

The significance of ATP concentration in cell-free and cell-based assays

ATP concentration greatly affects the determination of kinase inhibitory activity, given most kinase inhibitors bind to the ATP binding site of the kinase. Cell-free (biochemical) assays typically utilize ATP concentrations approximating the ATP K_m , however this is frequently far lower than the mM ATP concentrations found in cells.

In consideration of this difference, let's review the rationale of using ATP concentrations around the K_m in cell-free assays. K_m values approximate the affinity between ATP and the kinase, while K_i values describe the affinity between inhibitor and kinase. For ATP-competitive inhibitors, the dependencies between the half maximal inhibitory concentration (IC_{50}) and ATP concentrations are described by the Cheng-Prusoff equation ($IC_{50} = K_i + K_i/K_m \times [ATP]$). An inhibitor has different K_i values for every kinase, and every kinase has a different K_m for ATP. By representing the ATP concentration with its K_m value, the IC_{50} reflects $2 \times K_i$ value. Thus the IC_{50} value becomes a direct measure of affinity between the inhibitor and the investigated kinase¹⁾.

Table 1 is an example where the IC_{50} of Inhibitor Z is calculated using the Cheng-Prusoff equation. It hypothesizes that the ATP K_m values for Kinase A and Kinase B are $1\mu M$ and $10\mu M$, respectively, and

Enzyme	Kinase A	Kinase B
ATP K_m (μM)	1	10
Inhibitor Z K_i (μM)	0.1	0.2
<u>IC_{50} (μM) at each ATP conc. (μM)</u>		
1 (K_m for Kinase A)	0.2	0.2
10 (K_m for Kinase B)	1.1	0.4
1,000 (around the conc. in cells)	100	20

Table 1. IC_{50} calculation example ($IC_{50} = K_i + K_i/K_m \times [ATP]$)

the K_i values of Inhibitor Z for Kinase A and Kinase B are $0.1\mu M$ and $0.2\mu M$, respectively. The IC_{50} for Kinase A and Kinase B are calculated as $0.2\mu M$ and $0.4\mu M$ respectively when ATP concentration was assumed to be each K_m value. This result reflects $IC_{50} = 2 \times K_i$ value, thus establishing that the binding affinity of the inhibitor against various kinases can be

ranked on the basis of IC_{50} .

While this information is useful, prior to advancing a drug candidate, its cellular effects must be understood. Cellular ATP concentration is generally thought to be a few mM, and consistently higher than the ATP K_m of most kinases. Table 1 shows that the IC_{50} values for Kinase A and Kinase B are calculated to be $100\mu M$ and $20\mu M$ at $1mM$ ATP, based on the K_i values. This illustrates that at high ATP concentrations, the IC_{50} discrepancies are greater for kinases with lower K_m values compared to those with higher K_m values. Roberts M reported that the target occupancy of crizotinib at $1\mu M$ decreased drastically in a cell-based assay compared to the inhibition in a cell-free assay²⁾ (Fig.1).

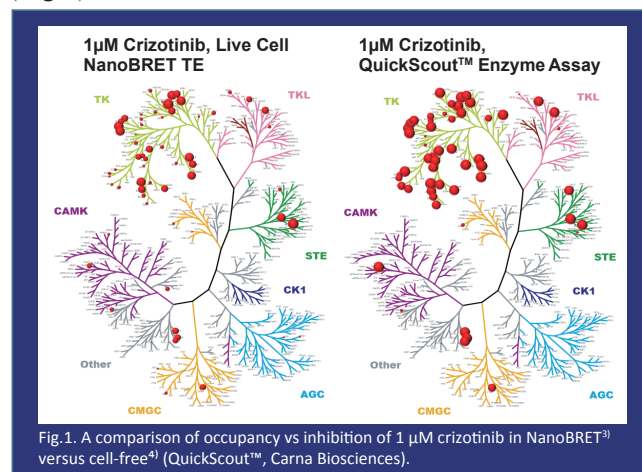


Fig.1. A comparison of occupancy vs inhibition of $1\mu M$ crizotinib in NanoBRET³⁾ versus cell-free⁴⁾ (QuickScout™, Carna Biosciences).

ATP concentrations in cell-free assays can be increased to more closely approximate cellular levels, however ATP concentration is not the only variable between these 2 assay systems. A multitude of factors affect compound binding affinity in cells. Given that cell membranes act as a barrier to binding, it is critical to confirm whether the inhibitor chosen by a cell-free assay maintains its inhibitory activity in a cell-based assay as well.

To address this need, Carna Biosciences offers a cell-based assay service known as the NanoBRET™ Target Engagement (TE) Intracellular Kinase Assay

System (Promega), in addition to cell-free recombinant kinase products and services (Profiling). In our cell-free assay service, we offer not only testing at ATP K_m values but also at 1mM ATP, while our cell-based assay services can examine compound affinity to the target kinase in a cell. The target kinase fused with NanoLuc[®] luciferase can be expressed in cultured cells and the binding of test compounds to the kinase can be detected by BRET (Bioluminescence Resonance Energy Transfer) (Fig.2).

Our cell-based assay makes it possible to evaluate the binding affinity of a compound to the target kinase inside intact cells, as it is performed without disrupting cell membranes. Additionally, it enables the determination of residence time, which is

especially important as a new pharmacodynamic parameter. Residence time data is emerging as a highly effective tool to differentiate new drug candidates from known drugs and other compounds which may have the same IC_{50} .

- Profiling Services
- NanoBRET[™] Target Engagement (TE) Intracellular Kinase Cell-Based Assay Services

References:

- 1) Protein Kinases as Drug Targets. 2011, ISBN: 978-3-527-31790-5, B. Klebl.
- 2) SLAS2020, Broad Kinome Selectivity and Residence Time Analysis in Live Cells with NanoBRET, Robers M.
- 3) Cell Chem Biol 2018; 25(2): 206-214.e11. Vasta JD.
- 4) PLoS One. 2014; 9(3): e92146. Uitdehaag JC.

