



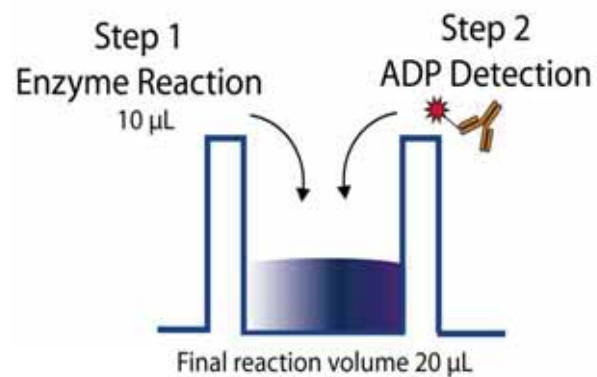
---

*Biology at Work*

# Analyzing Kinase Inhibitor Residence Times using the Transcreener ADP Assay

Meera Kumar  
BellBrook Labs

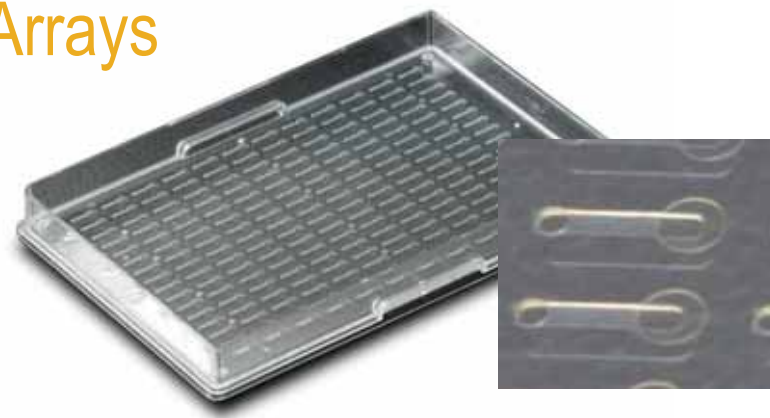
# T RANSCREENER<sup>®</sup> HTS Assays



 BellBrook  
Labs

*Biology at Work*

# I UVO<sup>™</sup> Microconduit Arrays



# T RANSCREENER® Assays

*Four assays, thousands of targets*



*Our Transcreener® assays enable detection of thousands of different nucleotide dependent enzymes.*

Transcreener Assay	Target Families
AMP/ GMP	Phosphodiesterases, Sialyltransferases
UDP	Glycosyltransferases, Ligases (DNA, Amino Acid, Protein)
GDP	GTPases (Small G Proteins), GAPs, Fucosyltransferases
ADP	Kinases (Protein, Lipid, CHO), ATPases, Carboxylases, Helicases
EPIGEN	Methyltransferases, Acetyltransferases

**Stuck on your screen?** Bypass HTS assay development roadblocks with BellBrook's [Transcreener® Assay Development Services](#).

## Webinar Presenters



### Roland Carrillo

- Director of Sales & Business Development
- PhD in Genetics from NC State University
- Email: [roland.carrillo@bellbrooklabs.com](mailto:roland.carrillo@bellbrooklabs.com)
- Phone: (608) 227-4504



*Biology at Work*



### Meera Kumar

- Senior Application Scientist
- MS in Molecular Biology from San Jose State University
- Email: [meera.kumar@bellbrooklabs.com](mailto:meera.kumar@bellbrooklabs.com)
- Phone: (608) 227-4506

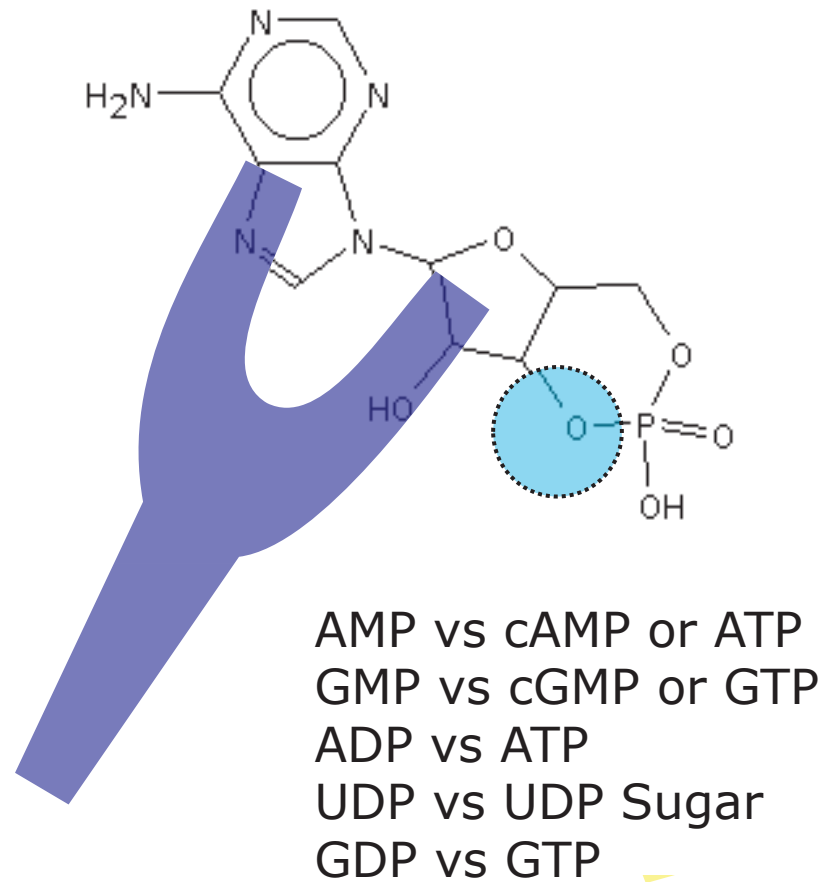
Website: [www.bellbrooklabs.com](http://www.bellbrooklabs.com)

- Overview of Transcreener technology.
- Overview of Drug Residence Times.
- Jump dilution Method using Transcreener assays
- Data Analysis
- Conclusions

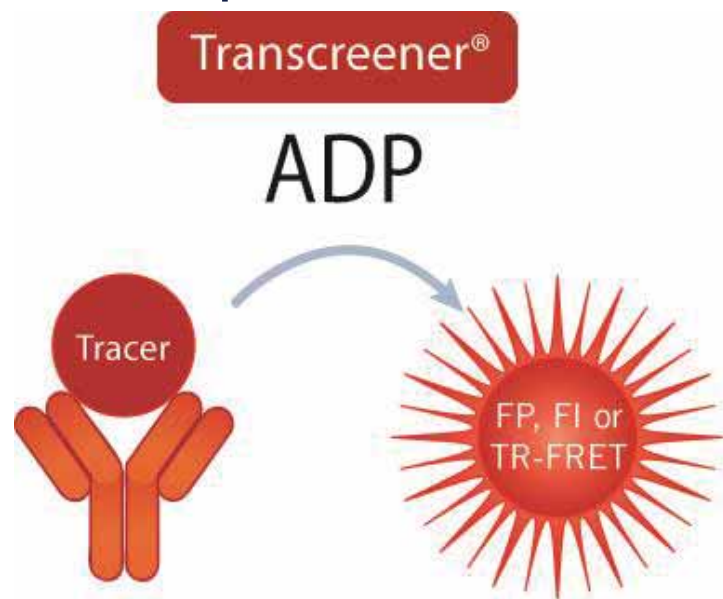
# Transcreener Technology

# Selective Nucleotide Detection

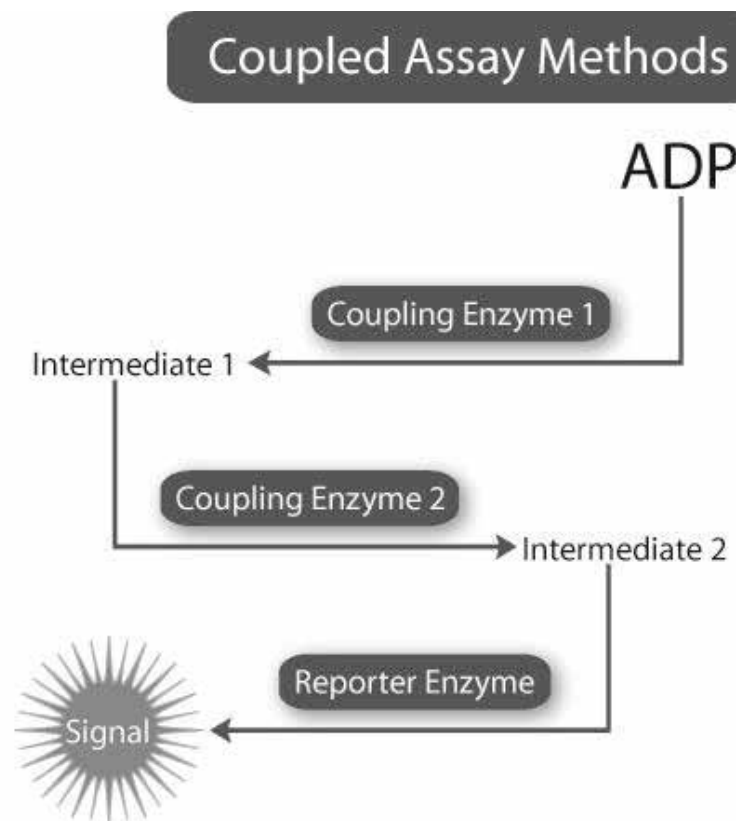
The Transcreener<sup>®</sup> assays rely on antibodies that are able to differentiate between nucleotides on the basis of subtle structural differences. This makes detection of the product nucleotide possible even in the presence of excess substrate nucleotide.



# Direct Detection of Nucleotides decreases potential for interference



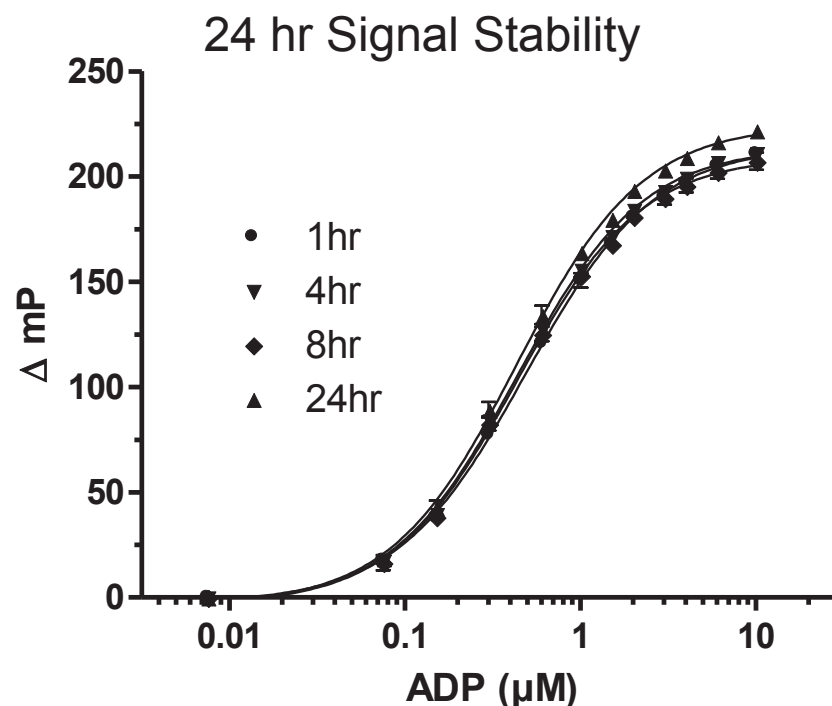
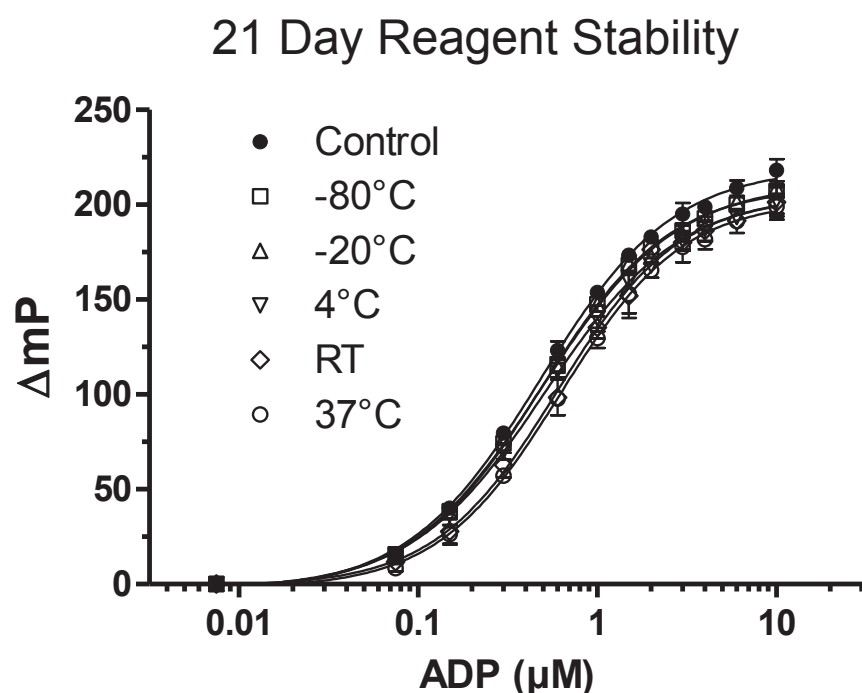
The Transcreener<sup>®</sup> assay relies on **direct detection** of ADP. Binding of tracer to antibody causes a change in fluorescence. There are just two components, and no intermediate steps.



All other ADP assays use **indirect detection** and are more complex. In a series of enzymatic steps, UDP is converted to a detectable product. Each step is subject to inhibition by library compounds.

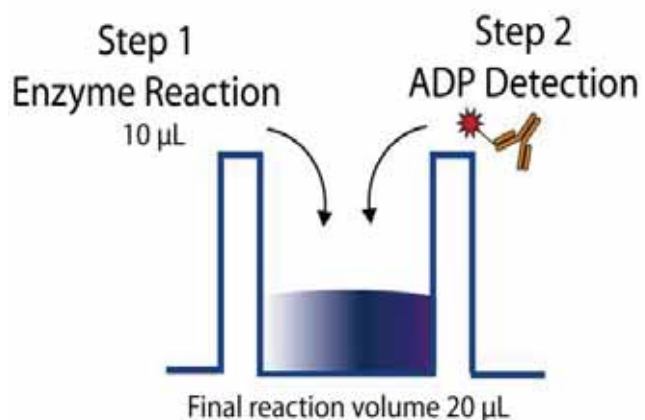


# Transcreener ADP<sup>2</sup> FP Assay: Overnight Reagent and Signal Stability

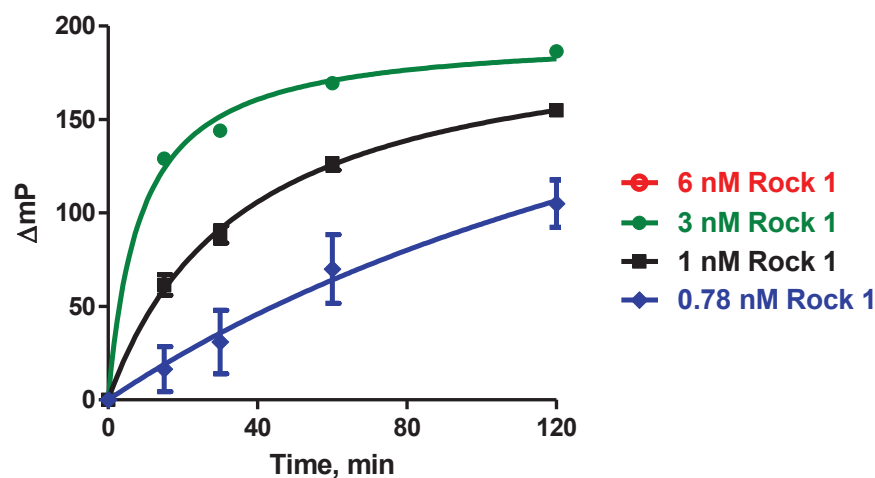


Standard curves for conversion of 10uM ATP to ADP demonstrate the outstanding stability of Transcreener detection reagents prior to addition to reaction and the stability of the signal following addition to kinase reactions. Data is for the FP assay, the FI and TR-FRET assays also have at least overnight reagent and signal stability. This provides outstanding flexibility for automated HTS platforms, especially with large numbers of plates, where there may be a lag between addition of detection reagent and plate-reading.

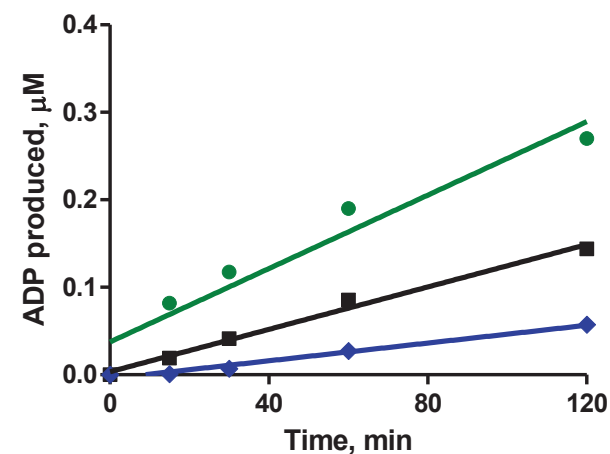
# True mix and read detection format, endpoint or continuous detection.



Transcreener assays are true mix-and-read format, with the enzyme quenching and ADP detection components added as a single reagent for endpoint assays. The assay can also be used in a continuous detection mode, which makes optimization of enzyme reactions simpler and reduces reagent consumption.



Standard Curve



Reaction is linear for 2 hrs with up to 1nM Rock 1; higher concentrations exceed initial velocity conditions.

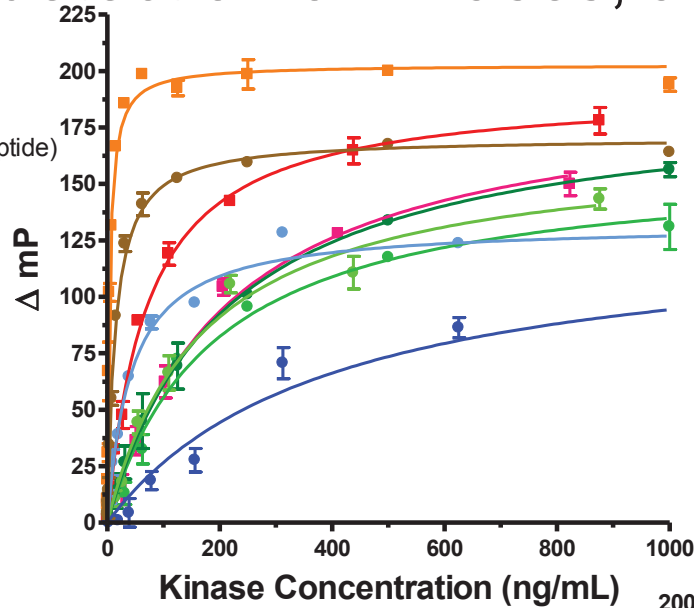
# Transcreener ADP<sup>2</sup> Assay: Universal detection of kinases, acceptors

## Peptide Substrates

- Abl1 (Abltide)
- AKT/PKB (Akt/SKG peptide)
- PKA (kemptide)

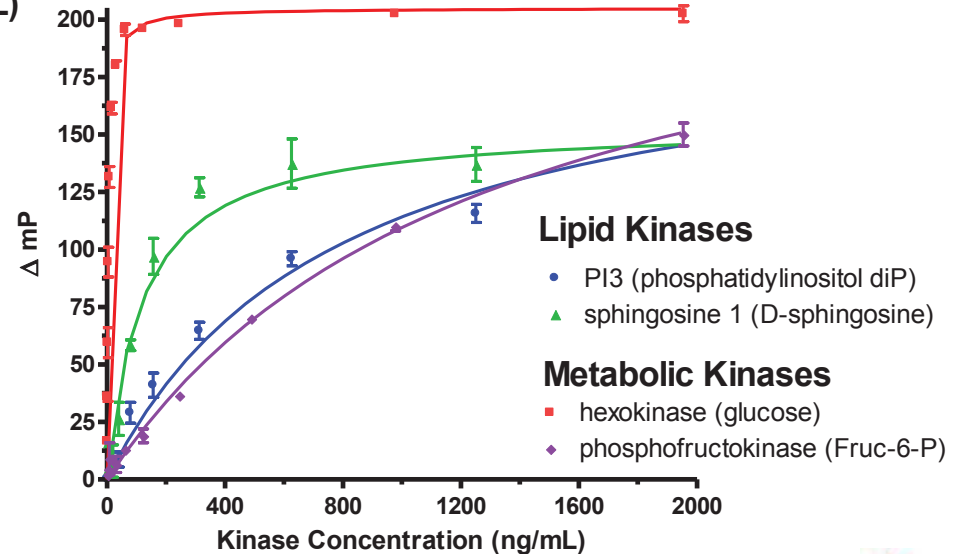
## Protein Substrates

- cdk5/p35 (Histone 1)
- COT (MEK1)
- p38alpha (MBP)
- PKA (histone H1)
- PKA (MBP)
- RAF1 (MEK1)



Universal detection means you can use any kinase and any acceptor substrate, including native proteins, which provide a more physiologically relevant measure of kinase activity.

Universal detection means straightforward detection of lipid and carbohydrate kinases in addition to protein kinases.



## Lipid Kinases

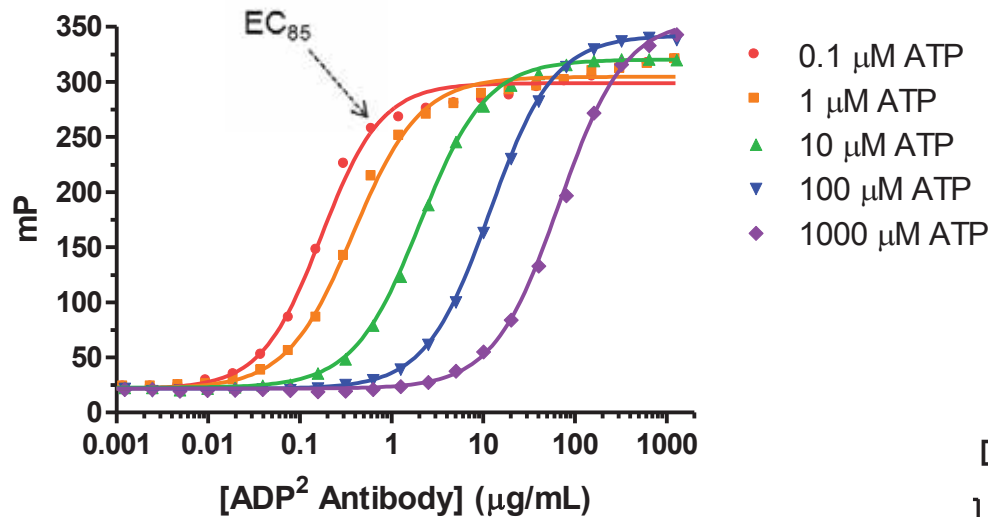
- PI3 (phosphatidylinositol diP)
- ▲ sphingosine 1 (D-sphingosine)

## Metabolic Kinases

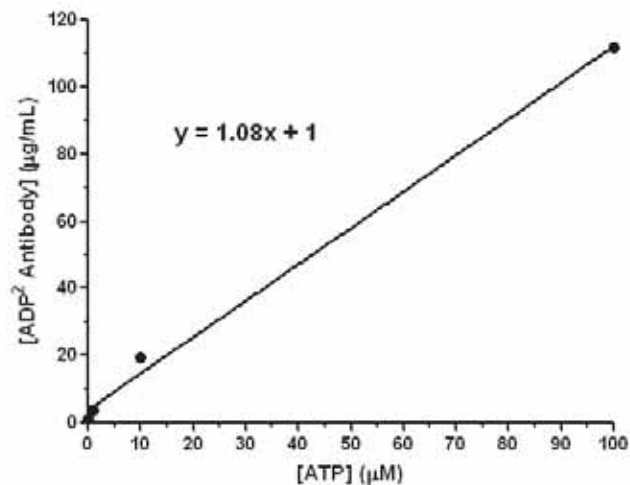
- hexokinase (glucose)
- ◆ phosphofruktokinase (Fruc-6-P)

# Dynamic Range can be Tuned for Robust Initial Velocity Detection at any ATP Concentration

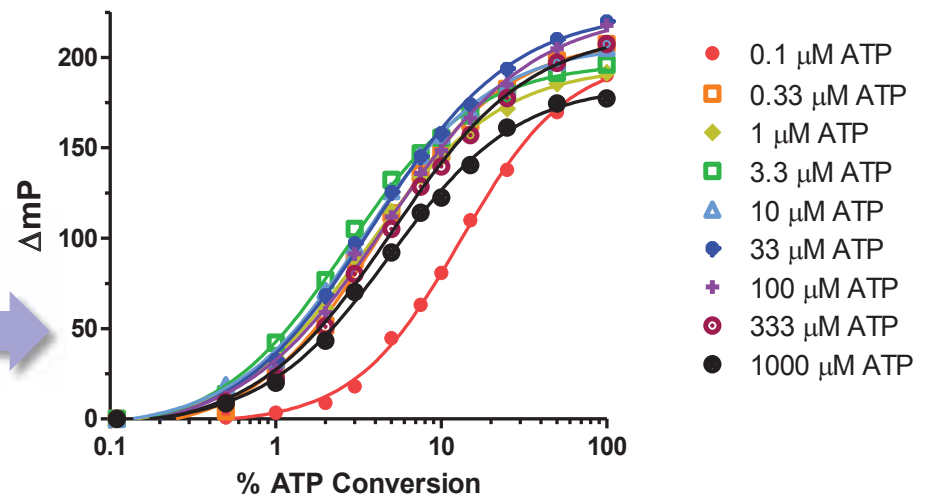
Ab titrations in the presence of ATP



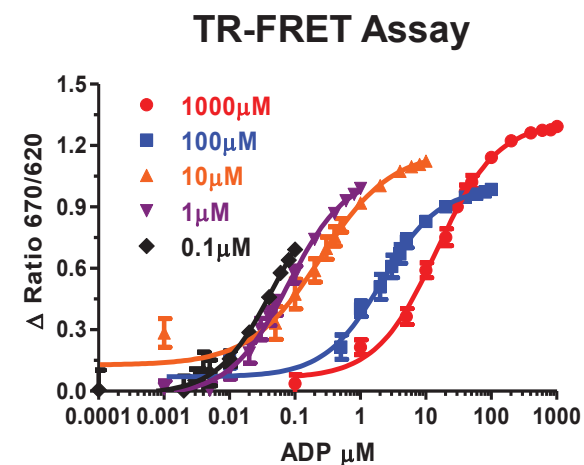
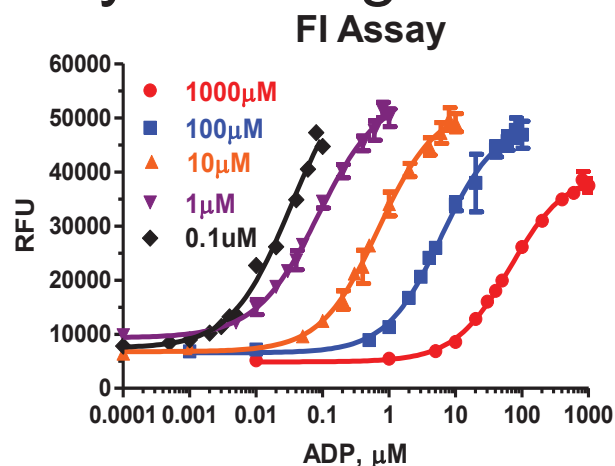
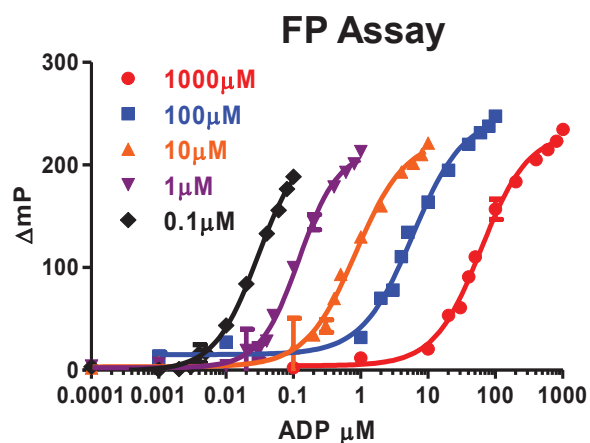
The ADP antibody concentration sets the dynamic range of the assay. EC<sub>85</sub> is optimal for initial velocity detection; ie, a good signal @ 10% conversion of ATP.



ATP/ADP Standard Curves  
[ADP<sup>2</sup> Ab] @ EC<sub>85</sub> for indicated [ATP]



# Transcreener ADP<sup>2</sup> Assay: Greater sensitivity reduces enzyme usage



	1 $\mu M$ ATP/ADP standard curve		10 $\mu M$ ATP/ADP standard curve		100 $\mu M$ ATP/ADP standard curve	
	Z' at 10% Conv	LLD ( $\mu M$ )	Z' at 10% Conv	LLD( $\mu M$ )	Z' at 10% Conv	LLD( $\mu M$ )
Transcreener FP	0.86	0.02 $\pm$ 0.07	0.85	0.01 $\pm$ 0.12	0.89	1.0 $\pm$ 0.3
Transcreener TR-FRET	0.71	0.10 $\pm$ 0.06	0.72	0.10 $\pm$ 0.09	0.72	1.0 $\pm$ 0.3
Transcreener FI	0.92	0.03 $\pm$ 0.01	0.88	0.05 $\pm$ 0.04	0.92	0.5 $\pm$ 0.4
Luc-ADP Detection Assay	ND	0.40 $\pm$ 0.87	0.30	0.50 $\pm$ 0.32	0.62	5.0 $\pm$ 0.7
Luc-ATP Depletion Assay	ND	0.25 $\pm$ 0.40	ND	1.50 $\pm$ 0.30	0.52	7.0 $\pm$ 0.6

Standard curves for conversion of ATP to ADP demonstrate robust detection at 10% conversion starting at ATP concentrations from 0.1 to 1,000  $\mu M$ . By comparison, Luciferase-based ADP and ATP detection methods require greater conversion of ATP, especially at lower ATP concentrations; this translates into higher enzyme consumption.

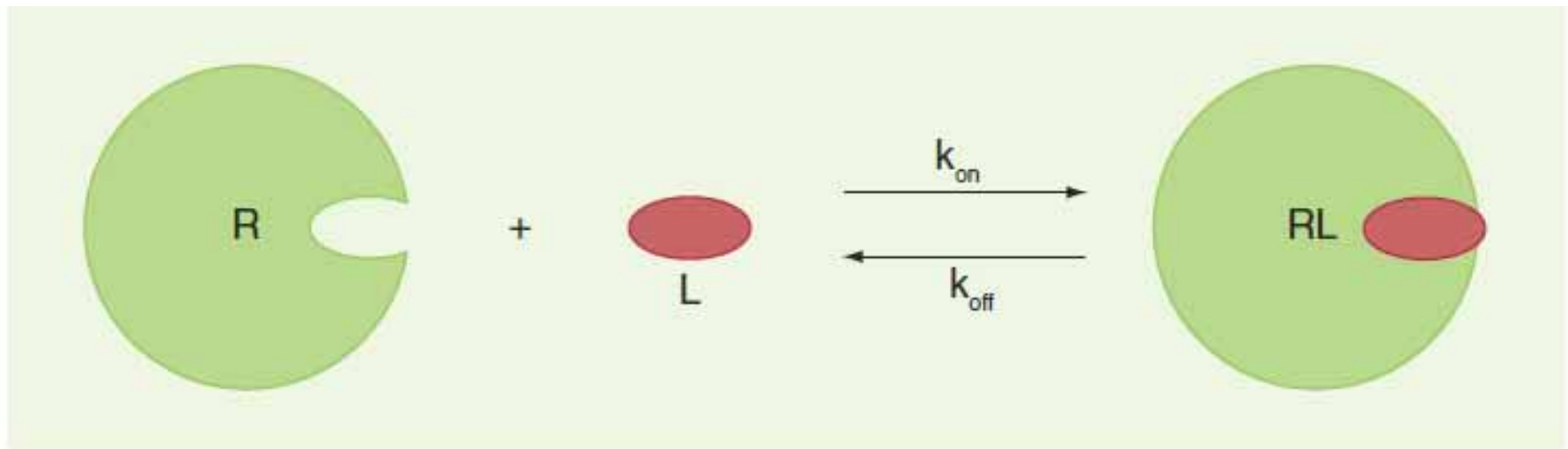
# Residence Time

# What is Drug Residence Time?

---

- **Drug–target residence time:** Lifetime of the binary complex between a drug molecule and a macromolecular target.
- Experimentally, the residence time is measured as the reciprocal of the rate constant for drug–target complex dissociation ( $1/k_{off}$ ).
- Drug potency is a function of drug occupancy on a target (residence time) as well as its affinity ( $IC_{50}$ ).

## The Static View of Drug-Target Interactions

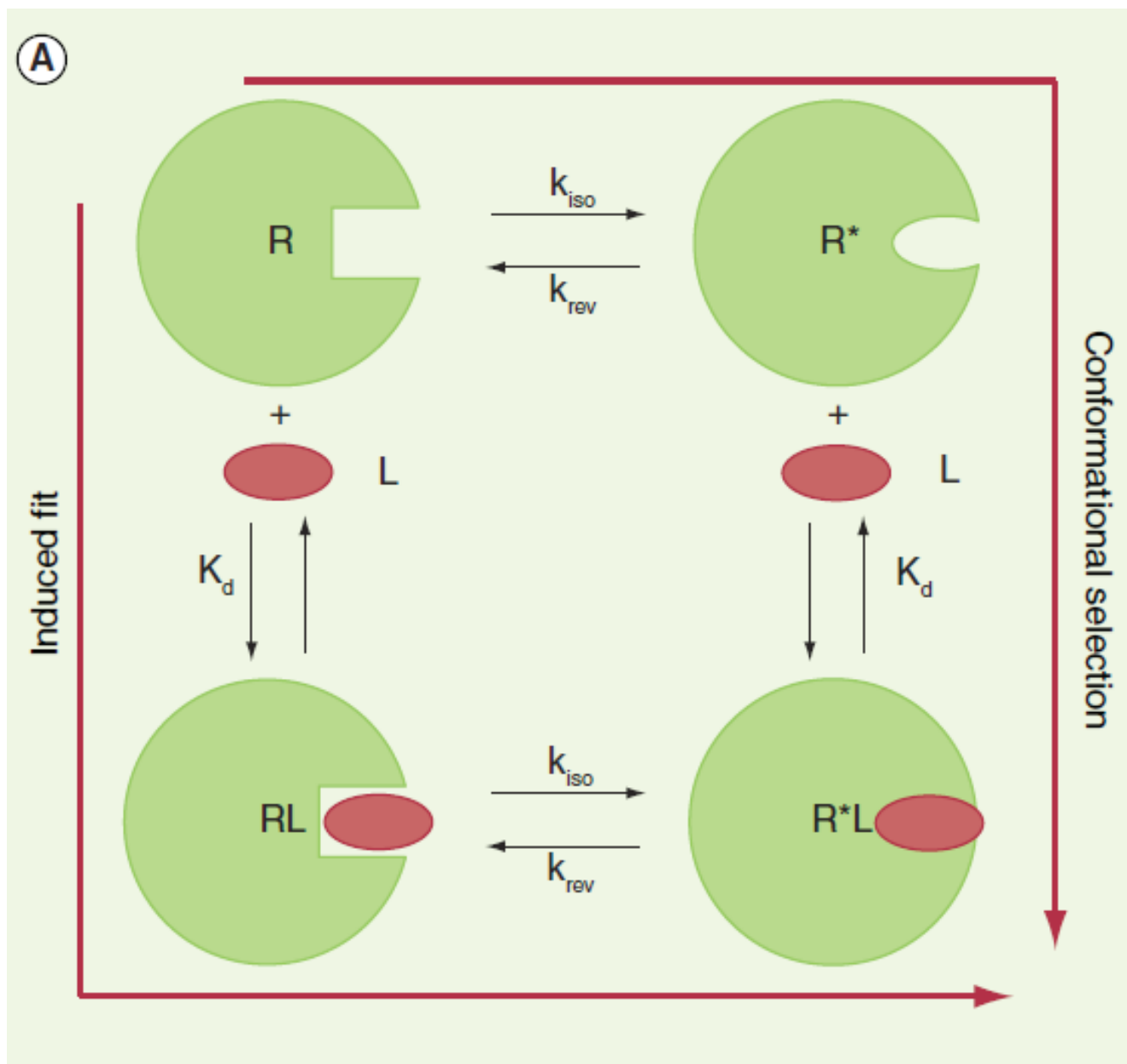


**Figure 1. Static lock-and-key model of receptor–ligand binding in which  $k_{on}$  and  $k_{off}$  each occur in a single kinetic step.**

$k_{on}$ : Complex association rate constant;  $k_{off}$ : Complex dissociation rate constant.



# EI Formation: Conformational Adaptation in Drug-Target Interactions



# Methods available to measure Drug residence times

---

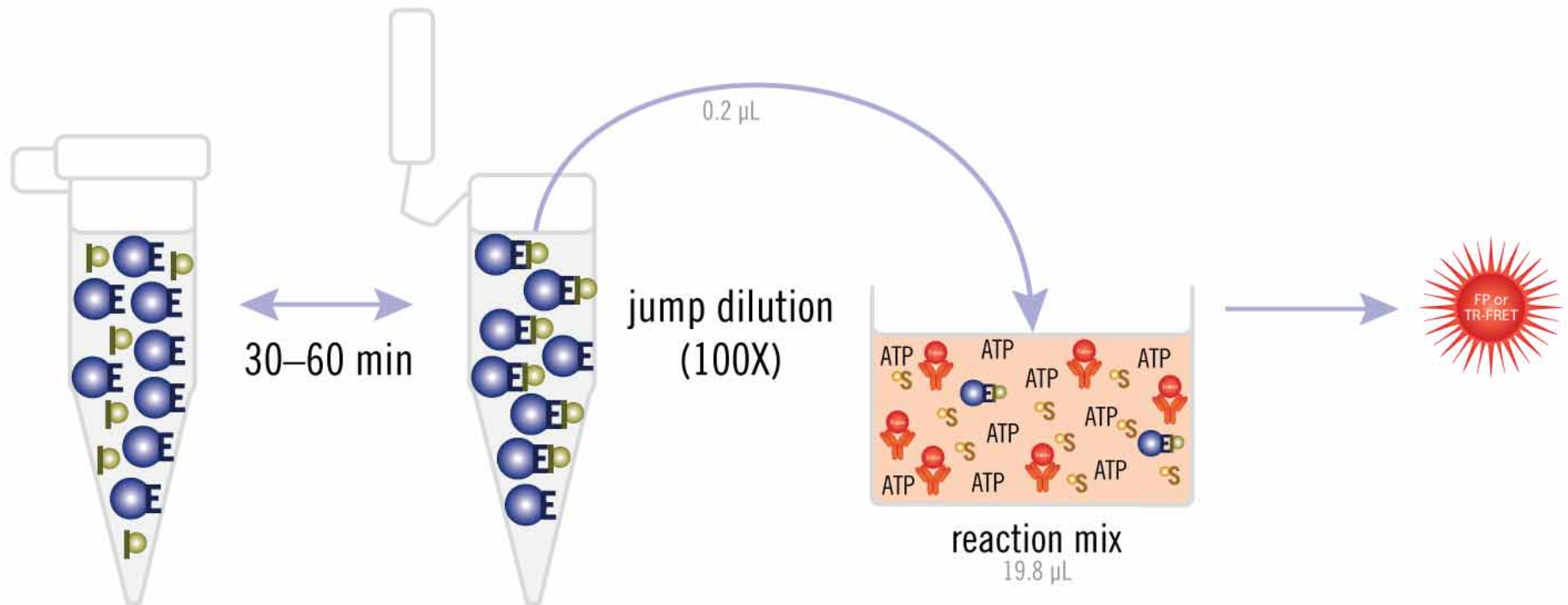
- Surface Plasmon Resonance
  - Requires that you immobilize proteins on a sensor surface.
  - Requires Expensive and specialized instruments, accessories and consumables to record measurements
  - Low Throughput
  - Only provides binding kinetics information
- Changes in intrinsic fluorescence
- Non Equilibrium reaction kinetics.
- Fluorescence Binding assays-Not activity based assays

# Streamlined protocol for residence time determination

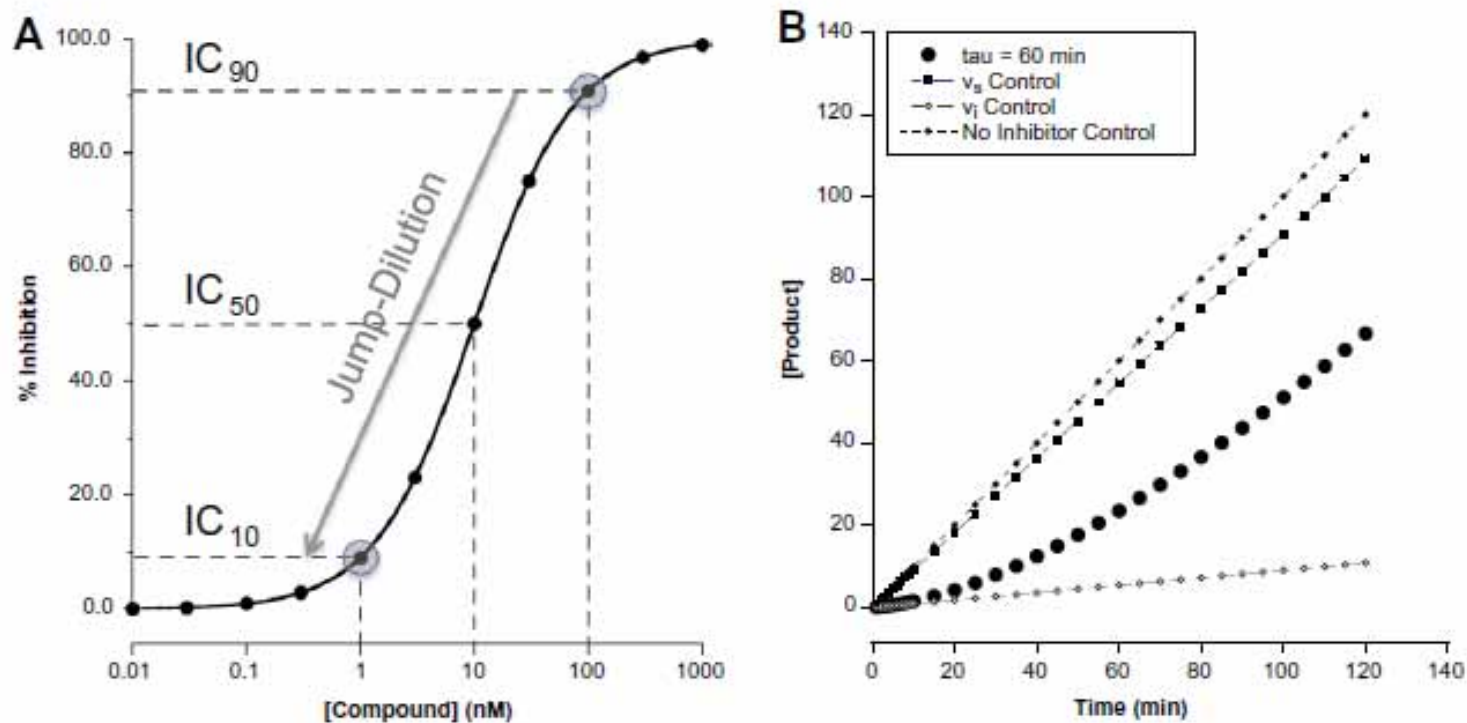
---

- Determine  $EC_{80}$  of target concentration.
- Determine  $IC_{50}$  values of the compounds.
- Preincubate  $10^* IC_{50}$  compound +  $100^*[EC_{80}]$
- Jump Dilution: Dilute 100X into enzyme reaction and Transcreener detection reagents.
- Continuously read over time and monitor the change in fluorescence.
- Data Analysis and calculate residence times.

# Jump dilution protocol for determining residence times



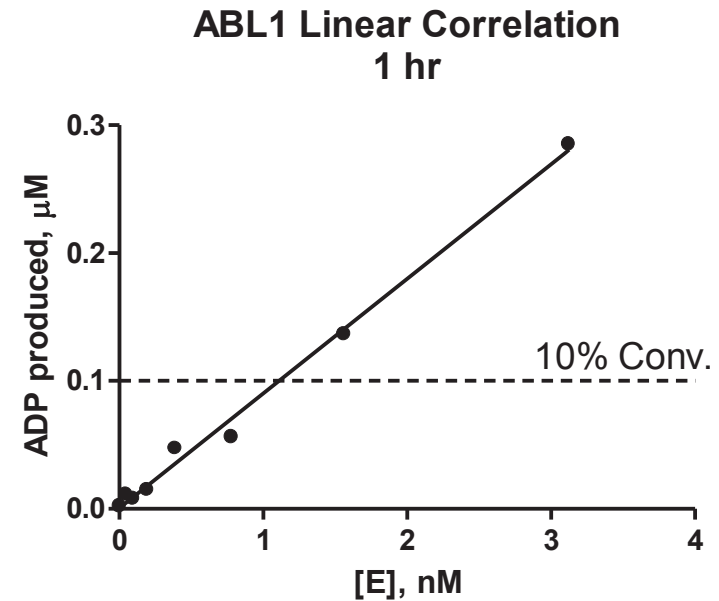
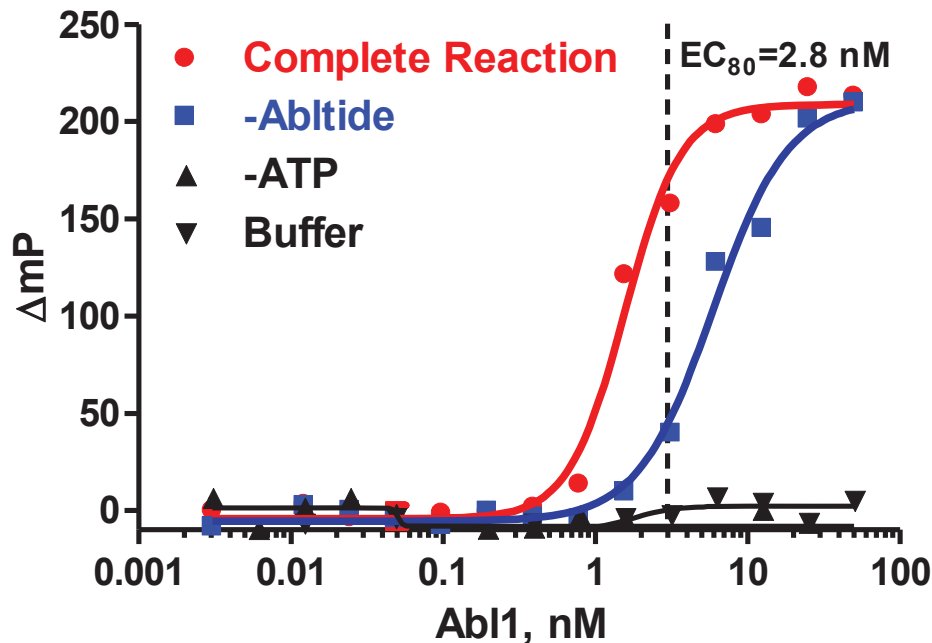
# Jump dilution



**Fig. 1.** (A) Typical concentration–response plot for target inhibition by a compound of interest illustrating the starting and ending concentrations of compound, relative to the IC<sub>50</sub>, used in a jump dilution experiment. (B) Example of a jump dilution progress curve for control samples and for a test compound displaying a residence time of 60 min. Both panels of this figure are adapted and redrawn from Copeland [2].

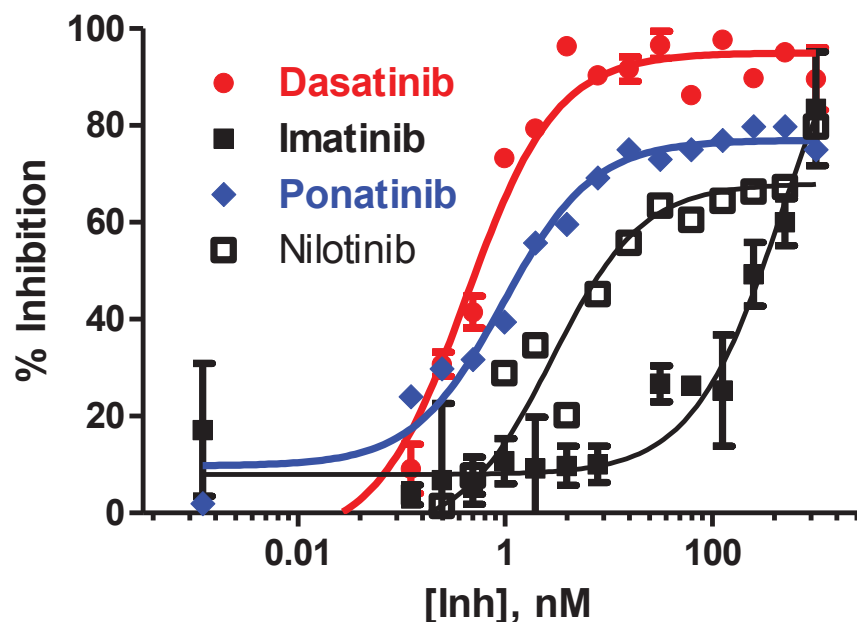
*Jump dilution analysis / R.A. Copeland et al. / Anal. Biochem. 416 (2011) 206–210*

# Case study 1: Abl1



**Figure 3. Determining the Optimal concentration of Abl1 and Drugs for the “Jump Dilution” Experiment.** **A.** Titration of Abl1 enzyme, performed in the presence of 1  $\mu M$  ATP and 10  $\mu M$  Abltide, determined an optimal  $EC_{80}$  concentration as 2.8 nM. **B.** Linear relationship between Abl1 concentration and ADP formation. The polarization values were converted into ADP (product formed) using a standard curve set up under similar conditions.

# Determining the optimal drug concentration for residence time experiments



	Dasatinib	Imatinib	Ponatinib	Nilotinib
IC <sub>50</sub> (nM)	0.4209	474.9	0.9547	2.847

A dose response curve for Abl1 enzyme (2.8 nM) in the presence of 1  $\mu$ M ATP and 10  $\mu$ M Abltide enabled determination of IC<sub>50</sub> values of 0.42 nM (dasatinib), 475 nM (imatinib), 2.8 nM (Nilotinib) and 0.95 nM (ponatinib).

# Preincubation for [EI] Complex formation

- Enzyme for [EI]
  - $100 \times EC_{80} = 2.8 \text{ nM} \times 100 = 280 \text{ nM}$
- Drug for [EI]
  - $10 \times IC_{50}$  (Dasatinib) = 45 nM
  - $10 \times IC_{50}$  (Imatinib) = 45  $\mu\text{M}$
  - $10 \times IC_{50}$  (Ponatinib) = 100 nM
  - $10 \times IC_{50}$  (Nilotinib) = 25  $\mu\text{M}$

**Preincubation Protocol.** The amount of Abl1 used for preincubation is at least  $50 \times EC_{80}$ , so that it still gives a robust signal after dilution. The inhibitor concentration is  $10 \times IC_{50}$ , such that the enzyme is saturated with the inhibitor. The mix was incubated for 1 hour at room temperature to insure formation of the EI complex.



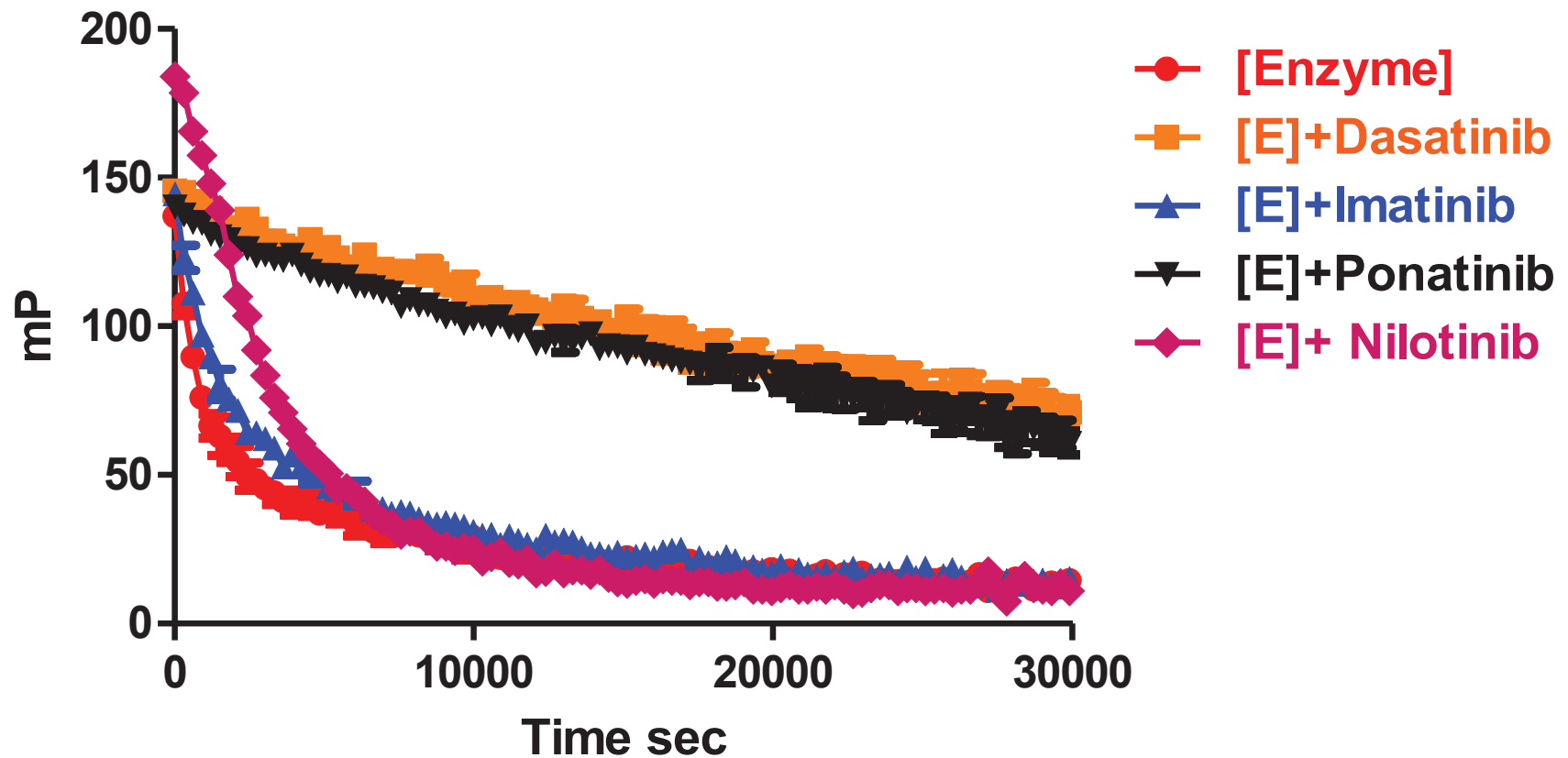
# Jump dilution

---

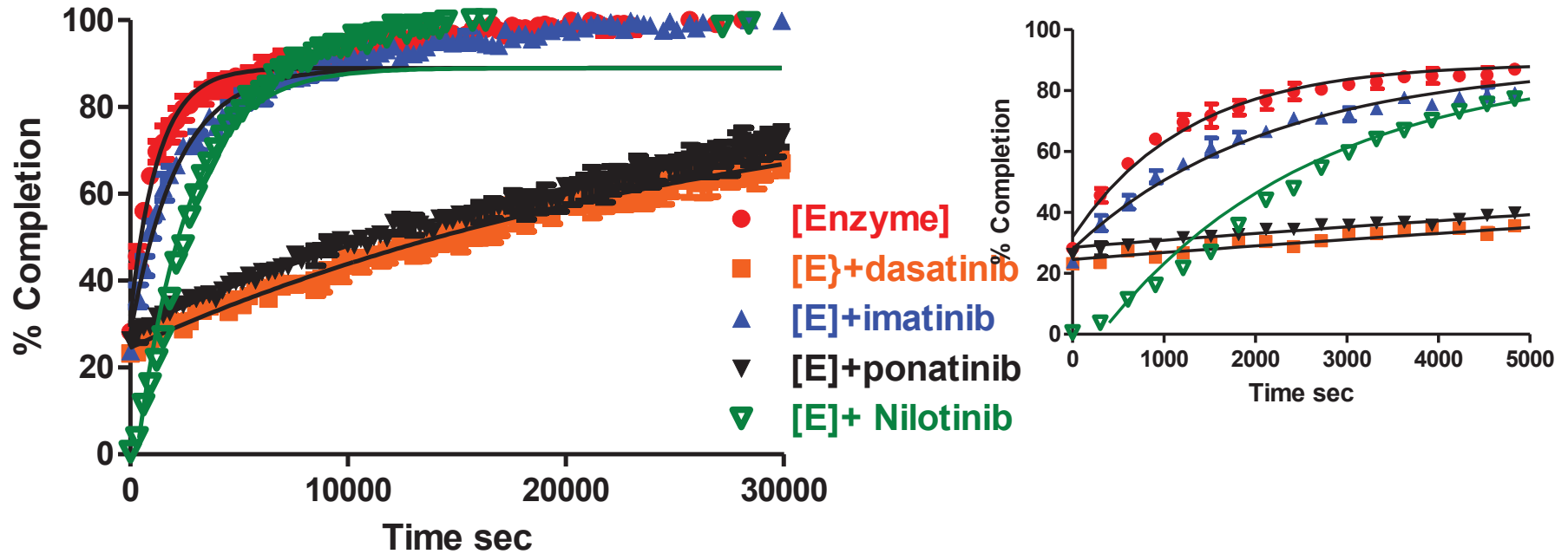
- The dilution of the EI complex is done such that the inhibitor concentration is at least 100-fold below its  $IC_{50}$  value.
- A 100-fold dilution (0.2  $\mu$ L of Abl/Inhibitor mixture into 19.8  $\mu$ L of Detection Mixture) was made in an LV-384 well plate.
- The Detection mixture comprised of 5  $\mu$ M ATP and 10  $\mu$ M Abltide in kinase buffer also contained 2 nM tracer and 3.2  $\mu$ g/mL of ADP<sup>2</sup> antibody.
- The plate was mixed well and read kinetically every 5 minutes for 4 hours in a Tecan Safire plate reader using the fluorescent polarization mode with EXC at 630 nm and EMS at 670 nm.

# Raw data showing mP values

Raw data: mP decreases as ADP is produced.



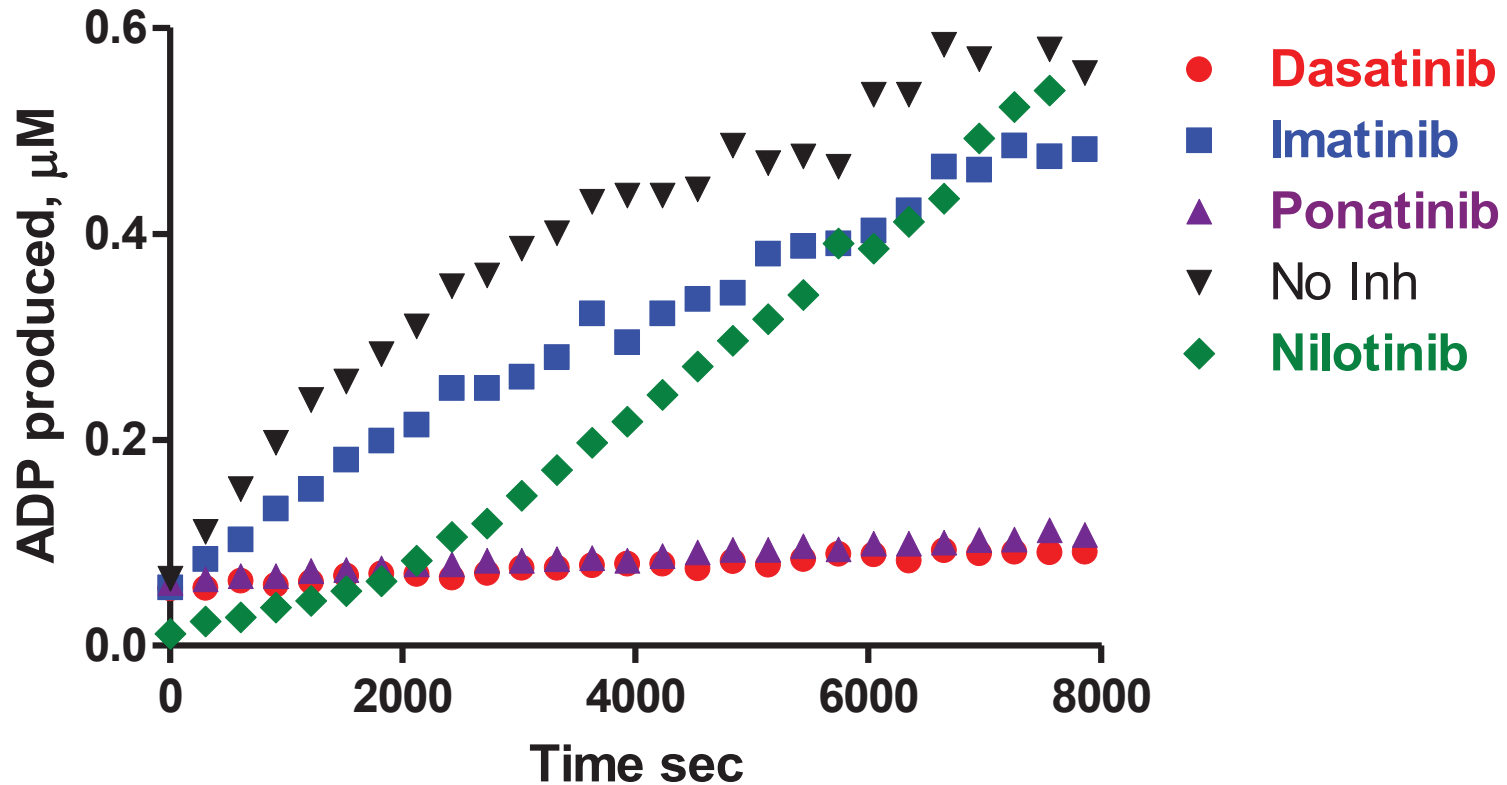
# Data analysis



	[Enzyme]	[E]+dasatinib	[E]+imatinib	[E]+ponatinib	[E]+ nilotinib
K	0.0007879	3.561e-005	0.0004612	3.899e-005	0.0004319
Half Life	879.8	19467	1503	17776	1605
Tau	1269	28085	2168	25645	2315

Data analysis: One phase decay using raw data.  
 Normalized data based on the following controls-  
 a) Inh without enzyme =0% activity.  
 b) Enzyme after completion= 100% activity.

# Converting mP values to ADP (product formed)

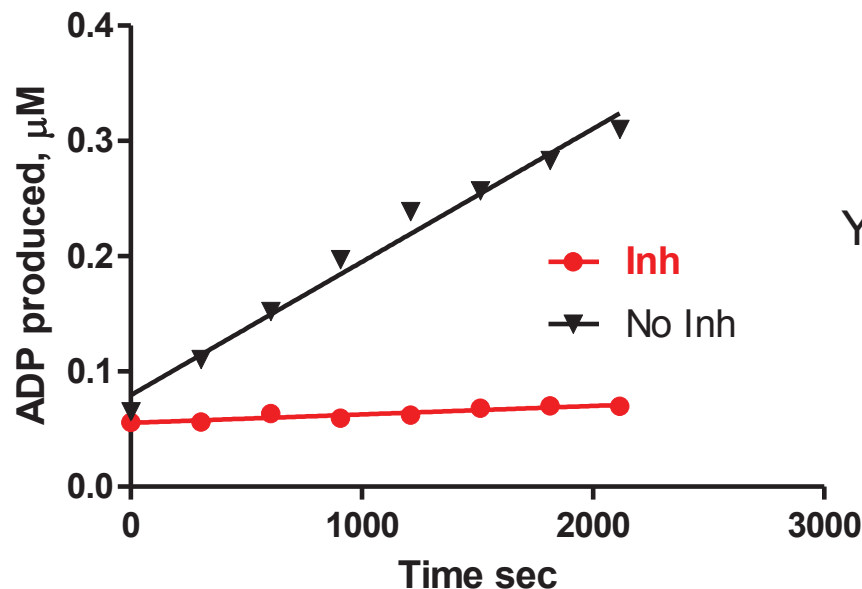


# How to determine residence time from the product formed

- Determine  $K_{off}$  values by fitting the curve to equation-

$$P = V_s t + (V_0 - V_s) \frac{(1 - e^{-kt})}{k}$$

- $V_0$  represents fully inhibited enzyme velocity,  $V_s$  represents the uninhibited enzyme velocity.

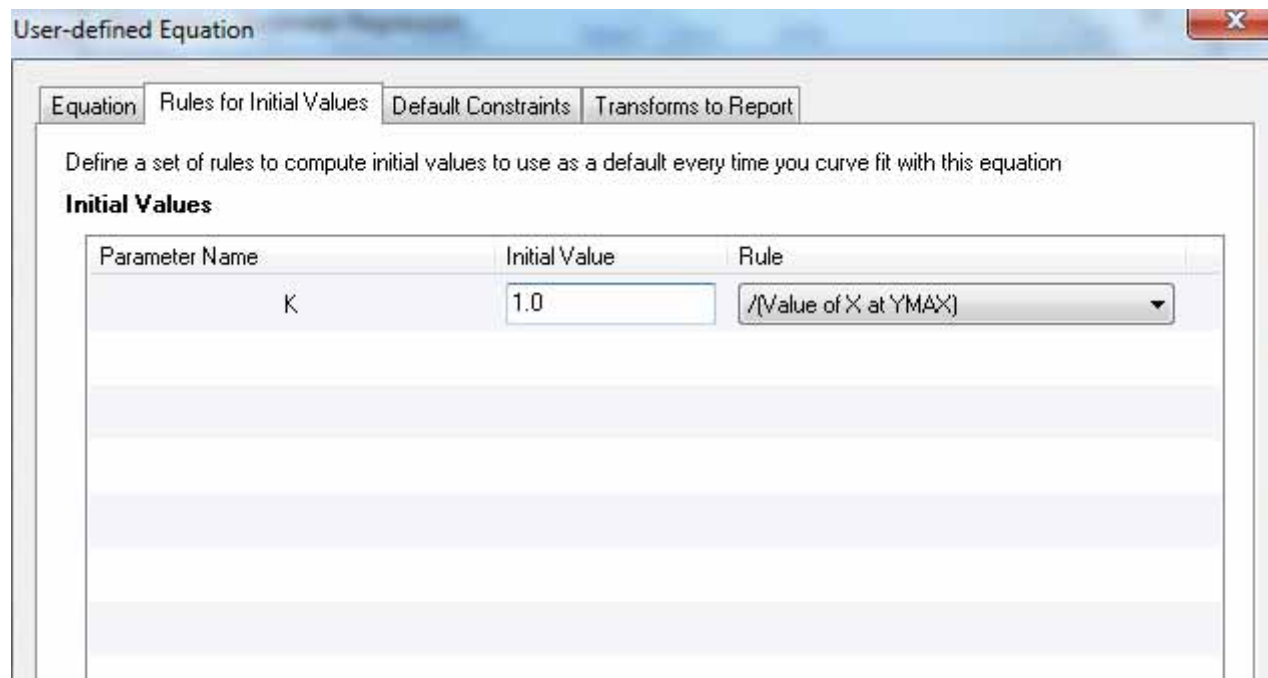


$$Y = 0.0001156X - 0.000108243/K(1 - \exp(-K*X))$$

$V_s$	$0.000007357 \pm 0.000001187$
$V_0$	$0.0001156 \pm 0.000006600$

# Curve fitting to this equation in Graph Pad Prism

- Open the analysis window in Graph Pad prism and select new equation.
- Enter the new equation.
- Set “rules”- Define K as 1/Value of X at Ymax



User-defined Equation

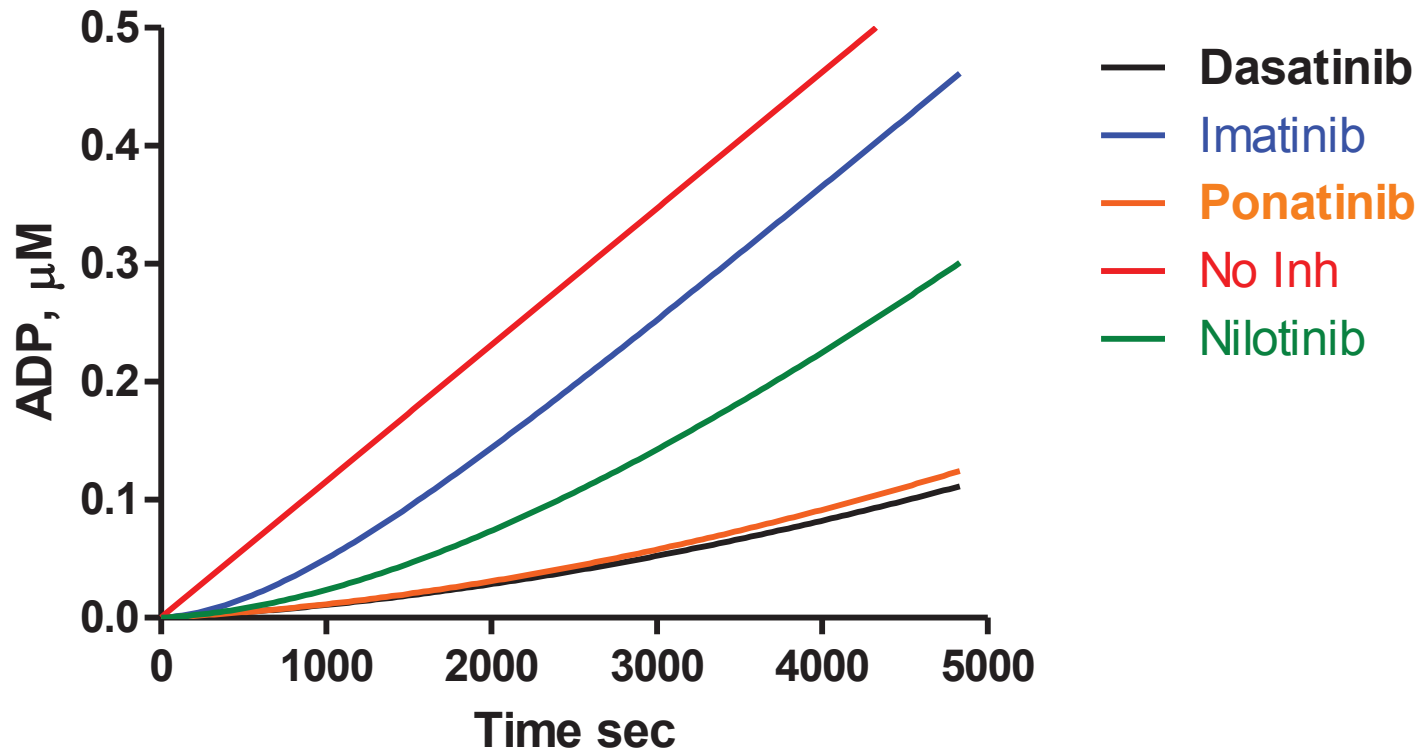
Equation Rules for Initial Values Default Constraints Transforms to Report

Define a set of rules to compute initial values to use as a default every time you curve fit with this equation

**Initial Values**

Parameter Name	Initial Value	Rule
K	1.0	/(Value of X at YMAX)

# Determining residence time



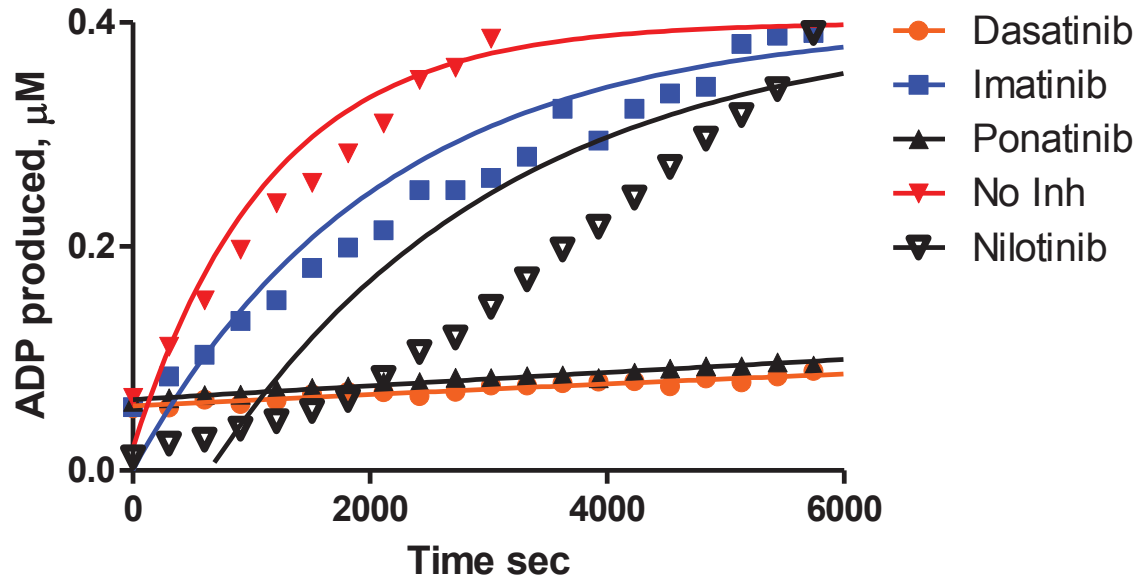
	Dasatinib	Imatinib	Ponatinib	Nilotinib
<b>K</b>	<b>0.00006669</b>	<b>0.001106</b>	<b>0.00007959</b>	<b>0.0003377</b>
<b>Tau, min</b>	<b>250</b>	<b>15</b>	<b>209</b>	<b>50</b>

Lit values

	Tau, min <sup>-1</sup>
Dasatinib	837
Ponatinib	500
Imatinib	14

# Determining residence time using one phase decay

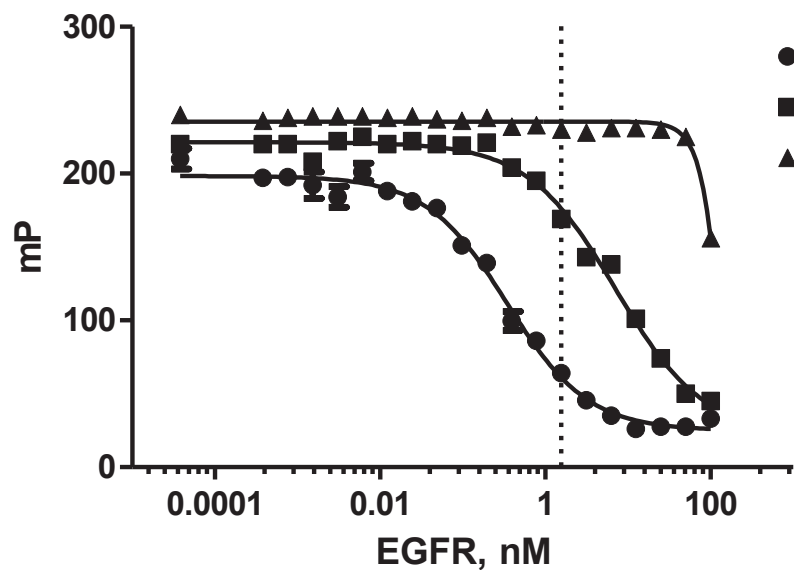
(product formation)



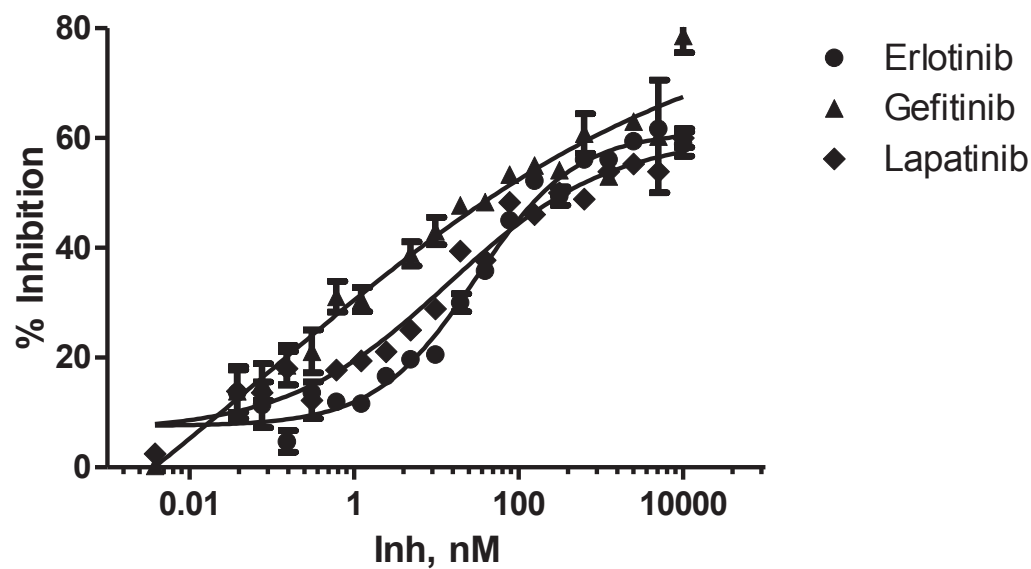
	Dasatinib	Imatinib	Ponatinib	Nilotinib
K	0.00003881	0.0004834	0.00001867	0.0004061
Tau, min	429	34	892	41



# Case Study 2: EGFR



	ATP+Srctide	ATP	Srctide
$EC_{80}$ (nM)	1.655	38.42	~ 1679



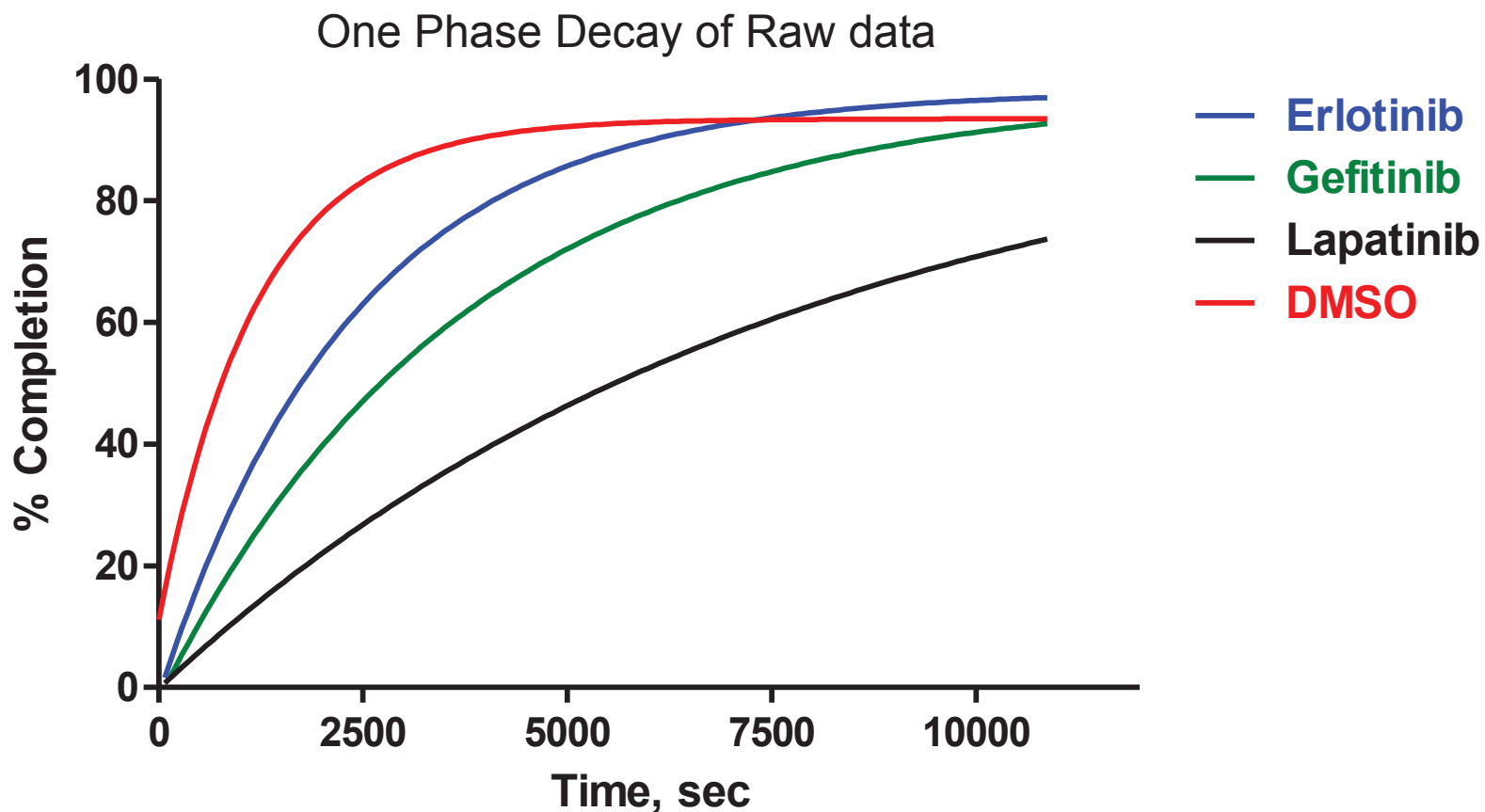
	Erlotinib	Gefitinib	Lapatinib
$IC_{50}$ (nM)	30.12	<1	13.75

# Preincubation for [EI] Complex formation

- Enzyme for [EI]
  - $100 \times EC_{80} = 1.5 \text{ nM} \times 100 = 150 \text{ nM}$
- Drug for [EI]
  - $10 \times IC_{50} \text{ (Erlotinib)} = 300 \text{ nM}$
  - $10 \times IC_{50} \text{ (Gefitinib)} = 100 \text{ } \mu\text{M}$
  - $10 \times IC_{50} \text{ (Lapitinib)} = 140 \text{ nM}$

**Preincubation Protocol.** The amount of EGFR used for preincubation is at least  $100 \times EC_{80}$ , so that it still gives a robust signal after dilution. The inhibitor concentration is  $10 \times IC_{50}$ , such that the enzyme is saturated with the inhibitor. The mix was incubated for 1 hour at room temperature to insure formation of the EI complex.

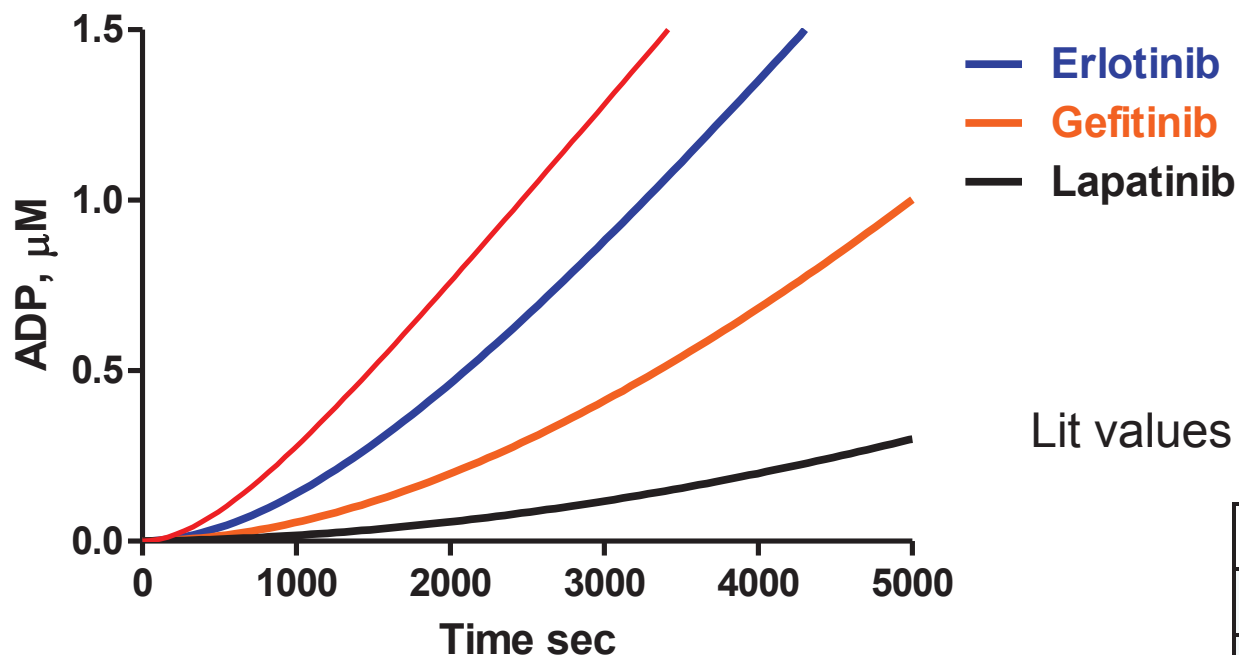
# Residence time for EGFR inhibitors using Transcreener assay (Normalization of raw data)



	Erlotinib	Gefitinib	Lapatinib
<b>K</b>	<b>0.0004181</b>	<b>0.00027</b>	<b>0.0001284</b>
<b>Tau(min)</b>	<b>40</b>	<b>62</b>	<b>130</b>

# Residence time for EGFR inhibitors using product formed

$$Y=0.0005288X-(0.000523/K)*(1-\exp(-K*X))$$



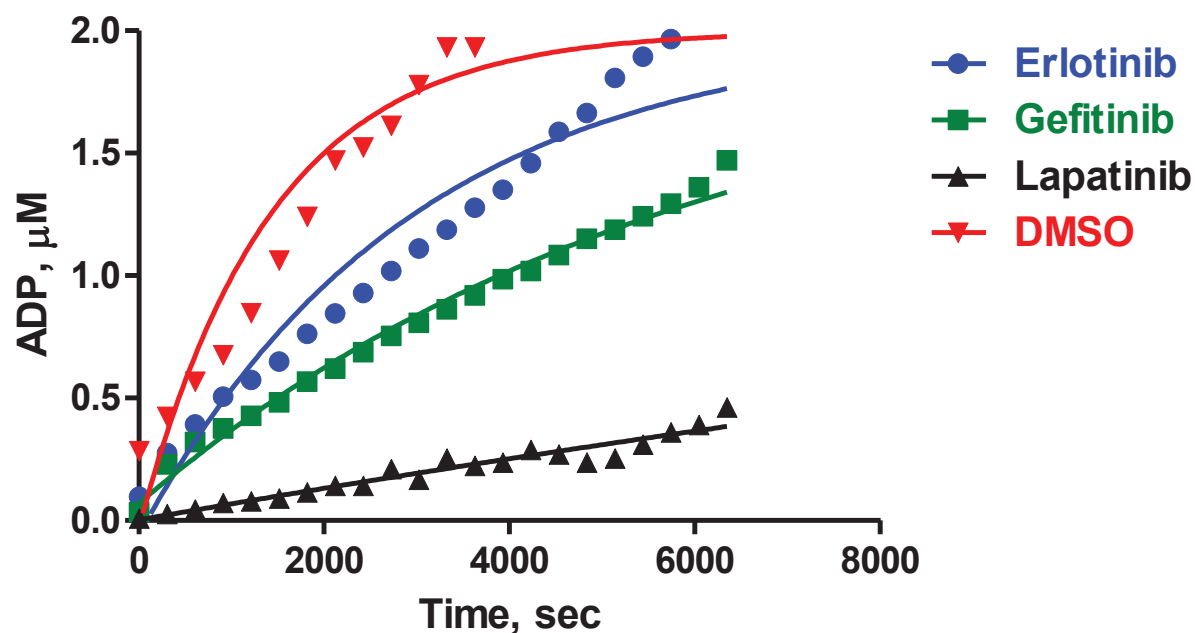
Lit values from Lanthascreen assay

	Tau, min
Gefitinib	14
Erlotinib	10
Lapatinib	52

	Erlotinib	Gefitinib	Lapatinib
K	0.00063	0.0002031	0.00004458
Tau(min)	27	82	374

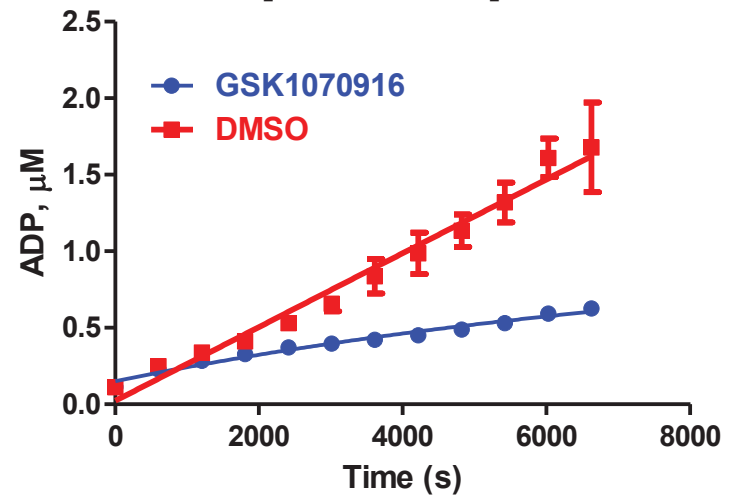
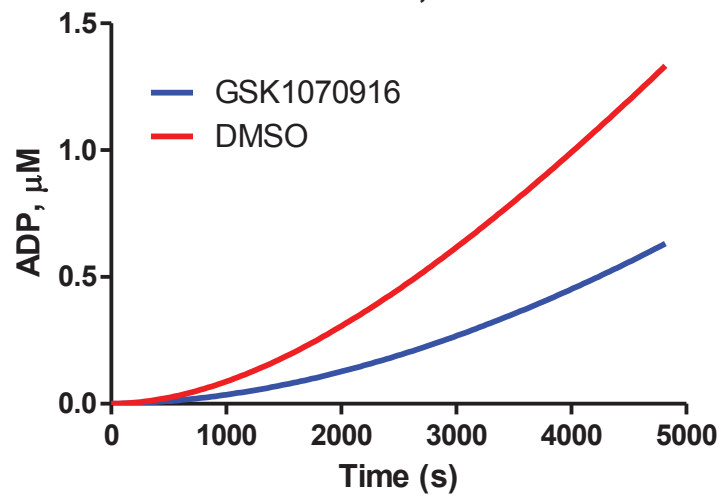
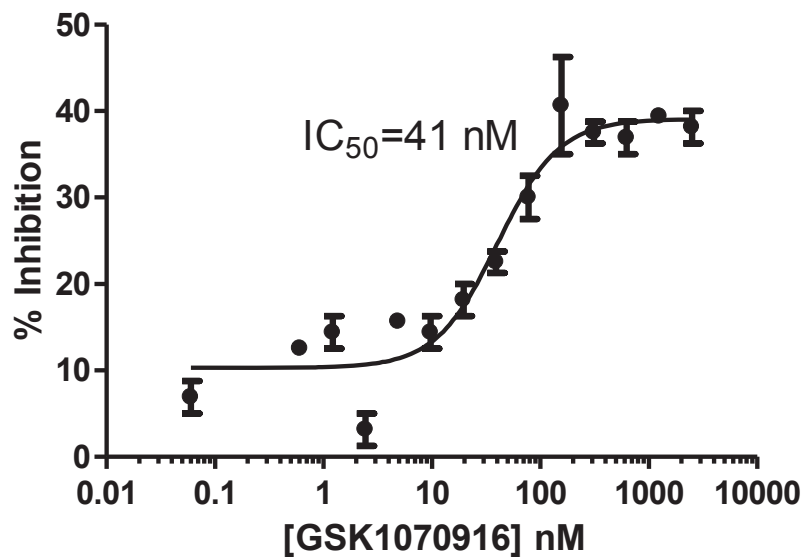
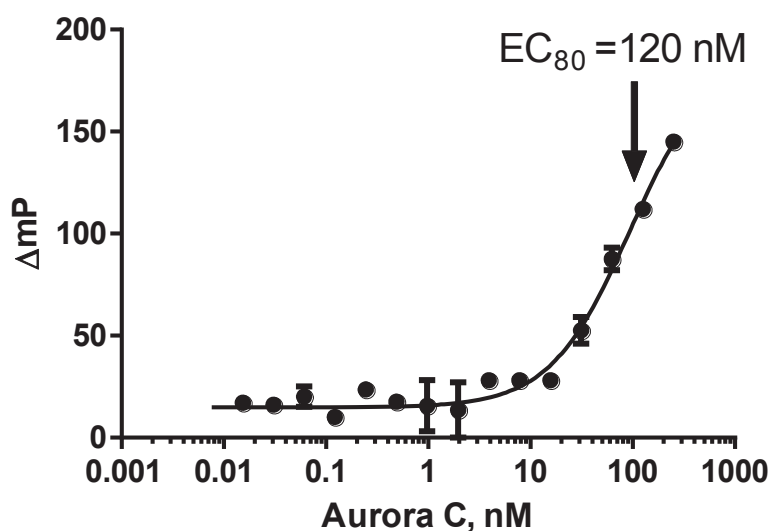
# Residence time for EGFR inhibitors using Transcreener assay

One phase Decay of product formed



	Erlotinib	Gefitinib	Lapatinib
K	0.0003412	0.0001693	0.00003341
Tau(min)	49	98	498

# Aurora C: Residence time for GSK-1070916



	GSK1070916
K	0.0001193
Tau (min)	140

	GSK1070916
K	0.0001
Tau (min)	154

Lit value-280 min  
 Biochem. J. (2009) **420**, 259–265

# Conclusions

---

- The Transcreener<sup>®</sup> ADP<sup>2</sup> Assay relies on highly selective immunodetection of ADP with FP, TR-FRET, and FI readouts. Direct detection has advantages over enzyme-coupled assays in terms of simplicity of use and resistance to compound interference.
- The assays can be tuned for detection of kinases or ATPases at any ATP concentration from 0.1 to 1000  $\mu\text{M}$ , making it well suited for profiling inhibitors with enzymes that have diverse ATP requirements.
- The ability to run reactions in kinetic mode enables determination of inhibitor residence times using a jump dilution method.
- The residence times determined for Abl1, EGFR and Aurora C inhibitors using the Transcreener ADP Assay in jump dilution experiments were consistent with literature values.

# References and Acknowledgements

- Morrison JF, Walsh CT. The behavior and significance of slow-binding enzyme inhibitors. *Adv Enzymol Relat Areas Mol Biol.* 1988;61:201-301.
- Impact of enzyme concentration and residence time on apparent activity recovery in jump dilution analysis. R.A. Copeland et al. / *Anal. Biochem.* 416 (2011) 206-210.
- Biochemical characterization of GSK1070916, a potent and selective inhibitor of Aurora B and Aurora C kinases with an extremely long residence time. K. Anderson et al. *Biochem. J.* (2009) 420, 259–265
- Evaluation of enzyme inhibitors in drug discovery. Robert A. Copeland.
- Acknowledgements:
  - **Robert A. Copeland-Epizyme**
  - **Tim Wigle-Warp Drive Bio**
  - **BellBrook Labs team.**



# Transcreener® Platform

*Four assays, thousands of targets*

