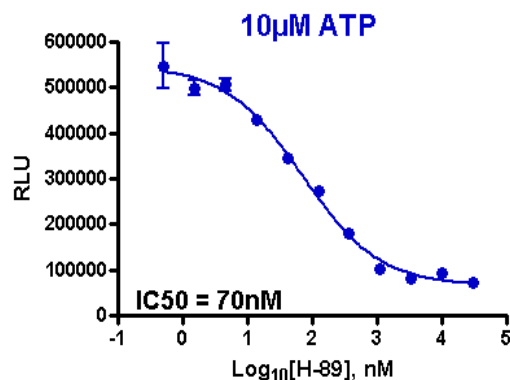
ADP-Glo™ Assay produces K_m values similar to literature

Determination of inhibitor's mechanism of action

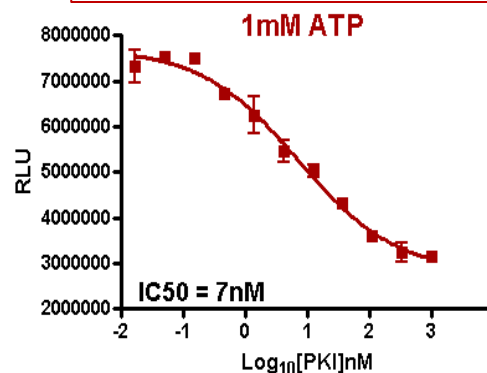
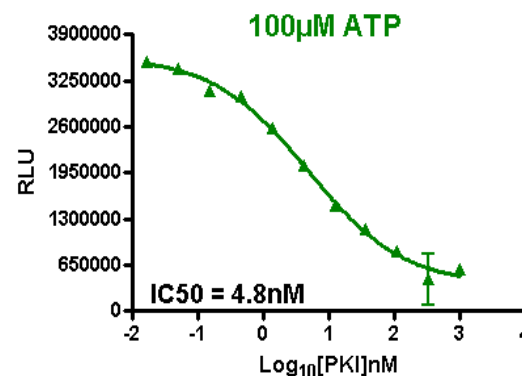
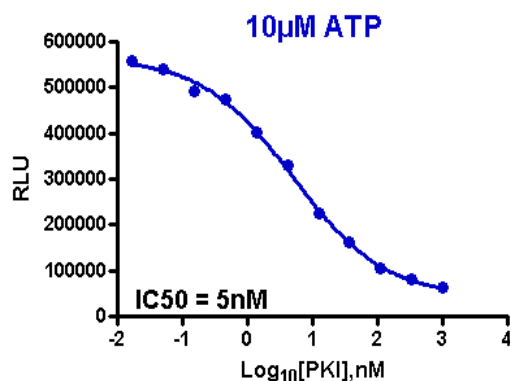
Mode of action studies



PKA ATP Competitive inhibitor H-89



PKA ATP non Competitive inhibitor PKI



Cellular Levels mM

ADP-Glo is a perfect assay to distinguish between ATP competitive and non competitive kinase inhibitors

ADP-Glo Assay in peer review

A subset of ADP-Glo use citations for kinases (a full list is available upon request)

A general framework for inhibitor resistance in protein kinases.

Balzano *et al.* Chemistry & Biology (2011) 18(8):966-975

Evidence that Aurora B is implicated in spindle checkpoint signaling independently of error correction.

Santaguida *et al.* The EMBO Journal (2011) 30, 1508-1519

Comparison of luminescence ADP production assay and radiometric scintillation proximity assay for cdc7 kinase.

Takagi *et al.* Combinatorial Chemistry & High Throughput Screening (2011) 14(8):669-687

A homogeneous and nonisotopic assay for phosphatidylinositol 4-kinases

Tai *et al.* Anal Biochem. (2011) 417(1):97-102

Deoxycytidine kinase regulates the G2/M checkpoint through interaction with cyclin-dependent kinase 1...

Yang *et al.* Nucleic Acid Research 2012, 40(19):9621-9632

Development and Validation of a High-Throughput Intrinsic ATPase Activity Assay for the Discovery of MEKK2...

Ahmad *et al.*, J Biomol Screen. 2012 Nov 7

Domain-Based Biosensor Assay to Screen for Epidermal Growth Factor Receptor Modulators in Live Cells

Antczak *et al.*, ASSAY and Drug Development Technologies 2012, 10(1): 24-36.

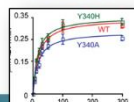
STK33 kinase inhibitor BRD-8899 has no effect on KRAS-dependent cancer cell viability

Luo *et al.*, PNAS, 2012 vol. 109 (8), 2860-2865 .

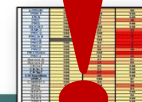
High-Throughput
Screening



Mode of action
studies



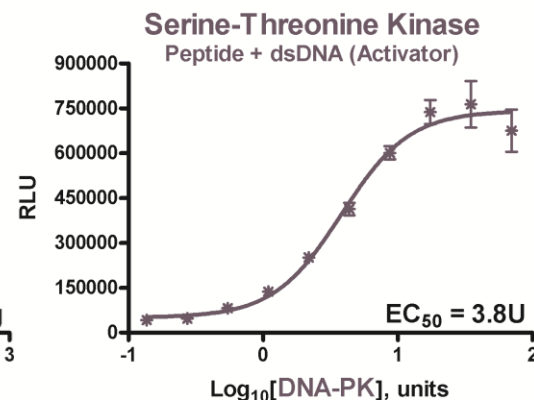
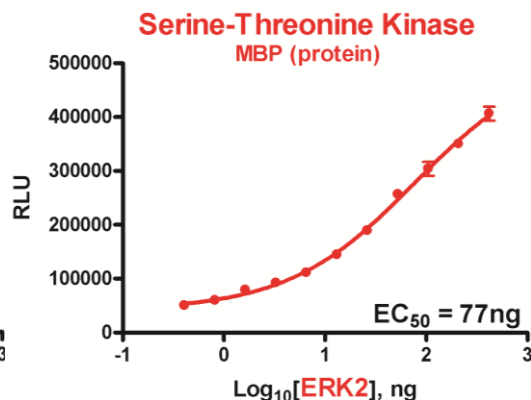
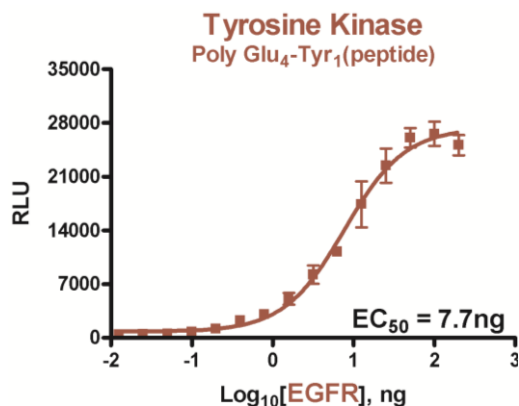
Kinase inhibitor
profiling



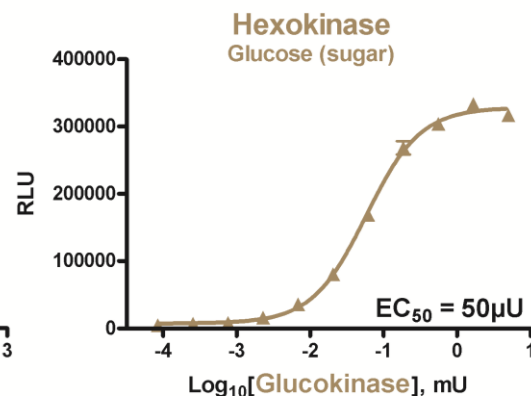
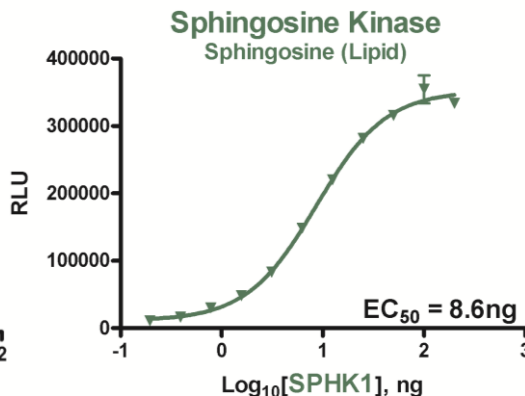
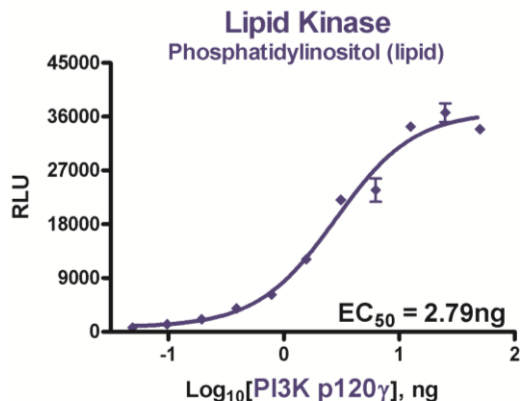
Diverse kinase-substrate combinations with ADP-Glo

Kinase Inhibitor Profiling

Protein Kinases

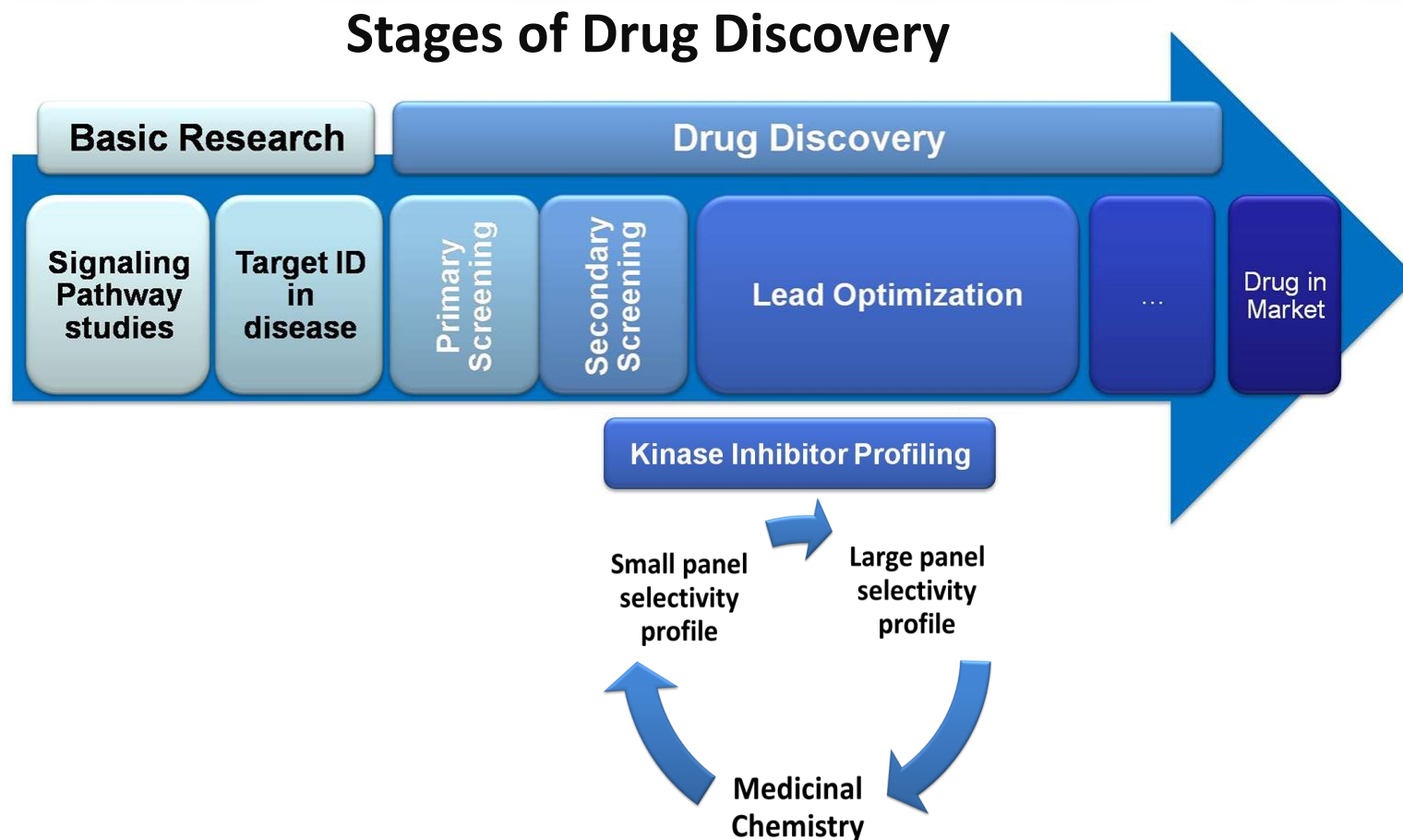


Non-Classical Kinases



ADP-Glo Kinase Assay detects the activity of any Kinase regardless of the substrate chemical structure

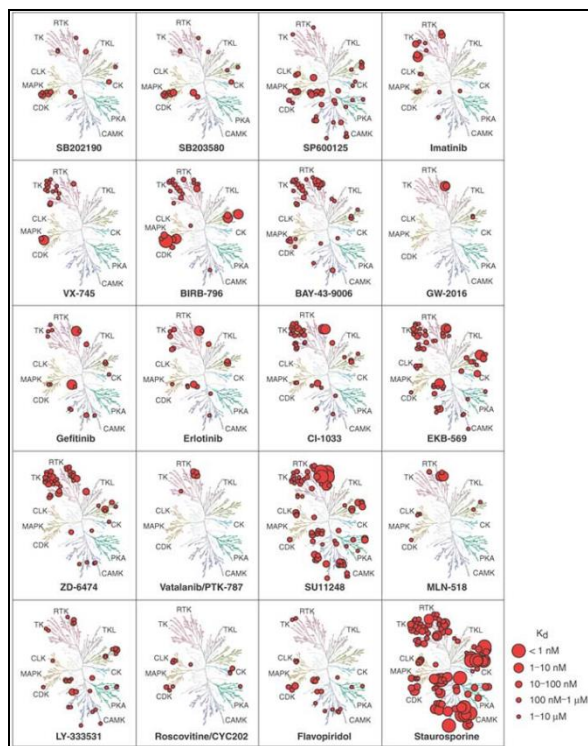
Kinase inhibitor Profiling is an important Step of Drug Discovery



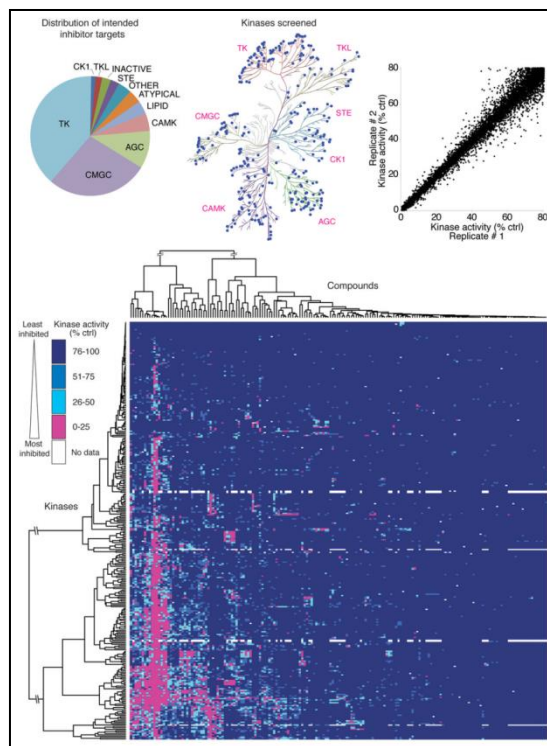
Kinase selectivity profiling allows for a balance between potency and selectivity of lead compounds during drug discovery

Kinase Profiling data generated in many ways

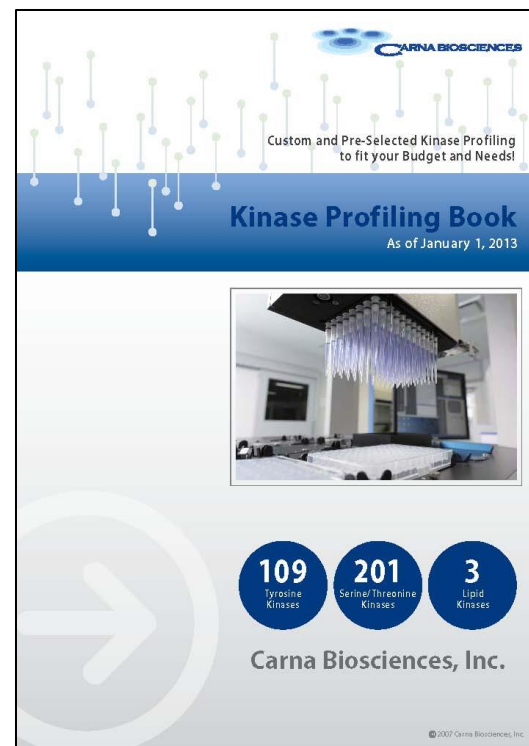
Many profiling platforms exist already



Binding



Radiometric



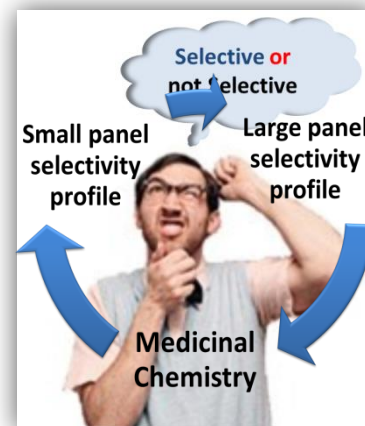
Mobility Shift

Large panel kinase profiles may not be performed routinely
Choosing the right approach for best results

Issues related to Kinase Profiling

Current Challenges in creating selectivity profiles for Kinase inhibitors

- ✓ Addressing confidentiality issues related to outsourcing.
- ✓ In house profiling requires kinase assay developments and rigorous optimization which can be time consuming.
- ✓ The lack of having enough enzymes available to do the profiling.
- ✓ Generating Selectivity profiles is costly and time consuming.
- ✓ Immediate accessibility of the data upon completion of Profiling.



There is a need in an **easy and flexible** solution that allows kinase selectivity profiles to be generated routinely without the time and cost challenges.

Kinase Enzyme System (KES) format



ADP-Glo Kinase Assay 0-1mM ATP

0.5ml	10mM UltraPure ATP
0.5ml	10mM ADP
5ml	ADP-Glo Reagent
10ml	Kinase Detection Buffer
1cake	Kinase Detection Substrate



Akt1 Kinase Enzyme System (Example)

0.1ml	AKT1 Kinase (10µg)
1ml	AKT (SGK) substrate (1mg)
1.5ml	Kinase Assay buffer
25µl	100mM DTT
25µl	2.5M MnCl ₂ *
500µl	Kinase Activator**

*For Tyrosine Kinases
**Lipids for PKC Kinases

KES: a complete kinase solution

Lipid Kinase Systems

Complete solution for measuring PI kinases without the use of radioactivity,
lipid substrate modification or lipid extraction

PI3K class I enzymes:

p110 α /p85 α
p110 α (E545K)/p85 α
p110 α (H1047R)/p85 α
P110 β / p85 α
P110 γ / p85 α
P110 δ / p85 α

+

PI:PS Lipid
Kinase
substrate
0.5ml x 1mM

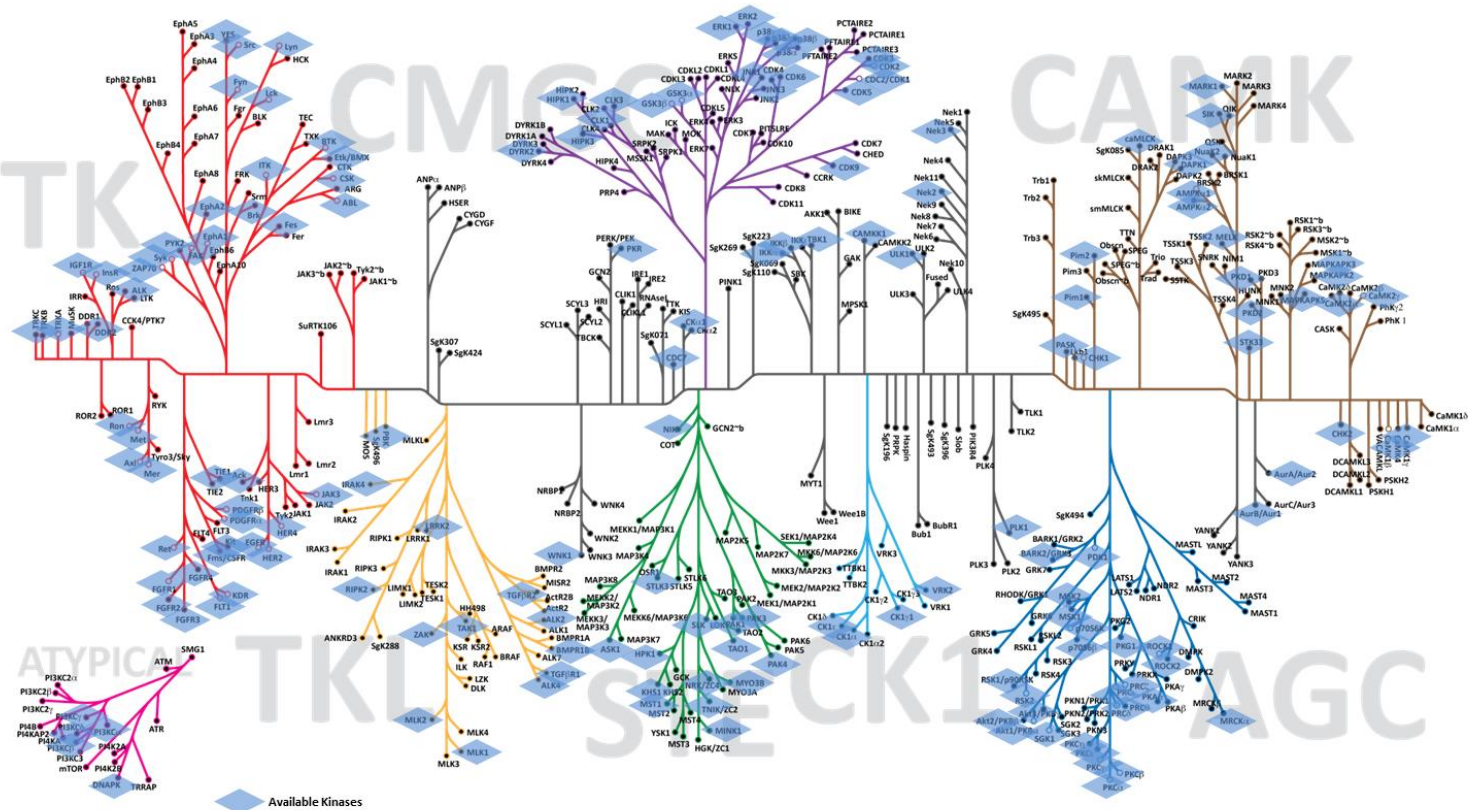
or

PIP2:PS Lipid
Kinase
substrate
0.25ml x 1mM

ADP-Glo Kinase Assay

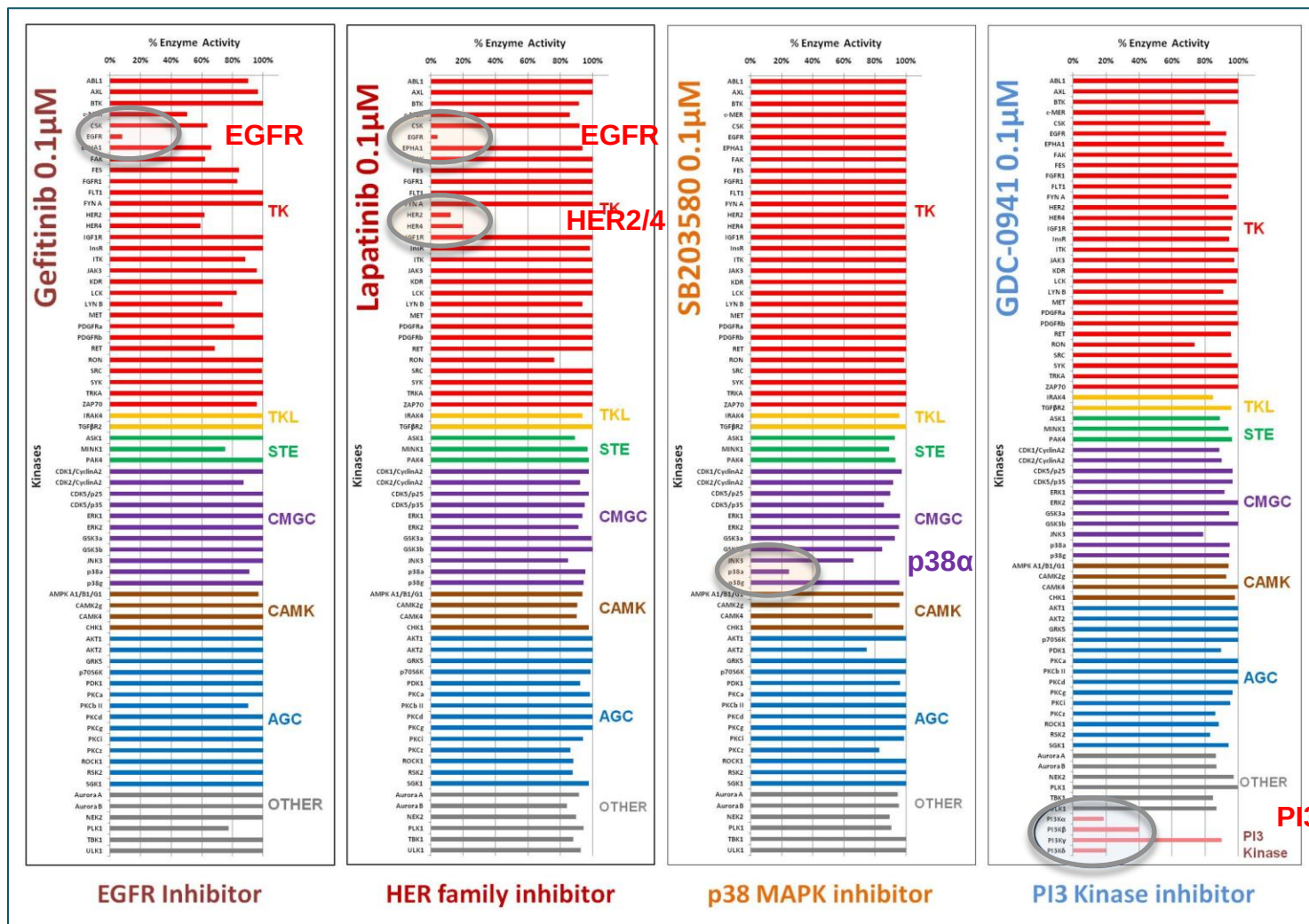
Addressing the kinome with large kinase panel

Profiling Panel should include close and distant kinases to assess compound selectivity



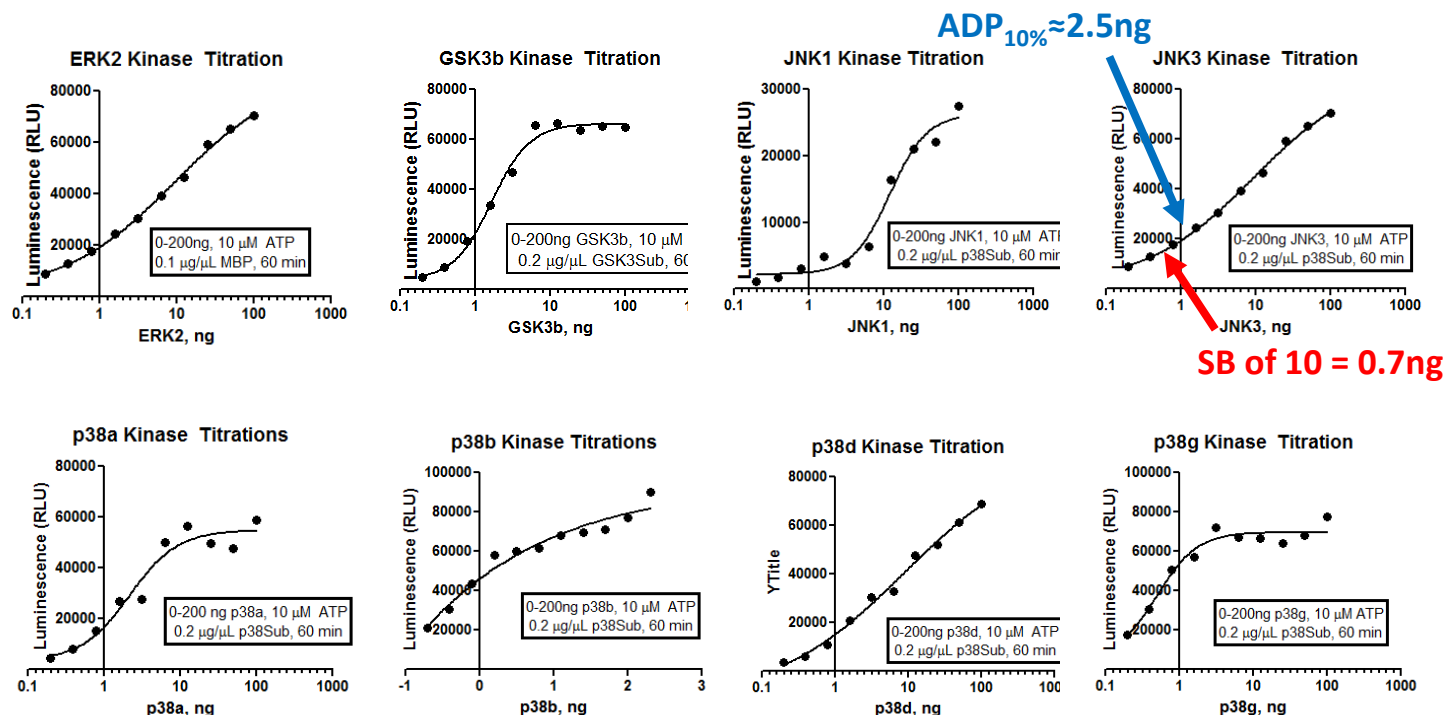
Broad Human Kinome coverage with >170 Kinase Enzyme Systems

Profiling inhibitors against a subset Kinase panel with ADP-Glo Kinase Assay



Optimizing enzyme concentration for kinase inhibitor studies

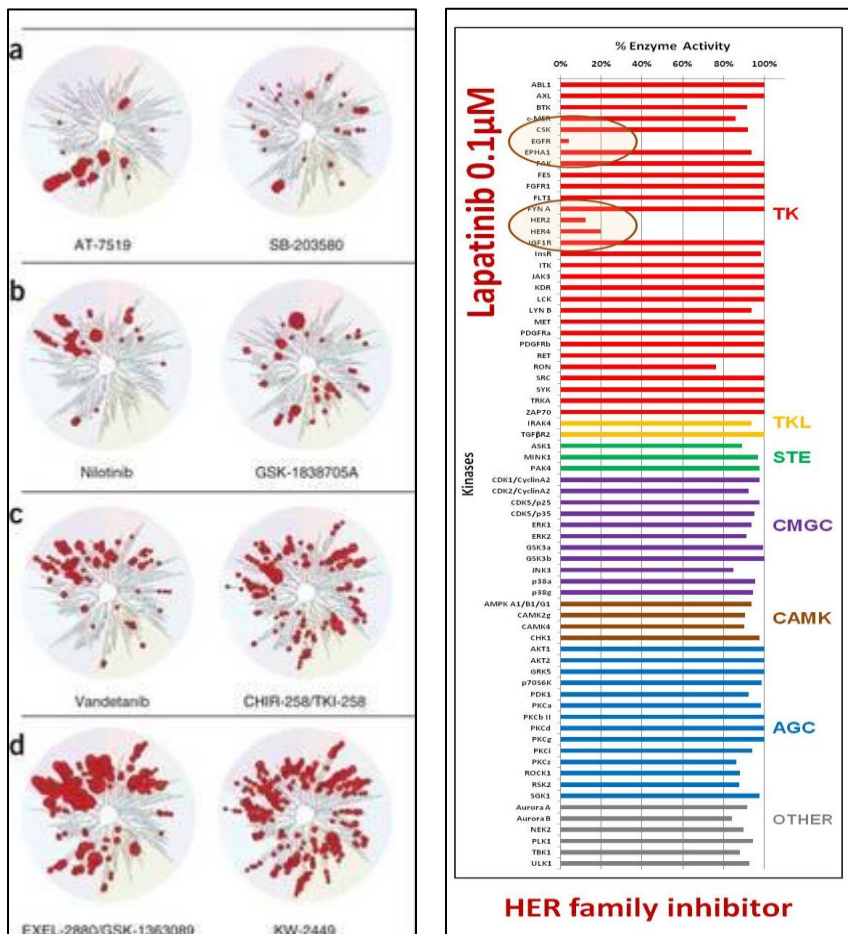
kinases need to be titrated to find the optimal amount to be used



➤ **Determination of SB and % conversion values:** kinase amount needed to generate a signal-to-background ratio of ≥ 10 and optimal % ATP conversion.

Large kinase profiling data vs. routine small selectivity panels

Did we achieve our goal of ease of use and flexibility?



- ✓ Costly and time consuming
- ✓ Extensive assay development required.

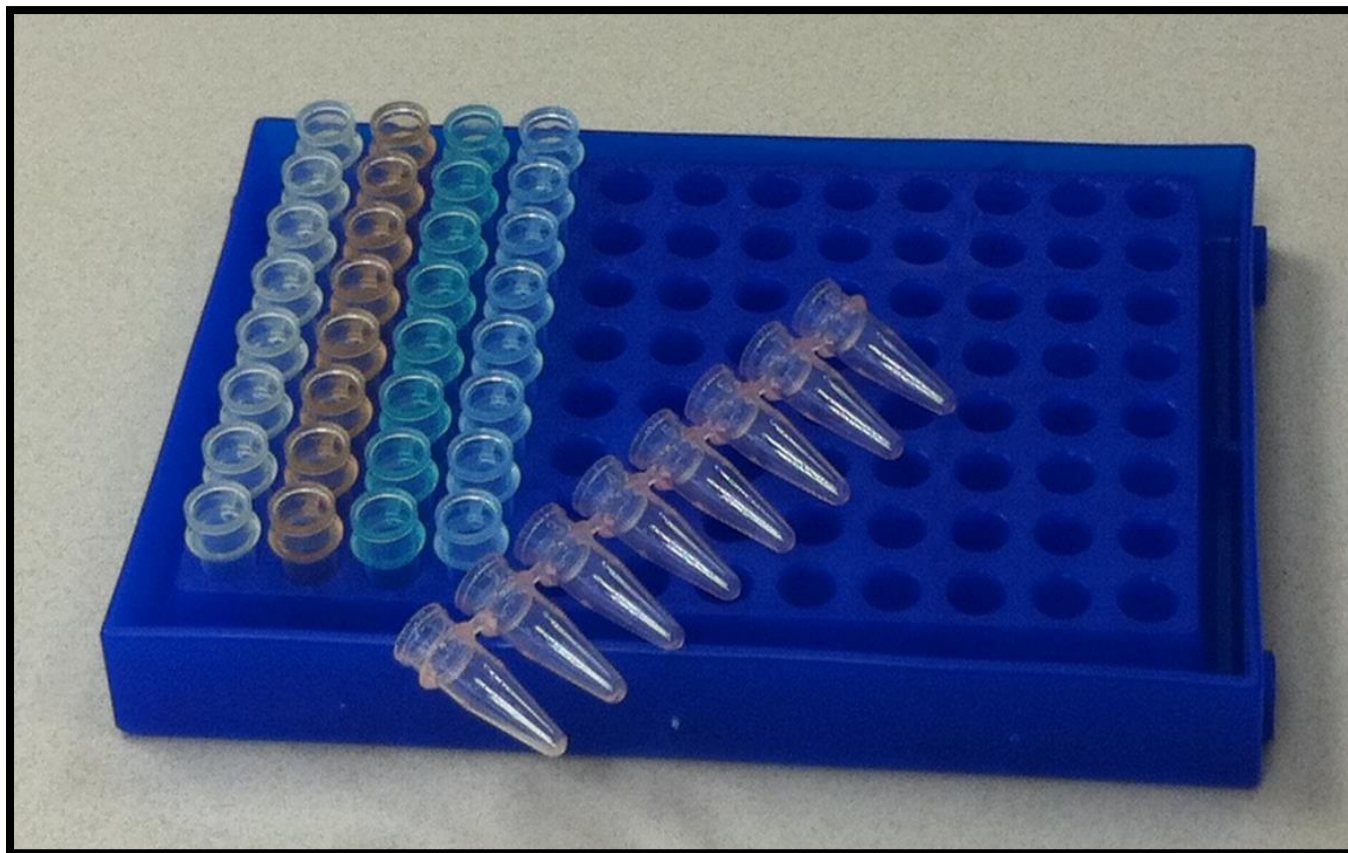
Developing a simple kinase profiling assay

Choosing the right kinases for flexible and targeted inhibitor profiling

Important kinase targets organized by kinase families in 8 well strips

**17 Kinase strips
(112* Kinases)**

***14 Family groups
and 3 mixed panels**



- Ready to profile with simple protocol (No development time required)
- Cost effective
- Quick results with no compound sending out

Kinase profiling Systems: 2 types

Family based strips

General panel

TK-1	TK-2	TK-3	TK-4	CMGC-1	CMGC-2	TKL-1	STE-1	CAMK-1	CAMK-2	AGC-1	AGC-2	OTHER/CK-1	OTHER-2	General Panel		
EGFR	ABL1	AXL1	c-MER	ERK2	CDK1/CyclinA2	ALK2	ASK1	CHK1	AMPK A1/B1/G1	AKT1	PKCa	Aurora A	CDC7/DBF4	FGFR1	CDK2/CyclinE1	AKT1
HER2	BRK	EPHA1	FGFR1	GSK3b	CDK2/CyclinE1	ALK4	HPK1	CHK2	AMPK A1/B1/G2	p70S6Kb	PKCb II	Aurora B	EIF2AK2	JAK3	GSK3b	PKCa
HER4	BTK	FAK	FGFR2	JNK1	CDK3/CyclinE1	IRAK4	MINK1	MAPKAPK2	AMPK A2/B1/G1	PDK1	PKCd	CK2a1	IKKb	LCK	p38a	ROCK1
IGF1R	CSK	ITK	FGFR4	JNK3	CDK5/p25	MLK2	MST1	MARK1	CAMK2a	PKA	PKCe	DNA-PK	NEK2	SYK	AMPK A1/B1/G2	Aurora A
InsR	FYN A	JAK3	FLT1	p38a	CDK5/p35	RIPK2	NIK	MELK	CAMK2g	PKC	PKCg	CK1α1	NEK3	MINK1	CAMK4	CK2a1
KDR	LCK	PYK2	FMS	p38b	CDK6/Cyclin D3	TAK1-TAB1	PAK1/CDC42	PASK	CAMK4	PRKG1	PKCi	CK1epsilon	PLK1	PAK1/CDC42	CHK1	IKKb
PDGFRa	LYN B	SYK	MET	p38d	CDK9/Cyclin K	TGFB2	PAK3	PIM1	DAPK1	ROCK1	PKCtheta	CK1g1	TBK1	IRAK4	DAPK1	CK1a1
PDGFRb	SRC	TRKA	RET	p38g	CLK3	ZAK	TNIK	PKCmu (PKD1)	STK33	RSK2	PKCz	VRK2	ULK1	TAK1-TAB1	MAPKAPK2	CK1g1

Relevant kinases organized in:

- Family based strips for intra-family inhibitor promiscuity analysis.
- General panel for quick selectivity assessment.
- Multiple strips are used for large selectivity profiling.

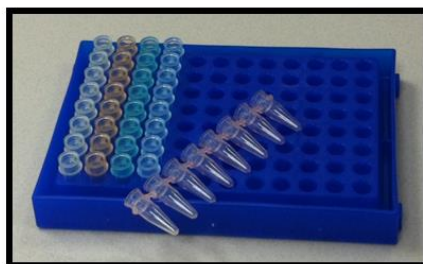
kinase profiling strip Concept

How does it work?

1	9	17	25	33	41	49	57	65	73	81	89
2	10	18	26	34	42	50	58	66	74	82	90
3	11	19	27	35	43	51	59	67	75	83	91
4	12	20	28	36	44	52	60	68	76	84	92
5	13	21	29	37	45	53	61	69	77	85	93
6	14	22	30	38	46	54	62	70	78	86	94
7	15	23	31	39	47	55	63	71	79	87	95
8	16	24	32	40	48	56	64	72	80	88	96

50x

Kinase Stocks (5 μ l)



S1	S9	S17	S25	S33	S41	S49	S57	S65	S73	S81	S89
S2	S10	S18	S26	S34	S42	S50	S58	S66	S74	S82	S90
S3	S11	S19	S27	S35	S43	S51	S59	S67	S75	S83	S91
S4	S12	S20	S28	S36	S44	S52	S60	S68	S76	S84	S92
S5	S13	S21	S29	S37	S45	S53	S61	S69	S77	S85	S93
S6	S14	S22	S30	S38	S46	S54	S62	S70	S78	S86	S94
S7	S15	S23	S31	S39	S47	S55	S63	S71	S79	S87	S95
S8	S16	S24	S32	S40	S48	S56	S64	S72	S80	S88	S96

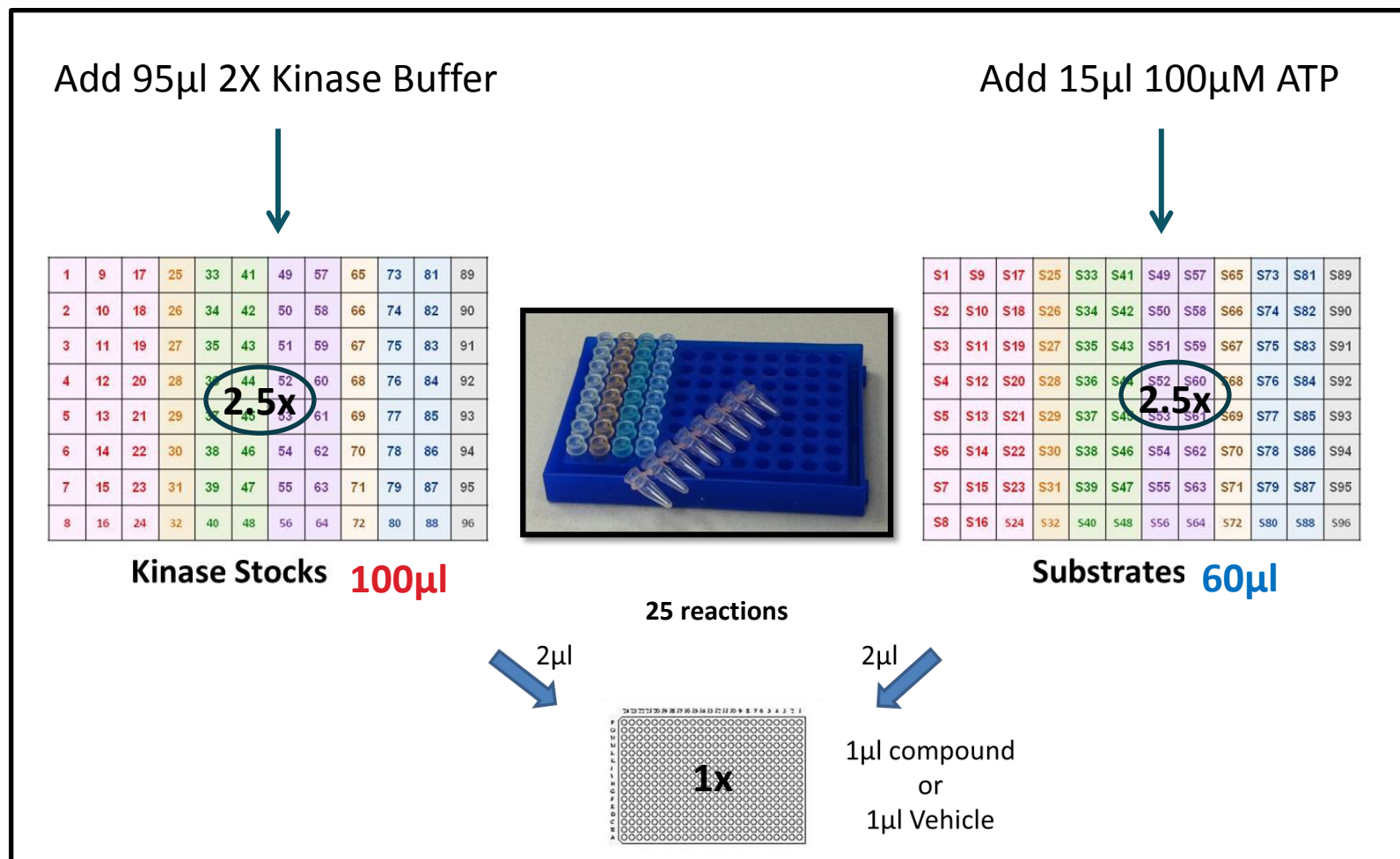
3.3x

Substrates (45 μ l)

Simple reaction assembly

kinase profiling strip Concept

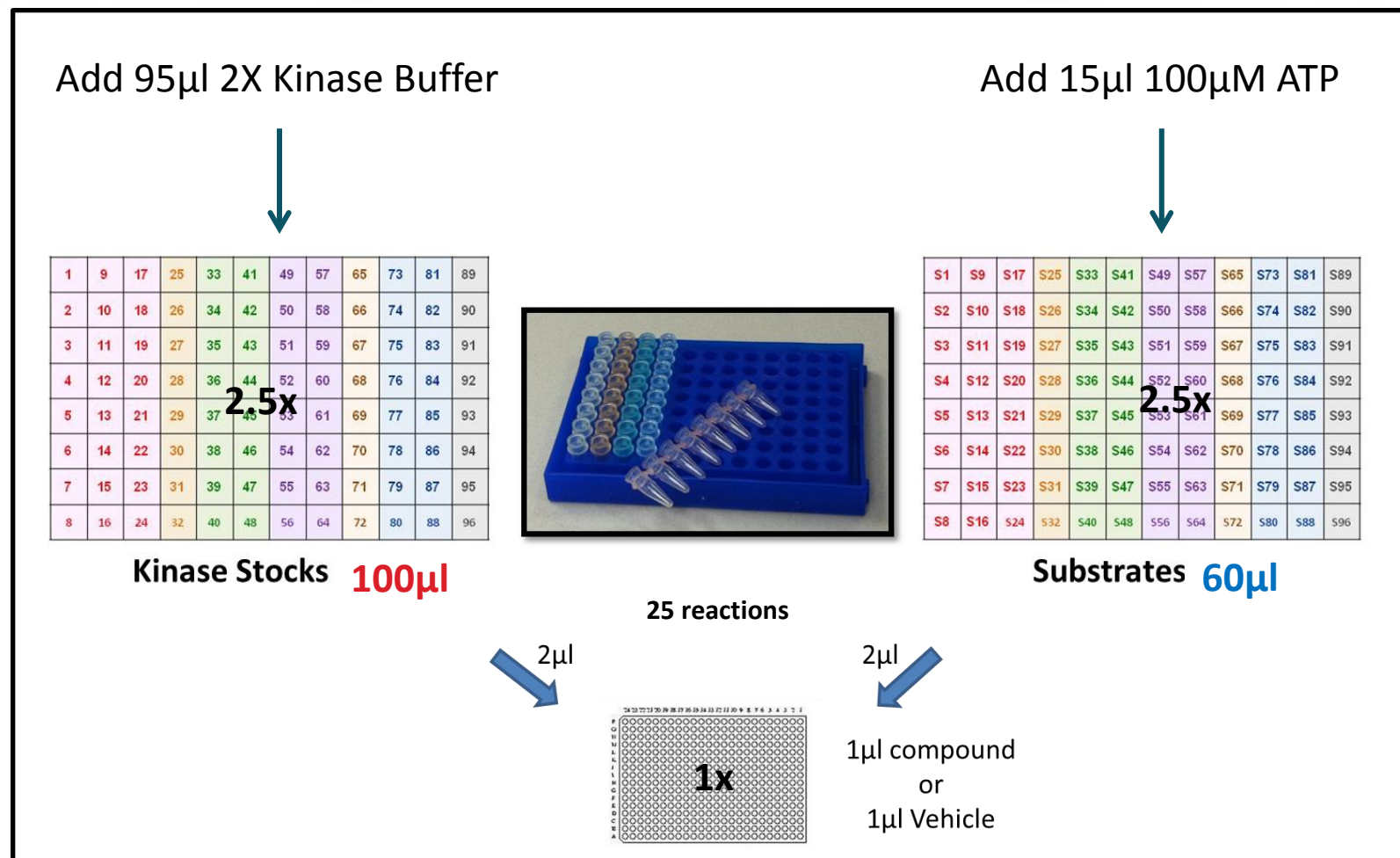
How does it work?



Simple reaction assembly

kinase profiling strip Concept

How does it work?

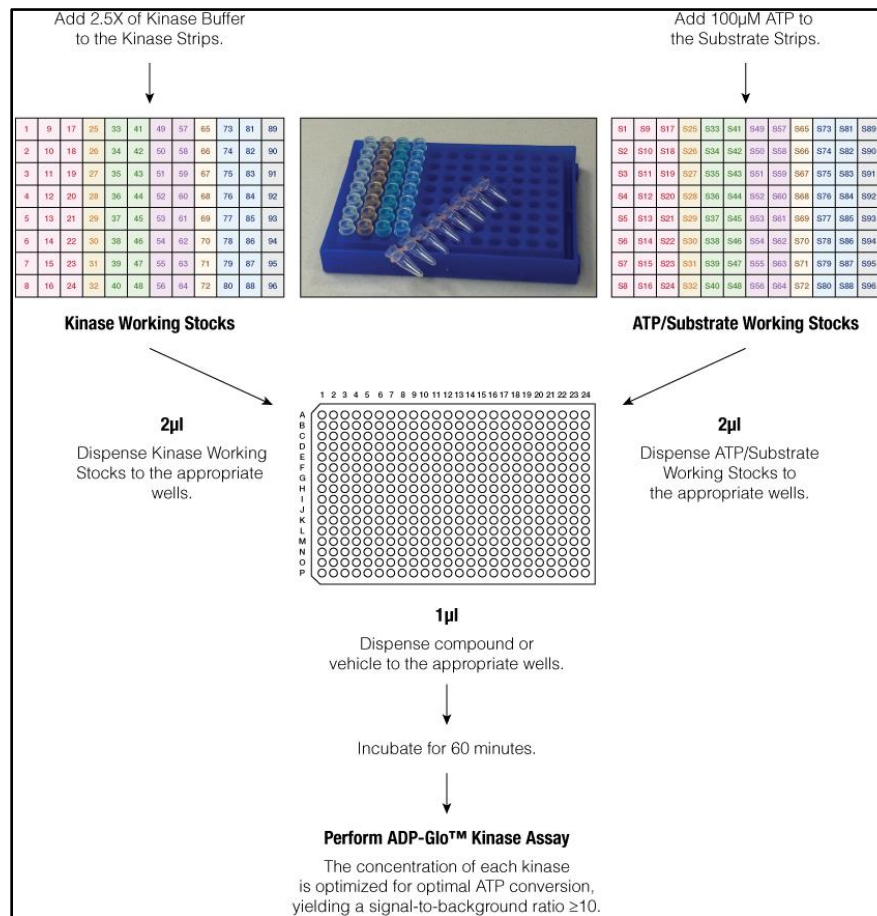


Each enzyme will generate an optimal % ATP conversion with an SB>10 fold

Simple profiling protocol for flexible and targeted inhibitor profiling

Kinase Profiling Strip Systems:

- ✓ **One-time use design:** Eliminating multiple freeze/thaw cycles ensures optimal kinase activity per experiment.
- ✓ **Flexible kinase inhibitor profiling:** Each strip has enough material to profile:
 - **Dose response for 1 compound against the 8 kinases at once.**
 - **10 compounds at a single dose**
- ✓ **Easy-to-automate inhibitor profiling.**



Profiling with ADP-Glo platform made simple

Streamline Your Kinase Profiling Workflow

Liquid Handling

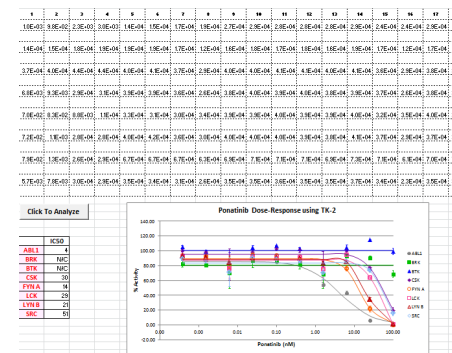


Detection



**GloMax[®]
Discover**

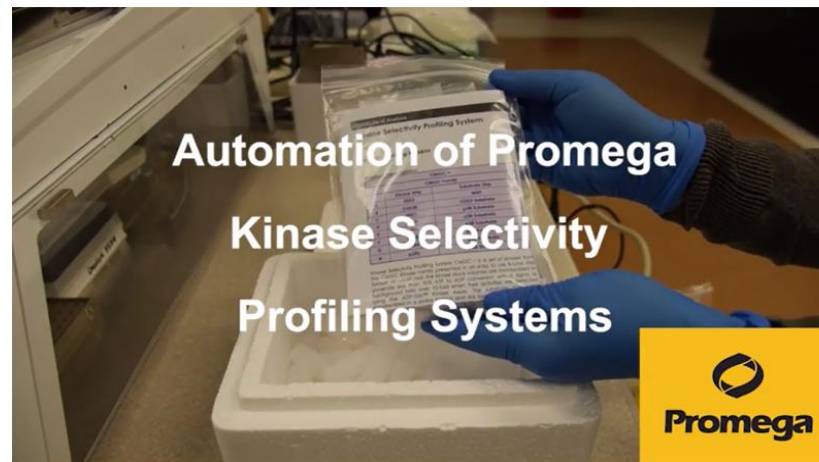
Data Analysis



**GloMax[®]
SMART Protocol**

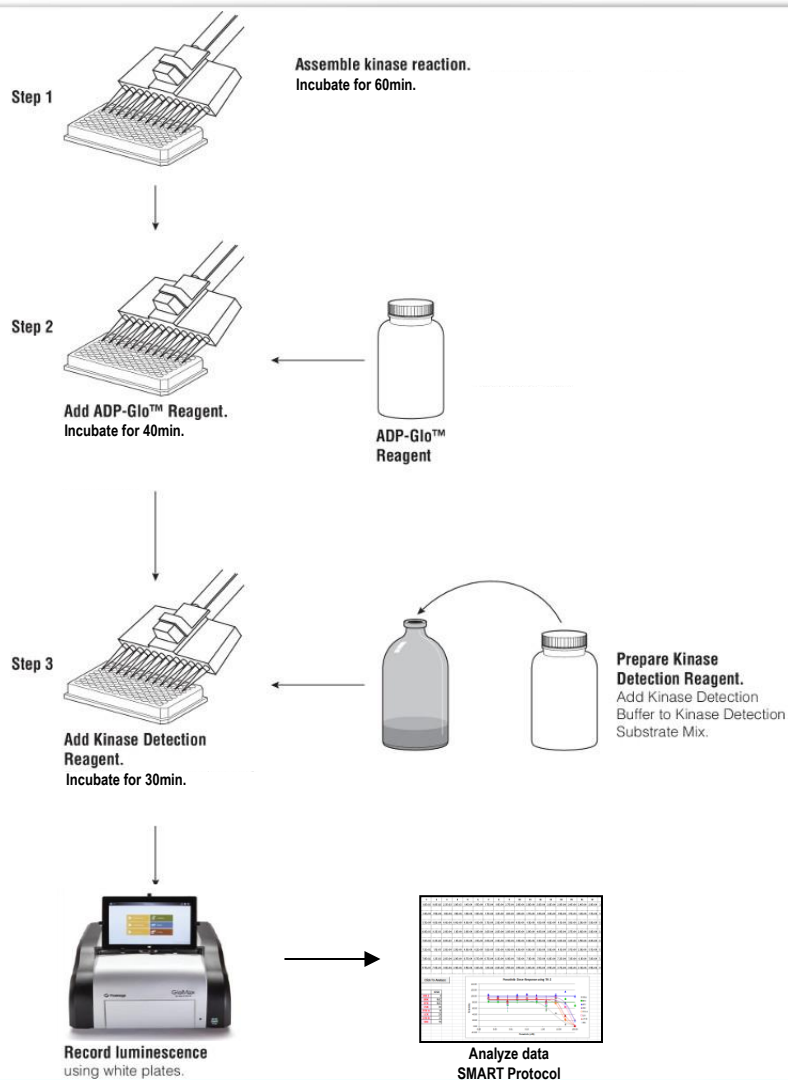
Automation of kinase profiling strip

How does it work?



[Click on the link below to watch the video](https://p.widencdn.net/pwhnbo)
<https://p.widencdn.net/pwhnbo>

ADP-Glo Assay Scheme



- Two simple additions after the Kinase reaction
- Record luminescence and export automatically data to SMART Protocol

Easily Collect Kinase Profiling Data and Quickly Analyze Results using GloMax[®] Systems

Select the built-in Kinase Selectivity Profiling SMART Protocol

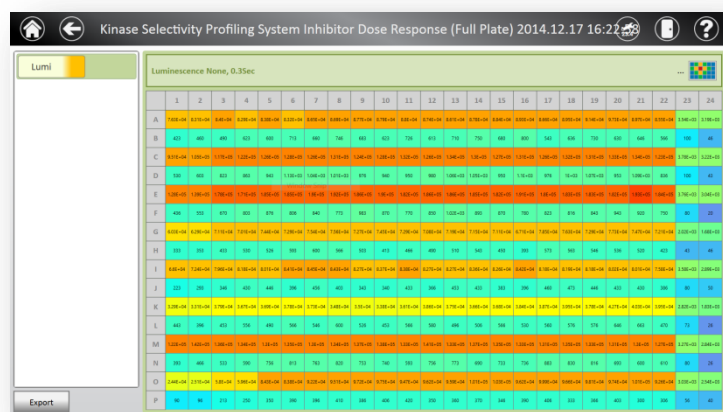
- Screen kinase inhibitors
- Automatically analyze results



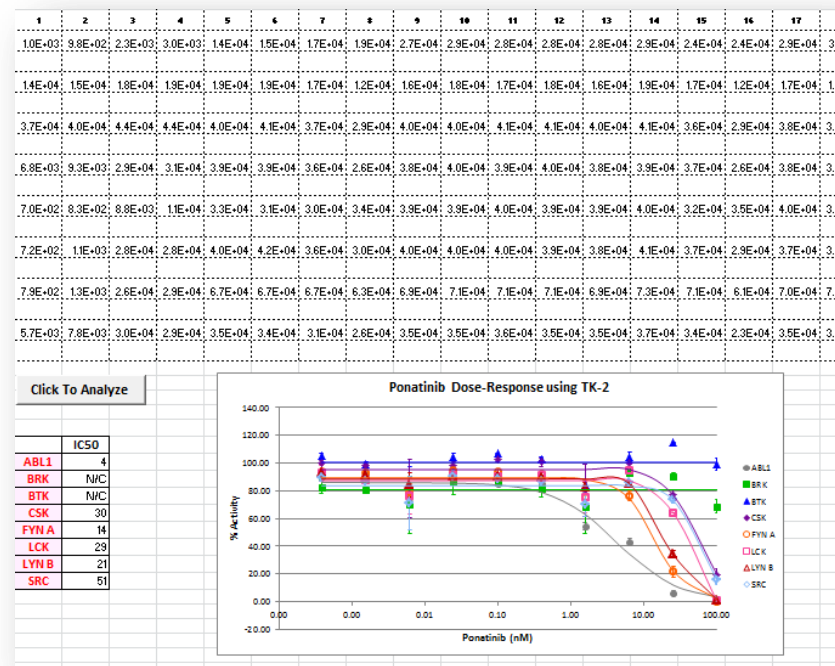
Over 50 Promega assay protocols are optimized and included with the GloMax[®] Discover System to streamline data collection and simplify your research

Easily Collect Kinase Profiling Data and Quickly Analyze Results using GloMax[®] Systems

View a heat map display to quickly assess the data



View analyzed dose response results after data collection



Kinase Selectivity Profiling Systems (KPS) *

Single Dose SMART Protocol

*Also available as a downloadable template on promega.com

Single Dose SMART Protocol

Raw data is imported automatically in the template. Click to analyze.

Single Dose SMART 2 Strip CV updating.xlsm [Read-Only] - Microsoft Excel

File Home Insert Page Layout Formulas Data Review View Add-Ins Acrobat

Clipboard Paste Copy Format Painter Font Calibri 11 A A Scientific Normal 2 Normal 3 Normal Bad Good Neutral Insert Delete Format AutoSum Fill Clear Sort & Filter Select

M17 99404

Kinase Selectivity Profiling Systems (KPS)
Single Dose SMART Protocol

Luminescence
Emission Filter: None
Integration Time(s): 0.3

Reading:	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	
A	5.34E+03	2.88E+04	1.64E+04	6.85E+03	3.05E+04	3.57E+04	3.54E+04	3.45E+04	4.05E+04	3.34E+04	2.76E+04	2.13E+04	5.34E+03	2.88E+04	1.64E+04	6.85E+03	3.05E+04	3.57E+04	3.54E+04	3.45E+04	4.05E+04	3.34E+04	2.76E+04	2.13E+04	5.34E+03
B	5.21E+03	3.19E+04	1.80E+04	6.69E+03	3.00E+04	3.46E+04	3.49E+04	3.32E+04	3.53E+04	3.53E+04	3.46E+04	2.59E+04	5.21E+03	3.19E+04	1.80E+04	6.69E+03	3.00E+04	3.46E+04	3.49E+04	3.32E+04	3.53E+04	3.53E+04	3.46E+04	2.59E+04	5.21E+03
C	5.98E+03	2.59E+04	1.09E+04	7.59E+03	3.73E+04	2.59E+04	2.66E+04	2.86E+04	2.69E+04	3.23E+04	2.90E+04	1.05E+04	5.98E+03	2.59E+04	1.09E+04	7.59E+03	3.73E+04	2.59E+04	2.66E+04	2.86E+04	2.69E+04	3.23E+04	2.90E+04	1.05E+04	5.98E+03
D	4.98E+03	2.37E+04	1.06E+04	7.89E+03	3.54E+04	2.49E+04	2.63E+04	2.63E+04	2.62E+04	3.28E+04	2.81E+04	1.06E+04	4.98E+03	2.37E+04	1.06E+04	7.89E+03	3.54E+04	2.49E+04	2.63E+04	2.63E+04	2.62E+04	3.28E+04	2.81E+04	1.06E+04	4.98E+03
E	5.95E+03	3.78E+04	3.55E+04	3.31E+04	2.11E+04	4.18E+04	4.26E+04	4.22E+04	5.94E+04	3.86E+04	4.01E+04	3.81E+04	5.95E+03	3.78E+04	3.55E+04	3.31E+04	2.11E+04	4.18E+04	4.26E+04	4.22E+04	5.94E+04	3.86E+04	4.01E+04	3.81E+04	5.95E+03
F	8.24E+03	3.96E+04	3.40E+04	2.77E+04	2.13E+04	4.22E+04	5.85E+04	4.32E+04	4.12E+04	3.92E+04	4.15E+04	3.54E+04	8.24E+03	3.96E+04	3.40E+04	2.77E+04	2.13E+04	4.22E+04	5.85E+04	4.32E+04	4.12E+04	3.92E+04	4.15E+04	3.54E+04	8.24E+03
G	5.10E+03	9.46E+04	3.73E+04	1.02E+05	9.78E+04	1.01E+05	9.41E+04	1.07E+05	1.00E+05	1.00E+05	1.02E+05	3.26E+04	5.10E+03	9.46E+04	3.73E+04	1.02E+05	9.78E+04	1.01E+05	9.41E+04	1.07E+05	1.00E+05	1.00E+05	1.02E+05	3.26E+04	5.10E+03
H	4.95E+03	9.07E+04	6.06E+04	9.55E+04	9.25E+04	9.69E+04	9.04E+04	9.94E+04	9.74E+04	8.82E+04	1.08E+05	3.27E+04	4.95E+03	9.07E+04	6.06E+04	9.55E+04	9.25E+04	9.69E+04	9.04E+04	9.94E+04	9.74E+04	8.82E+04	1.08E+05	3.27E+04	4.95E+03
I	5.42E+03	7.81E+03	5.41E+03	7.95E+03	7.65E+03	7.65E+03	7.45E+03	7.82E+03	7.82E+03	7.82E+03	8.04E+03	5.04E+03	5.42E+03	7.81E+03	5.41E+03	7.95E+03	7.65E+03	7.65E+03	7.45E+03	7.82E+03	7.82E+03	7.82E+03	8.04E+03	5.04E+03	5.42E+03
J	5.12E+03	7.66E+03	5.09E+03	7.47E+03	7.70E+03	7.87E+03	7.50E+03	7.54E+03	7.59E+03	7.29E+03	7.78E+03	4.71E+03	5.12E+03	7.66E+03	5.09E+03	7.47E+03	7.70E+03	7.87E+03	7.50E+03	7.54E+03	7.59E+03	7.29E+03	7.78E+03	4.71E+03	5.12E+03
K	6.05E+03	1.37E+05	6.34E+03	1.12E+05	1.18E+05	1.17E+05	7.20E+04	1.17E+05	1.19E+05	1.41E+05	1.16E+05	5.87E+03	6.05E+03	1.37E+05	6.34E+03	1.12E+05	1.18E+05	1.17E+05	7.20E+04	1.17E+05	1.19E+05	1.41E+05	1.16E+05	5.87E+03	6.05E+03
L	5.62E+03	1.17E+05	5.88E+03	1.00E+05	1.08E+05	1.19E+05	7.27E+04	1.17E+05	1.19E+05	1.25E+04	1.18E+05	5.67E+03	5.62E+03	1.17E+05	5.88E+03	1.00E+05	1.08E+05	1.19E+05	7.27E+04	1.17E+05	1.19E+05	1.25E+04	1.18E+05	5.67E+03	5.62E+03
M	5.84E+03	6.99E+04	5.05E+03	3.61E+04	7.10E+03	6.12E+04	4.21E+04	5.43E+04	2.44E+04	5.72E+03	6.32E+03	4.66E+03	5.84E+03	6.99E+04	5.05E+03	3.61E+04	7.10E+03	6.12E+04	4.21E+04	5.43E+04	2.44E+04	5.72E+03	6.32E+03	4.66E+03	5.84E+03
N	5.23E+03	6.99E+04	5.12E+03	4.61E+04	6.72E+03	6.02E+04	4.59E+04	5.77E+04	4.64E+04	5.30E+03	6.12E+03	5.23E+03	5.23E+03	6.99E+04	5.12E+03	4.61E+04	6.72E+03	6.02E+04	4.59E+04	5.77E+04	4.64E+04	5.30E+03	6.12E+03	5.23E+03	5.23E+03
O	6.10E+03	8.91E+04	4.62E+03	3.42E+04	5.51E+03	7.15E+04	4.41E+04	6.18E+04	2.77E+04	5.47E+03	6.70E+04	4.32E+03	6.10E+03	8.91E+04	4.62E+03	3.42E+04	5.51E+03	7.15E+04	4.41E+04	6.18E+04	2.77E+04	5.47E+03	6.70E+04	4.32E+03	4.32E+03
P	5.62E+03	8.46E+04	4.81E+03	5.19E+04	5.82E+03	7.73E+04	6.82E+04	5.84E+04	5.14E+04	5.73E+03	7.67E+04	4.56E+03	5.62E+03	8.46E+04	4.81E+03	5.19E+04	5.82E+03	7.73E+04	6.82E+04	5.84E+04	5.14E+04	5.73E+03	7.67E+04	4.56E+03	4.56E+03

Click to Analyze Data Click to Clear Results

Kit Analysis
of Inhibitors
KPS

Ready KPS Single Dilution Results Statistical Analysis

70% 10:20 AM 5/8/2014

Single Dose SMART Protocol

Answer simple questions about the experiment: Strip kind, inhibitors, etc...

Single Dose SMART 2 Strip CV updating.xlsxm [Read-Only] - Microsoft Excel

File Home Insert Page Layout Formulas Data Review View Add-Ins Acrobat

Clipboard Font Alignment Number Styles Cells Editing

M17 99404

Kinase Selectivity Profiling Systems (KPS)
Single Dose SMART Protocol

Luminescence
Emission Filter: None
Integration Time(s): 0.3

Reading: 1

1 2 3 4 5 6 7 16 17 18 19 20 21 22 23 24

Single Dose Inhibitor

Columns 1-12
Enter the # of inhibitors tested in the box below.
Select the Kinase Profiling Strip used for the assay.

Columns 13-24
Enter the # of inhibitors tested in the box below.
Select the Kinase Profiling Strip used for the assay.

☐ Inhibitors are the same for both assays

Ok

Click to Analyze Data
Click to Clear Results

Kit Analysis
of Inhibitors
KPS

KPS Single Dilution Results Statistical Analysis

Ready 70% 10:22 AM 5/8/2014

Single Dose SMART Protocol

Data generated in a format that can be even published

TK-3

	Dasatinib	Tofacitinib	Erlotinib	Roscovitine	Compound	Compound	Compound	Compound	Compound	Compound
AXL1	47.63	8.35	99.78	119.21	119.28	114.05	120.41	116.15	103.28	73.24
EPHA1	27.50	11.83	161.43	104.16	109.66	114.32	109.07	141.73	120.61	26.69
FAK	87.50	73.65	44.68	110.49	137.48	112.72	136.74	100.65	106.58	93.72
ITK	50.28	106.74	102.87	107.36	99.50	118.84	107.04	107.31	114.03	31.73
JAK3	-0.69	98.93	97.53	100.81	89.28	97.79	98.70	92.65	106.92	-8.76
PYK2	8.23	82.79	88.58	90.62	54.89	91.75	93.37	6.07	91.80	-0.08
SYK	-0.69	55.28	2.14	85.77	59.84	78.48	46.42	-0.01	88.07	-0.80
TRKA	-1.81	45.88	-0.24	84.60	62.11	66.96	41.59	-0.32	81.46	-1.75

TK-1

	Dasatinib	Tofacitinib	Erlotinib	Roscovitine	Compound	Compound	Compound	Compound	Compound	Compound
EGFR	47.63	8.35	99.78	119.21	119.28	114.05	120.41	116.15	103.28	73.24
HER2	27.50	11.83	161.43	104.16	109.66	114.32	109.07	141.73	120.61	26.69
HER4	87.50	73.65	44.68	110.49	137.48	112.72	136.74	100.65	106.58	93.72
IGF1R	50.28	106.74	102.87	107.36	99.50	118.84	107.04	107.31	114.03	31.73
InsR	-0.69	98.93	97.53	100.81	89.28	97.79	98.70	92.65	106.92	-8.76
KDR	8.23	82.79	88.58	90.62	54.89	91.75	93.37	6.07	91.80	-0.08
PDGFRa	-0.69	55.28	2.14	85.77	59.84	78.48	46.42	-0.01	88.07	-0.80
PDGFRb	-1.81	45.88	-0.24	84.60	62.11	66.96	41.59	-0.32	81.46	-1.75

Legend:

- >60% Active
- 20-60% Active
- <20% Active

Dose Response SMART Protocol

Raw data is imported into the template. Click to analyze.

Answer simple questions about the experiment: Strip kind, inhibitors concentration...

Kinase Selectivity Profiling Systems (KPS)
Dose-Response SM

Full Plate Dose-Response

Strip 1

What compound was used for the dilution series?
Tofacitinib

Enter the final concentration of compound in Column 1 (in nM).
10000

Enter the dilution factor used. Ex. 20nM to 5nM would be a dilution factor of 4
4

Select the Kinase Profiling Strip used for the assay.
TK-3

Strip 2

What compound was used for the dilution series?
PF-477736

Enter the final concentration of compound in Column 1 (in nM).
10000

Enter the dilution factor used. Ex. 20nM to 5nM would be a dilution factor of 4
4

Select the Kinase Profiling Strip used for the assay.
TK-1
TK-2
TK-3
CMGC-1
CMGC-2

Ok

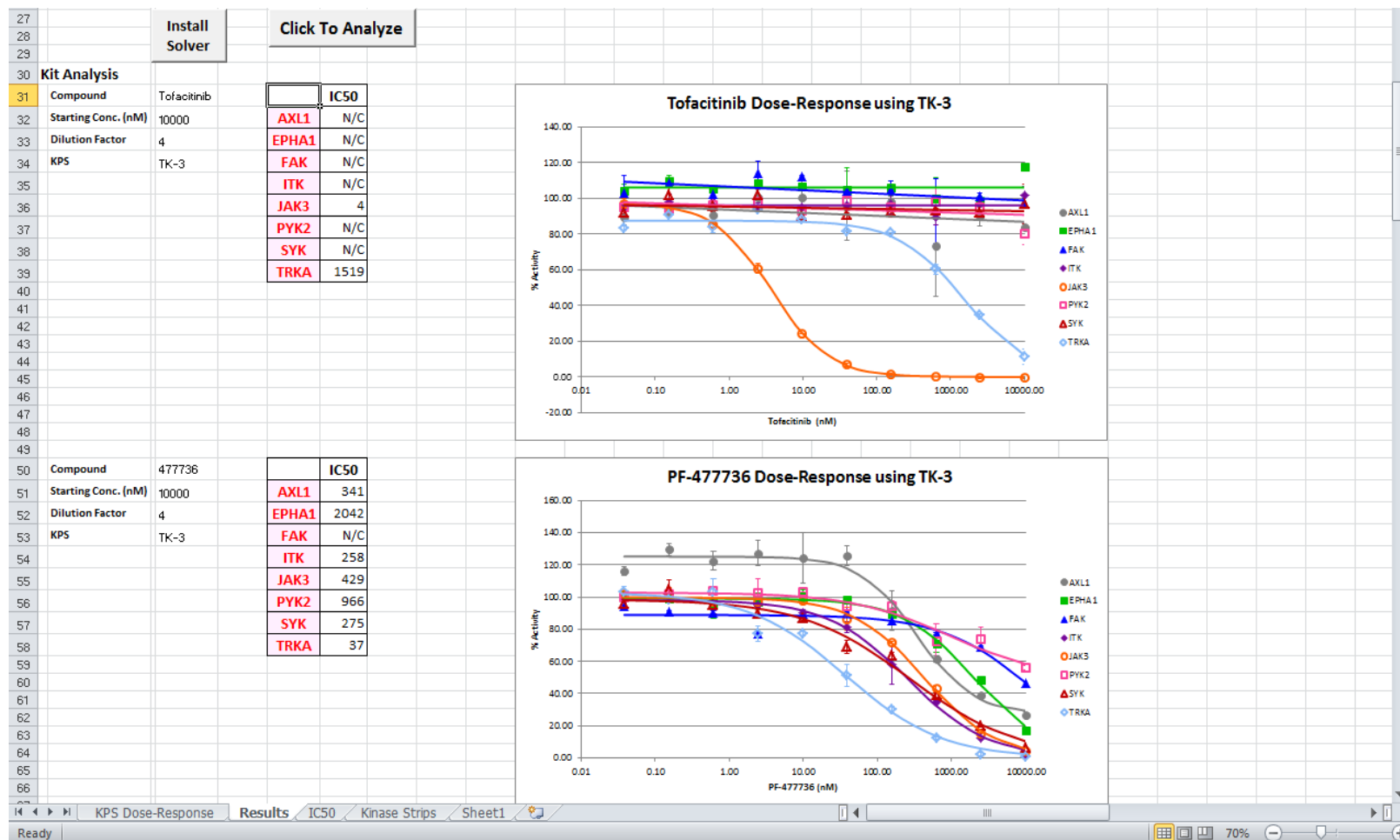
Install Solver

Kit Analysis
Compound Tofacitinib

Ready 10:41 AM 5/8/2014

Dose Response SMART Protocol

Data generated in a format that can be even published with accurate IC50 determination



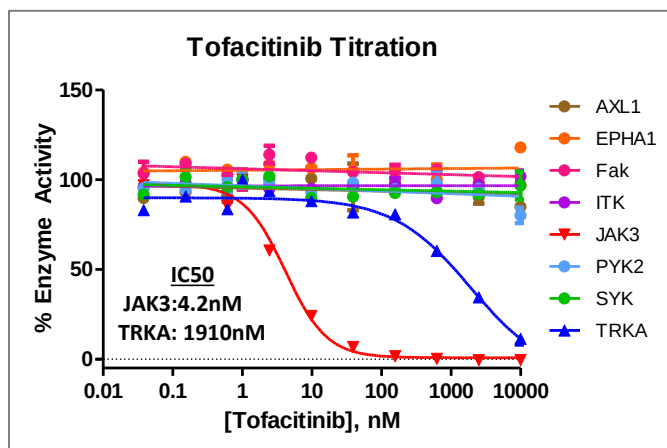
Flexible profiling: One Dose or Dose response

Single dose
Profiling
(Up to 8/strip)

	Kinase	Gefitinib	Dasatinib	Tofacitinib	SB203580	Roscovitine	PF-477736	Tozasertib	Enzastaurin
ST-3 TK family	AXL1	108	92	105	94	102	54	107	101
	EPHA1	84	1	102	79	97	63	77	100
	Fak	95	84	118	106	109	86	91	115
	ITK	102	101	99	94	100	33	24	99
	JAK3	99	113	0	87	94	36	88	103
	PYK2	133	134	122	106	134	120	80	116
	SYK	104	72	84	87	94	34	77	103
	TRKa	95	82	60	89	87	9	1	89

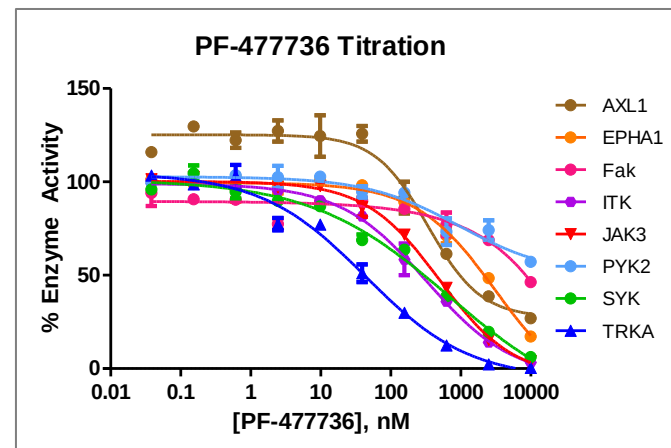
Inhibitor Concentration = 1 μ M

Selectivity



Dose response
Profiling
(Up to 2/strip)

Promiscuity



- Enabling flexible Kinase Profiling with the Strip Systems.
- Verified selectivity profiles against 8 kinases at once.

Large selectivity panels created with the profiling strips

64 Kinases with 8 compounds at a single dose profiling

ADP-Glo profiling systems

Strip #	Kinase	SB203580	Dasatinib	PF-477736	Gefitinib	Dasatinib	Tofacitinib	Roscovitine	Tozasertib
ST-1 TK Family	EGFR	99	104	2	2	94	102	94	103
	HER2	87	106	2	2	89	89	97	101
	HER4	96	104	2	2	101	101	101	96
	IGF1R	102	105	2	2	100	102	96	97
	InsR	98	100	95	97	94	97	101	88
	KDR	84	93	4	86	83	83	91	34
	PDGFRa	94	84	1	62	2	90	93	36
	PDGFRb	85	88	2	85	2	77	52	40
ST-2 TK Family	ABL1	98	102	27	97	106	107	98	102
	BRK	54	105	36	73	4	115	105	52
	BTX	99	99	33	83	3	101	111	46
	CSK	95	105	41	97	3	99	102	81
	FYN A	90	93	2	74	0	97	97	30
	LCK	90	88	22	74	0	89	82	27
	LYN B	95	91	28	75	0	97	96	78
	SRC	89	96	10	90	0	98	91	52
ST-3 TK Family	AXL1	94	101	54	108	97	105	102	107
	EPHA1	79	100	63	84	1	102	97	77
	Fak	106	115	88	95	84	106	109	91
	ITK	94	99	33	102	10	99	104	24
	JAK3	87	103	35	99	1	9	8	88
	PYK2	106	116	120	133	133	122	134	80
	SYK	87	103	34	104	72	81	94	77
	TRKa	89	89	9	95	82	60	87	1
ST-4 CMGC Family	CDK1/Cyclin A2	111	109	89	95	85	112	102	102
	CDK2/Cyclin A2	99	101	64	111	99	77	24	96
	CDK3/Cyclin E1	89	90	92	99	116	99	44	106
	CDK5/p25	97	98	75	89	104	10	26	99
	CDK5/p35	101	105	75	107	99	9	25	100
	CDK6/Cyclin D3	104	96	79	106	111	95	88	102
	CDK9/Cyclin K	110	72	83	100	91	149	41	110
	CLK3	90	98	87	106	92	98	98	92
ST-5 CMGC Family	ERK2	97	105	26	103	99	109	99	102
	GSK3b	55	8	86	86	103	105	96	102
	JNK1	85	107	95	89	99	104	107	97
	JNK3	30	106	88	81	99	93	106	88
	p38a	99	54	2	99	64	105	101	103
	p38b	10	103	103	94	54	100	101	105
	p38d	105	105	95	105	101	95	107	109
	p38g	94	104	98	112	101	103	104	107
ST-7 AGC Family	PKCa	96	4	77	88	107	106	103	102
	PKCb II	101	78	91	101	100	105	96	96
	PKCd	97	31	77	71	106	90	97	92
	PKCe	94	9	66	99	97	97	106	100
	PKCg	100	34	37	89	96	78	101	89
	PKCi	93	97	84	101	97	99	98	99
	PKCtheta	102	3	66	97	95	87	100	102
	PKCa	101	98	95	108	98	95	95	116
ST-8 CAMK Family	CHK1	97	97	0	95	96	105	100	101
	CHK2	100	106	8	80	98	107	109	90
	MAPKAPK2	101	106	100	91	98	101	106	98
	MARK1	93	96	2	97	97	78	103	97
	MELK	98	109	43	101	104	86	94	96
	PASK	98	99	90	101	111	93	96	92
	PIM1	108	70	65	96	112	99	96	104
	PKCmu (PKD1)	97	98	76	82	102	98	97	97
ST-9 Other/CK Family	Aurora A	99	108	4	102	95	80	105	4
	Aurora B	93	111	2	95	94	94	105	2
	CK2a1	94	115	86	103	108	105	105	99
	DNA-PK	102	91	85	106	104	99	113	115
	CK1a1	86	103	93	104	98	102	95	107
	CK1epsilon	58	95	96	100	105	96	86	107
	CK1g1	92	99	97	102	105	94	88	104
	VRK2	95	105	96	105	107	99	94	109

- Selective and promiscuous compounds are identified.
- Inhibitor potency compared to different members of the same family or different families in the kinome.

Inhibitors tested at 1µM

Inhibitors	Gefitinib	Dasatinib	Tofacitinib	SB203580	Roscovitine	PF-477736	Tozasertib	Enzastaurin
Kinase	EGFR	Abl/Src family	JAK	p38	CDK	CHK	Aurora	PKC

Validation of data created with ADP-Glo Profiling Systems

Comparison of ADP-Glo™ profiling to published radiometric data

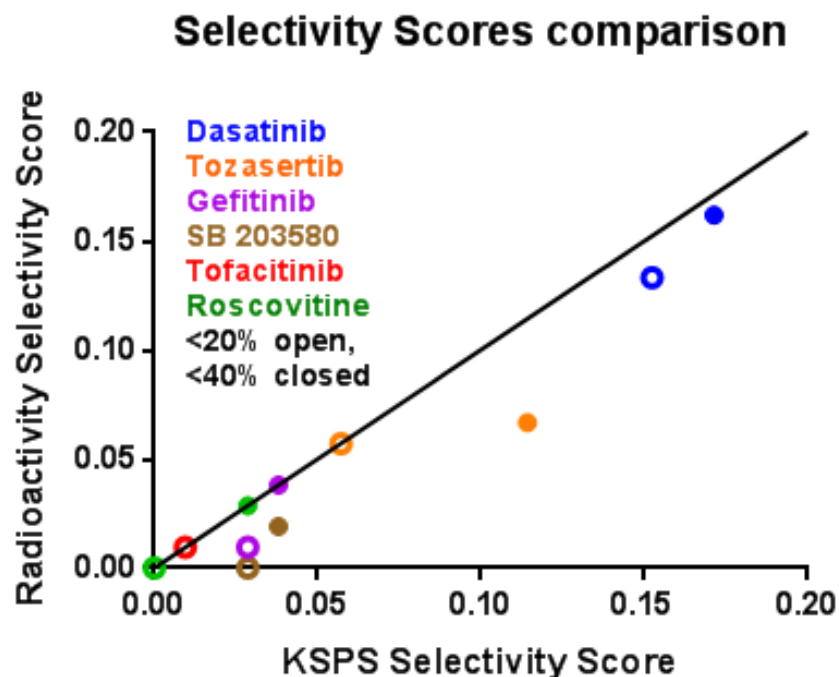
ADP-Glo™ profiling systems

Radiometric data¹

Strip #	Kinase	S8203580	Dasatinib	PF-47736	Gefitinib	Dasatinib	Tofacitinib	Resorcinol	Tozasertib	Dasatinib	Tofacitinib	Resorcinol	Tozasertib	Kinase
ST-1 TK Family	EGFR	99	104	43	2	64	102	94	103	21	96	97	78	EGFR
	HER2	87	106	102	88	106	89	97	101	80	109	103	102	HER2
	HER4	96	104	64	28	98	103	101	98	98	103	98	98	HER4
	IGF1R	102	105	98	103	102	102	96	97	97	101	98	44	IGF1R
	InsR	98	100	95	97	94	97	101	88	97	97	94	52	InsR
	KDR	84	93	4	86	83	83	91	34	103	106	97	85	KDR
	PDGFRa	94	84	1	62	2	90	93	36	2	104	98	71	PDGFRa
	PDGFRb	85	88	27	97	1	77	92	43	2	99	97	92	PDGFRb
ST-2 TK Family	ABL1	96	102	27	97	4	106	107	20	3	103	102	94	ABL1
	BRK	54	105	36	73	4	115	105	92	5	101	103	95	BRK
	BTk	99	99	33	83	5	101	111	46	2	99	83	99	BTk
	CSK	95	105	41	97	3	99	102	83	7	101	89	78	CSK
	FYN A	90	93	74	74	2	67	97	30	4	106	4	106	FYN A
	LCK	90	88	22	74	2	89	92	27	0	80	81	52	LCK
	LYN B	95	91	75	75	9	97	96	78	2	99	105	63	LYN B
	SRC	89	96	90	90	0	98	91	52	4	100	107	27	SRC
ST-3 TK Family	AXL1	94	101	54	108	92	105	102	107	61	98	100	44	AXL1
	EPHA1	79	100	63	84	102	97	77	98	103	103	96	100	EPHA1
	Fak	106	115	86	95	94	118	109	93	102	99	105	99	Fak
	ITK	94	99	33	102	101	99	100	24	102	99	105	99	ITK
	JAK3	87	103	35	99	113	4	94	88	93	2	97	92	JAK3
	PKY2	106	116	120	133	134	122	134	80	68	100	100	95	PKY2
	SYK	87	103	34	104	72	84	94	72	95	107	103	97	SYK
	TRKa	89	89	9	95	82	60	87	1	97	94	98	7	TRKa
ST-4 CMGC Family	CDK1/Cyclin A2	111	109	89	95	85	112	64	102	96	96	83	99	CDK1/Cyclin A2
	CDK2/Cyclin A2	99	101	64	111	99	77	24	96	92	86	24	88	CDK2/Cyclin A2
	CDK3/Cyclin E1	89	90	92	99	116	99	44	106	92	92	58	95	CDK3/Cyclin E1
	CDK5/p25	97	98	75	97	104	100	26	104	100	100	101	91	CDK5/p25
	CDK5/p35	101	105	75	107	99	96	25	100	100	99	91	100	CDK5/p35
	CDK6/Cyclin D3	104	96	79	106	111	93	84	102	103	111	96	121	CDK6/Cyclin D3
	CDK9/Cyclin K	110	72	83	100	83	104	41	110	84	100	79	93	CDK9/Cyclin K
	CLK3	90	98	87	106	92	98	98	92	106	105	105	110	CLK3
	ERK2	97	105	101	108	99	109	99	102	93	91	98	91	ERK2
	GSK3b	55	4	86	86	103	105	96	102	97	100	96	104	GSK3b
	JNK1	84	107	95	89	99	104	107	97	106	108	98	95	JNK1
	JNK3	30	106	86	81	95	93	106	88	100	100	101	107	JNK3
ST-5 CMGC Family	p38a	64	102	99	64	24	105	101	103	37	93	96	97	p38a
	p38b	10	103	103	103	101	105	103	103	77	103	103	90	p38b
	p38d	105	105	95	105	105	105	107	109	99	94	97	80	p38d
	p38g	94	104	98	112	101	103	104	107	100	97	95	96	p38g
	PKCa	96	5	77	88	107	106	103	102	103	79	87	82	PKCa
	PKCb II	105	4	75	91	106	100	105	95	89	88	96	90	PKCb II
	PKCd	97	31	77	73	106	90	97	92	100	75	102	98	PKCd
	PKCe	94	9	66	99	97	97	106	100	99	98	98	98	PKCe
	PKCg	100	14	37	89	96	78	101	89	86	91	96	94	PKCg
	PKCi	93	97	84	101	97	99	98	99	101	103	92	99	PKCi
	PKCtheta	102	8	66	97	95	87	100	102	97	89	90	86	PKCtheta
	PKCz	102	92	95	108	95	96	105	108	101	92	100	76	PKCz
ST-6 CAMK Family	CHK1	97	97	97	99	96	105	103	94	96	94	94	101	CHK1
	CHK2	100	106	8	80	98	107	109	90	103	104	97	100	CHK2
	MAPKAPK2	103	106	100	91	98	101	106	98	100	97	101	101	MAPKAPK2
	MARK1	93	96	6	97	97	78	103	97	107	119	115	114	MARK1
	MELK	98	109	43	101	106	86	94	92	81	79	85	58	MELK
	PASK	98	99	90	101	112	105	91	96	110	105	93	101	PASK
	PIM1	108	70	65	96	112	99	96	104	102	102	105	99	PIM1
	PKCmu (PKD1)	97	98	76	82	102	98	98	97	97	97	99	102	PKCmu (PKD1)
	Aurora A	99	108	4	102	95	80	105	4	93	84	89	4	Aurora A
	Aurora B	93	111	9	95	94	80	105	4	93	99	95	9	Aurora B
	CK2a1	115	115	100	108	105	104	105	95	N/A	N/A	N/A	114	CK2a1
	ST-9 Other/CK Family	DNA-PK	102	91	85	106	104	99	113	115	N/A	N/A	N/A	N/A
CK1a1		86	103	93	104	98	102	95	107	N/A	N/A	91	97	CK1a1
CK1epsilon		58	95	96	100	105	96	86	107	90	101	81	99	CK1epsilon
CK1g1		92	99	97	102	105	96	88	104	97	101	94	99	CK1g1

Validation of data created with ADP-Glo Profiling Systems

Comparison of Selectivity profiling Scores* between ADP-Glo™ and radiometric data
(Over 100 kinases compared)

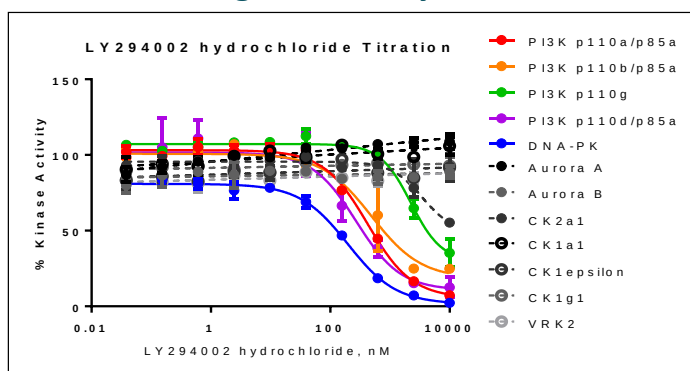


Selectivity Data generated with ADP-Glo™ platform correlates perfectly with radioactivity-based selectivity data.

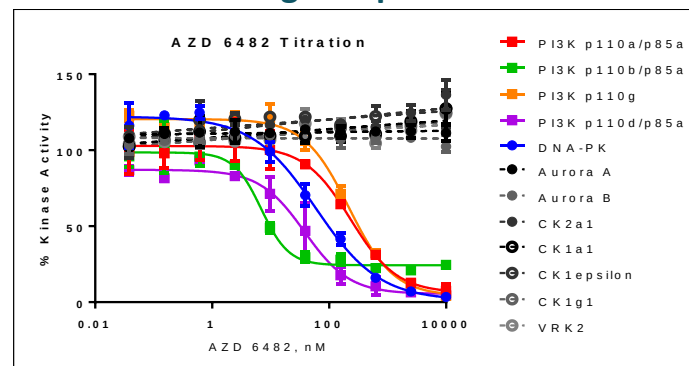
*Selectivity scores: # of kinases with % activity left < 20 or 40% / # of kinases tested

Simultaneous profiling compounds against protein and lipid kinases.

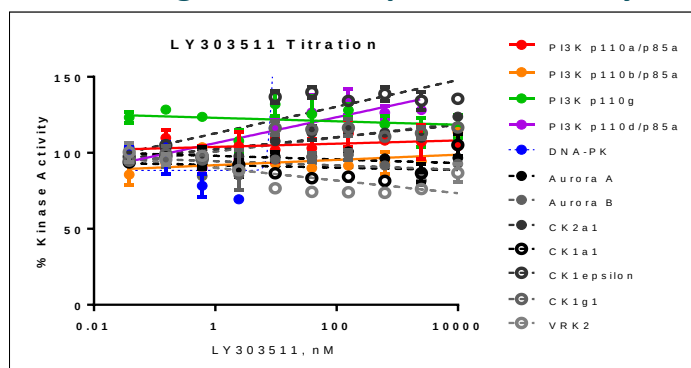
Profiling PI3K family inhibitor



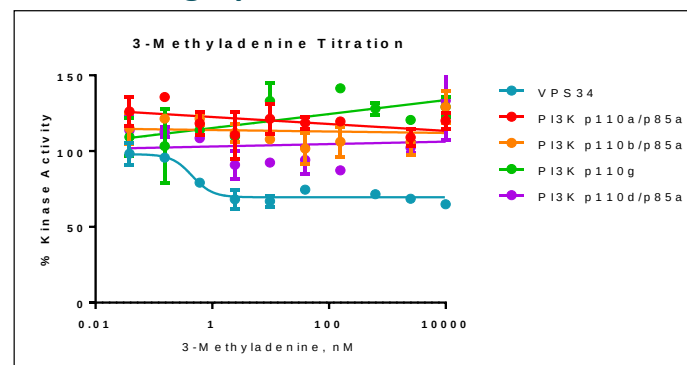
Profiling PI3Kβ inhibitor



Testing control compound inactivity



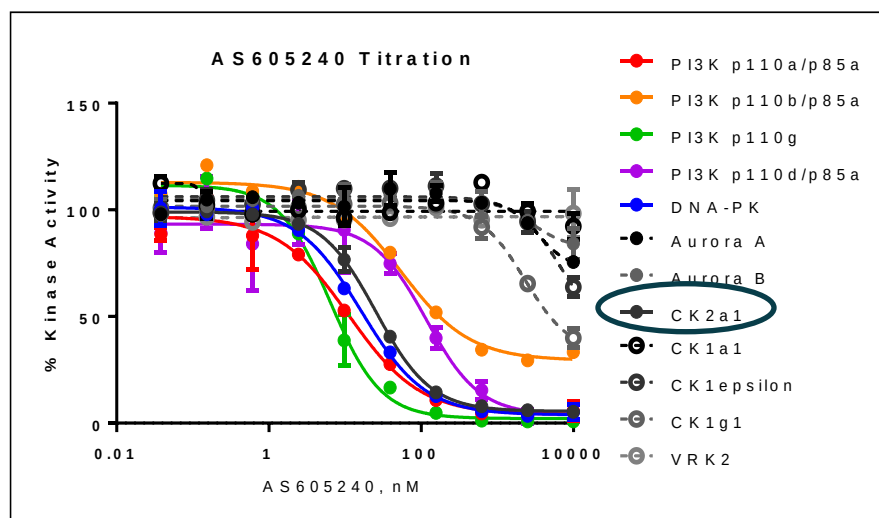
Profiling Lipid Kinase Class III inhibitor



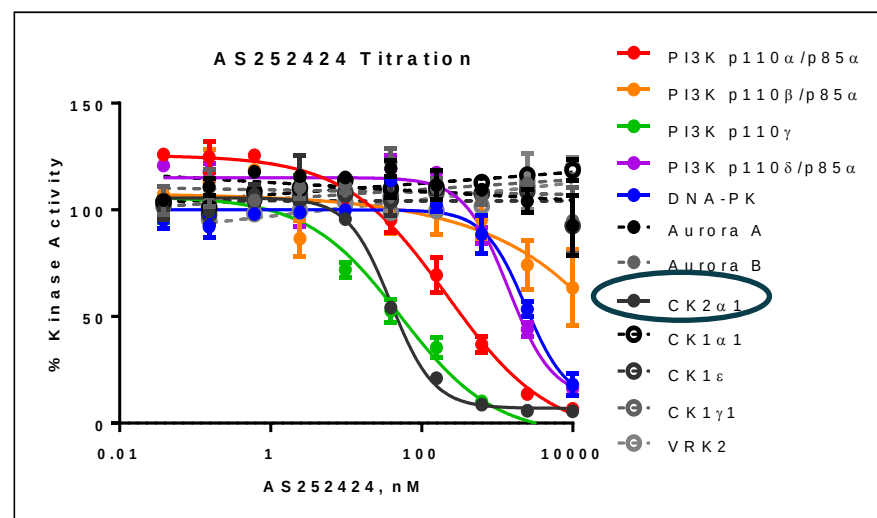
Profiling kinases in strips to identify compounds' selectivity towards members of PI3 Kinase family.

Simultaneous profiling compounds against protein and lipid kinases.

Profiling PI3Ky inhibitor



Profiling PI3Ky inhibitor



Simultaneous Profiling of Protein and Lipid kinases identifies promiscuity of PI3K specific inhibitor towards other protein kinases such as CK2α1.

Advantages of Kinase profiling strips

Kinase Selectivity Profiling Strip Systems have the following advantages:

- ✓ **Fast and simple reaction assembly:** Two quick dilutions provide working stocks of kinases and substrate/co-factor solutions sufficient for 25 kinase reactions.
- ✓ **One-time use design:** Eliminating multiple freeze/thaw cycles ensures optimal kinase activity per experiment.
- ✓ **Optimized kinase activity for inhibitor profiling:** All kinases have been optimized to provide optimal ADP production when assayed at 10 μ M ATP.
- ✓ **Formatted strips provide access to eight kinases at a time:** Kinase from singular kinase families are grouped together for a more relevant selectivity profiles.
- ✓ **Flexible kinase inhibitor profiling:** Each strip has enough material to profile 4 compounds at a single dose or create a dose response for 1 compound against the 8 kinases at once.

Advantages of Kinase profiling strips

ADP-Glo™ Kinase Selectivity Profiling Systems generate data similar to the “gold standard” Radiometric assay

“Now even chemists can do kinase profiling?”



General features of ADP-Glo™ Kinase platform

ADP-Glo™ platform offers so many positive attributes that make it ideal for all stages of drug discovery

Basic Research, Primary and secondary screenings and for profiling of lead compounds.



Summary

ADP-Glo™ Kinase Platform



1. ADP-Glo Kinase Assay

Kinase reactions up to 1mM ATP



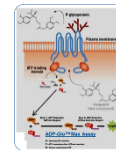
2. Kinase Enzyme Systems

174 complete kinase assays

Kinase
ERK2
GSK3b
JNK1
JNK3
p38a
p38b
p38d
p38g

3. Kinase profiling Systems

8 well kinase strips for easy profiling



4. ADP-Glo Max Assay

*ATPase reactions up to 5mM ATP
(ABC transporters,...)*



Streamlining Kinase Profiling Workflow by Integrating Bioluminescent Kinase Strips with Multimode Detection System

Hicham Zegzouti, Ph.D.

Assay Design Group, Promega, Madison WI

SLAS 2015 Tutorial

February 10, 2015