

Cell Proliferation MTS Assay

HT-29 Cells: Human colorectal cancer cell line from ATCC.

Assay format: Cell proliferation MTS assay.

Assay Protocol: 5000 HT-29 cells were seeded into each well of the 384-well plate. Cells were treated with or without compound Saha or Staurosporine for 72 hours. 5 µl of CellTiter 96® AQueous One Solution Reagent was added to each well and incubated for 3 hours. The colored formazan product was used to determine the viable cells present in each well. Absorbance at 492 nm was recorded by Envision 2104 Multilabel Reader. The IC50 (EC50) values were calculated and the IC50 curves were plotted using the GraphPad Prism 4 program based on a sigmoidal dose-response equation.

