

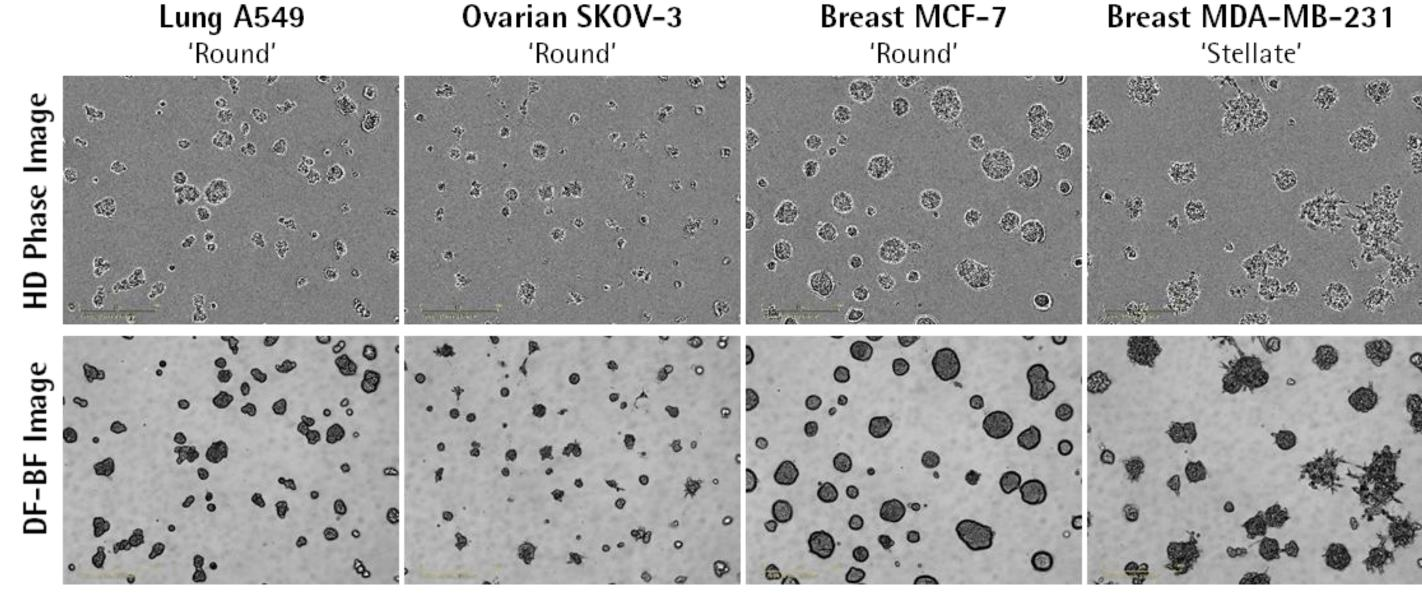
Development and optimization of matrigel-based multi-spheroid 3D tumor assays using real-time live-cell analysis

K. Patel, M. Oliver, G. Lovell, N. Holtz, T. Jackson, N. Dana, T. Dale, D. Trezise Essen BioScience, Welwyn Garden City, AL7 3AX UK or Ann Arbor, MI, 48108, USA

Summary & Impact

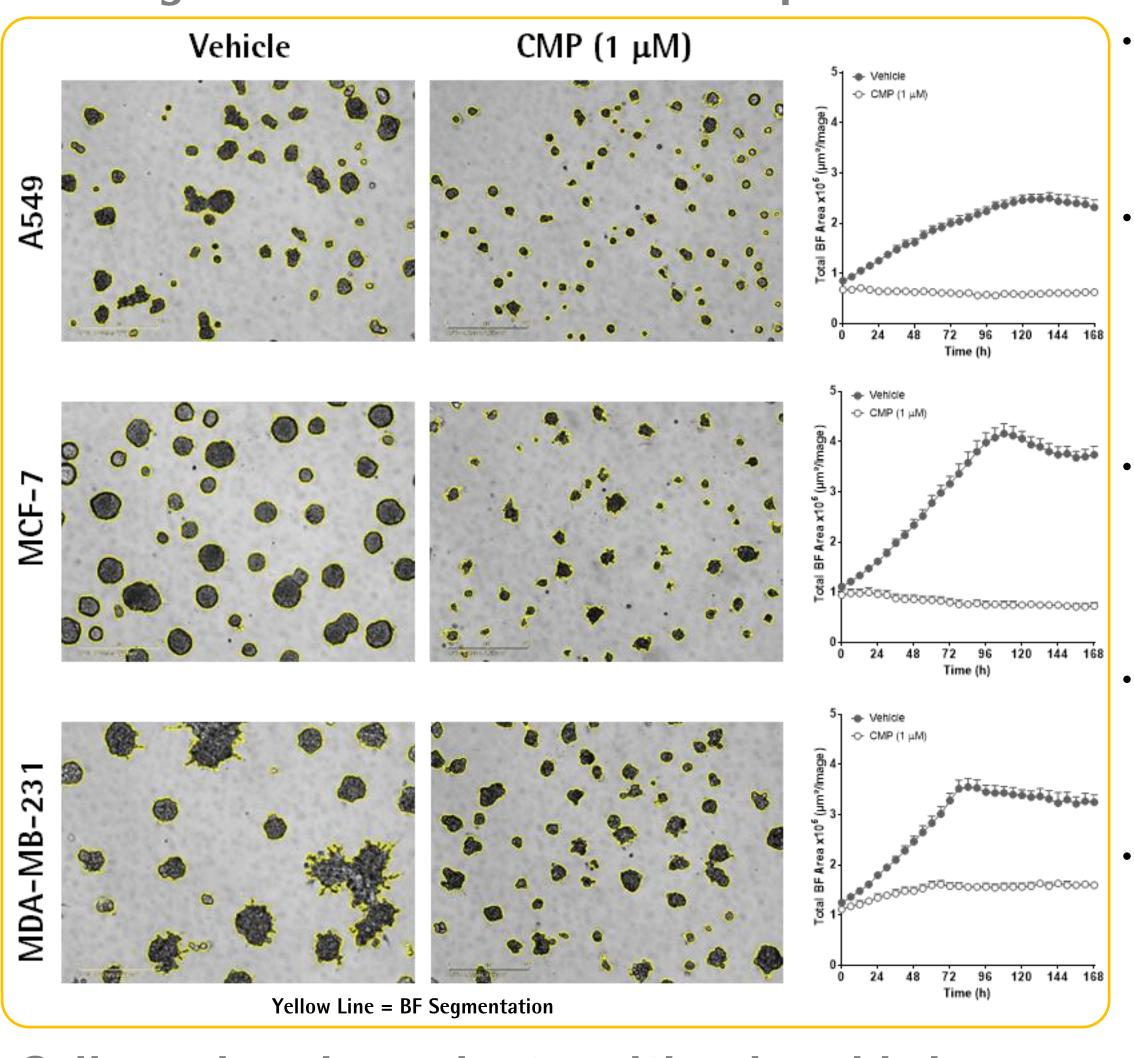
- The tumour-associated extracellular matrix (ECM) microenvironment provides critical biochemical cues as well as an essential structural scaffold for solid tumours to survive and grow.
- Here we describe a robust 3D ECM-based technique for culturing multiple tumour spheroids formed of lung, ovarian or breast cancer cell lines in a 96-well format.
- IncuCyte's DF Brightfield image acquisition tool enables the ability to monitor and quantify changes in spheroid size
- and morphology (brightfield) as well as viability (fluorescence) using real-time live-cell analysis.
- The use of IncuCyte® Cell Health reagents such as Annexin V to label apoptotic cells elucidates mechanism of action of compound treatments.
- Furthermore, this approach should facilitate more translational investigation of primary- and patient-derived organoid tumours.

Novel DF-Brightfield image capture



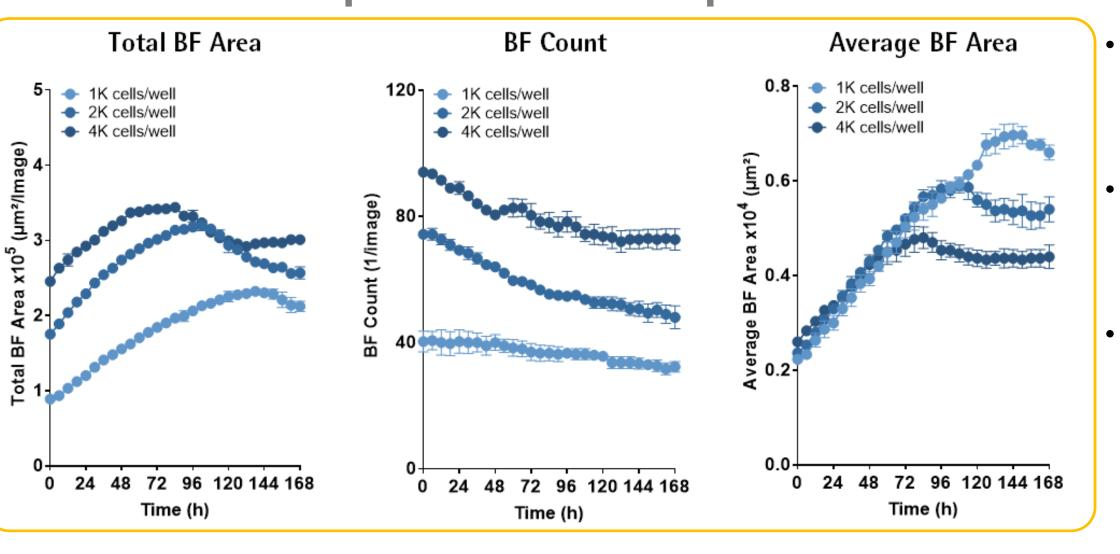
- High quality HD phase and DF-Brightfield (BF) images of multi-spheroids (MS) formed from a range of tumour cell lines (5 days post seeding) on a Matrigel® base.
- IncuCyte's proprietary image acquisition technique, DF-BF for 3D cultures, generates high contrast, extended depth of focus images.
- 3 days post seeding, A549, SKOV-3 and MCF-7 cells formed round aggregates, while MDA-MB-231 MS exhibited stellate branching distinctive of an invasive morphology.

DF-Brightfield enables label free quantification



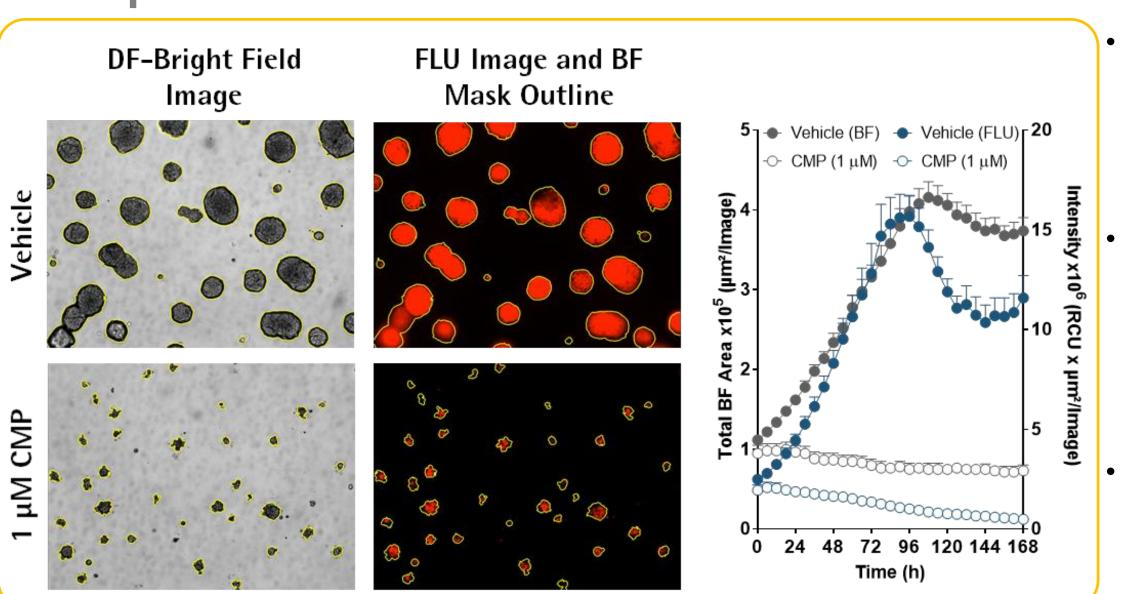
- A549, MCF7, and MDA-MB-231 cells were seeded (2K cells per well) and MS allowed to form for 3 days.
- MS were subsequently treated with the cytotoxic agent camptothecin (CMP, 1 μM) or vehicle control (0.1% DMSO) and images (DF-BF) collected every 6h for 7 days.
- IncuCyte multi-spheroid software enables kinetic quantification of growing and shrinking MS via size measurements (Total BF Area).
- A549, MCF-7 and MDA-MB-231 MS increased 1.9-, 2.0and 1.8-fold in size over 3 days, respectively.
- Treatment with CMP (1 μ M) inhibited growth of all MS.

Cell number dependent multi-spheroid size



- A549 MS were seeded at various densities (1K 4K cells per well).
- MS size and rate of growth, was proportional to the number of cells seeded.
- The decrease in number of MS over time reflects the merging of larger neighbouring MS.

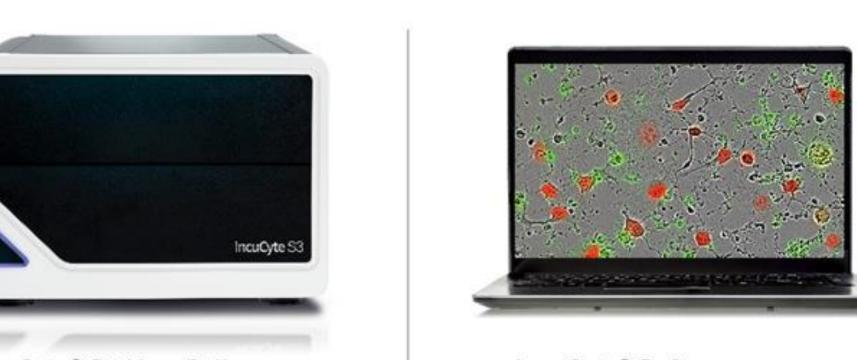
FP expression as an alternative measure for cell viability



- MCF-7 MS stably expressing nuclear restricted RFP (MCF-7-NucLight Red) were treated with CMP (1 μM) or vehicle (0.1% DMSO) for 7 days.
- IncuCyte analysis software reports both MS size (BF Area) and viability (Fluorescence intensity within BF Area) without the need to mask the fluorescent object
- fluorescent object.

 Fluorescence (RFP) intensity measurements provide a potential surrogate for MS health.

IncuCyte® System for Continuous live-cell analysis: Methodology



IncuCyte® S3 Live-Cell

Analysis System

A fully automated phase contrast and two-color fluorescence imager that

IncuCyte® Software

Fast, flexible and powerful control hub for continuous live-cell analysis comprising image acquisition,

New IncuCyte S3 Spheroid software module

processing and data visualization.

SSEN Service Annual Service State of Service S

IncuCyte® Reagents and Consumables

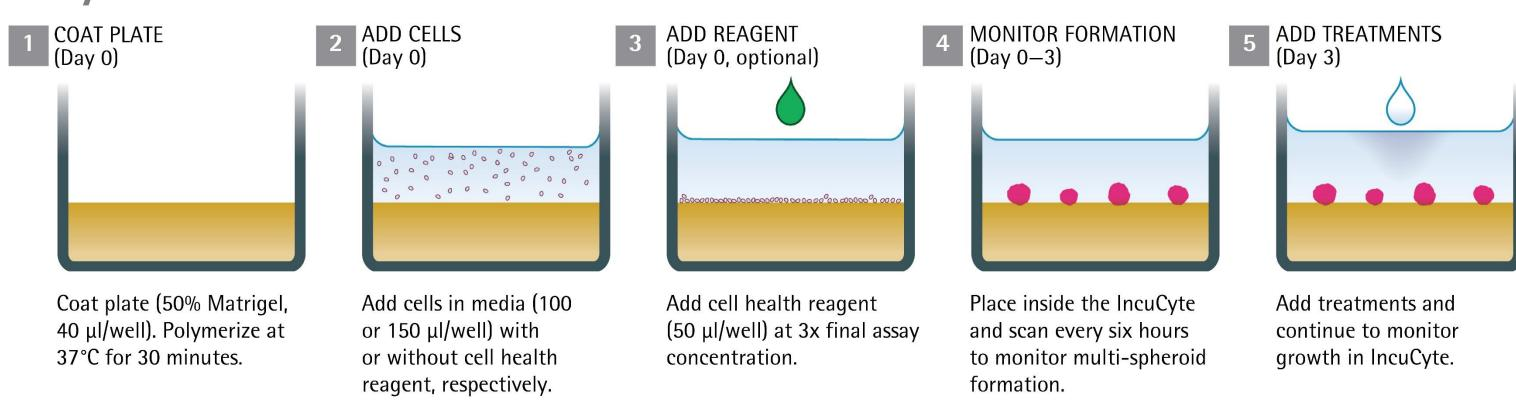
A suite of non-perturbing cell labeling and reporter reagents. Includes nuclear-targeted GFP and RFPs for cell counting plus no-wash cell health reagents for apoptosis and cytotoxicity.

Assay Workflow

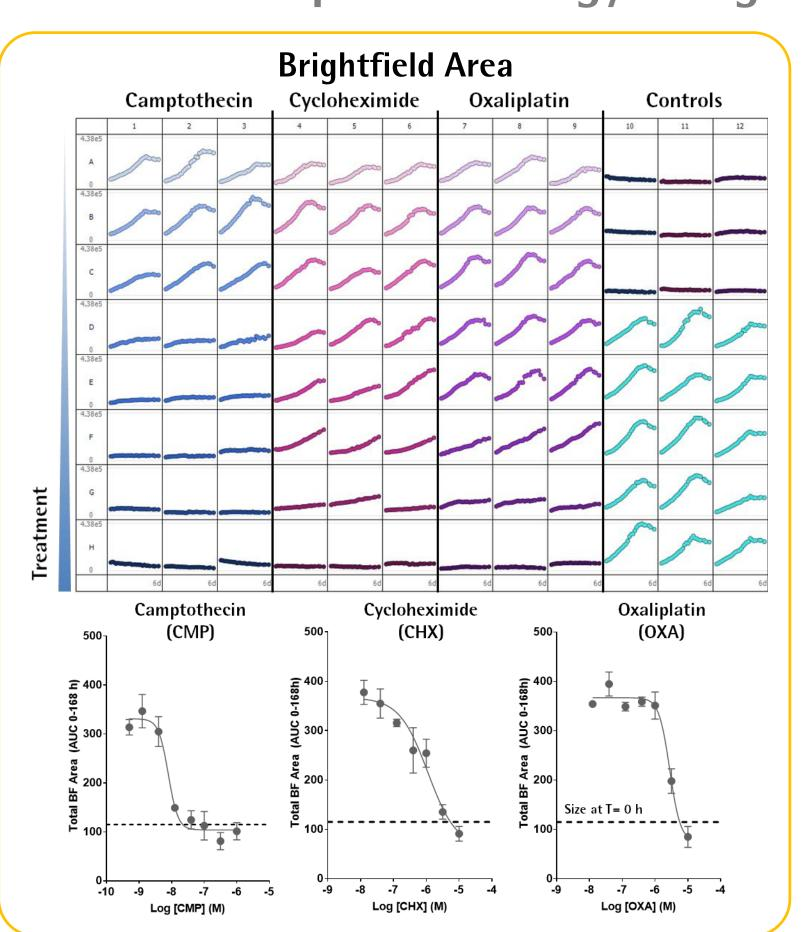
resides within a standard cell incubator

plates and flasks repeatedly over time.

for optimal cell viability. Designed to scan



Quantitative pharmacology using label free and fluorescent readouts



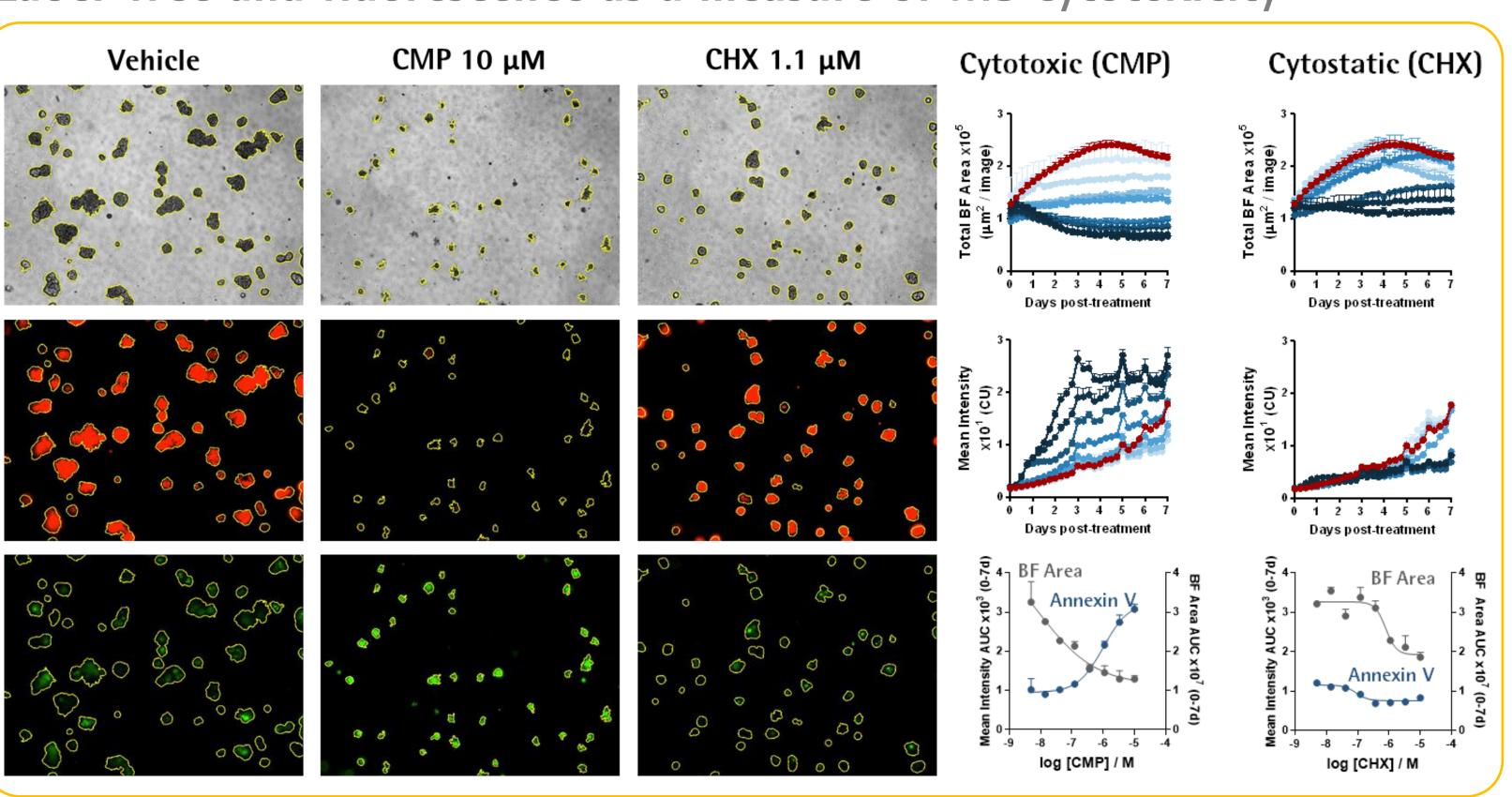
Tuesting of the control of the contr

Integrated Intensity

Controls

- MCF-7-NucLight Red MS were allowed to form for 3 days prior to treatment (7 days) with known cytotoxic compounds.
- Time-course plate-views enable rapid visualisation of treatment effects on both MS size (Total BF Area) and viability (FLU Intensity within BF Boundary).
- Concentration response curves represent area under curve (AUC) analysis of the time-course data.
- All compounds caused a concentration dependent inhibition of growth and viability with rank order of potency CMP > CHX ≥ OXA.

Label-free and fluorescence as a measure of MS cytotoxicity



- A549-NR cells (2K cells per well) were seeded in the presence of IncuCyte Annexin V Green reagent (1%) and MS allowed to form for 3 days.
- MS were treated with CMP, CHX or vehicle control and images (DF-BF, red and green fluorescence) acquired every 6h for 7 days.
- Both CMP (cytotoxic) and CHX (cytostatic) caused a concentration-dependent inhibition of MS growth (Total BF time courses).
- A loss of RFP signal, lack of growth and a simultaneous increase in Annexin V green fluorescence intensity (apoptosis) was observed in CMP treated MS.
- Despite CHX inhibiting MS growth, RFP expression remained high, whilst little or no increase in Annexin V fluorescence was observed, suggesting minimal cell death. These observations are consistent with the cytostatic properties of CHX.
- CRCs compare the cytotoxic vs cytostatic mechanisms of CMP and CHX respectively.