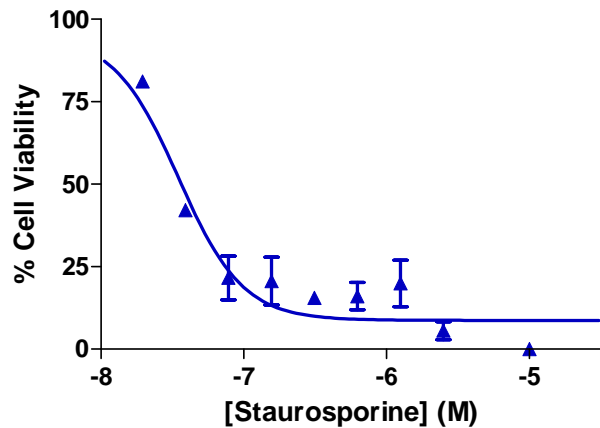


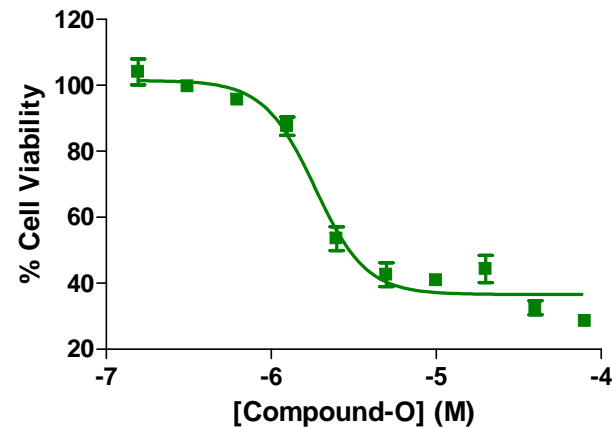
MTT Assay

HCT116 cells, a human colorectal carcinoma cell line, were treated with Staurosporine for 72hours and the cell viability was measured by MTT assay.



	Staurosporine
HILLSLOPE	-1.931
EC50	3.491e-008

HCT116 cells, a human colorectal carcinoma cell line, were treated with Compound-O for 72hours and the cell viability was measured by MTT assay.



	Compound-O
HILLSLOPE	-2.800
EC50	1.842e-006