

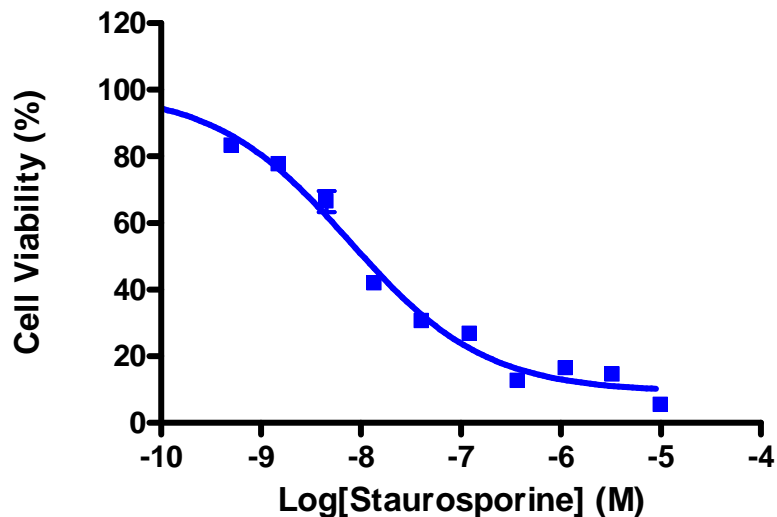
# CyQUANT Direct Cell Proliferation Assay

**Cell line:** HT-29 human colon cancer cell line, and Jurkat human T cell leukemia cell line from ATCC.

**Assay format:** 384-well format.

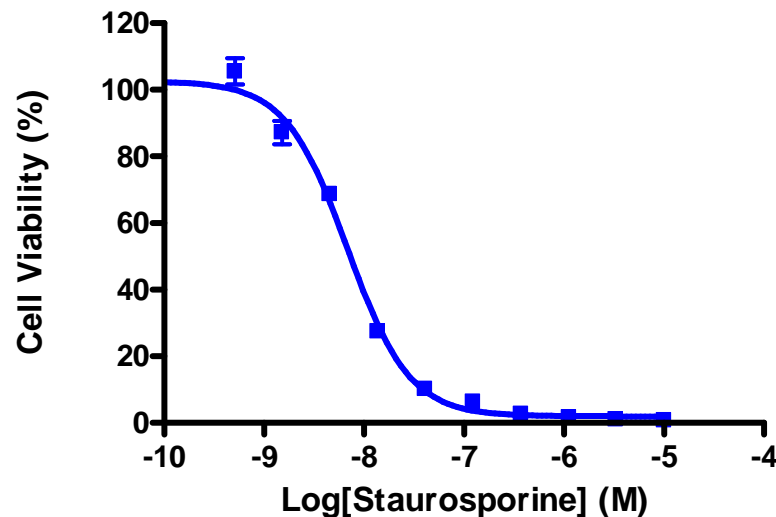
**Assay Protocol:** 2000 HT-29 cells or Jurkat cells were seeded into each well of a 384-well plate. Cells were treated with or without compound staurosporine for 48 hours. CyQUANT Assay Detection Reagent was added to each well and incubated for 60 minutes at 37°C. Fluorescence was measured by BioTek Synergy4 Microplate Reader. The IC50 (EC50) values were calculated and IC50 curves were plotted using the GraphPad Prism 4 program based on a sigmoidal dose-response equation.

**HT-29**



	Staurosporine (M)
HILLSLOPE	-0.6425
EC50	7.745e-009

**Jurkat**



	Staurosporine (M)
HILLSLOPE	-1.403
EC50	6.838e-009