

## ➤ The Service in Brief

The biochemical ATP-competition assay service includes the biochemical determination of the effect of different ATP concentration on the  $IC_{50}$  of a compound against a given kinase. The resulting data give information about a possible ATP-competitive mode of action of the test compound, since ATP-competitive compounds show increasing  $IC_{50}$  values if tested at increasing ATP-concentrations.

## ➤ Background

The  $K_i$ -values of purely ATP competitive inhibitors can be derived from their  $IC_{50}$  values using the Cheng-Prusoff equation (Cheng Y, Prusoff WH (1973) *Biochem Pharmacol* 22: 3099-3108).

Non-ATP-competitive inhibitors (e.g. allosteric inhibitors) do not show this behaviour, i.e. their  $IC_{50}$  values do not change under different ATP-concentrations in a given assay setup. Mixed forms of inhibitors are also possible, which may show an intermediate dependency of the  $IC_{50}$  value on the ATP-concentration.

$$K_i = IC_{50} / (1 + \frac{[ATP]}{K_m})$$

### Figure 1: Cheng-Prusoff equation

$K_i$  = dissociation constant for binding of the inhibitor to the kinase, "true  $IC_{50}$ ";  $IC_{50}$  = measured  $IC_{50}$  value for the inhibitor at a distinct assay ATP-concentration;  $K_m$  = apparent  $ATP-K_m$  of the particular kinase in a given assay setup

## ➤ Example

$IC_{50}$  values of Staurosporine and the covalent MEK1 inhibitor Selumetinib were determined at different ATP-concentrations. The  $IC_{50}$  value of Staurosporine increased 14fold with increasing ATP-concentrations indicating an ATP-competitive mode of action. In contrast increasing ATP-concentration had no effect on the  $IC_{50}$  values of Selumetinib indicating that ATP does not compete with this compound.

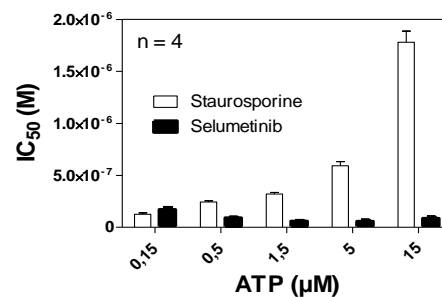


Figure 2: Dependency of the  $IC_{50}$  values of MEK1 inhibitor Selumetinib compared to Staurosporine

## ➤ Description of Service

The Service includes three steps:

- Determination of the linearity of the assay at the ATP-concentration that is 10fold higher than the apparent  $ATP-K_m$  of the kinase
- $IC_{50}$  determination of the compound at the apparent  $ATP-K_m$  of the kinase to determine the optimal compound concentration range for  $IC_{50}$  measurement.
- Determination of  $IC_{50}$  values at five different ATP-concentrations. Routinely, ATP-concentrations of 0.1/0.3/1/3/10fold of the  $ATP-K_m$  are used for quadruplicate  $IC_{50}$  determination, spanning a maximal ATP-concentration range of factor 100.

All assays are performed as a radiometric, filter-plate based protein kinase activity assay.  $IC$  values are determined using 10 different compound concentrations.

Based on the Cheng-Prusoff equation, the  $IC_{50}$  values of an ATP-competitive compound would theoretically change under these conditions as follows:

- 0.1 x  $ATP-K_m$  ➤  $IC_{50} = K_i \times 1.1$
- 0.3 x  $ATP-K_m$  ➤  $IC_{50} = K_i \times 1.3$
- 1 x  $ATP-K_m$  ➤  $IC_{50} = K_i \times 2$
- 3 x  $ATP-K_m$  ➤  $IC_{50} = K_i \times 4$
- 10 x  $ATP-K_m$  ➤  $IC_{50} = K_i \times 11$