



Blended fluorescent and phase-contrast IncuCyte images of HT1080 cells after 24 hour exposure to Staurosporine (SSP) , Camptothecin (CMP) and Cycloheximide (CHX). Green pixels denote DNA labeling of cells which have lost membrane integrity.