

Cell Viability Assay (Measurement of ATP Amount)

Cat #: Lum_ATP

Cells: Primary Cells or Cell lines

Description of Sample Data:

3T3 Cells were treated with Staurosporine for overnight in the presence of 0.1% Fetal Bovine Serum. Lysates of the 3T3 cells were analyzed for the amount of ATP by luminescence detection.

Reference Compound IC50:

Assay	Staurosporine, IC50 (nM)
Luminescence_ATP Assay	0.9

Reference Compound IC50 for Viability of 3T3 Cells

