

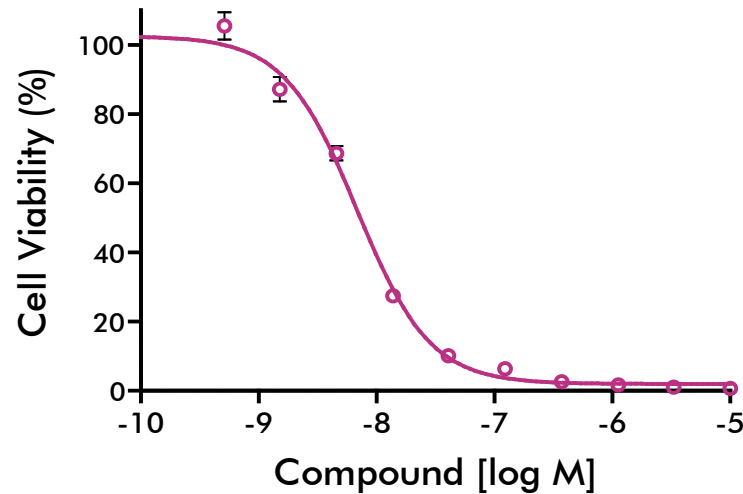
# Cell Proliferation & Viability Assays — CyQUANT Direct Cell Proliferation Assay

**Cell line:** Jurkat human T cell leukemia cell line

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

**Assay Protocol:** 2000 Jurkat 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. 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 value was calculated and IC50 curve was plotted using the GraphPad Prism 4 program based on a sigmoidal dose-response equation.



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