

➤ The Target

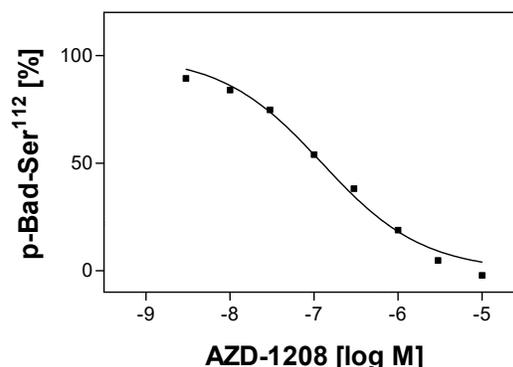
Pim kinases are constitutively active serine/threonine kinases that promote growth factor-independent proliferation by phosphorylation and subsequent inhibition of a range of cellular proteins. Activity of Pim kinases is physiologically regulated via protein stability. Especially Pim1 and Pim2 kinase have been found to be overexpressed or mutated in a variety of human tumors including B cell lymphomas, leukemias, prostate cancer and oral cancer.

➤ Cellular Phosphorylation Assay

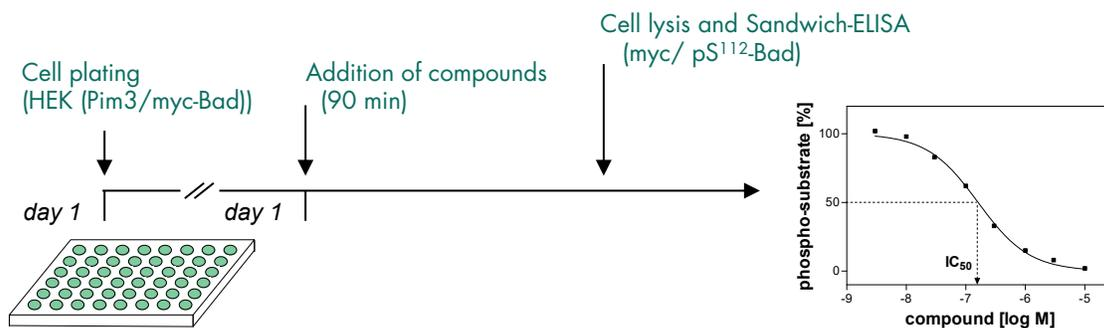
Transiently transfected Human embryonic kidney cells (HEK) express full length Pim3 kinase as well as Pim-substrate Bad labeled with a myc-tag. Constitutive activity of exogenous Pim3 results in phosphorylation of exogenous Bad at Ser¹¹². Upon compound incubation, cells are lysed and Bad phosphorylation is quantified by Sandwich-ELISA technique. The assay is validated based on the cognate inhibitor of Pim kinases AZD1208 (see Fig.1).

Figure 1: Assay validation.

Cognate Pim inhibitor AZD1208 blocks Pim3 and inhibits phosphorylation of Bad at Ser¹¹² with highly reproducible IC₅₀ values. The graph shows a representative result.



➤ You ship your compounds – Reaction Biology performs the testing



- IC₅₀ values are determined by testing 8 compound concentrations in semi-logarithmic steps (each concentration in duplicates).
- Quality assurance is provided by calculation of Z' factors for Low/High controls on each assay plate and by including a full IC₅₀ curve for a reference inhibitor to monitor adequate dose/response relation in your assay run.