

# Promega Corporation

***ADP Detection Platform for  
kinase inhibitor screening, mode  
of action studies and profiling***

Hicham Zegzouti, Ph.D.

May 2013

***Hicham.zegzouti@promega.com***

# ***Kinases orchestrate complex biological processes***

Kinases play a critical role in Human biology

- Important components of Cell Signal transduction
- Regulation of many cellular processes through phosphorylation of diverse substrates (proteins (S/T, Y), Lipids, Sugars...)



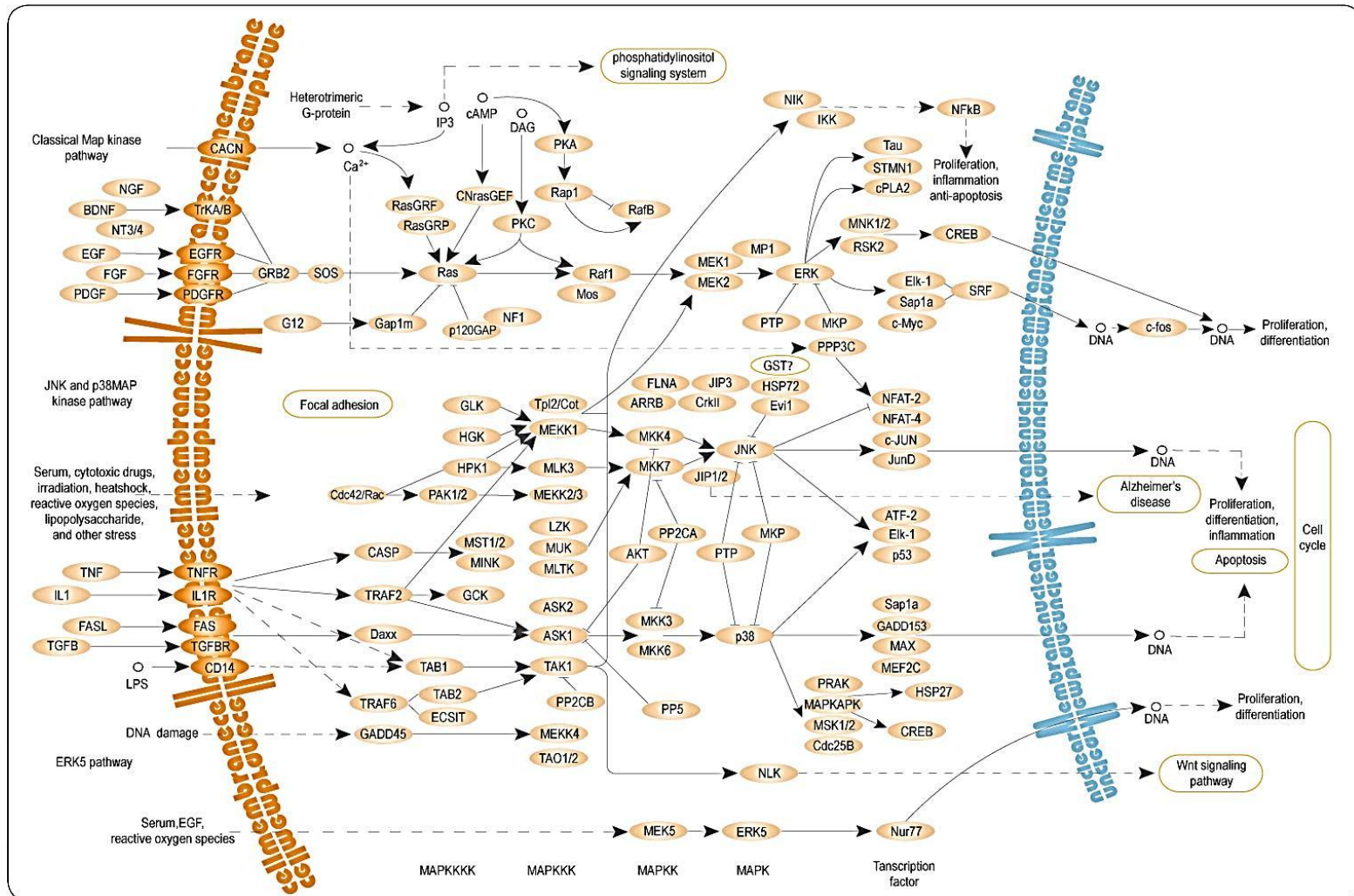
- ~518 protein Kinase in human genome (388 S/T, 90 Tyr, 40 atypical) & more
- More than third of all human proteins are phosphorylated





# MAP Kinase pathway example

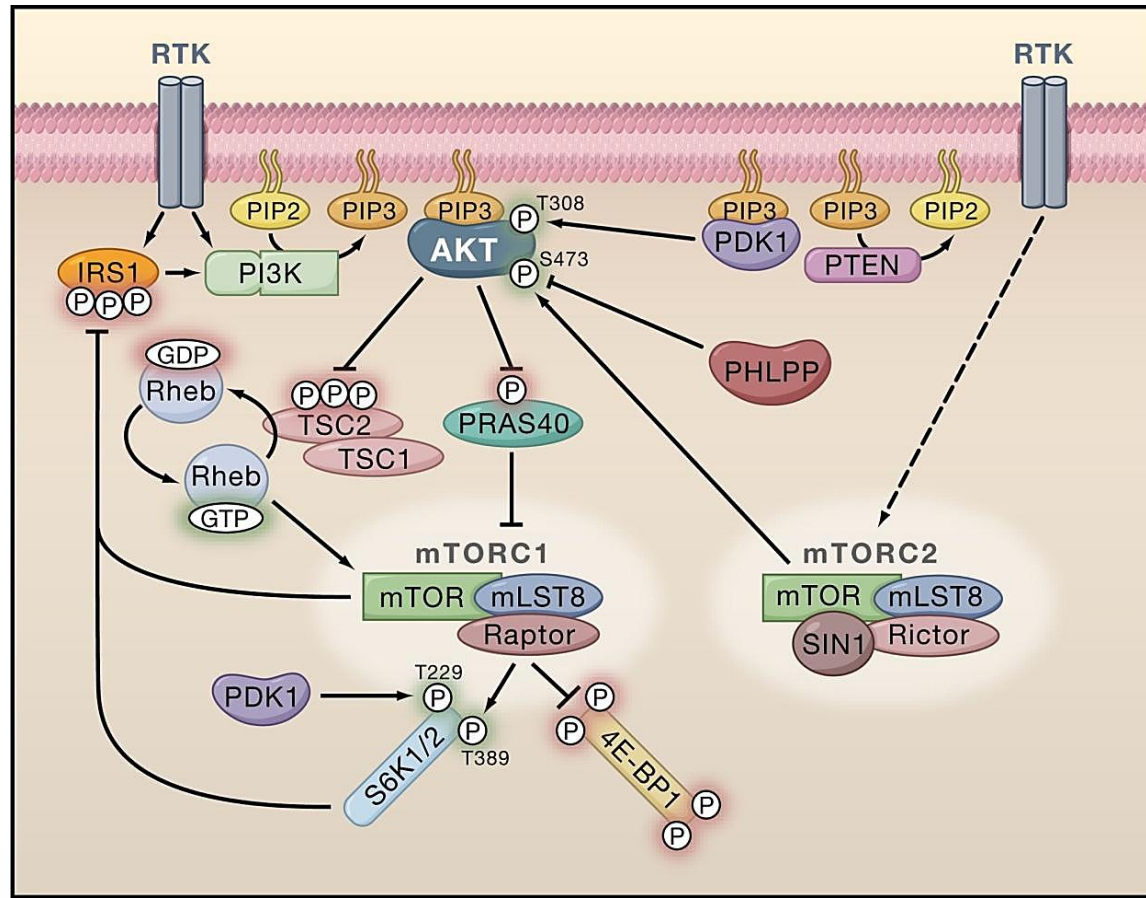
Signal Transduction (Phosphorylation cascades) from the receptors to the transcription factors





# Akt-Lipid Kinase Signaling example

## Cross talk between Signal Transduction pathways



Regulation of:

- cell survival
- Proliferation
- insulin-dependent metabolic responses.

Manning and Cantley (2007) *Cell* **129**, 1261-1274.

# ***Kinases orchestrate complex biological processes***

**Kinases play a critical role in Human biology**

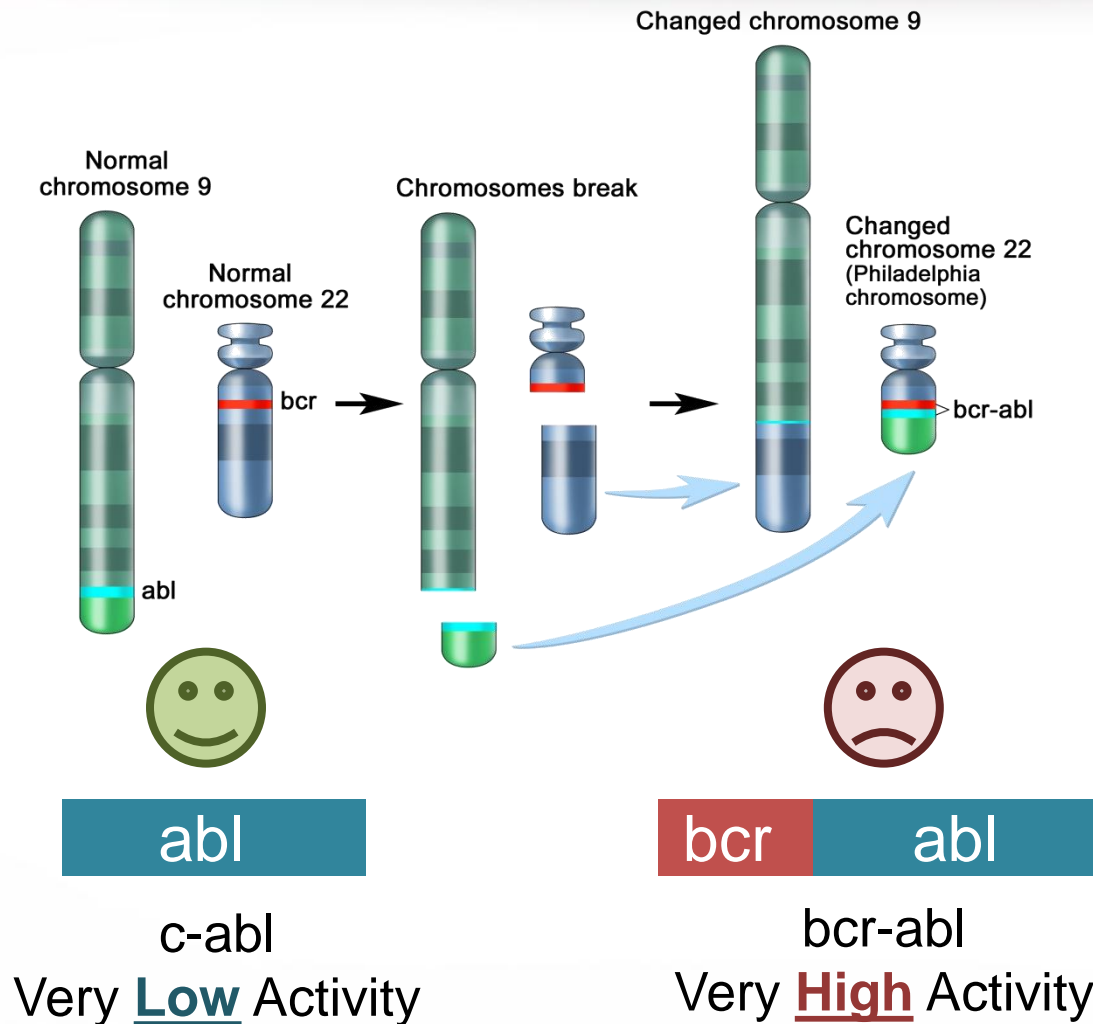
**Kinases are directly involved in many diseases**

- **Under pathological conditions Kinases can be deregulated/mutated**
- **Alteration in phosphorylation states results in abnormalities**
- **Over 400 Human Diseases are Linked to Defects in Kinases- (and Ppases) Dependent Signaling Pathways**



**Kinases became the largest target in the drug discovery market**

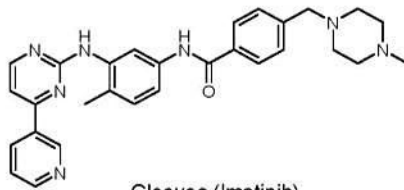
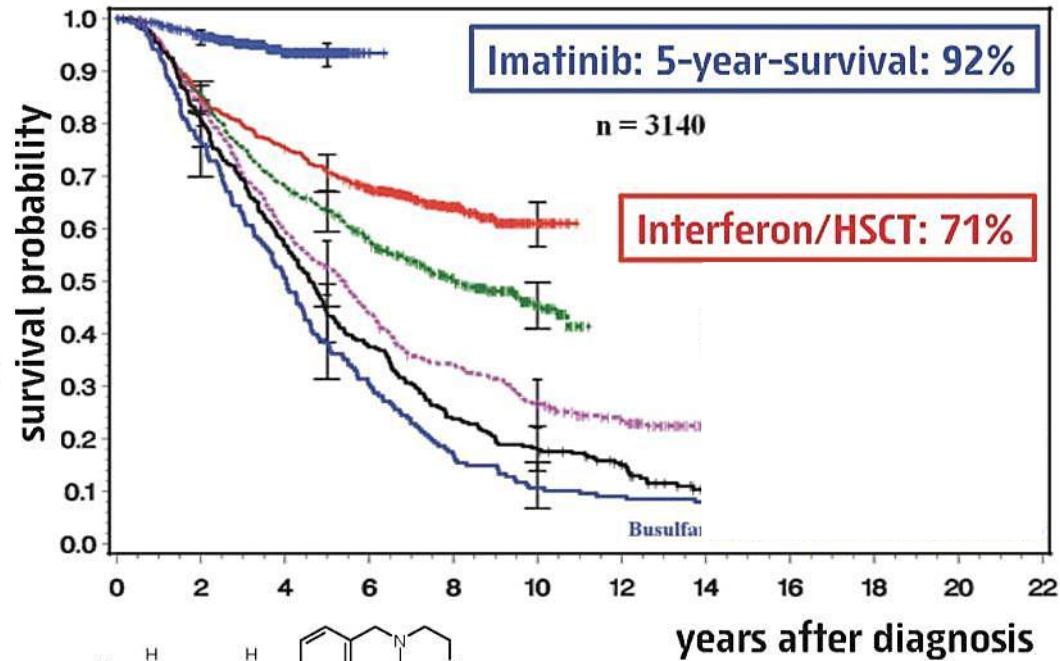
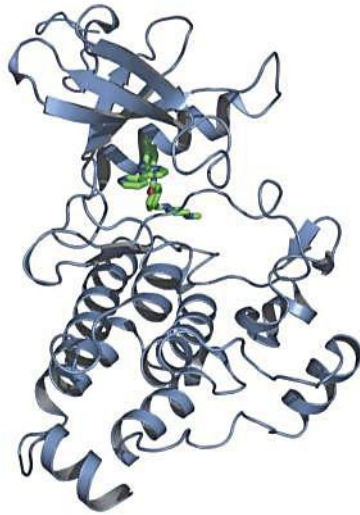
# ***bcr-abl: An oncogene that started it all***



- First chromosomal translocation identified
- Deregulated constitutively active Tyrosine Kinase
- Expressed in Chronic Myeloid Leukemia (CML)
- Expression sufficient to cause CML



# Imatinib (Gleevec): Improved Long-term survival



Gleevec (Imatinib)

Gleevec inhibits the DFG out state of the Kinase (Inactive form)

DFG (Asp-Phe-Gly): Mg-ATP binding site

# Importance of Kinases in Drug discovery

## US Marketed Small Molecule Kinase Inhibitors

**Table 1. Targeted Therapeutics in Cancer.\***

Gene	Genetic Alteration	Tumor Type	Therapeutic Agent
<b>Receptor tyrosine kinase</b>			
<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
<b>Non-receptor tyrosine kinase</b>			
<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
<b>Serine-threonine-lipid kinase</b>			
<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

• Kinases

• Disease  
Neuro

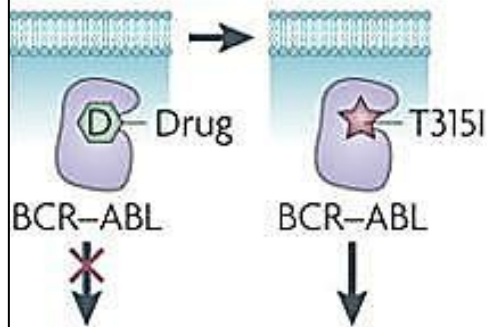
market

ascular,

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

# ***Tumors become resistant to Kinase inhibitors***

**a** Mutation that blocks drug binding



**Resistance mechanisms:**

- 1. Mutation in kinase target**
- 2. Induction of Bypass mechanism**
- 3. Activation in up/down-stream effectors**

➤ **Patient relapse requires second-line therapies:**

- **Second generation of inhibitors needed**
- **Therapeutic combinations of multiple kinase targets**



# ***Combinatorial therapy as a new trend for recalcitrant cancer treatment***



## **HIV treatment example:**

**Elvitegravir  
(integrase inhibitor)**

**Cobicistat  
(drug metabolism inhibitor)**

**Emtricitabine and Tenofovir  
(nucleoside RT inhibitors)**

**vs.**

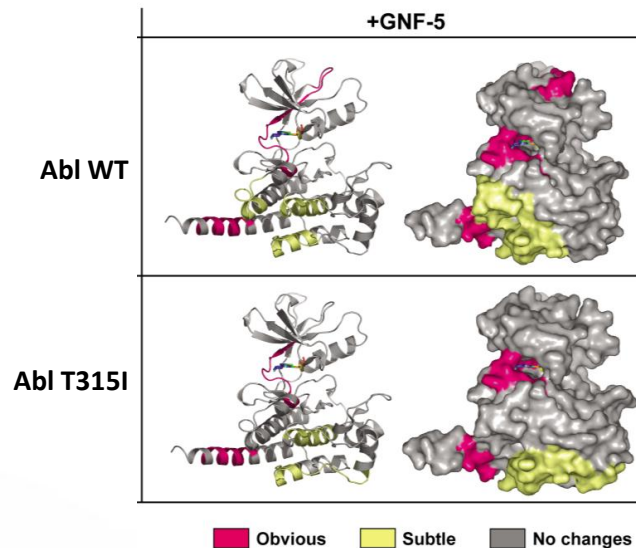
## **Cancer treatment example:**

**Kinase inhibitors are used  
as a single therapy**

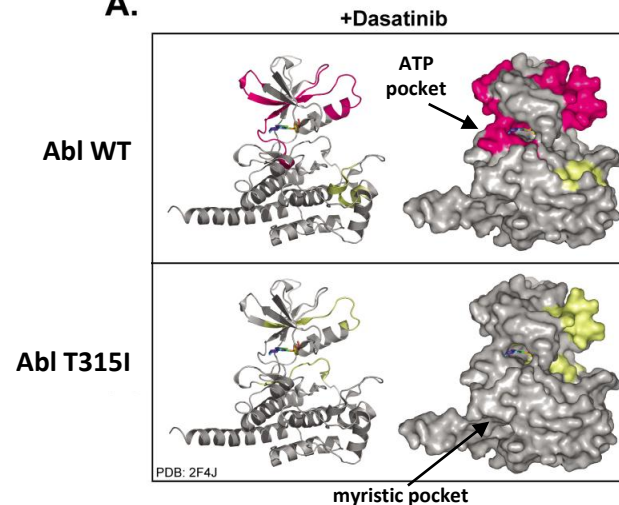
# Combinatorial therapy as a new trend for recalcitrant cancer treatment (Abl example)

**in combination with:** Allosteric inhibitors  
binding to non ATP binding sites:

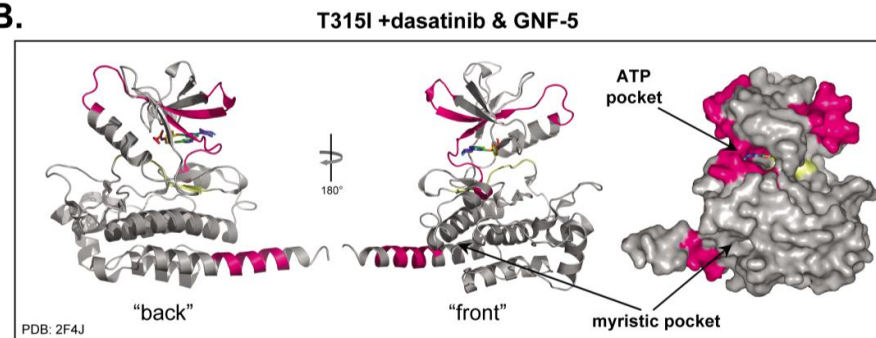
- PIF pocket
- Substrate pocket
- **Myristic pocket**



## A. ATP site inhibitors



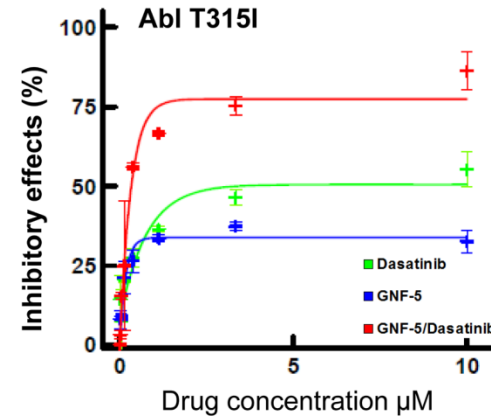
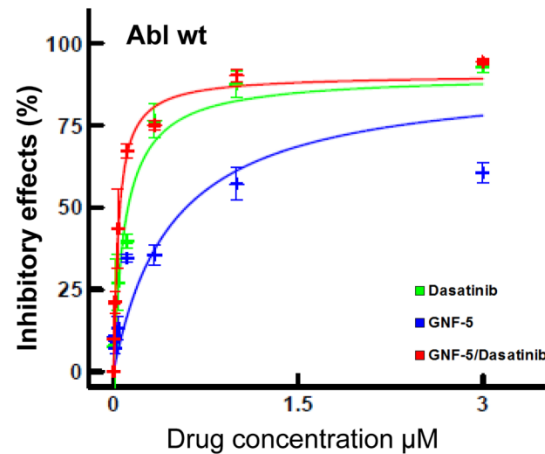
## B.



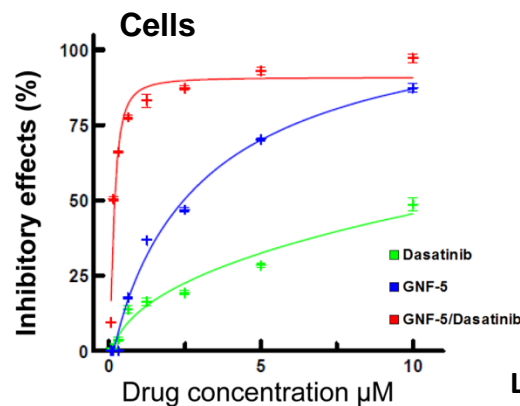
Simultaneous binding of Dasatinib and GNF-5 to T315I caused conformational changes in Abl such that dasatinib binding to T315I became similar as when it is bound to WT Abl.

# Combinatorial therapy as a new trend for cancer treatment

**A.**



**B.**



(Similar effect with imatinib and nilotinib)

Lacob et al., PLoS One. 2011; 6(1): e15929

**Additive interaction between GNF-5 and the ATP-competitive inhibitor Dasatinib to inhibit proliferation of Bcr-Abl T315I mutant**



# *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

Two way paradigm

Selective compounds



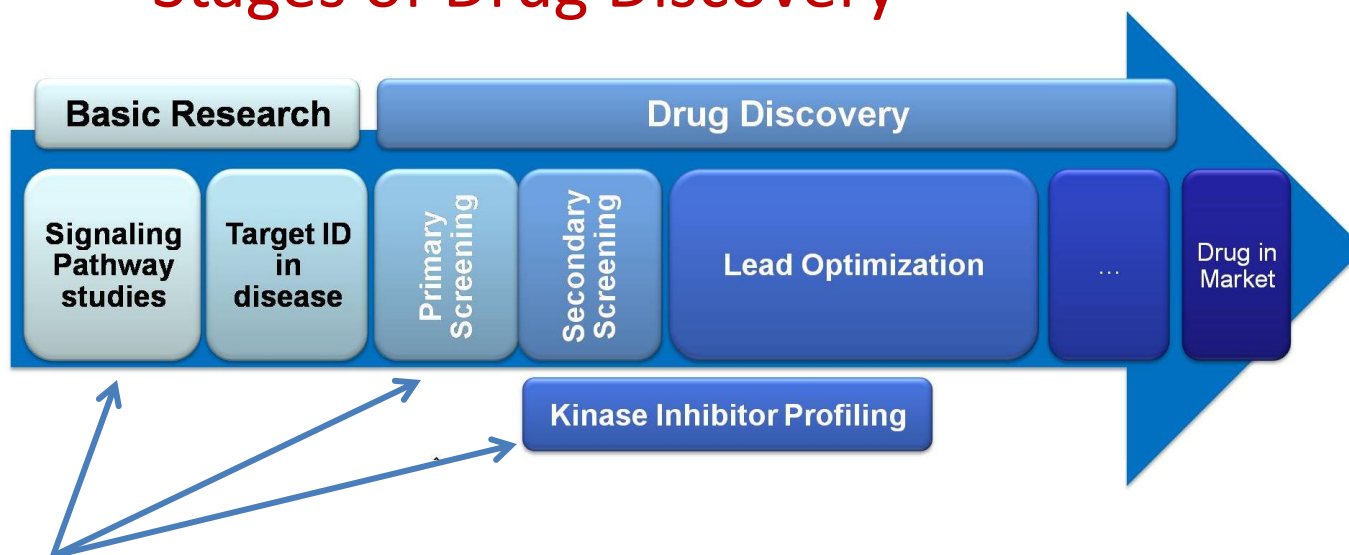
Potency

Choosing the right assay for success

# ***Kinase Studies in Basic Research and Drug Discovery***

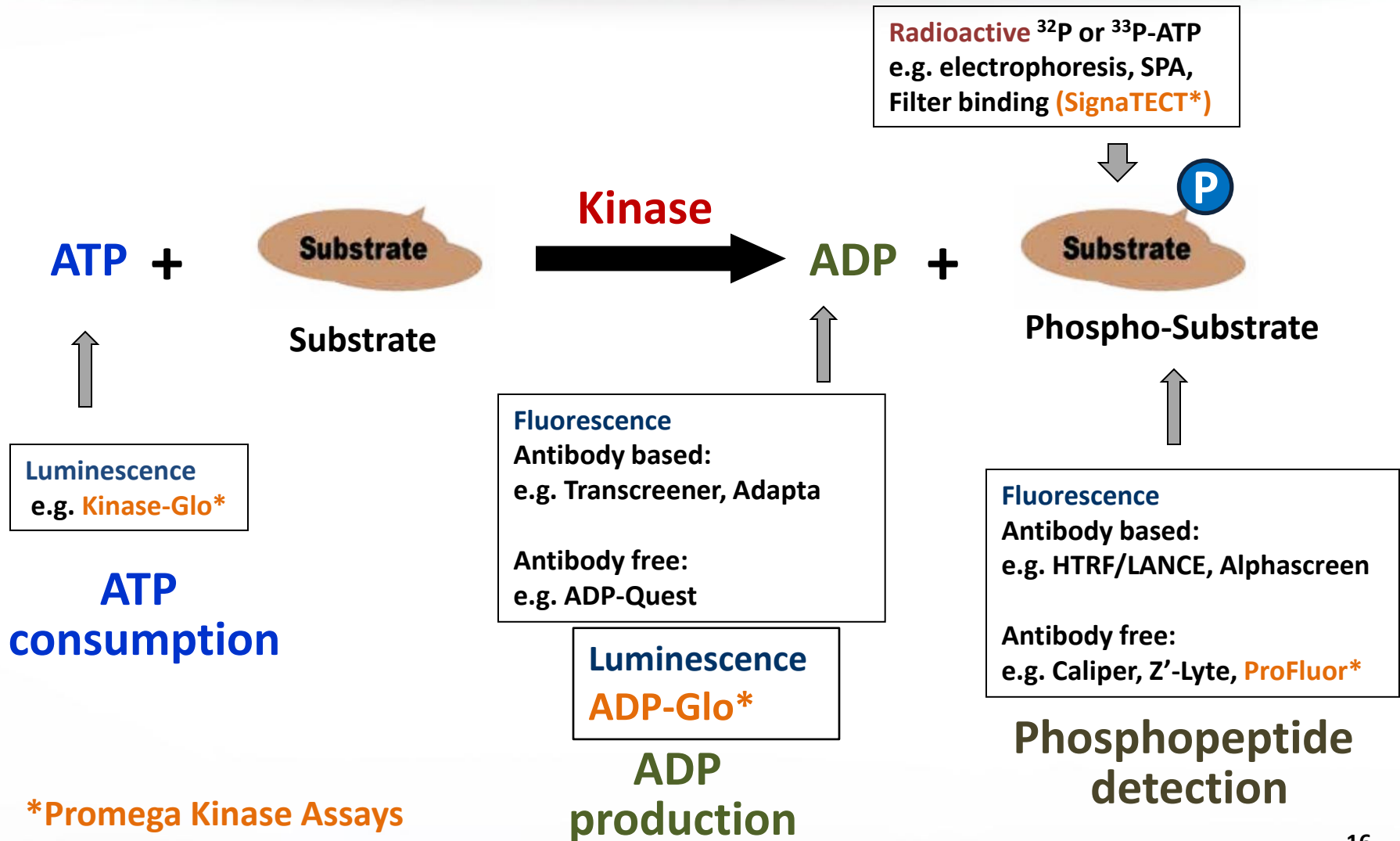
## **Biochemical Kinase Assays**

### **Stages of Drug Discovery**



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

# Detection of Kinase activity



# ***Current Kinase Assays features and drawbacks***

## **Drawbacks of Current Assays:**

- If not radioactive they require specific antibodies
- Require fluorescently-labeled peptides
- False Hits with Fluorescence based assays
- Very expensive
- Require special detection technology
- Not tolerant to High concentrations of ATP

## **Features of an Ideal Assay:**

- ✓ Homogeneous, Nonradioactive, Robust
- ✓ Universal (any enzyme substrate combinations)
- ✓ Applicable for different kinases and diverse substrates
- ✓ Use multiphosphorylated substrates
- ✓ Minimal False Hits
- ✓ Distinguish between ATP competitive and noncompetitive inhibitors

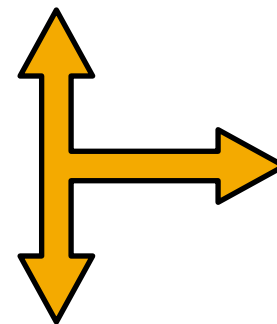


## *Kinase Research: parts of the equation*

➤ Kinase families (targets)

➤ Kinase Assays (Glo)

➤ ~~Pharmaceuticals~~ Kinase Inhibitors (Drugs)

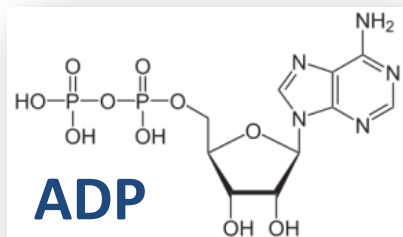
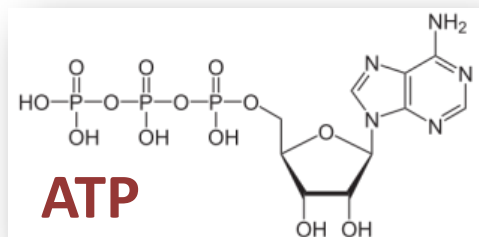


# ***Bioluminescent Kinase Assay Platform***

## **Kinase Glo<sup>®</sup> Assay**

### **Monitoring ATP Depletion**

- Kinase Glo<sup>®</sup> (10μM ATP)
- Kinase Glo<sup>®</sup> Plus (100μM ATP)
- Kinase Glo<sup>®</sup> Max (500μM ATP)

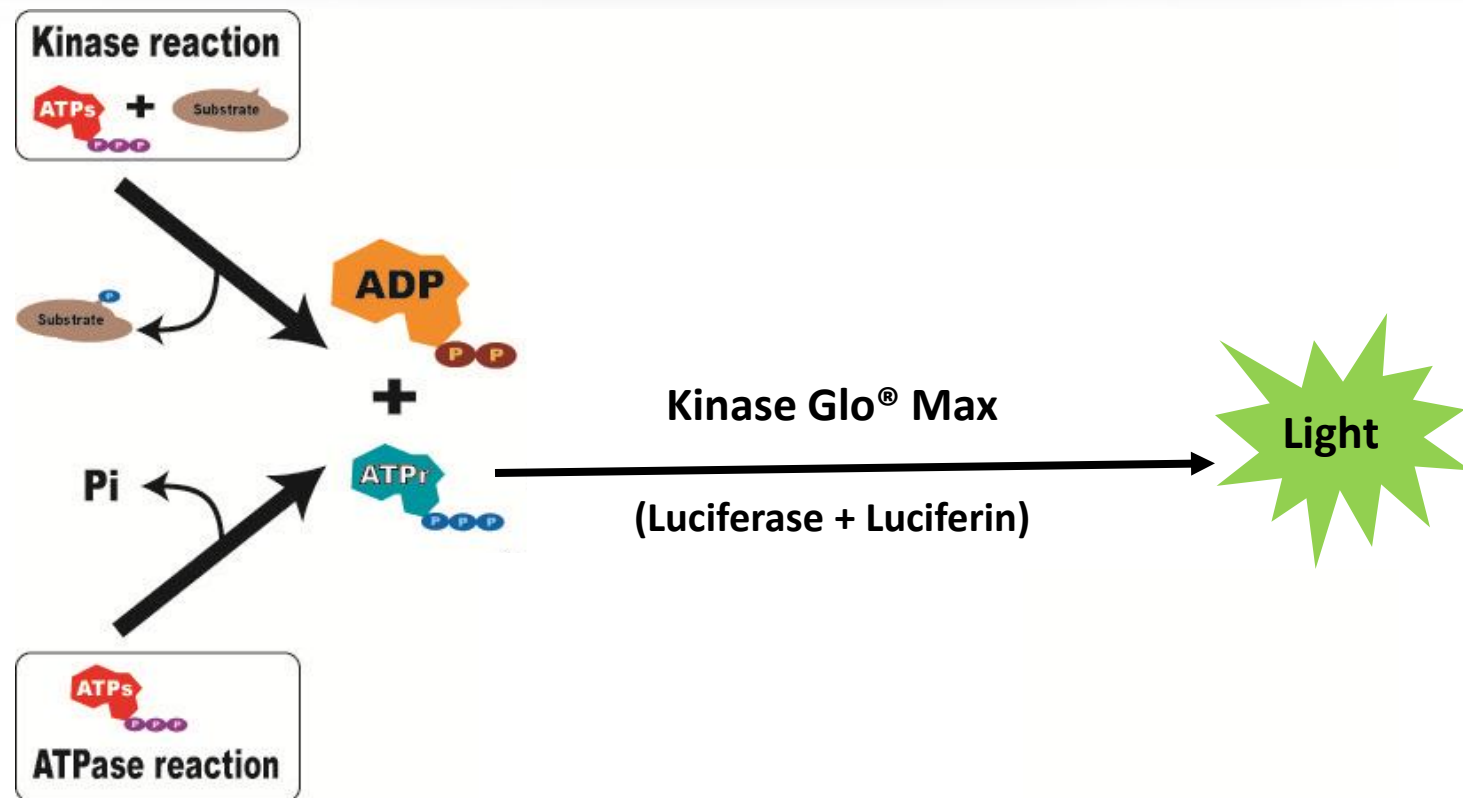


## **ADP Glo<sup>™</sup> Assay**

### **Monitoring ADP Production**

**Micro to Millimolar  
ATP Concentration**

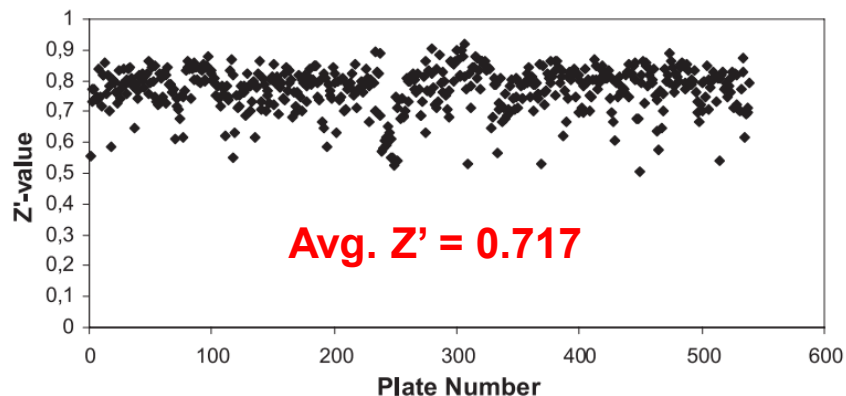
# ATP detection: Kinase-Glo<sup>®</sup> Platform



Light output is correlated with the amount of **ATP** remaining and is inversely correlated with the amount of kinase activity

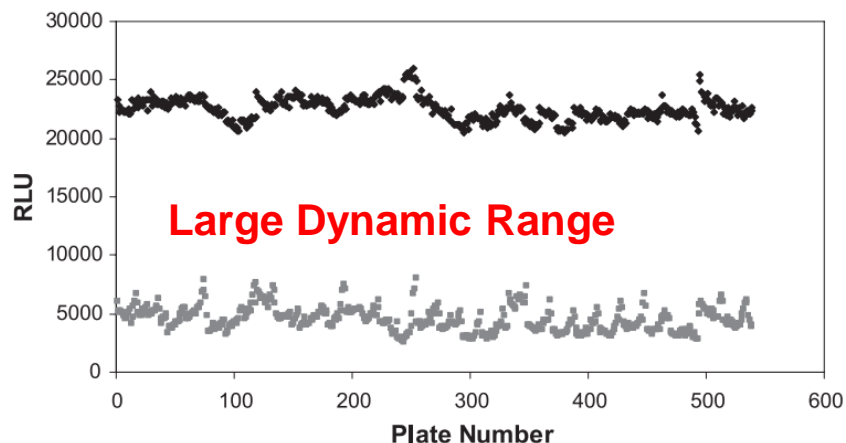
# Sensitivity and Robustness of Kinase-Glo<sup>®</sup>

## Z' Values calculated for Test Plates



- ✓ Kinase-Glo<sup>®</sup> Assay generates a high signal to background ratios
- ✓ Kinase-Glo<sup>®</sup> Assay is a robust assay ideal for HTS

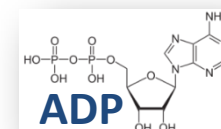
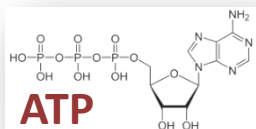
## Averages for controls (+ and -)



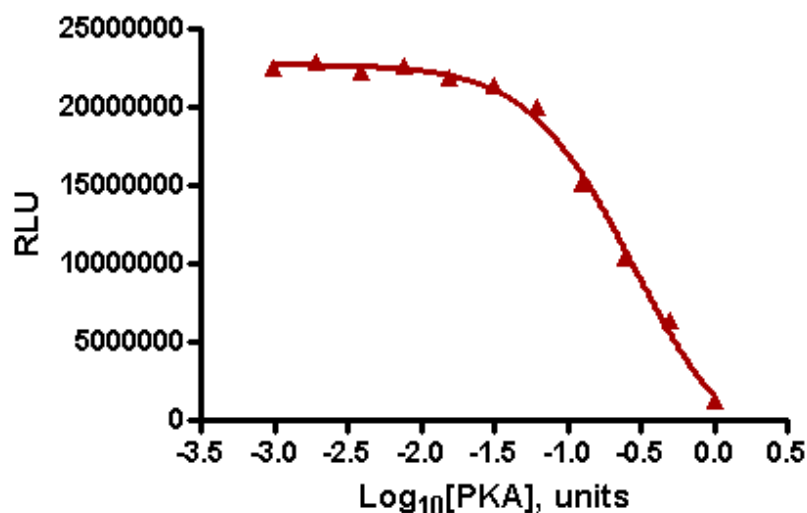
Baki A, *et al* (2007)  
*Assay and Drug Dev. Technol.* **5**: 75-83



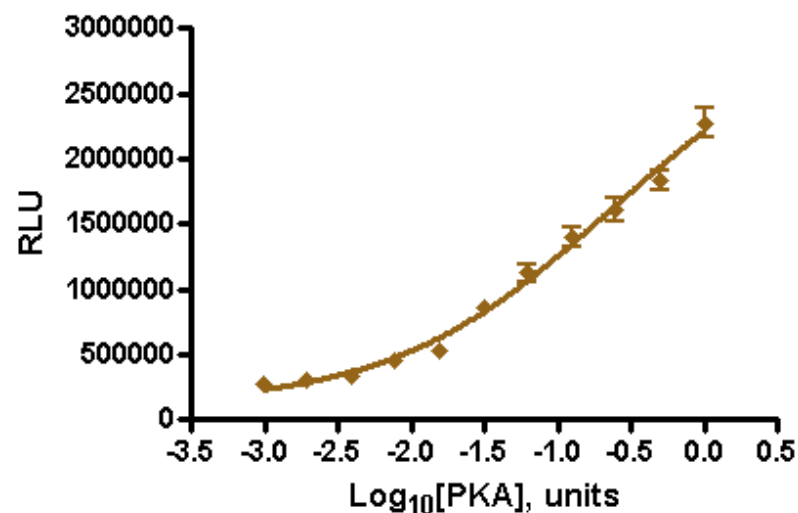
# Luminescent Kinase assays: Two Different Approaches



▲ Kinase Glo<sup>®</sup> Max

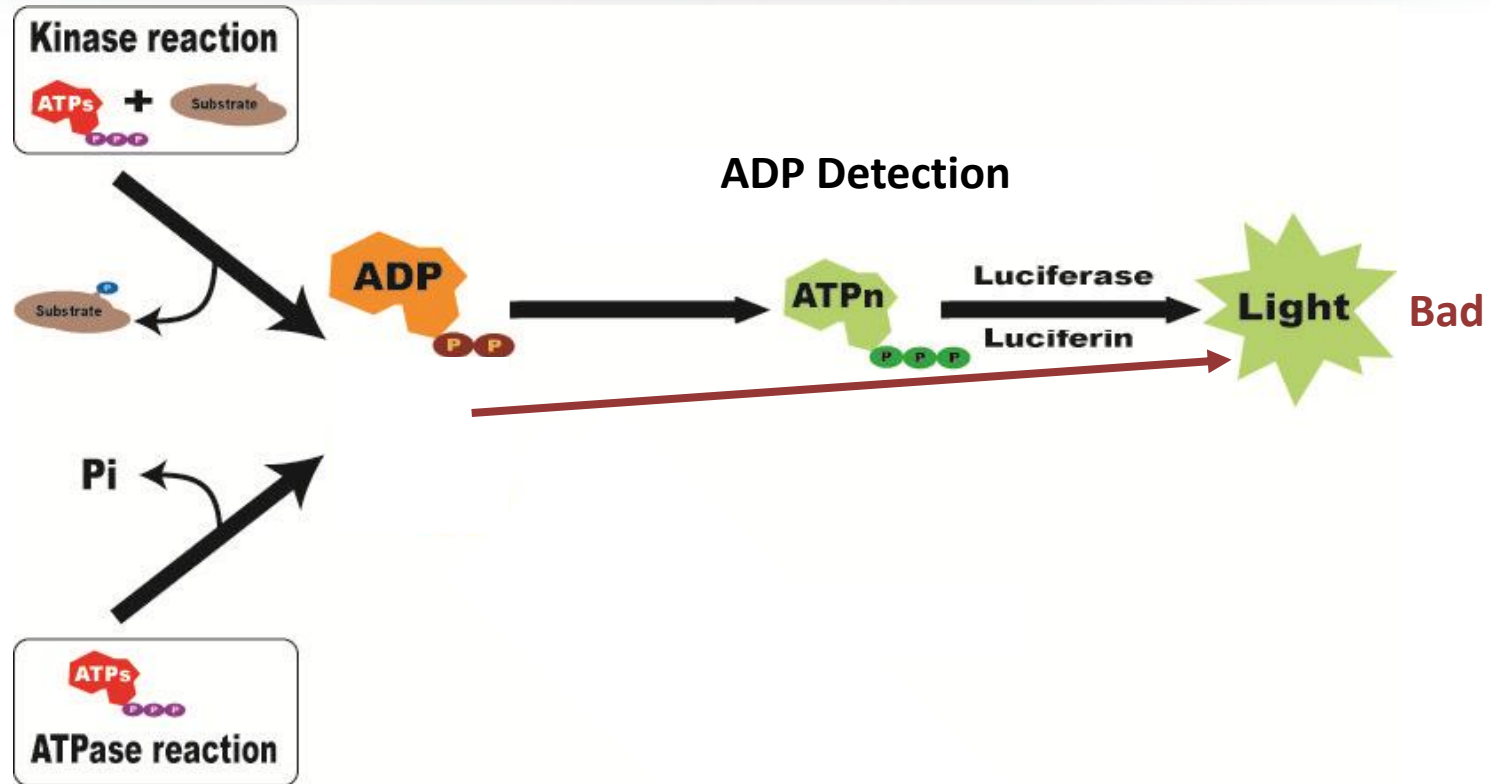





◆ ADP-Glo<sup>™</sup>



- Kinase-Glo<sup>®</sup> measures the remaining **ATP** after a kinase Reaction
- ADP-Glo<sup>™</sup> measures **ADP** produced in a kinase Reaction

# Luminescent ADP detection assay requirements



 : Starting ATP amount  
 : ATP remaining after kinase reaction  
 : Newly synthesized ATP

# ADP-Glo™ Assay Format- 1:1:2

**384-well plate**

**5µl kinase reaction**

+

**5µl ADP-Glo™ Reagent**

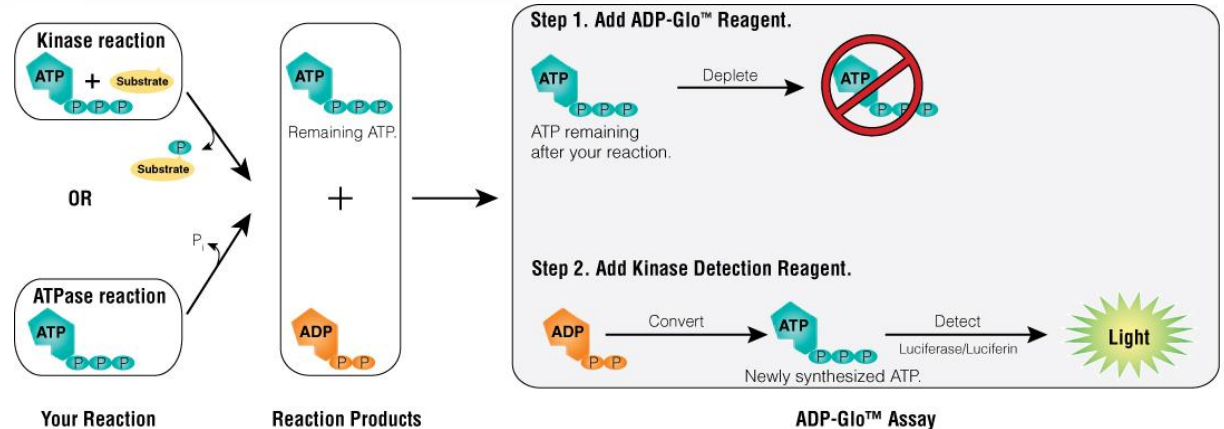
**40 min. Incubation**

+

**10µl Kinase Detection Reagent**

**30-60 min. Incubation**

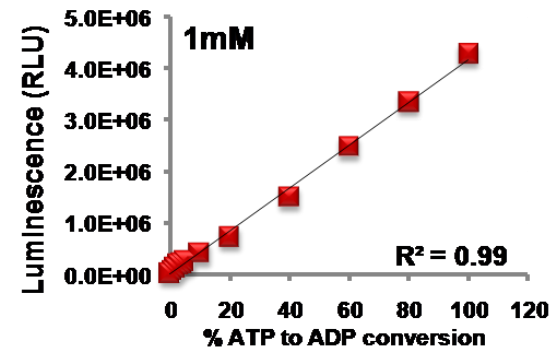
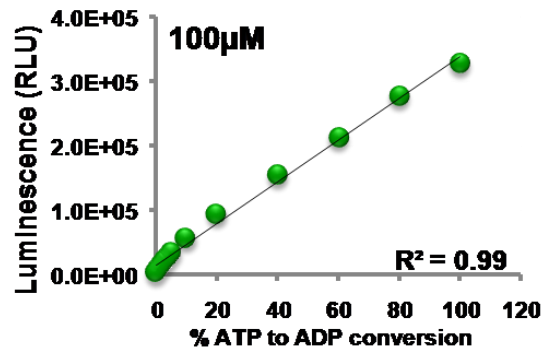
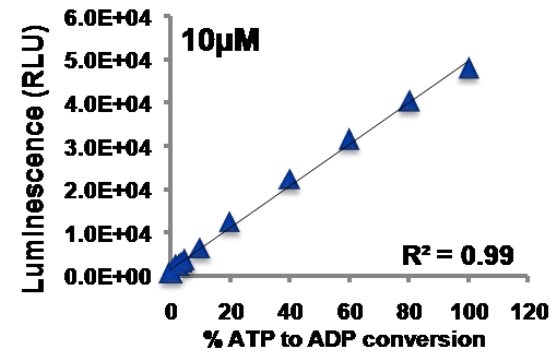
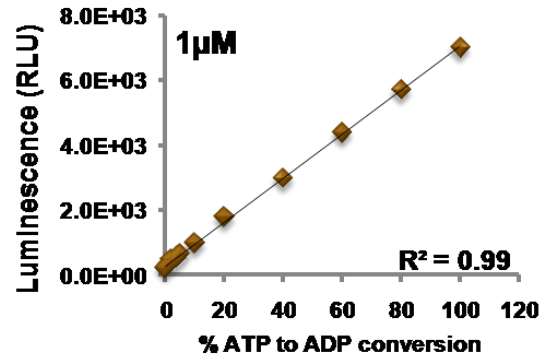
**Record Luminescence**



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

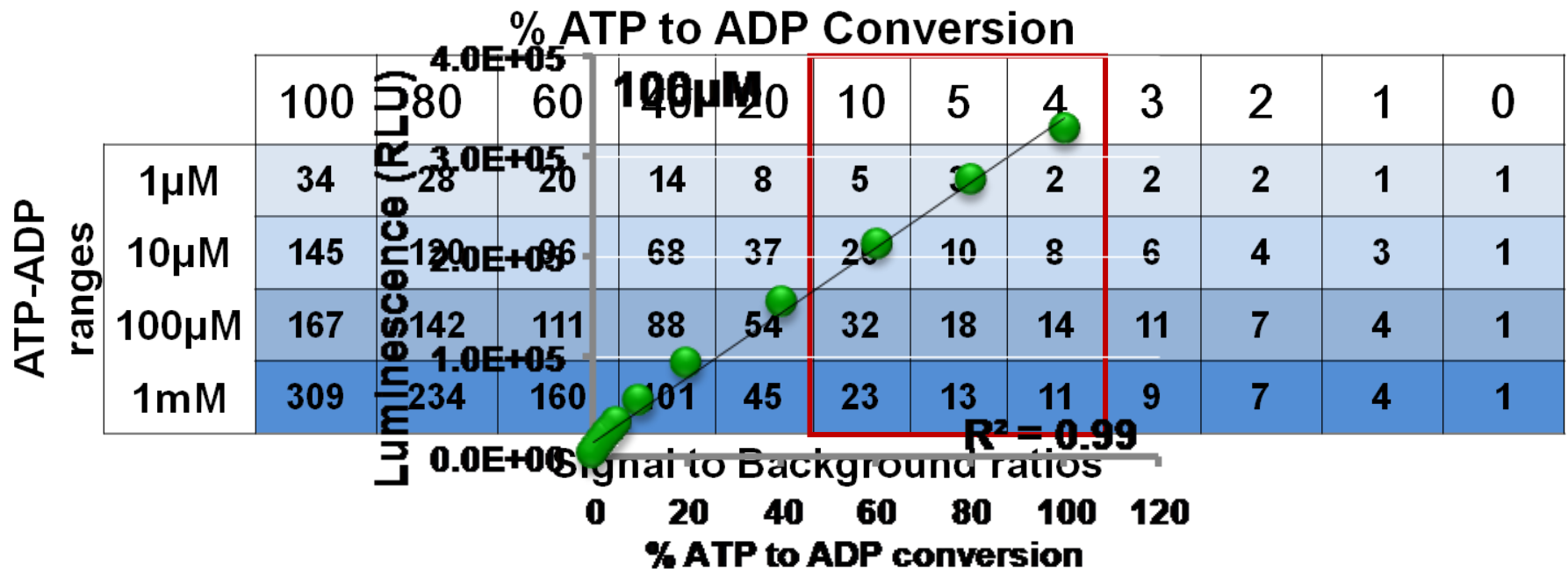


**ADP-Glo™ can be used virtually with any ATP concentration**



# Sensitivity of the ADP-Glo™ assay

Signal to Background ratios produced at different  
ATP to ADP conversion %



ADP-Glo™ can detect as low as 20nM ADP in  
5μl (0.1pmole) with a high Z' value

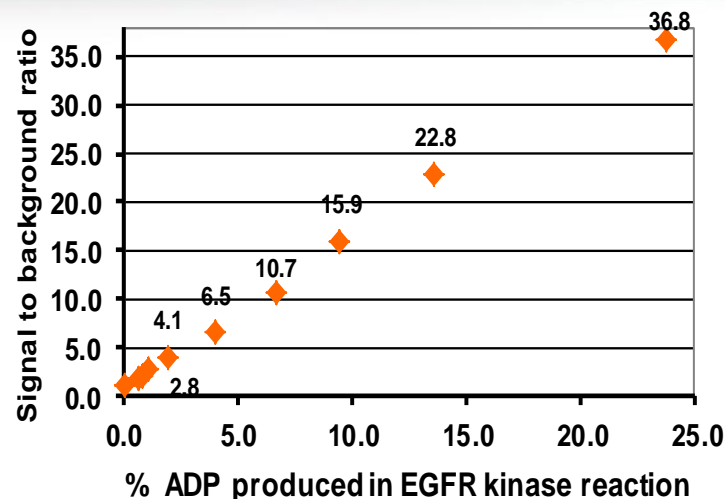
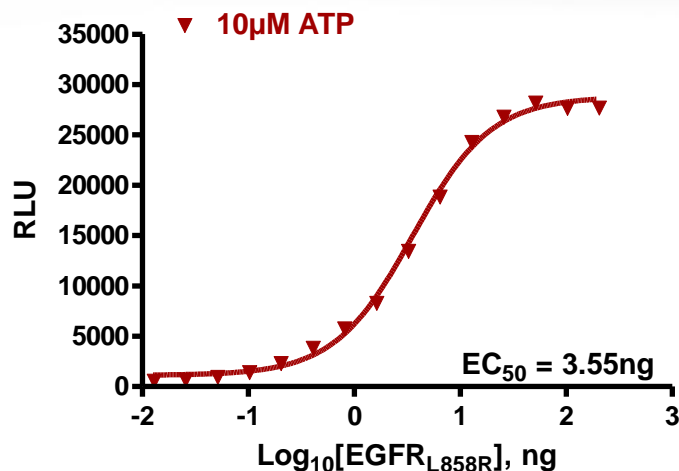
# MORE Sensitivity of the ADP-Glo™ assay

ADP-Glo™ Assay Performance Using Four Different Sources of ATP.

	ATP Concentrations								
	1mM			10μM					
% ADP in ATP + ADP									
Pro						5%	20%	10%	5%
Pro				54	32	18	37	20	10
Comp		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 highly improved with the new Ultra Pure ATP**

# ADP-Glo™ detection of Tyrosine kinase activity



EC<sub>50</sub>

EGFR <sub>L858R</sub> (ng/reaction)	100	50	25	12.5	6.25	3.13	1.56	0.78	0.39	0.20	0.1	0.05	0
RLU (Average)	27753	28278	26871	24358	18944	13551	8384	5849	3923	2404	1502	1045	368
% ADP produced	87.4	89.1	84.6	76.7	39.8	23.8	13.6	9.5	6.7	4.0	1.9	1.1	0
S/B	75.4	76.8	73.0	66.2	51.5	36.8	22.8	15.9	10.7	6.5	4.1	2.8	1

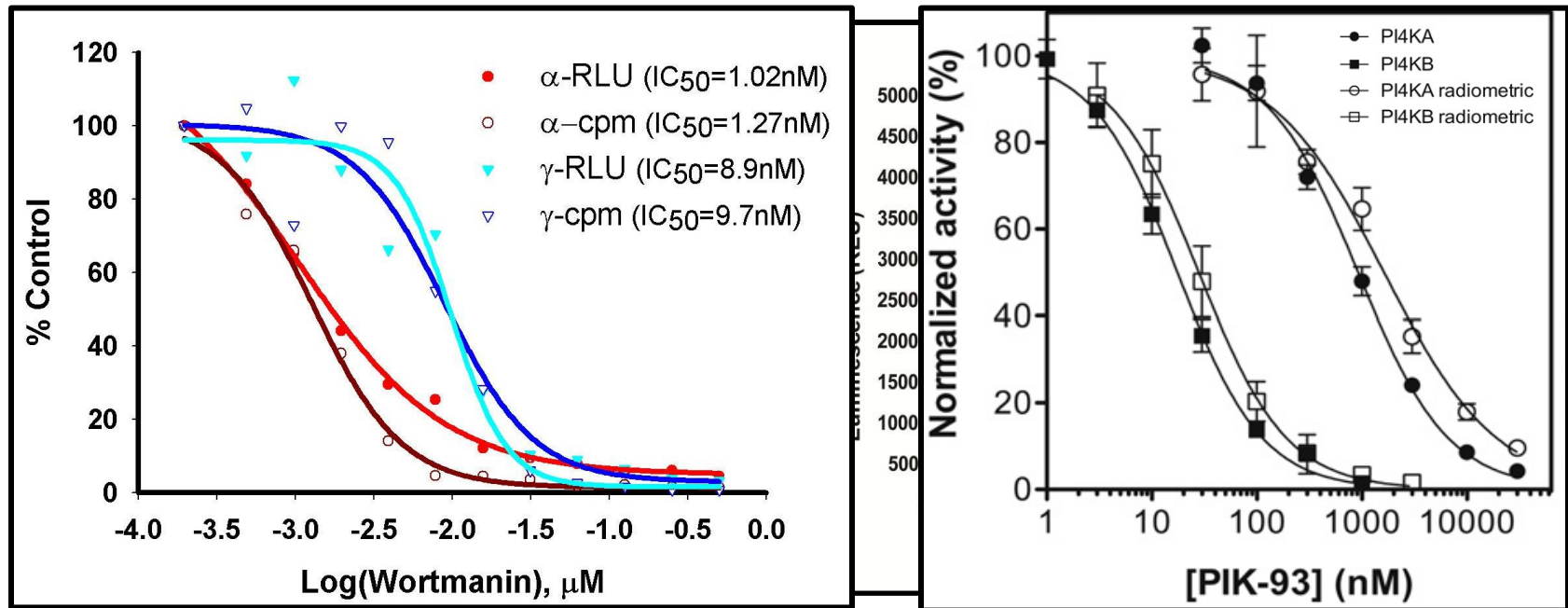
SB<sub>10</sub>

EGFR (ng)

0.36

Amount of enzyme to generate  
5-10% conversion

# 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.

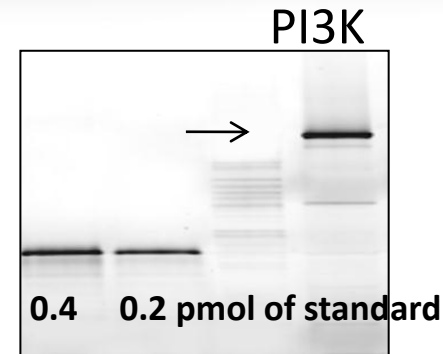
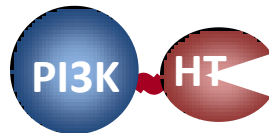
Tai AW, Bojjireddy N, Balla T. (2011) A homogeneous and nonisotopic assay for phosphatidylinositol 4-kinases. [Anal Biochem.](#) 417(1):97-102.

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



# Validation of ADP-Glo with immobilized Kinase

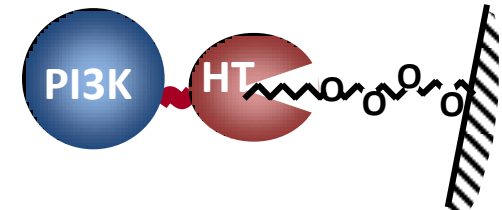
Expression as HaloTag fusion (HT)



In gel Visualization

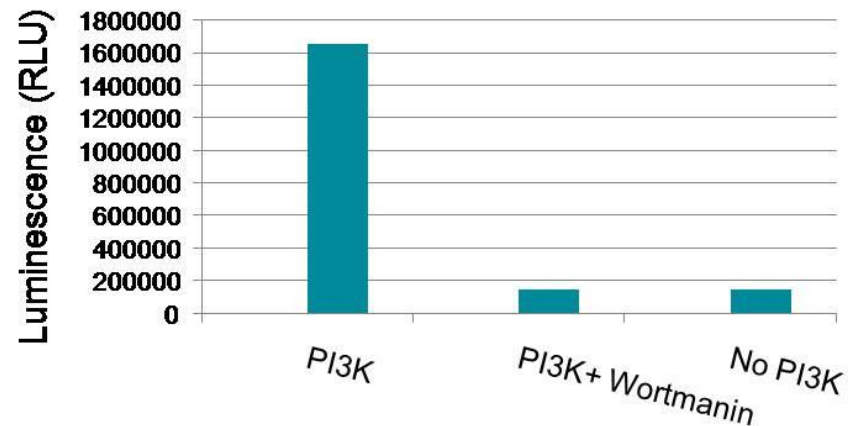


Immobilization on HaloLink beads



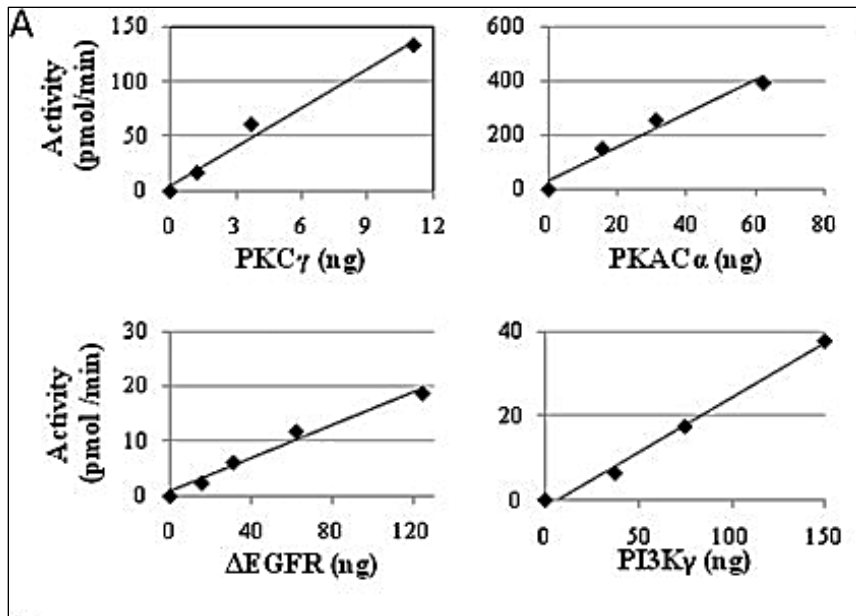
Activity measurement

ADP-Glo



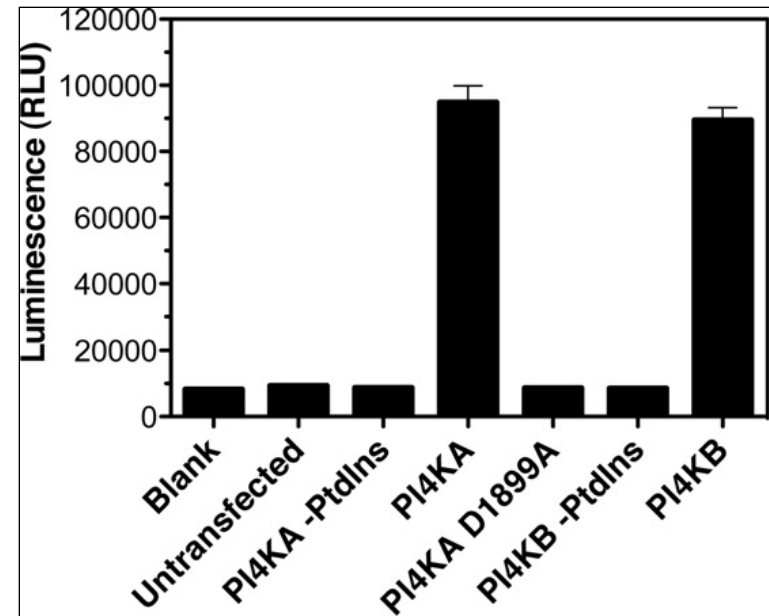
# Validation of ADP-Glo with mammalian expressed Kinases

## HaloTag



[Ohana RF](#), et al (2011) HaloTag-based purification of functional human kinases from mammalian cells. [Protein Expr Purif.](#) 76(2):154-64.

## FLAG Tag



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

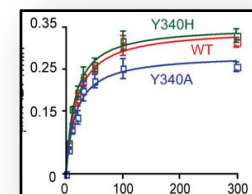
# One assay platform - many applications

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

**High-Throughput Screening**

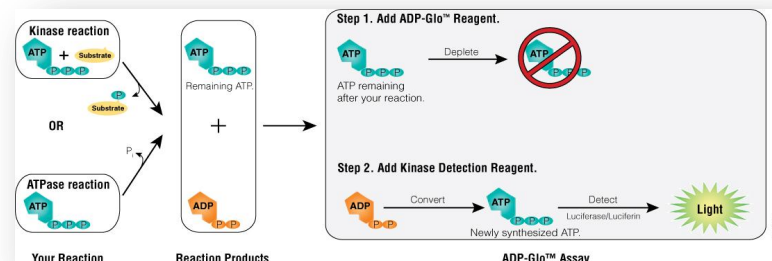


**Mode of action studies**



**Kinase inhibitor profiling**

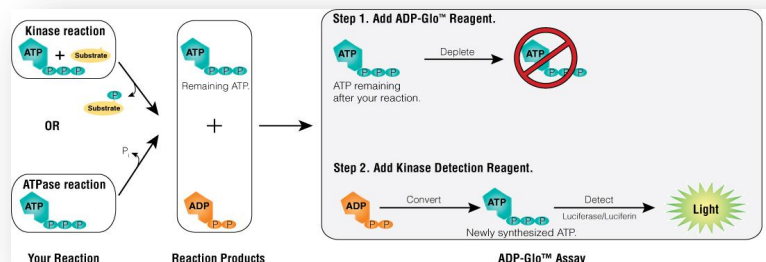
Protein	IC50	IC90	IC95
PKA	100	100	100
PKC	100	100	100
PKD	100	100	100
PKF	100	100	100
PKG	100	100	100
PKL	100	100	100
PKM	100	100	100
PKN	100	100	100
PKO	100	100	100
PKP	100	100	100
PKQ	100	100	100
PKR	100	100	100
PKS	100	100	100
PKT	100	100	100
PKU	100	100	100
PKV	100	100	100
PKW	100	100	100
PKX	100	100	100
PKY	100	100	100
PKZ	100	100	100
PKAA	100	100	100
PKAB	100	100	100
PKAC	100	100	100
PKAD	100	100	100
PKAE	100	100	100
PKAF	100	100	100
PKAG	100	100	100
PKAH	100	100	100
PKAI	100	100	100
PKAJ	100	100	100
PKAK	100	100	100
PKAL	100	100	100
PKAM	100	100	100
PKAN	100	100	100
PKAO	100	100	100
PKAP	100	100	100
PKAQ	100	100	100
PKAR	100	100	100
PKAS	100	100	100
PKAT	100	100	100
PKAU	100	100	100
PKAV	100	100	100
PKAW	100	100	100
PKAX	100	100	100
PKAY	100	100	100
PKAZ	100	100	100



**ADP-Glo™ Assay Platform**

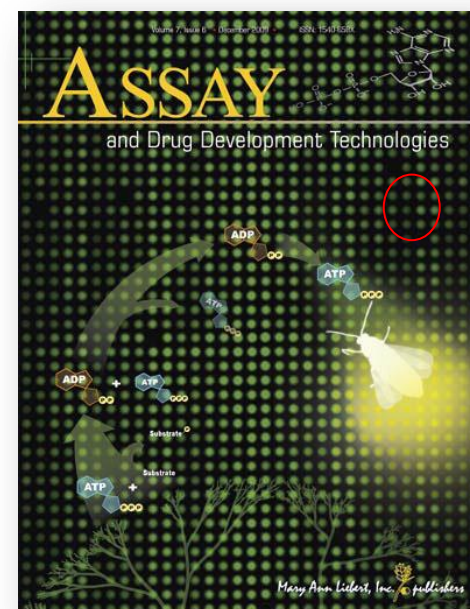
# One assay platform - many applications

## Screening Kinase Inhibitors with ADP-Glo™ Kinase assay



## ADP-Glo™ Assay Platform

## High-Throughput Screening

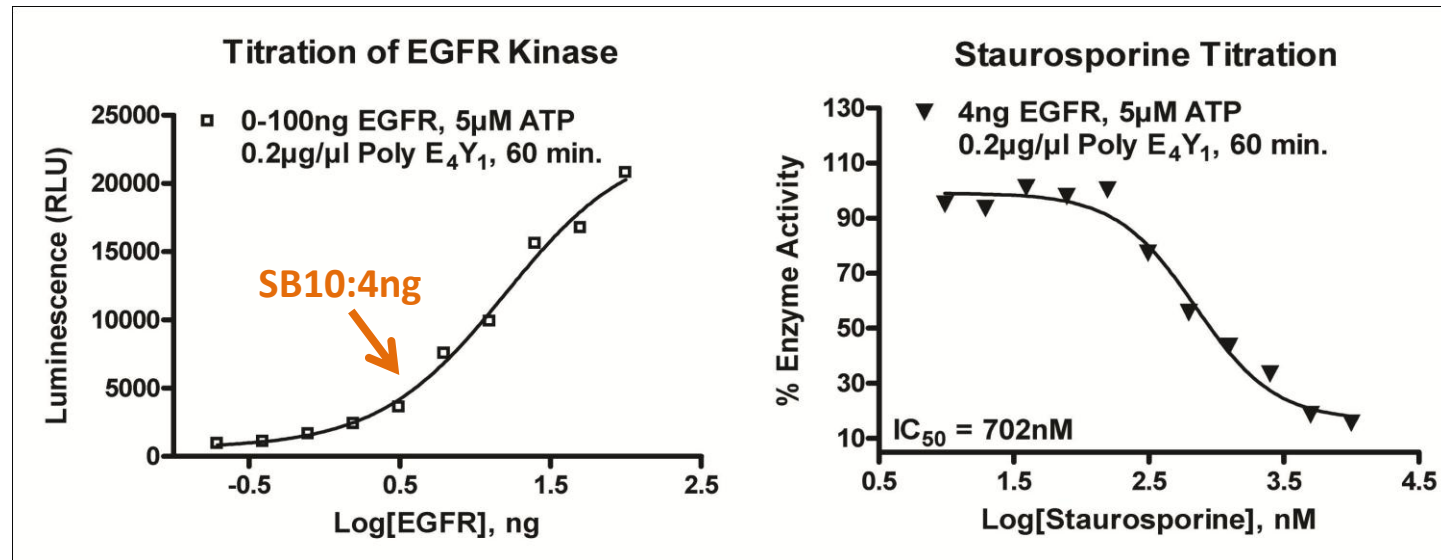


Free Access here:

<http://www.liebertonline.com/toc/adt/7/6>

# Developing kinase assay for screening and profiling

**SB 10 value** allows a high dynamic range with less variability and more accuracy in an inhibitor screening or Profiling (High Z' factor)



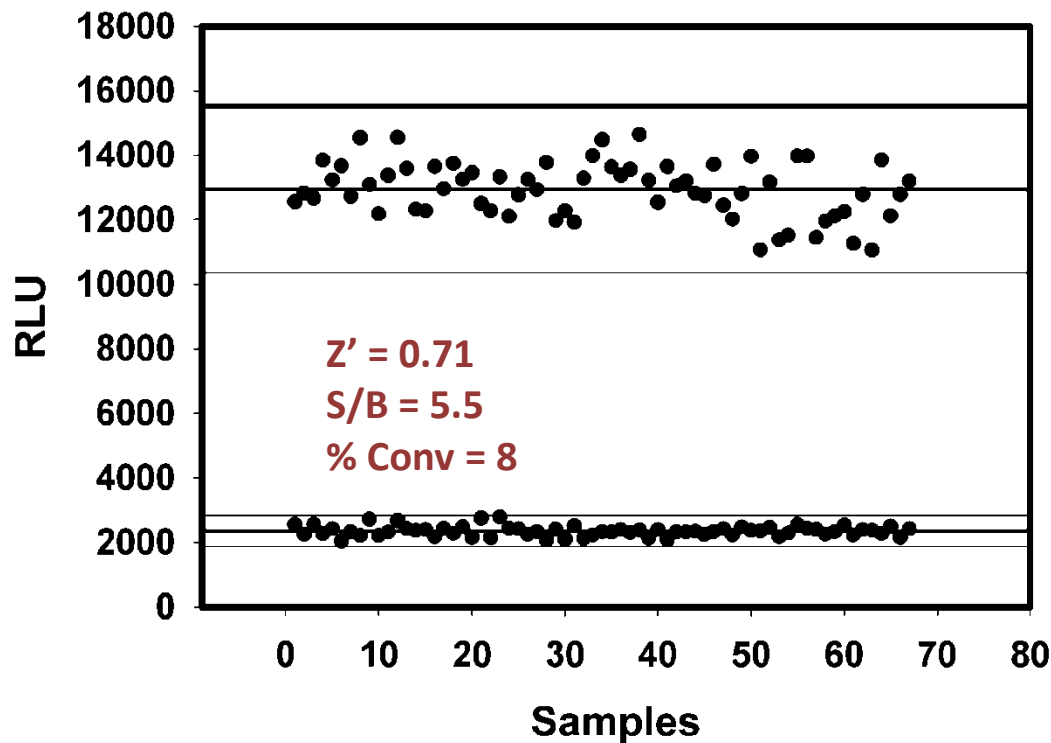
**Kinase Titration:** Using the defined ATP concentrations, perform ½ serial dilution of Enzyme.

**SB10 value:** Corresponds to the amount of Kinase needed to generate a 5-10% substrate conversion.



# Sensitivity, Robustness of the ADP-Glo™ Assay

During HTS, a Z' values >0.5 shows a robust assay



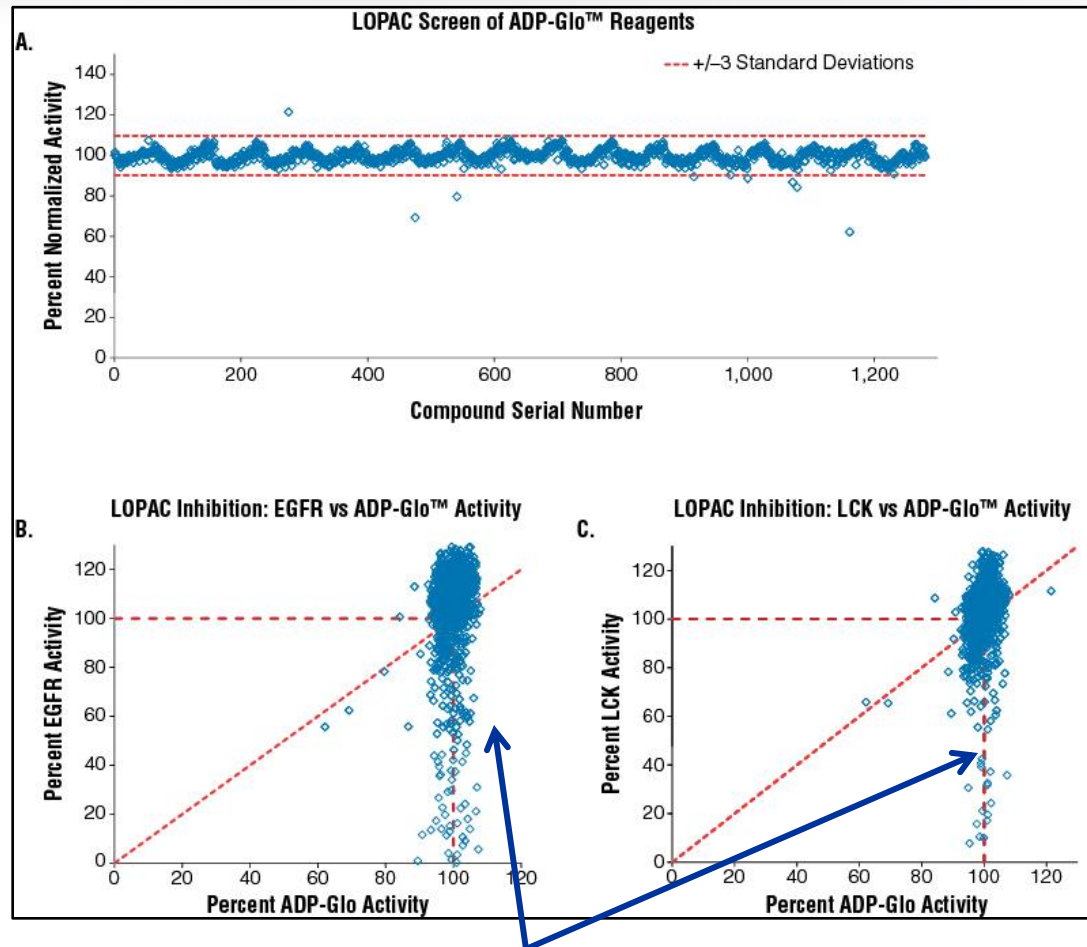
$$Z\text{-factor} = 1 - \frac{3(\sigma_p + \sigma_n)}{|\mu_p - \mu_n|}$$

$\sigma_p$ : STDEV of sample    $\sigma_n$ : STDEV of control  
 $\mu_p$ : Mean of sample    $\mu_n$ : Mean of control

Zhang JH, Chung TD, Oldenburg KR (1999) *J Biomol Screen.* **4**, 67–73

**ADP-Glo™ can detect low Enzyme activity with a high Z' value**

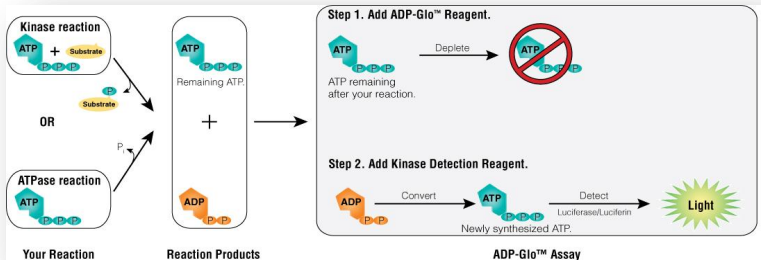
# Screening for Kinase inhibitors using luminescent ADP detection



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

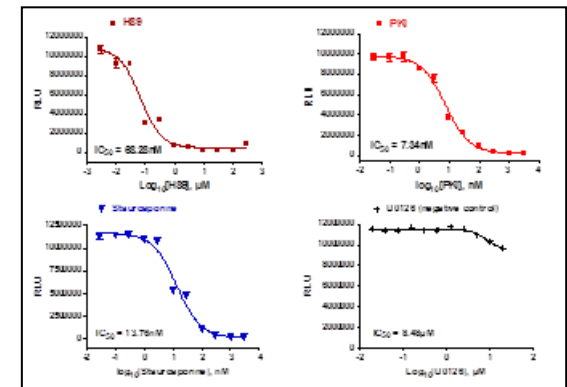
# One assay platform - many applications

## Mode of Action Studies with ADP-Glo™ Kinase Assay



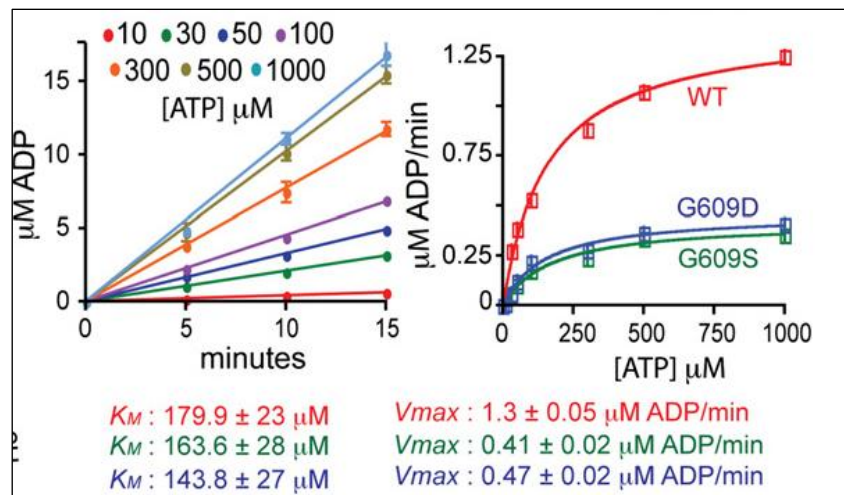
## ADP-Glo™ Assay Platform

## Mode of action studies

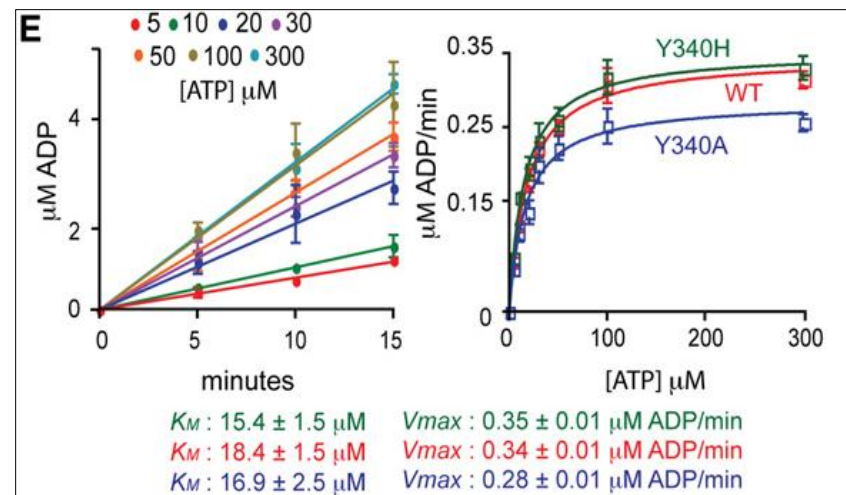


# Determination of Biochemical values using ADP-Glo™

## C-Src



## Haspin

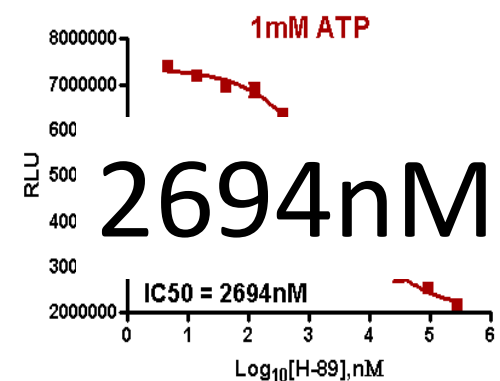
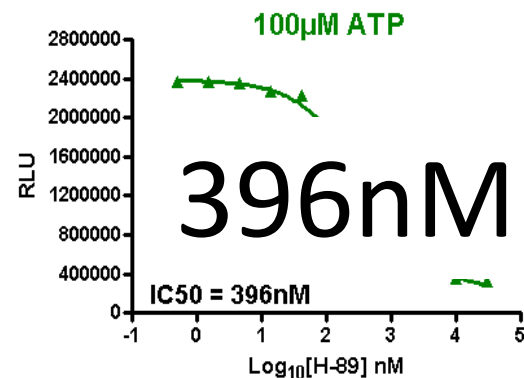
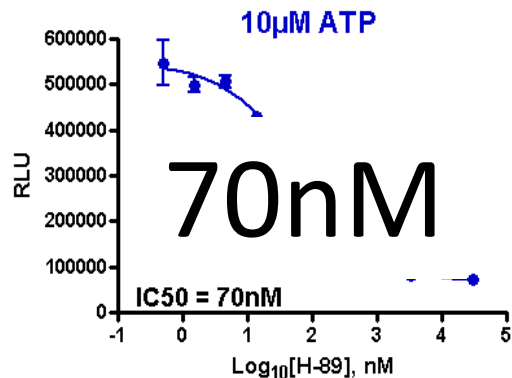


Balzano D et al., (2011). A general framework for inhibitor resistance in protein kinases. *Chem Biol.* 18(8):966-75.

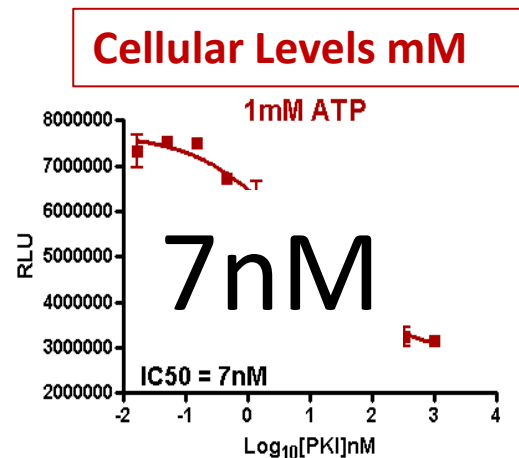
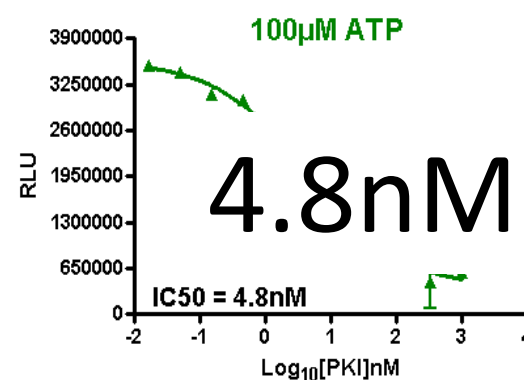
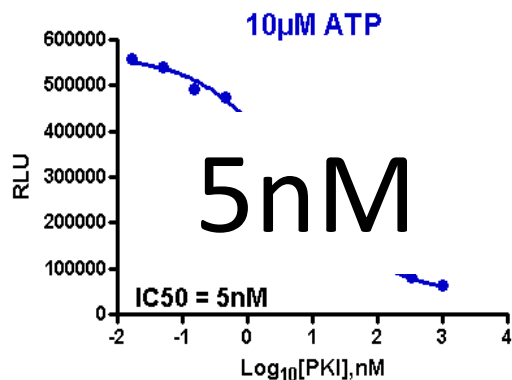
**ADP-Glo™ produces  $K_M$  values similar to literature**

# Determination of inhibitor's mechanism of action

## PKA ATP Competitive inhibitor H-89



## PKA ATP non Competitive inhibitor PKI

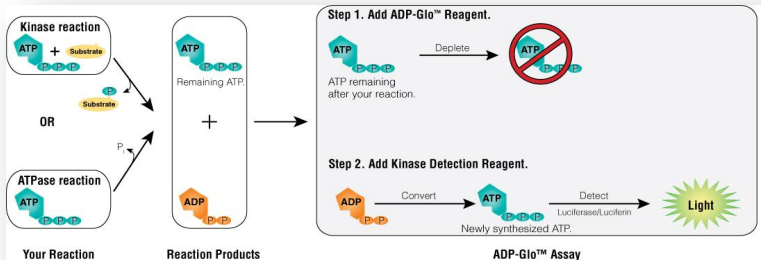


ADP-Glo™ is a perfect assay to distinguish between ATP competitive and non competitive kinase inhibitors



# One assay platform - many applications

## Profiling Kinase Inhibitors with ADP-Glo™ Kinase Assay



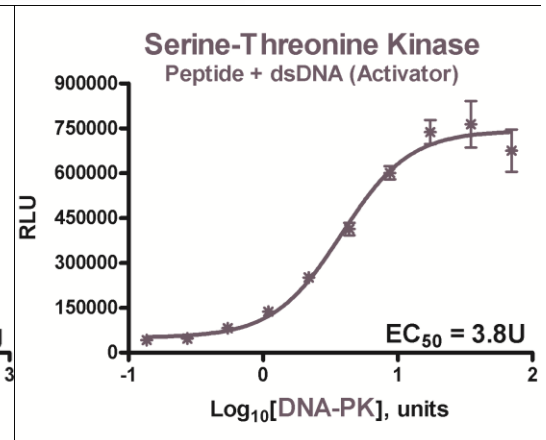
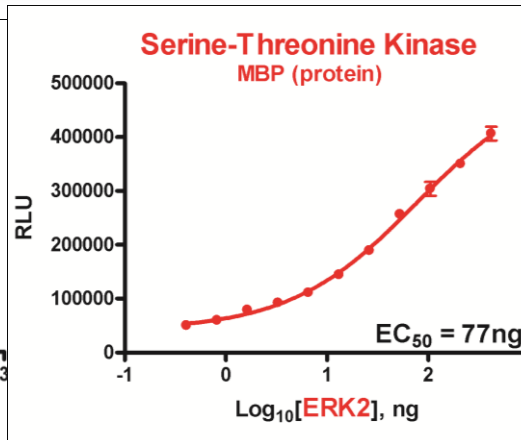
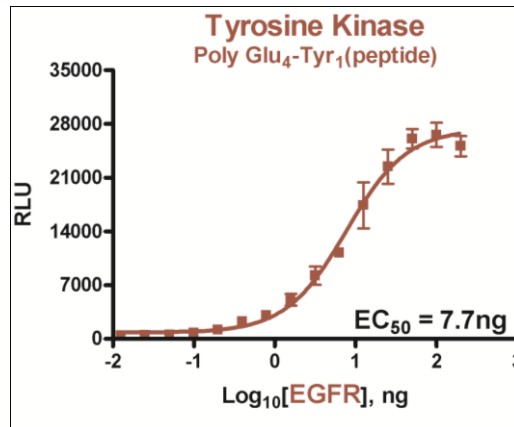
## ADP-Glo™ Assay Platform

## Kinase inhibitor profiling

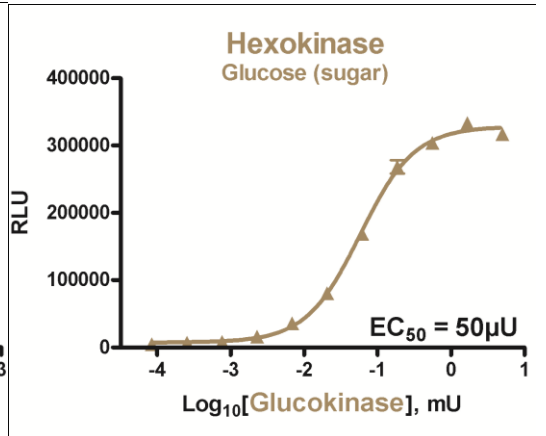
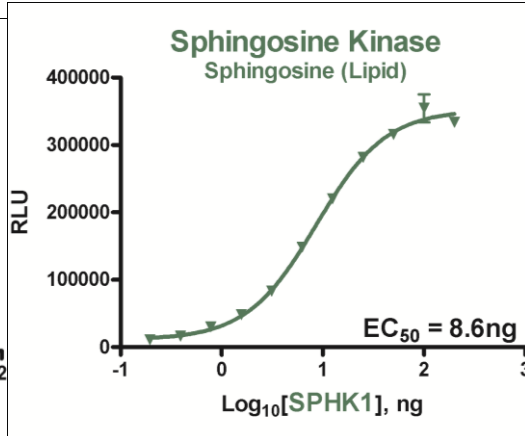
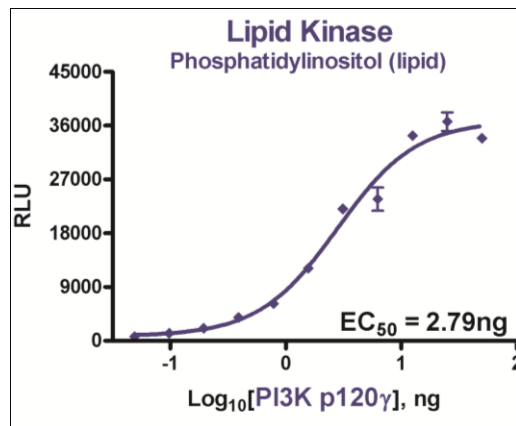
Kinase	LCK inhibitor	CDK/CRK inhibitor	Ro-32-0432
ASK1	100	83	99
HPK1	100	24	92
KHS1	96	77	100
MEK1	100	89	99
MEKK1	64	82	100
MINK1	100	96	81
NIK	100	92	100
MEKK2	60	95	95
MST1	92	63	98
MYO3b	100	91	95
PAK1/CDC42	66	100	100
PAK3	100	90	100
SLK	91	85	93
TAOK1	92	25	100
TNIK	88	100	94
ERK2	81	100	85
GSK3b	68	20	25
JNK1	92	100	100
JNK3	91	98	100
p38a	100	100	100
p38g	100	100	100
p38d	98	100	100
CDK2/A2	12	1	73
CDK3E1	45	17	80
CDK5/p25	76	2	98
CDK5/p35	73	1	100
CDK6/D3	58	100	60
CDK9/K	37	43	63
CLK3	57	100	94
AMPK A1/B1/G1	100	68	71
AMPK A1/B1/G2	100	67	100
AMPK A2/B1/G1	100	46	100
CAMK2a	100	81	100
CAMK2g	100	72	100
CAMK4	100	67	100
DAPK1	100	100	100
STK33	99	100	100
CHK2	100	100	100
MAPKAPK2	100	100	100
MARK1	100	73	100
PASK	100	87	100
PKCmu	100	93	100

# Diverse kinase-substrate combinations with ADP-Glo™

## Protein Kinases

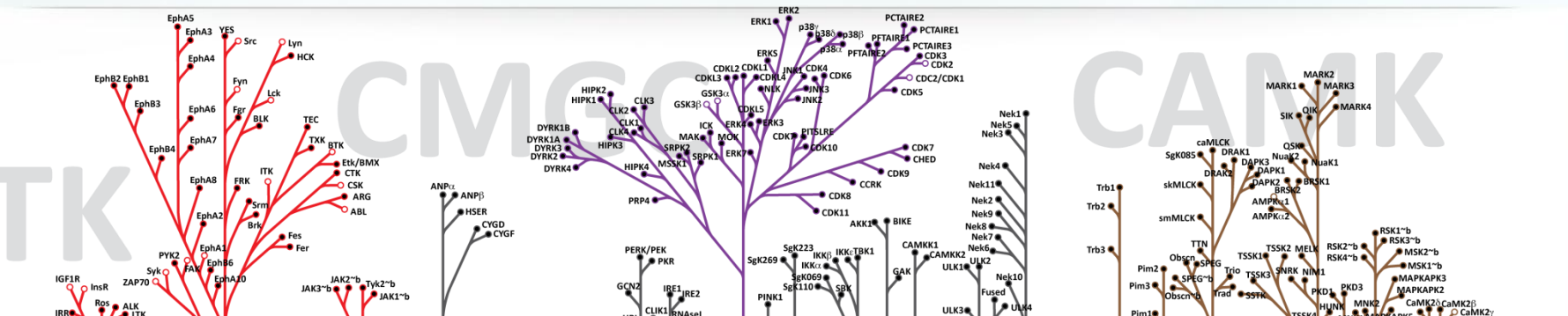


## Non-Classical Kinases

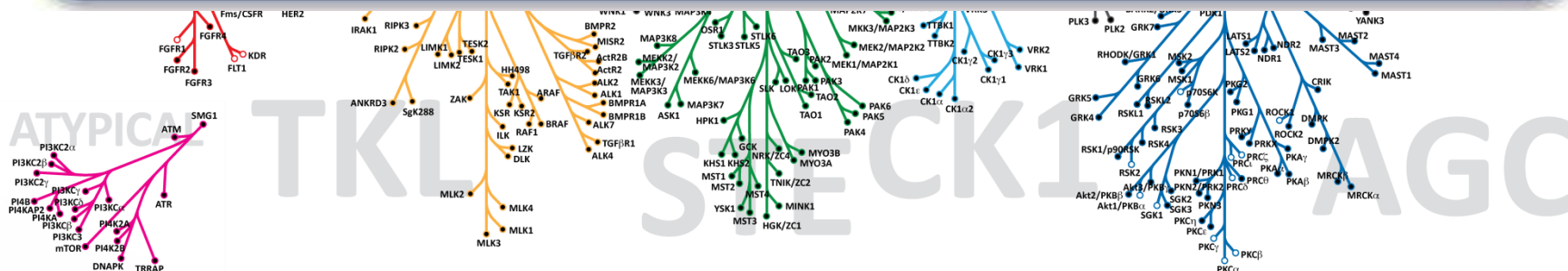


**ADP-Glo™ Kinase Assay detects the activity of any Kinase regardless of the substrate chemical structure**

# Profiling the Human Kinome



## Kinase Inhibitor Profiling



**ADP-Glo™ is a universal Kinase Assay that detects the activity of any Kinase regardless of the substrate chemical structure**

# ***Profiling Kinase inhibitors is a critical step in Drug Development***

## **Assessing selectivity and cross reactivity of Kinase inhibitors**

Knowing the exact kinase inhibition profile of a compound is critical for the understanding of its mode of action and to assess any side effects that may result from its extra activities.

**Gleevec (BCR-ABL), PD 98059 (p38), U0126 (MEK1), Rapamycin (mTOR), were shown to be very selective inhibitors, while PP1 and PP2 (Src family) inhibit other kinases in different families, CSK, SAPK2a/p38, CK1, KIT and BCR-ABL.**

Identifying new targets during a compound profiling could lead to novel therapeutic applications.

**Gleevec (BCR-ABL inhibitor) was marketed by Novartis initially for the treatment of chronic myeloid leukemia (CML), then its indication was expanded to gastrointestinal stromal tumors (GIST) because of its effect on c-Kit kinase.**

Profiling already in development or clinically approved compounds against drug resistant Kinase targets could lead to new treatment of relapsed patients with ineffective first-line targeted therapies.

**MK-0457 (Aurora Kinase inhibitor) in clinical development for the treatment of solid tumors is shown to inhibit an ABL kinase mutant (T315I), that it could be potentially used to treat Gleevec-resistant cancer.**

# *Providing Complete solutions for the kinase field:*

*Promega + Signalchem collaboration*

## **Kinase Enzyme Systems (KES)**

Kinase Enzyme System Manufactured By



Exclusively Distributed By



---

**Promega Corporation**

---

2800 Woods Hollow Road

---

Madison WI 53711 USA

---

Telephone: 608-274-4330

---

Toll-Free: 800-356-9526

---

Fax: 608-277-2516

---

Internet: [www.promega.com](http://www.promega.com)

---



# Kinase Enzyme System (KES) offering

2 catalog numbers per enzyme system

  
**Promega**  
 Catalog # V9101

## 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

+

 **SignalChem**  
*Specialists in Signaling Proteins*

Promega Catalog #

## 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 <sub>2</sub> *
500µl	Kinase Activator**

\*For Tyrosine Kinases  
 \*\*Lipids for PKC Kinases

=

  
**Promega**  
 Catalog # V9061

# Promega kinase panel

## (Kinase Enzyme Systems)

Kinase Enzyme Systems (Example)

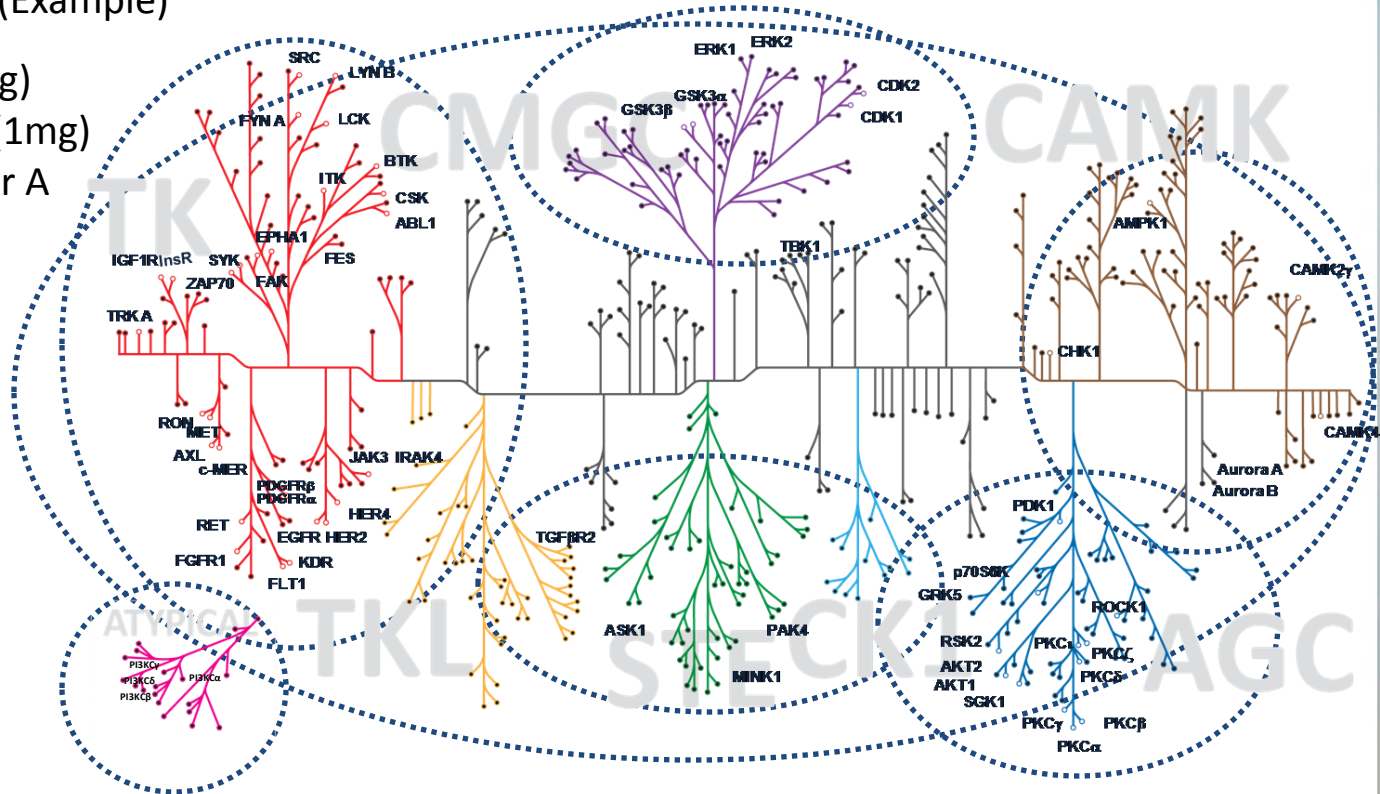
Akt1 Kinase (10 $\mu$ g)

Akt (PKB) substrate (1mg)

5 x Reaction Buffer A



**ADP-Glo™**  
**Kinase Assay**



**Promega Kinase Panel contains initially 70 Kinases that  
Cover the Human kinome**

# Validating Kinase Enzyme systems with ADP-Glo™

## Kinases were tested with ADP-Glo™ Kinase Assay

Kinase	Substrate	ATP Concentration (μM)	SB10 (ng)	IC50 Staurosporine (nM)
<b>CMGC</b>				
CDK1/CyclinA2	Histone H1	50	1	1.9
CDK2/CyclinA2	Histone H1	50	0.6	1.8
CDK5/p25	Histone H1	10	2	1
CDK5/p35	Histone H1	10	3	1.2
ERK1	MBP	50	2.8	>1000
ERK2	MBP	50	1.6	>1000
GSK3a	GSK3 Substrate	25	1	6.4
GSK3b	GSK3 Substrate	25	1	0.3
LYN B	SRC substrate	25	3	4.7
MET	Poly (Glu <sub>4</sub> , Tyr <sub>1</sub> )	10	4	280
PDGFRa	Poly (Glu <sub>4</sub> , Tyr <sub>1</sub> )	25	7	0.48
PDGFRb	Poly (Glu <sub>4</sub> , Tyr <sub>1</sub> )	25	6.5	0.3
RET	IGF1tide	25	2	2.3
RON	Axltide	25	2	76.5
SRC	SRC substrate	50	1.8	250.8
SYK	Poly (Glu <sub>4</sub> , Tyr <sub>1</sub> )	10	2.8	0.17
TRKA	Poly (Glu <sub>4</sub> , Tyr <sub>1</sub> )	50	0.6	0.3
ZAP70	Poly (Glu <sub>4</sub> , Tyr <sub>1</sub> )	3	2.9	103.5
<b>TKL</b>				
IRAK4	MBP	25	5.0	0.75
TGFβR2	MBP	50	30	>1000
<b>STE</b>				
ASK1	MBP	25	4.4	21.4
MINK1	MBP	50	0.6	2.8
PAK4	Akt substrate II	5	10	0.2
p70S6K	S6K substrate	25	20	82
PDK1	PDKtide	5	8	0.75
PKCa	CREBtide	25	0.2	5.94
PKCb II	CREBtide	50	0.1	6.18
PKCd	CREBtide	50	0.8	1.87
PKCg	PKCtide	50	0.3	6.2
PKCi	CREBtide	25	0.5	518
PKCz	CREBtide	5	0.2	>1000
ROCK1	S6K substrate	5	10	6.7
RSK2	RSK Substrate	25	1	11.3
SGK1	Akt (PKB) substrate	50	3.5	12.2
<b>OTHER</b>				
Aurora A	MBP	25	7	0.47
Aurora B	MBP	25	6.4	1.5
NEK2	MBP	50	5	>1000
PLK1	Dephospho Casein	5	40 (SB4)	>1000
TBK1	MBP	25	2.2	0.05
ULK1	MBP	10	5.0	18.6

- ADP-Glo
- To gene

ted

# 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 $\alpha$ /p85 $\alpha$   
 p110 $\alpha$  (E545K)/p85 $\alpha$   
 p110 $\alpha$  (H1047R)/p85 $\alpha$   
 P110 $\beta$ / p85 $\alpha$   
 P110 $\gamma$ / p85 $\alpha$   
 P110 $\delta$ / p85 $\alpha$

+

PI:PS Lipid  
Kinase substrate  
0.5ml x 1mM

or

PIP2:PS Lipid Kinase  
substrate  
0.25ml x 1mM



**Promega**

Catalog #XXXX

Enzymes are offered separately  
&  
together in a complete Class I  
profiling kit

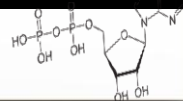
ADP-Glo™ Kinase Assay

# Promega kinase panel

## (Kinase Enzyme Systems)

Promega Kinase Panel only expanding by 70 Kinases that covers 9 families human kinome

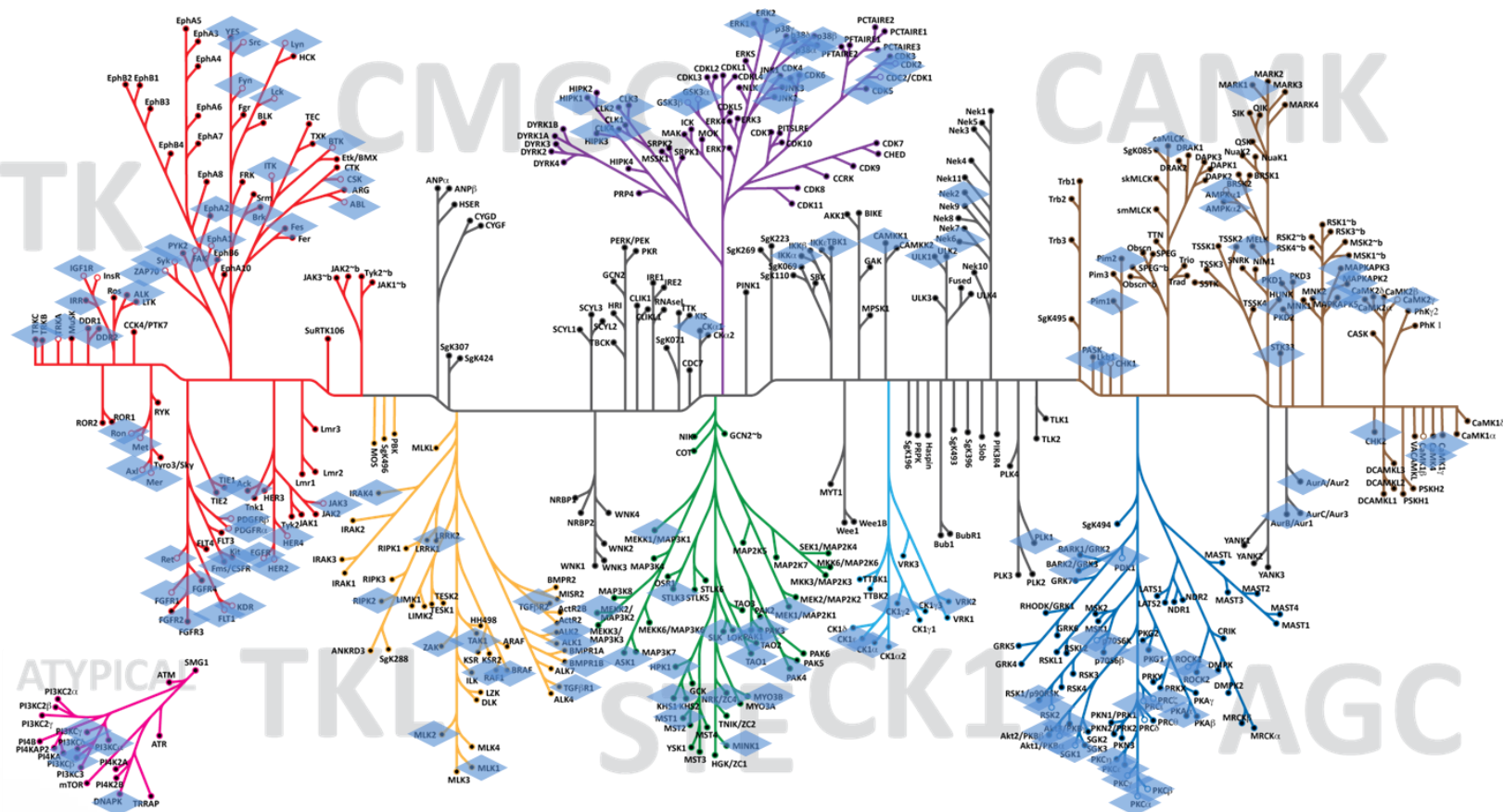
TK			TKL	CAMK	STE	CMGC	AGC	OTHER/CK1
ABL1	RET	KIT T670I	IRAK4	AMPK A1/B1/G1	ASK1	CDK1/CyclinA2	AKT1	Aurora A
AXL	RON	PDGFRA D842V	TGFβR2	CAMK2g	MINK1	CDK2/CyclinA2	AKT2	Aurora B
BTK	SRC	PDGFRA T674I	TGFβR1	CAMK4	PAK4	CDK5/p25	GRK5	NEK2
c-MER	SYK	RET Y791F	BRAF	CHK1	MEK1	CDK5/p35	p70S6K	PLK1
CSK	TRKA	RET V804L	BRAF V600E	AMPK A2/B1/G1	PAK3	ERK1	PDK1	TBK1
EGFR	ZAP70	MET M1250T	RAF1(EE)	MAPKAPK5	MST1	ERK2	PKCa	ULK1
EPHA1	FLT3	FGFR3 K650E	MLK2	PIM2	SLK	GSK3a	PKCb II	NEK6
FAK	FMS	FLT3 D835Y	TOPK	PKD2	HPK1	GSK3b	PKCd	DNA-PK
FES	FGFR2		ALK1	PIM1	PAK1/CDC42	JNK3	PKCg	IKKa
FGFR1	FGFR4		MLK1	CAMK1g	MEKK1	p38a	PKCi	CK2a1
FLT1	ACK		RIPK2	MLCK/MYLK	MEKK2	p38g	PKCz	CAMKK1
FYN A	DDR2		BMPR1b (ALK6)	CAMK2a	MYO3b	CDK9/Cyclin K	ROCK1	CK1α1
HER2	PYK2		TAK1-TAB1	MAPKAPK2	STK39/STLK3	JNK1	SGK1	VRK2
HER4	c-Kit		ALK2	MAPKAPK3	KHS1	CDK2/Cyclin E1	RSK2	CK1g2
IGF1R	TRKB		ZAK	DAPK1	NIK	p38b	AKT3	CK1epsilon
InsR	BRK		LRRK2	SIK	TAOK1	p38d	PKCe	CK2
ITK	EGFR L858R			MYLK3/caMLCK		CDK7	RSK1	
JAK3	EGFR L861Q			STK33		CDK4/Cyclin D3	PKCtheta	
KDR	EGFR T790M			CHK2		CDK6/Cyclin D3	ROCK2	
LCK	EGFR T790M L858R			PKCmu (PKD1)		CDK3/CyclinE1	p70S6Kb	
LYN B	ABL1 E255K			PASK		CLK1	GRK2	
MET	ABL1 G250E			MELK		CLK3	PKG	
PDGFRa	ABL1 T315I			MARK1		HIPK1	PKA	
PDGFRb	ABL1 Y253F			AMPK A1/B1/G2		HIPK3	PKC	





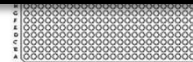
# Promega kinase panel

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



Broad Human Kinome coverage with 174 Kinase Enzyme Systems

## Streamlined profiling protocol



## 16 Kinase strips (128 Kinases)

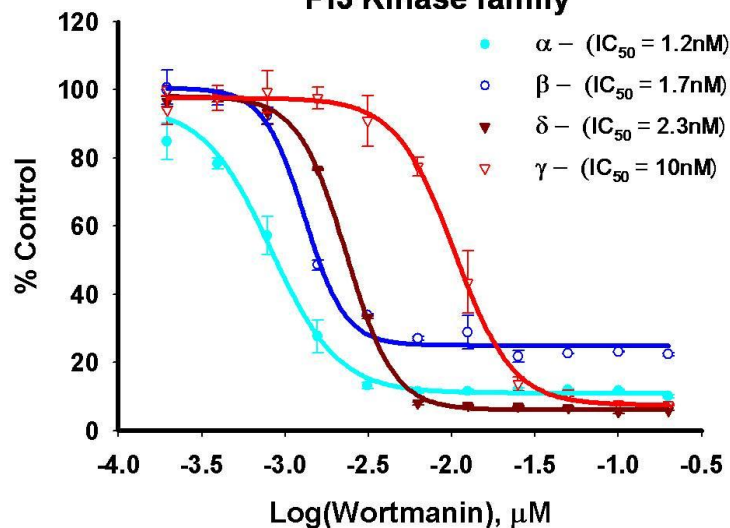
## Profiling with ADP-Glo platform made simple



# Profiling kinase inhibitors with ADP-Glo™ kinase platform

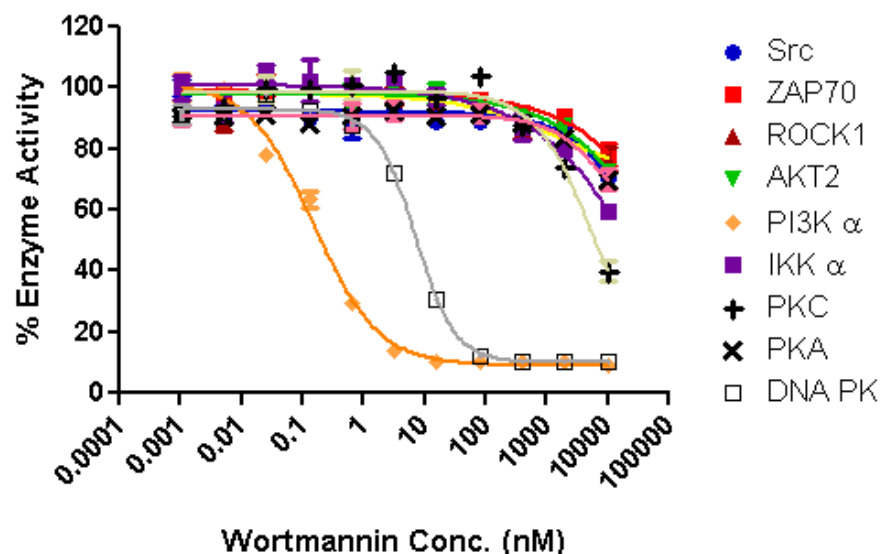
## Against a kinase family

**Selectivity Profile of Wortmanin towards PI3 Kinase family**



## Against a Panel of Different families

**Selectivity profile for Wortmannin against 9 kinase panel using ADP-Glo™**

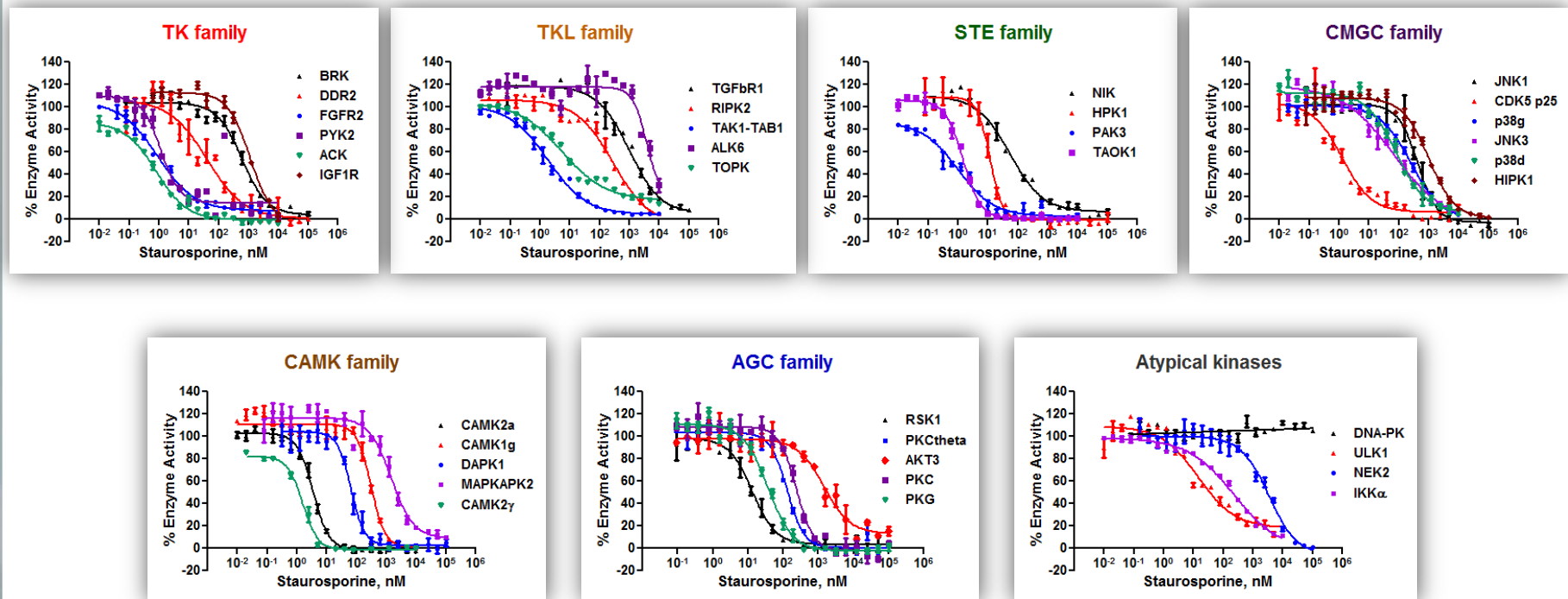


**As a “One Assay for all”, ADP-Glo™ is an ideal assay for Profiling inhibitors**



# Profiling kinase inhibitors with ADP-Glo™ kinase platform

## 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**

# Determination of inhibitor profiles of different kinase families

LCK inhibitor  
(Src family specific)

Kinase	LCK inhibitor	CDK/CRK inhibitor	Ro-32-0432
ER2	72	78	100
HR4	16	83	100
IGF1R	100	100	100
InsR	95	96	100
KDR	56	98	96
PDGFR	25	77	97
PDGFRβ	36	63	63
ABL1	100	100	59
BRK	1	100	72
BTX	3	87	88
CSK	11	100	89
FYN A	4	66	90
LCK	20	89	81
LYN B	12	96	71
SRC	15	92	100
AXL	100	94	100
EPHA1	68	85	98
FAK	100	100	100
ITK	84	86	90
JAK3	100	92	92
PK2	100	90	100
SYK	89	96	95
TRKA	36	100	50
FGFR1	38	27	98
FGFR2	76	96	98
FGFR4	81	93	100
FLT1	84	91	100
FLT3	13	93	75
FMS	62	100	100
MET	87	71	100
RET	17	99	100
BMPRIb	100	87	100
BRAP	98	98	100
IRAK4	100	92	100
LRRK2	71	84	100
MLK1	96	72	100
RAF1(EE)	92	92	100
TAK1-TAB1	100	72	100
TGFR1	100	76	100
ALK1	95	99	86
BRAP-V500E	97	100	97
MLK2	87	90	100
RIPK2	15	95	87
TGFR2	82	100	100
TOPK	99	100	100
ZAK	87	100	100

Cdk/CRK inhibitor  
(CDK family specific)

Kinase	LCK inhibitor	CDK/CRK inhibitor	Ro-32-0432
ASK1	100	83	99
HPK1	100	21	92
KHS1	96	77	100
MEK1	100	89	99
MEKK1	64	82	100
MINK1	100	96	81
NIK	100	92	100
MEKK2	60	95	95
MST1	92	63	98
MYO3b	100	91	95
PAK1/CDC42	66	100	100
PAK3	100	90	100
SLK	91	85	93
TAK1	92	25	100
TNIK	88	100	94
ERK2	81	100	85
GSK3β	68	20	25
JNK1	92	100	100
JNK3	91	98	100
p38α	100	100	100
p38β	100	100	100
p38γ	98	100	100
CDK2/A2	12	1	73
CDK3E1	46	17	80
CDK5/p25	76	2	98
CDK5/p35	73	3	100
CDK6/D3	58	100	60
CDK9/K	37	33	63
CLK3	57	100	94
AMPK A1/B1/G1	100	68	71
AMPK A1/B1/G2	100	67	100
AMPK A2/B1/G1	100	46	100
CAMK2a	100	81	100
CAMK2γ	100	72	100
CAMK4	100	67	100
DAPK1	100	100	100
STK33	99	100	100
CHK2	100	100	100
MAPKAPK2	100	100	100
MARK1	100	73	100
PASK	100	87	100
PKCmu	100	93	100

Ro-32-0432  
(PKC family specific)

Kinase	LCK inhibitor	CDK/CRK inhibitor	Ro-32-0432
p70S6K	100	82	90
PKA	100	100	100
PKC	100	89	100
PKCβ	100	91	100
PKG	100	87	100
ROCK1	100	100	100
RSK2	100	87	33
PKCa	100	100	5
PKCb II	100	94	0
PKCγ	100	100	7
PKCd	100	94	4
PKCe	100	87	8
PKCi	100	88	49
PKCtheta	100	100	2
PKCz	100	94	41
Aurora A	100	91	91
Aurora B	100	33	89
CK2a1	100	95	100
CK1α	100	83	100
CK1γ2	100	17	100
CK1epsilon	100	60	100
VRK2	100	100	97
CAMKK1	100	84	95
IKKα	87	100	44
IKKβ	100	100	84
NEK2	100	100	100
NEK6	100	88	98
PLK1	100	89	100
TBK1	100	82	100
ULK1	100	75	91

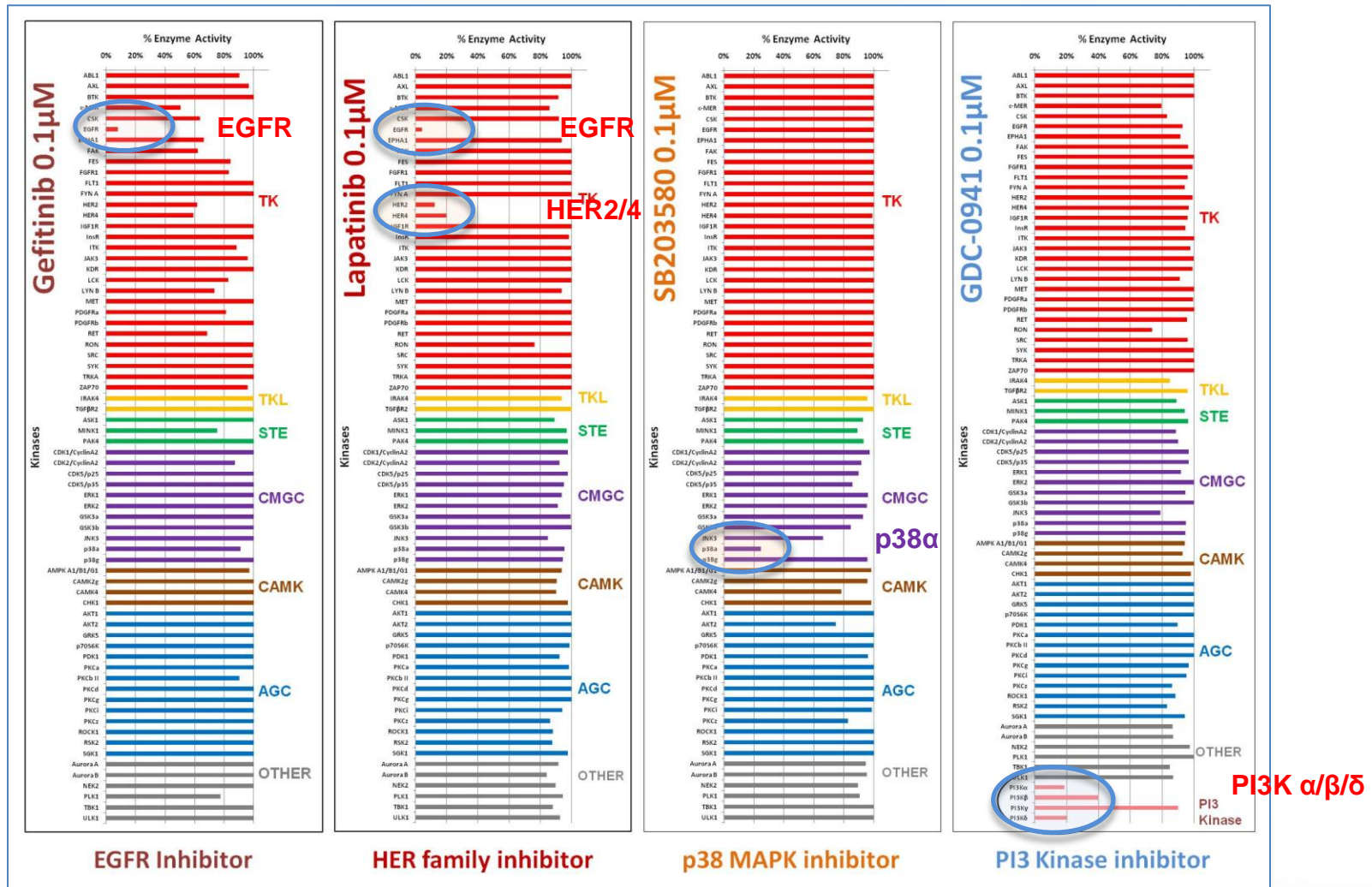
More inhibited

Less inhibited

ADP-Glo™ Kinase platform for convenient and meaningful selectivity profiles creation



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



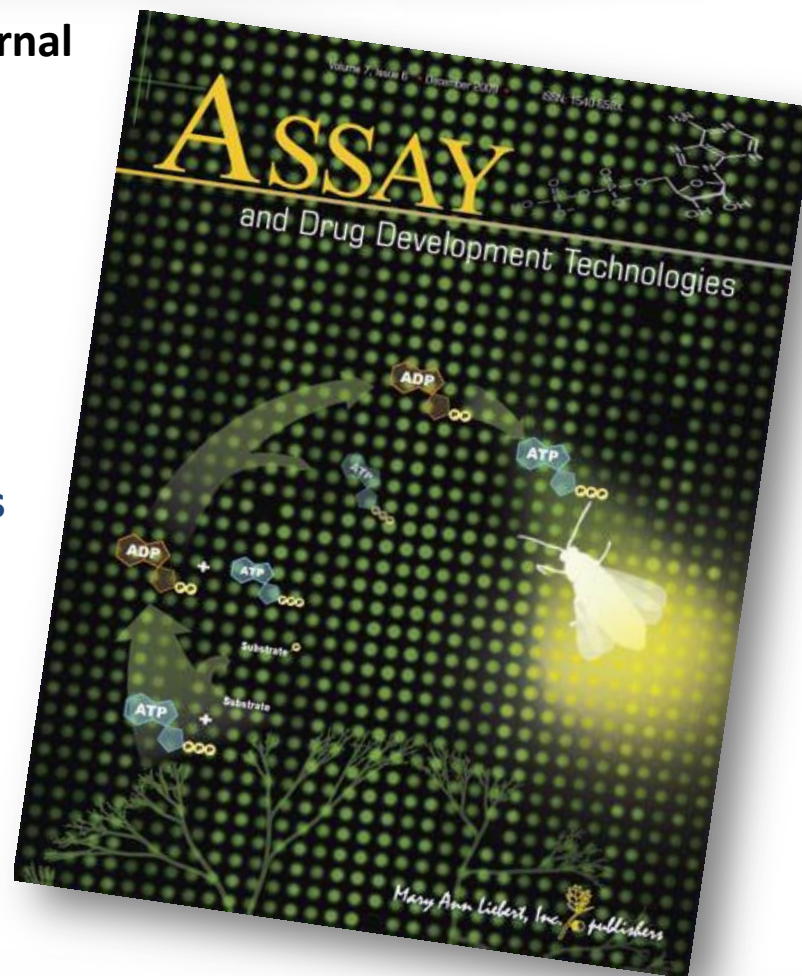
# *ADP-Glo™ Assay in peer review*

Assay and Drug development Technologies Journal  
(December 09 issue) contains  
**6 publications** related to ADP-Glo:

1. ADP-Glo Technology description
2. ADP-Glo for Lipid Kinases
3. uHTS screening with ADP-Glo
4. Comparison of ADP-Glo with other assays
5. Radioactivity vs. ADP-Glo
6. Profiling with ADP-Glo

Free Access here:

<http://www.liebertonline.com/toc/adt/7/6>



# ***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 .



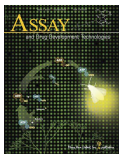
# General features of ADP-Glo™ assay

- ✓ Homogenous, non radioactive and Antibody Free
- ✓ Luminescent assay: Less Compound interference
- ✓ **ADP-Glo™ assay offers so many positive attributes that make it ideal for all stages of drug discovery**
- ✓ Universal: Any kinase/any substrate, ideal for profiling
- ✓ **Basic Research, Primary and secondary screenings and for profiling of lead compounds.**
- ✓ **High dynamic range: High Signal to Background at low % ATP to ADP conversion allows use of lower amount of enzyme during HTS**
- ✓ **Broad range of ATP conc. (linear from 10<sup>-10</sup> to 10<sup>-11</sup> M range) allows distinction between ATP competitive and non competitive inhibitors**
- ✓ High sensitivity: 20nM ADP detected with more than 2.5 fold difference

**Features and Applications: [www.promega.com/kinase](http://www.promega.com/kinase)**

# Summary

## ADP-Glo™ Platform



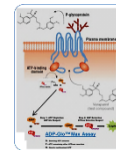
### **1. ADP-Glo™ Kinase Assay**

*Kinase reactions up to 1mM ATP*



### **2. Kinase Enzyme Systems**

*174 complete kinase assays*



### **3. ADP-Glo™ Max Assay**

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

**Next**



### **4. Complete solutions for Profiling**

*Large number of kinases ready to assay  
in a flexible “do it yourself” profiling kits*





**Thank you**

**Questions?**

[hicham.zegzouti@promega.com](mailto:hicham.zegzouti@promega.com)