

96-Well Live-Cell Assays for Immune Cell Killing of 3D Tumor Spheroids

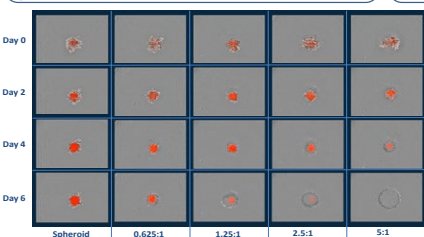
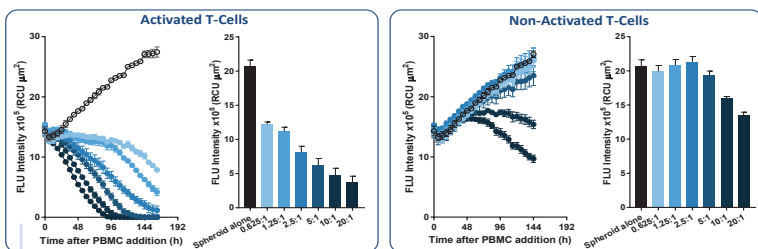
M. Oliver¹, K. Patel¹, N. Holtz², E. Endsley², T. Dale¹, and D. Trezise¹

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

Summary and Impact

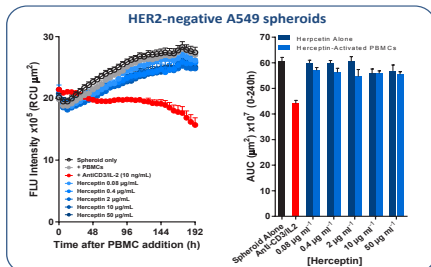
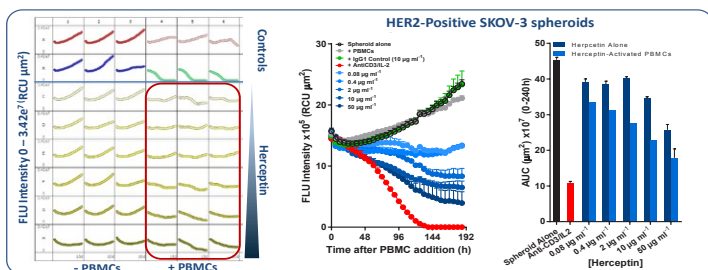
- Immunotherapies such as checkpoint inhibitors, CAR-Ts and immune-targeting Abs have great promise for cancer treatment. Translational cell-based assays are required to optimise these approaches.
- Here we describe image-based, immune cell-killing assays of 3D tumour spheroids, geared for assessing the efficacy of novel immune-modulators.
- Human tumour cell lines expressing RFP were used to form spheroids in 96-well ULA plates. Immune cells were then added and activated to kill. Spheroid viability was assessed over time (up to 10 days) by measuring the loss of RFP fluorescence using IncuCyte live-cell analysis.
- This method is exemplified with a range of immune cell types (PBMCs, T-cells, NK-cells) and activators, including anti-CD3 & IL-2
- In an ADCC format, Herceptin induced a concentration-dependent specific killing of Her-2 expressing tumours. Higher concentrations of Herceptin were required in 3D vs 2D ADCC assays.
- These data demonstrate how immune-cell killing and ADCC assays can be extended from traditional 2D mono-cultures to 3D spheroid assays, providing the potential for greater translational relevance. These assays will be highly valuable in the search for novel immune-modulators.

Effector-to-Target Ratio Dependent Cytotoxicity

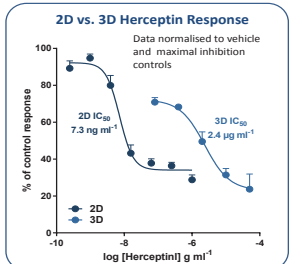


- Blended phase and fluorescent images of A549 NuLight Red™ spheroids in the presence and absence of immune cells.
- A549 cells (2.5K/well) seeded with PBMCs activated with anti-CD3 (10 ng/ml) and IL-2 (10 ng/ml). Cytotoxicity was quantified based on the red fluorescent intensity.
- Data demonstrates an E:T ratio-dependent destruction of tumour spheroids by the activated T-cell population.
- Note that E:T ratio optimisation is required as non targeted cell death was observed at E:T ratios >5:1.

Herceptin Induced ADCC in HER2-Positive SKOV-3 Cells



- A similar assay was conducted in a 2D culture model. SKOV-3 cells (1.6K/well) were seeded overnight prior to the addition of PBMCs (8K/well) and subsequent treatment with Herceptin.
- SKOV-3 tumour spheroids appear to exhibit ~300-fold lower Herceptin sensitivity in comparison to 2D.
- Note the apparent 34% inhibition of the 3D spheroid at the lowest test concentration (0.08 µg ml⁻¹). This suggests that a biphasic concentration response curve may exist, where the outermost cells behave as in the 2D model, whereas the spheroid centre has lower sensitivity.
- Additional experimentation is required to further understand the differential effects of Herceptin in 2D vs 3D models.



Continuous Live-Cell Analysis: Methodology



IncuCyte® S3 Live-Cell Analysis System

A flexible assay platform that sits inside a standard tissue culture incubator. IncuCyte automatically and continuously acquires and analyzes HD phase and fluorescent images of living cells cultured in microplates, dishes, or flasks.



IncuCyte® Software

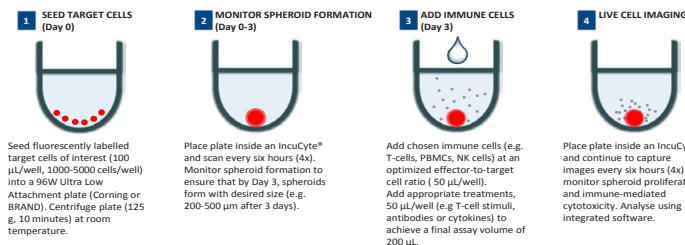
Fast, flexible, and powerful control hub for continuous live-cell analysis comprising image acquisition, processing, and date visualization.



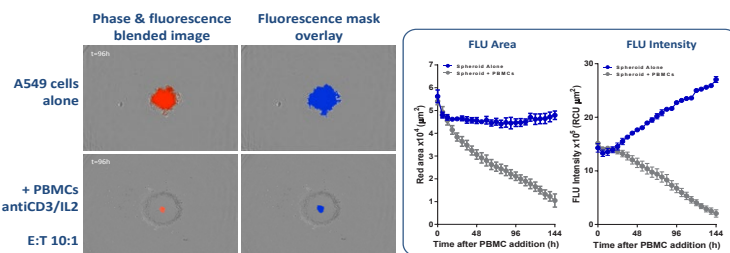
IncuCyte® Reagents & Consumables

A suite of non-perturbing cell labeling and reporter reagents. Includes nuclear-targeted GFP and RFPs for cell counting, no-wash caspase 3/7 substrate for apoptosis, and cell kits for angiogenesis.

96-Well 3D Cell Killing Assay Workflow

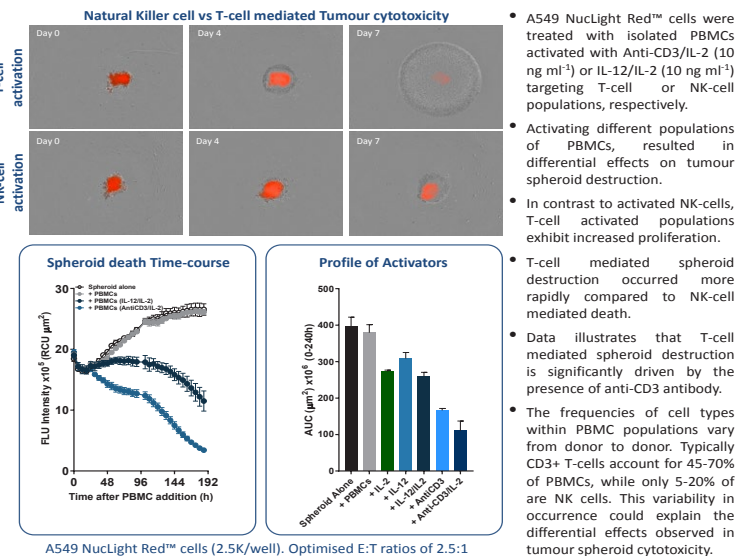


Fluorescence as a Measure of Spheroid Cytotoxicity



- Blended phase and fluorescent images, with corresponding masks, of A549 human lung epithelial carcinoma cells stably expressing RFP (A549 NuLight Red™, Essen BioScience).
- Note the increase in fluorescence intensity of the spheroid alone and the decline of fluorescence in the presence of immune cells.
- Spheroid proliferation and immune cell-mediated cytotoxicity can be quantified kinetically using the IncuCyte® size metrics (fluorescence intensity and fluorescence area) which require masking of the fluorescent spheroid.

Activator-Dependent Tumor Cytotoxicity



A549 NuLight Red™ cells (2.5K/well). Optimised E:T ratios of 2.5:1