

Simplifying high throughput 3D spheroid growth and shrinkage assays using live content imaging

T. Dale¹, K. Patel¹, B. O'Clair², T. O'Callaghan¹, D. Appledorn², D. Trezise¹

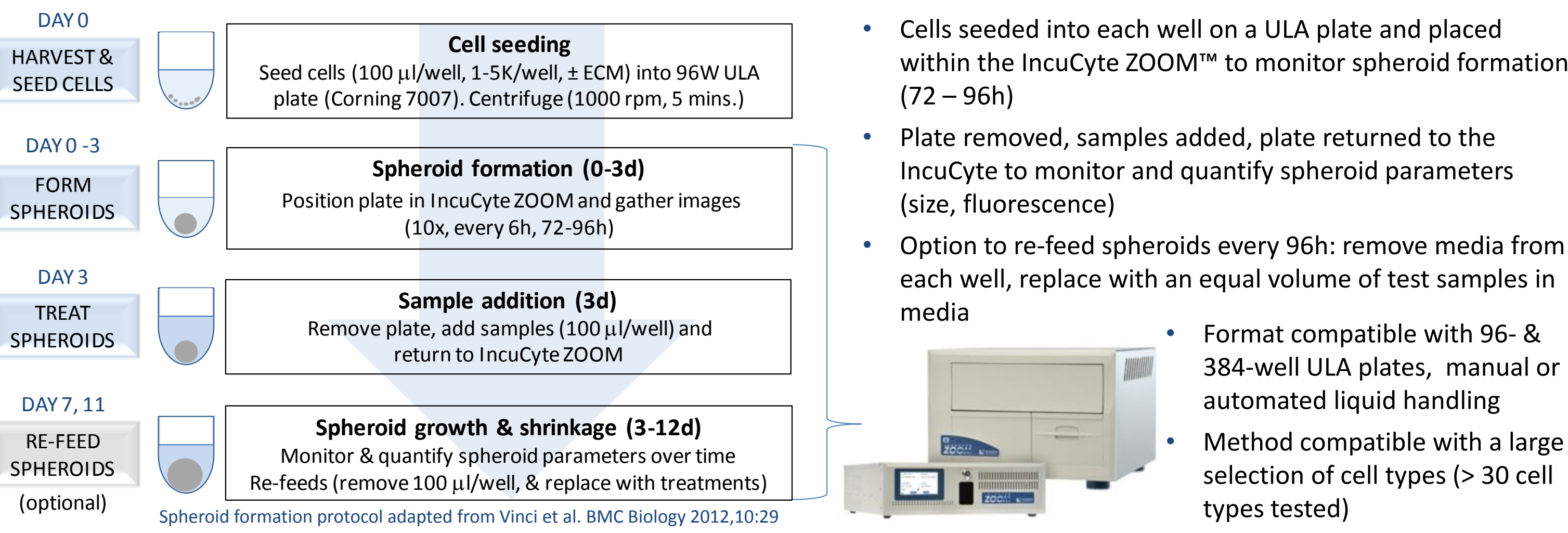
¹Essen BioScience Ltd, Welwyn Garden City, AL7 3AX UK & ²Essen BioScience Inc, Ann Arbor, Michigan, 48108 USA



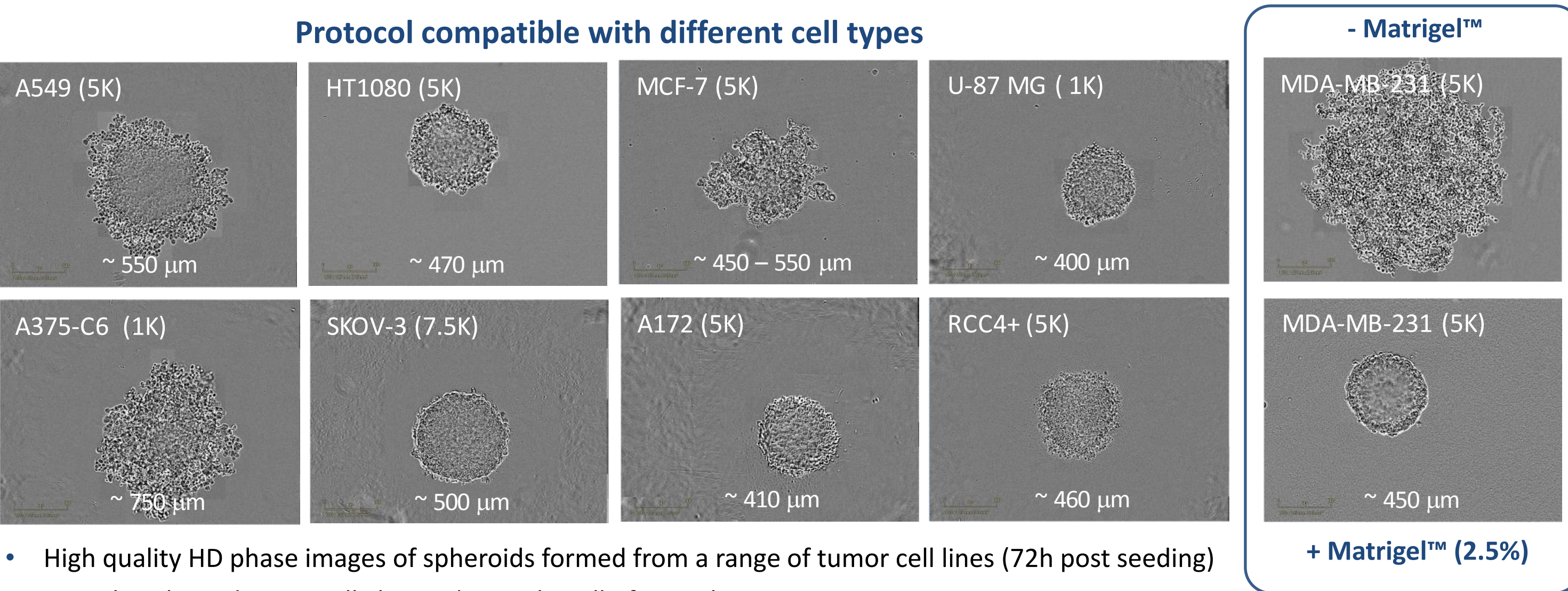
Summary & Impact

- Creating spheroids using Corning® Ultra Low Attachment plates provides a technically simple, robust and economical approach to creating 3D spheroids in 96-well plates (equivalent to 2D).
- Using fluorescently labelled cells and IncuCyte ZOOM™, we have assembled and validated fully-kinetic, spheroid growth and shrinkage assays.
- IncuCyte ZOOM™ metrics (spheroid size, (area/confluence) and fluorescence intensity (Mean Image Fluorescence, MIF)) are informative, and can be readily gathered to monitor spheroid growth and shrinkage over time.
- Spheroid fluorescence intensity can be measured without masking, and may be a useful indicator of spheroid 3D nature and/or cell health.
- The images and time-lapse movies provide important information and validation regarding spheroid morphology and the effects of test samples.
- These assays should prove useful for medium throughput, quantitative pharmacology of test samples for effects on the growth and shrinkage of tumor cells in 3D.
- Cell proliferation/ spheroid growth can be directly compared in 2D and 3D.

96-well spheroid assay workflow – an integrated solution

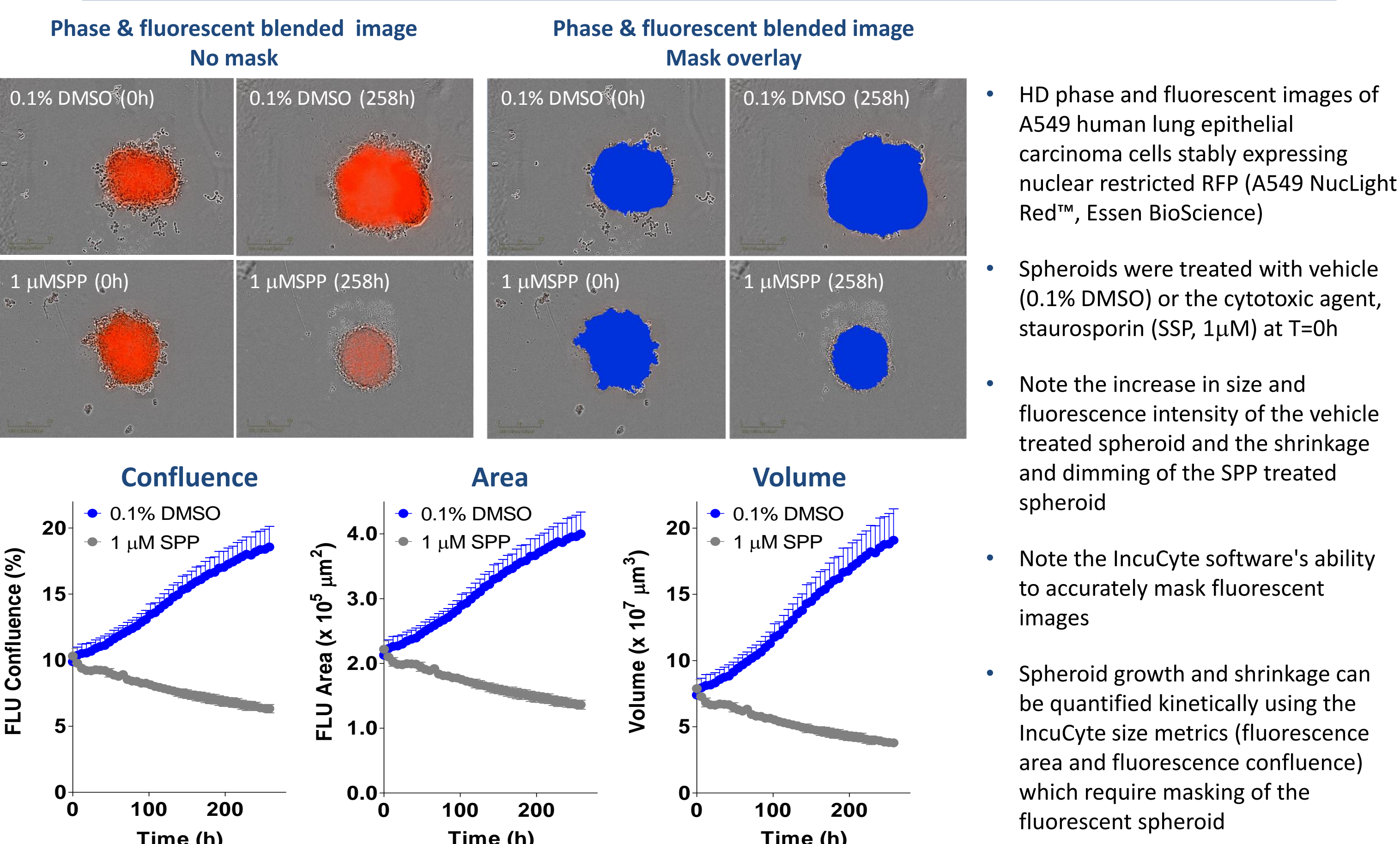


Successful spheroid formation using multiple cell types

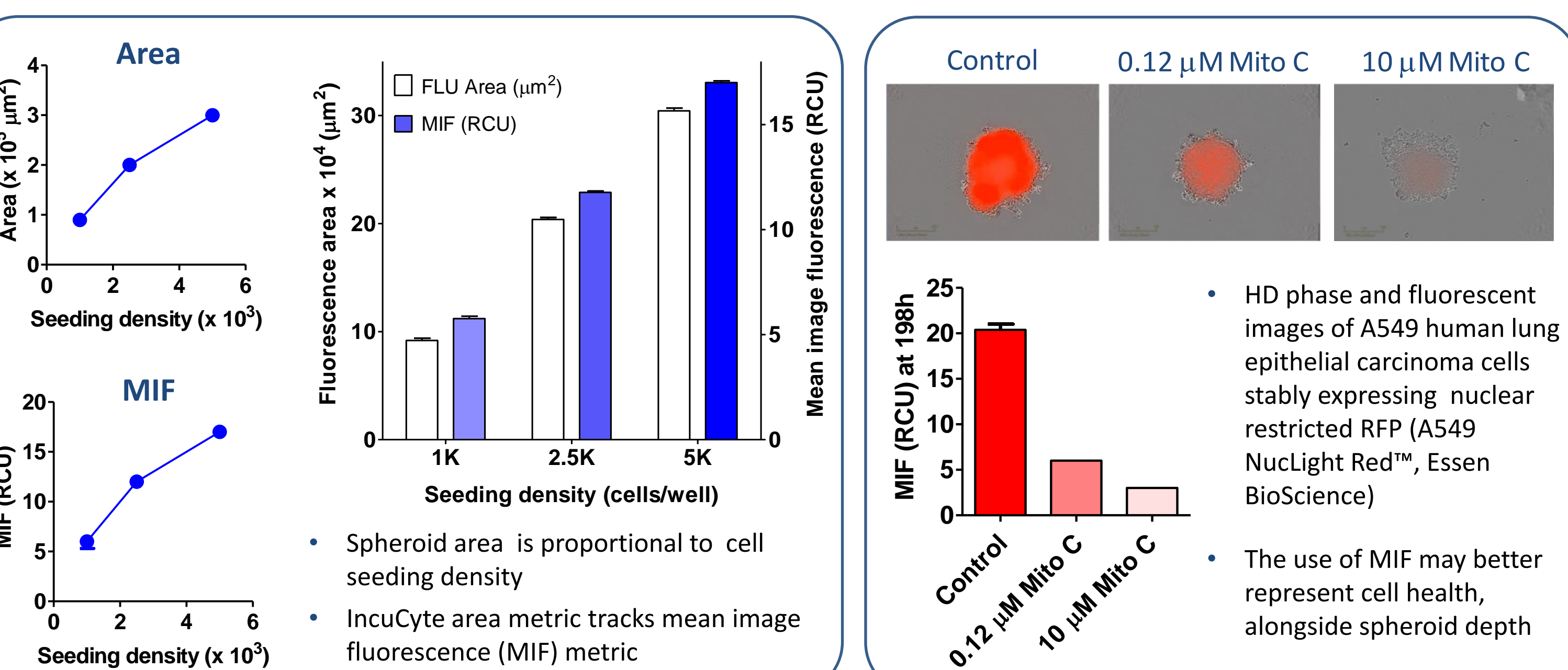


- High quality HD phase images of spheroids formed from a range of tumor cell lines (72h post seeding)
- A single spheroid is centrally located in each well of ULA plate
- Different 3D morphologies are observed (tight spheroids, compact and loose aggregates)
- Note, the 3D structure of loose aggregates can be improved by the inclusion of a basement membrane extract (e.g. 2.5% Matrigel™) at the stage of spheroid initiation

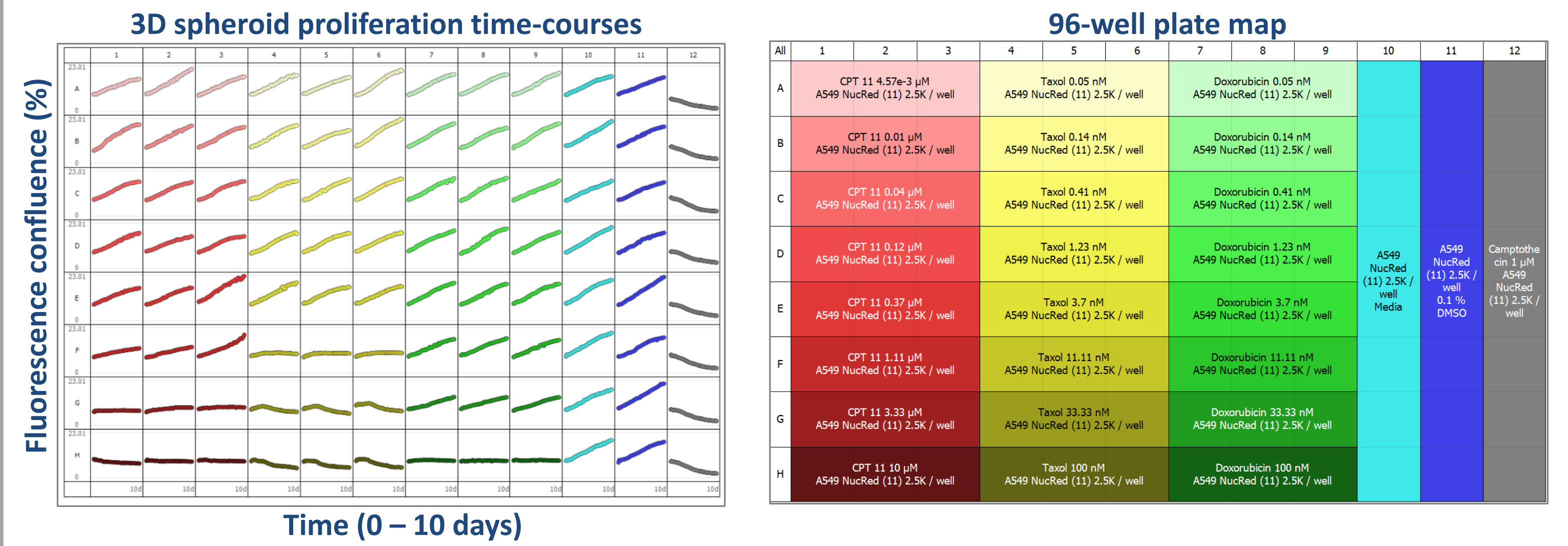
Quantifying spheroid growth and shrinkage over time



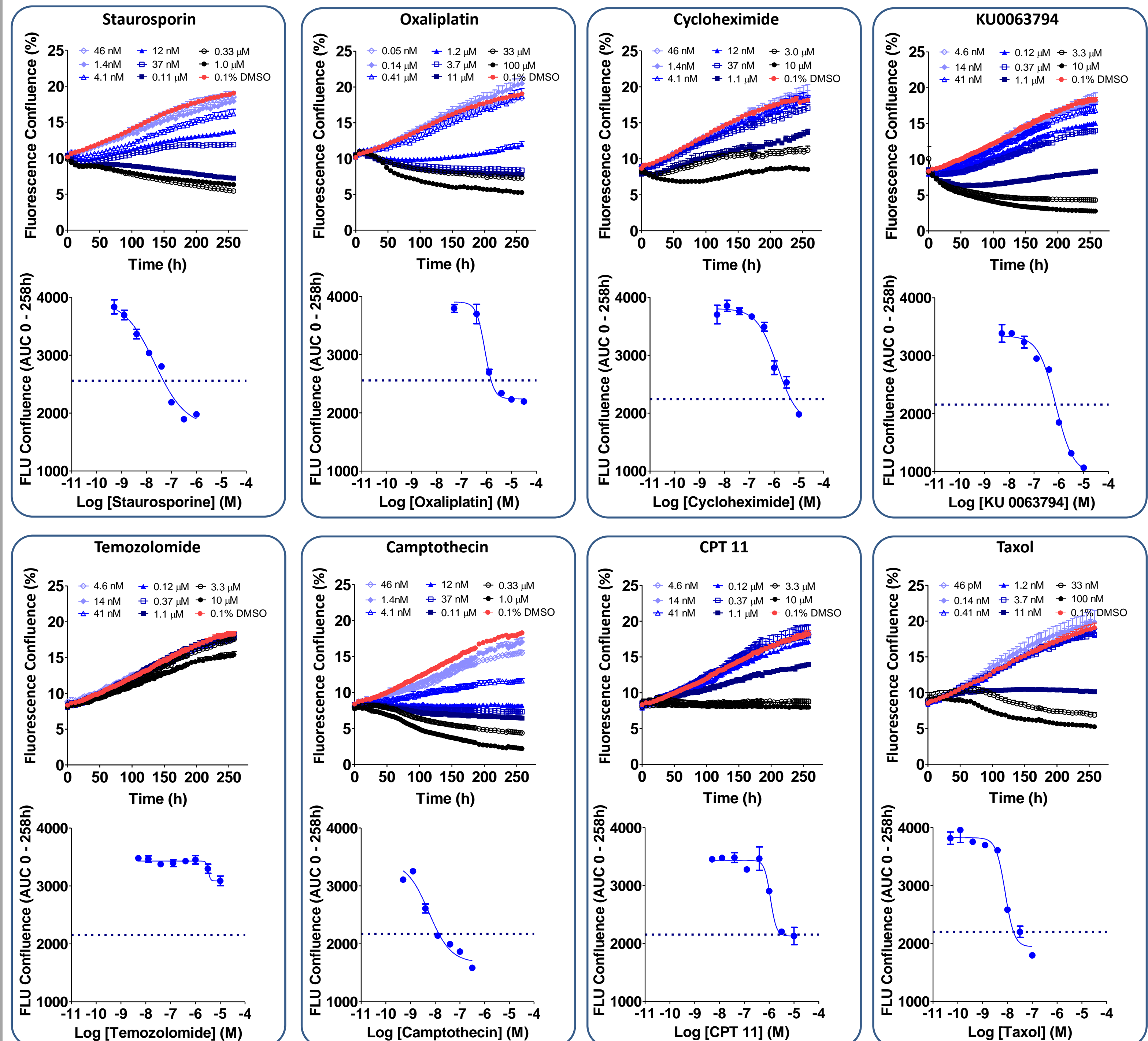
Validation of size metric



3D Pharmacology



- Plate views represent a facile visualisation the kinetic temporal data for spheroid growth and shrinkage (fluorescence confluence %) collected from a single 96-well microtitre plate



- A panel of compounds with different mechanisms of action were profiled in a 3D spheroid culture of human A549 cells stably expressing a red fluorescent label (A549 NuLight Red™)
- Time-course plots with the corresponding concentration response curves derived from the calculated area under the curve (AUC)
- The dotted line on the CRC plots represents the FLU confluence at time = 0 h, thus compounds yielding inhibition below this level (e.g. KU063794) are likely to be inducing spheroid shrinkage and cell death

Preliminary colony-forming assays in soft agar

