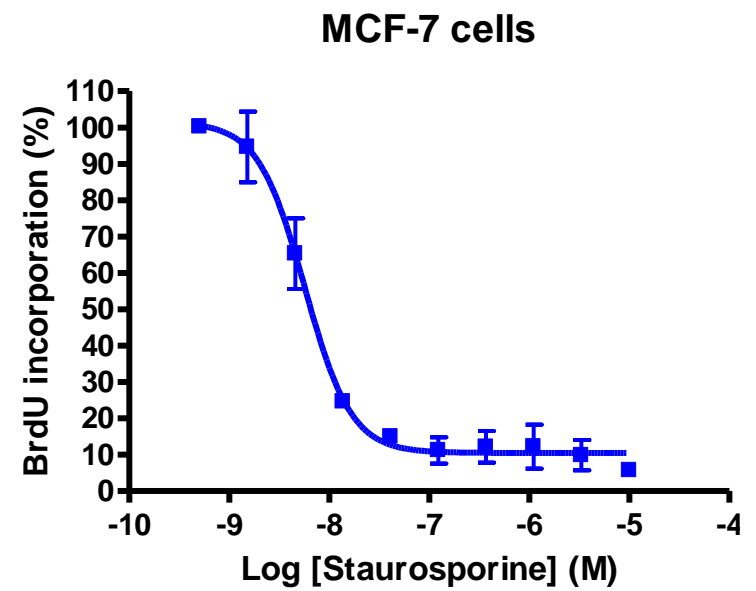


# BrdU Cell Proliferation ELISA Assay

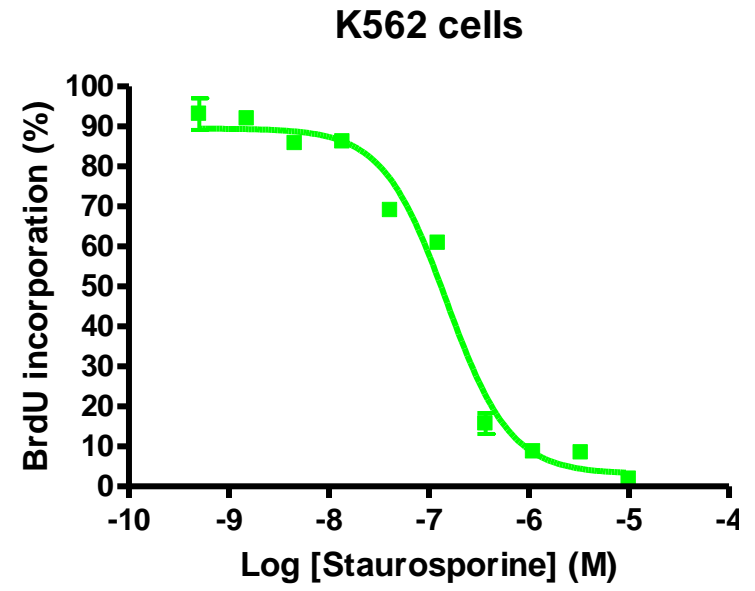
**Cell lines:** MCF-7 human breast cancer and K562 human chronic myelogenous leukemia (CML) cell lines from ATCC.

**Assay format:** BrdU cell proliferation ELISA assay.

**Assay Protocol:** MCF-7 and K562 cells were seeded into each well of a 96-well plate. The cells were treated with various concentrations of reference compound staurosporine for 72 hours. The BrdU incorporation into cellular DNA during cell proliferation was measured using BrdU ELISA assay. Absorbance signal was recorded by Envision 2104 Multilabel Reader. The IC50 (EC50) value was calculated and the IC50 curve was plotted using the GraphPad Prism 4 program based on a sigmoidal dose-response equation.



	Staurosporine
HILLSLOPE	-1.867
EC50	5.701e-009



	Staurosporine
HILLSLOPE	-1.362
EC50	1.504e-007