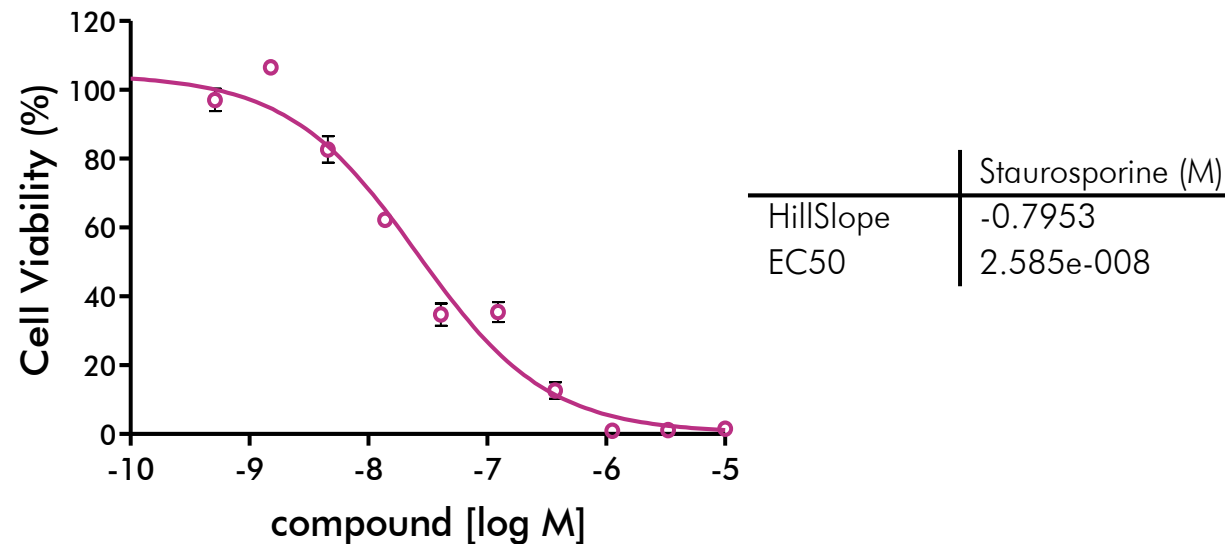


Cell Proliferation & Viability Assays — MTT Cell Proliferation Assay

Cell line: KMS-11 human multiple myeloma cell line.

Assay format: 384-well format.

Assay Protocol: 2000 KMS-11 cells were seeded into each well of a 384-well plate. Cells were treated with the indicated concentrations of reference compound staurosporine for 48 hours. MTT Dye Solution was added to each well and incubated for 6 hours at 37°C until formazan was produced in the cells. Solubilization/Stop Solution was added to each well to dissolve the formazan crystals. The 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.