

Determination of Drug Residence Time in a HTS Format using Kinetic Analysis with the Transcreener® ADP² Assay

Meera Kumar, Tom Zielinski, Roland Carrillo, and Robert G. Lowery
BellBrook Labs, Madison, WI, USA



Overview

Drug efficacy is influenced by the affinity of the drug to the target and longevity of the drug's interaction with the target, or residence time. In general, longer residence time results in improved efficacy, as the extended contact between drug and enzyme results in extended inhibition of target activity. This in turn allows pharmacological effect to be achieved over protracted time at lower doses, reducing off-target effects. Residence time can be determined by equilibrium binding measurements using methods such as surface plasmon resonance (SPR), which, while highly quantitative data, is low-throughput and requires costly instrumentation. Furthermore, SPR requires that protein targets or drugs to be covalently conjugated to surfaces, which can modify target function and/or drug binding. Here we describe, a "jump dilution" kinetic approach to determine the residence time for protein kinase inhibitor drugs using the Transcreener ADP² Assay, which relies on homogenous immunodetection of ADP. ADP displaces a tracer from a highly specific monoclonal antibody resulting in a change in fluorescence. Three homogenous mix-and-read formats that use a red-shifted tracer to minimize compound interference are available: fluorescence polarization (FP), time resolved FRET (TR-FRET), and fluorescence intensity (FI).

Transcreener ADP² Assays: Direct, Homogenous ADP Detection with FP, TR-FRET, or FI Readouts

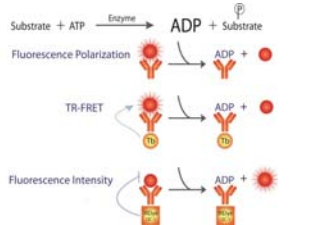


Figure 1. Transcreener® ADP² Assays: the Only Direct Detection Method Available for ADP. ADP displaces a tracer from a highly specific monoclonal antibody resulting in a change in fluorescence. Three homogenous mix-and-read formats that use a red-shifted tracer to minimize compound interference are available: fluorescence polarization (FP), time resolved FRET (TR-FRET), and fluorescence intensity (FI).

Jump Dilution Method using Transcreener ADP² Assay

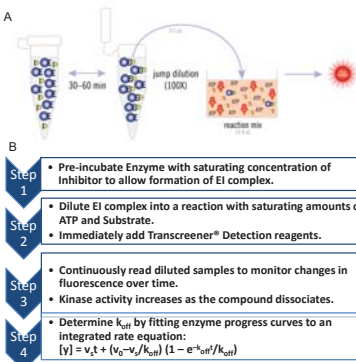


Figure 2. Using the Transcreener® ADP² Assay for "Jump Dilution". The reversibility of inhibition is determined by measuring the recovery of enzymatic activity after a rapid dilution of the EI complex into a reaction mix that contains ATP, acceptor substrate and Transcreener® reagents. This method, known as the "Jump Dilution" method, monitors the recovery of kinase activity over time. **A**, Shows the principle of Jump Dilution experiment and **B** shows a flowchart for running a Jump Dilution experiment using Transcreener assays.

Experimental Example: Determining the Dissociation Rate of Abl1 Inhibitors

Determining Optimal ABL1 and Drug Concentrations for the "Jump Dilution" Experiments

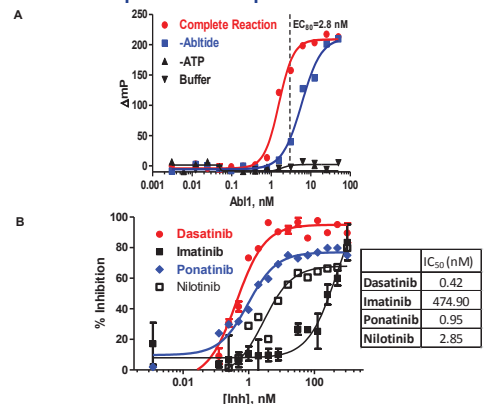


Figure 3. Determining the Optimal Concentration of Abl1 and Drugs for the "Jump Dilution" Experiment. A. Titration of Abl1 enzyme, performed in the presence of 5 μM ATP and 10 μM Abilite, determined an optimal EC₅₀ concentration as 2.8 nM. B. A dose response curve for Abl1 enzyme (2.8 nM) in the presence of 5 μM ATP and 10 μM Abilite. The reaction was run in a typical kinase buffer (50 mM Tris (pH 7.5), 5 mM MgCl₂, 0.01% Brij).

Determining Enzyme Velocities-Initial and Steady state velocities to create the Integrated Rate Equation.

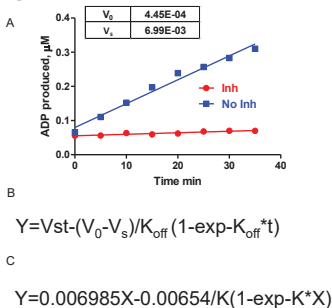


Figure 4. Determining Enzyme velocity V₀ and V_{ss}. A. V₀ was determined for fully inhibited reactions using Abl1 enzyme in the presence of 400 nM Dasatinib. B. Integrated rate equation used to determine the residence time values. C. Substituting the values of V₀ and V_{ss} to create an integrated equation

Jump Dilution Step: Recovery of Kinase Activity and Determining the K_{off} rates.

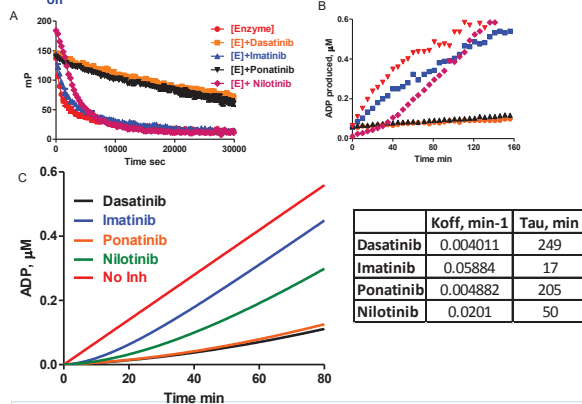


Figure 5. Determination of residence time for inhibitors. A. EI complexes were formed by pre-incubating with Abl1 (280 nM) with saturating inhibitor concentrations (10 x IC₅₀): 45 nM (Dasatinib), 45 μM (Imatinib), 25 μM (Nilotinib) or 100 nM (Ponatinib). EI mixtures were diluted 100-fold by adding 0.2 μL into 19.8 μL containing 10 μM Abilite and 5 μM ATP and Transcreener ADP² detection reagents; plates were read at 5 min intervals. B. Raw polarization was converted into product formed (ADP) using a standard curve. C. Progress curves were fit to the integrated rate equation shown in Figure 4C and K_{off} and Tau values were determined using Graph Pad Prism. Residence times (Tau values) are reciprocal of K_{off} values.

Dissociation Rate and residence times for EGFR inhibitors

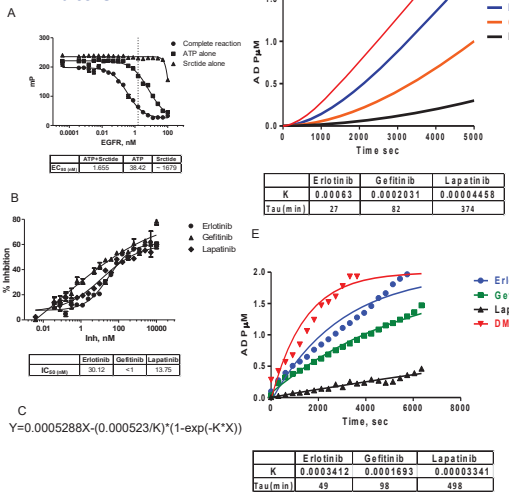


Figure 6. Residence Times for EGFR Inhibitors. A. An optimal (EC₅₀) concentration of 1.5 nM EGFR was determined in the presence of 5 μM ATP and 10 μM SrcTide. B. Dose response curve for EGFR inhibitors. C. An integrated rate equation for determining residence time for inhibitors. D. Progress curves after jump-dilution of EI complexes into reaction containing ATP, SrcTide and Transcreener ADP² detection reagents. Inhibitors were pre-incubated with EGFR at 10 x IC₅₀ to allow EI formation. E. An approximate determination of time occupied at the target can also be determined by fitting the converted data to one phase decay. This can be used to rank-order compounds residence times.

Dissociation Rate and residence time for GSK1070916 with Aurora C

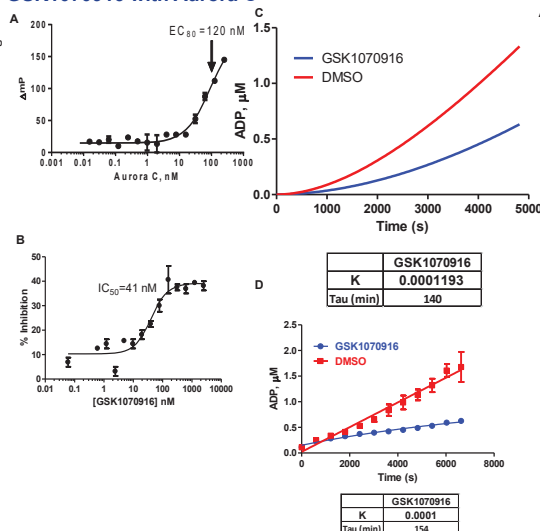


Figure 7. Residence Times for GSK1070916 with Aurora C. A. Aurora C was titrated in the presence of 20 μM ATP and 10 μM Kemptide to determine an optimal EC₅₀ concentration of 120 nM. B. A dose response curve for Aurora C enzyme yielded an IC₅₀ value of 42 nM. C. Residence time for GSK1070916 was determined for Aurora C in a "jump dilution" experiment using the Transcreener® ADP² Assay. Incubations were performed with Aurora C concentration of 1200 nM and inhibitor concentrations of 400 nM. A 40-fold jump dilution was made into 10 μM Kemptide, 20 μM ATP and Transcreener detection reagents. D. An approximate determination of time occupied at the target can also be determined by fitting the converted data to one phase decay.

Dissociation Rate and residence time with Aurora B

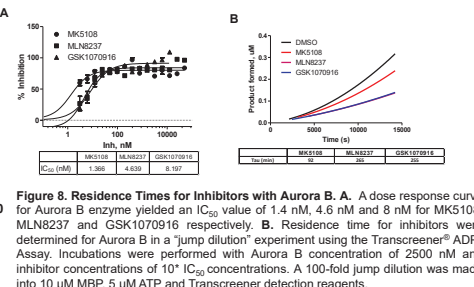


Figure 8. Residence Times for Inhibitors with Aurora B. A. A dose response curve for Aurora B enzyme yielded an IC₅₀ value of 1.4 nM, 4.6 nM and 8 nM for MK5108, MLN8237 and GSK1070916 respectively. B. Residence time for inhibitors were determined for Aurora B in a "jump dilution" experiment using the Transcreener® ADP² Assay. Incubations were performed with Aurora B concentration of 2500 nM and inhibitor concentrations of 10 x IC₅₀ concentrations. A 100-fold jump dilution was made into 10 μM MBP, 5 μM ATP and Transcreener detection reagents.

Conclusions

- The Transcreener® ADP² 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 assay can be tuned for detection of kinases or ATPases at any ATP concentration from 0.1 to 1000 μM, 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 ADP² Assay in jump dilution experiments were consistent with literature values.