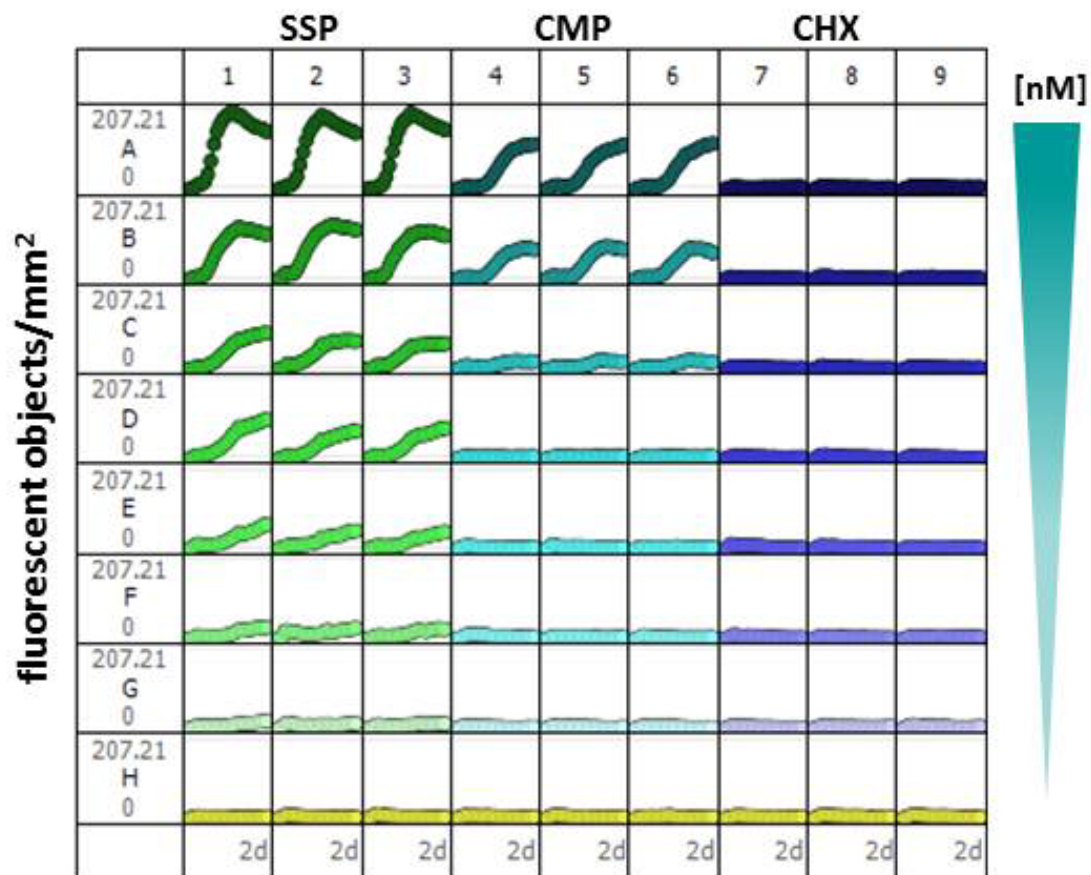


## HT 1080



96-well Cellplayer™ Cytotoxicity data. IncuCyte microplate graph showing kinetic measurement of cell death in HT1080 cells as determined by YOYO®-1 (Life Technologies) staining in response to increasing concentrations of Staurosporine (SSP), Camptothecin (CMP) and Cycloheximide (CHX).