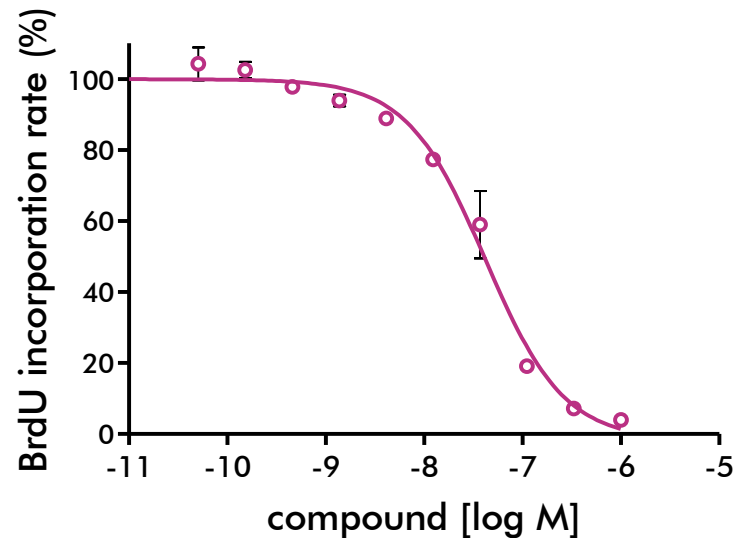


# Cell Proliferation & Viability Assays — BrdU Cell Proliferation ELISA Assay

**Cell lines:** B-474 breast cancer cell line.

**Assay format:** BrdU cell proliferation ELISA assay.

**Assay Protocol:** BT-474 cells were seeded into each well of a 96-well plate. The cells were treated with the indicated concentrations of reference compound staurosporine for 72 hours. The BrdU incorporation into cellular DNA during cell proliferation was measured using BrdU ELISA assay. Absorbance signal was recorded 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
HillSlope	-1.085
EC50	4.198e-008