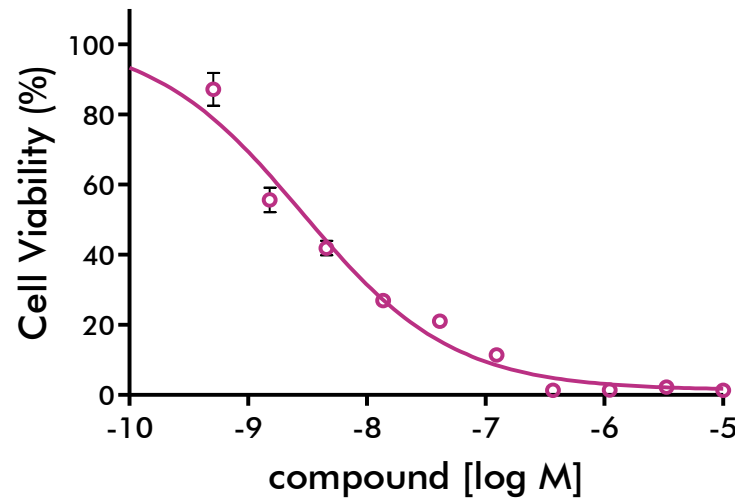


Cell Proliferation & Viability Assays — ATPlite Cell Proliferation Assay

Cell line: HeLa human cervical cancer cell line.

Assay format: 384-well format.

Assay Protocol: 2000 HeLa 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. ATPlite Assay Reagent was added to each well and incubated for 15 minutes. Luminescence was measured by 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.



	Staurosporine (M)
HillSlope	-0.6843
EC50	2.855e-009