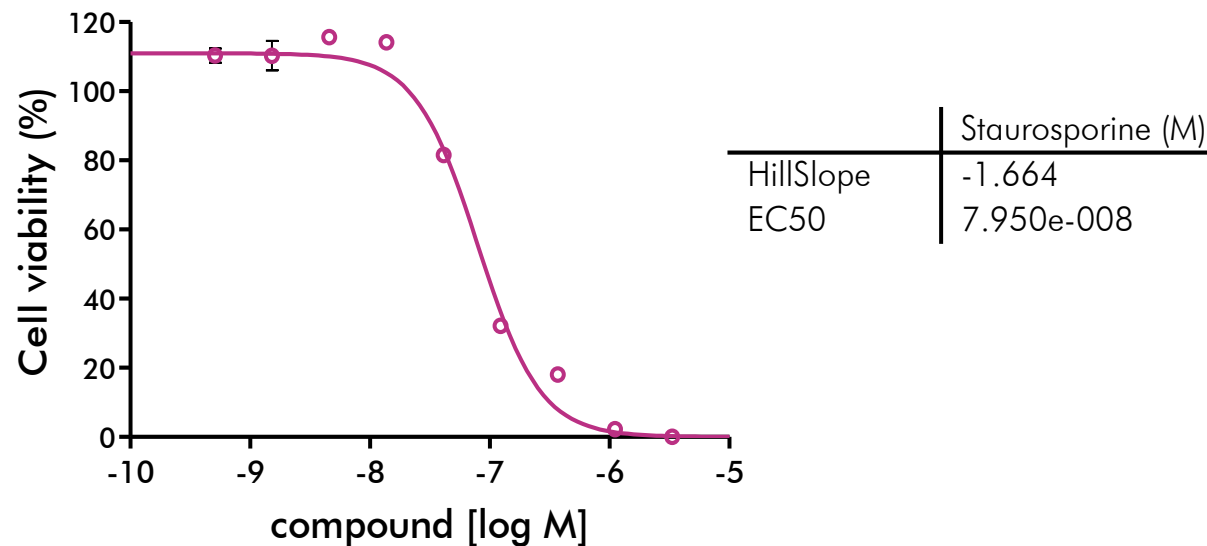


Cell Proliferation & Viability Assays — WST-8 Cell Proliferation Assay

Cell line: HepG2 human hepatocellular cancer cell line.

Assay format: 384-well format.

Assay Protocol: 5000 HepG2 cells were seeded into each well of a 384-well plate. Cells were treated with the indicated concentrations of reference compound staurosporine for 72 hours. CCK-8 Solution (product from Dojindo) was added to each well and incubated at 37 C for 4 hours. Absorbance of each sample was measured using Envision 2104 Multilabel Reader. The IC50 value was calculated and IC50 curve was plotted using the GraphPad Prism 4 program based on a sigmoidal dose-response equation.



© USA, June-2020. Material may not be reproduced or distributed without written permission from Reaction Biology Corp.