

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

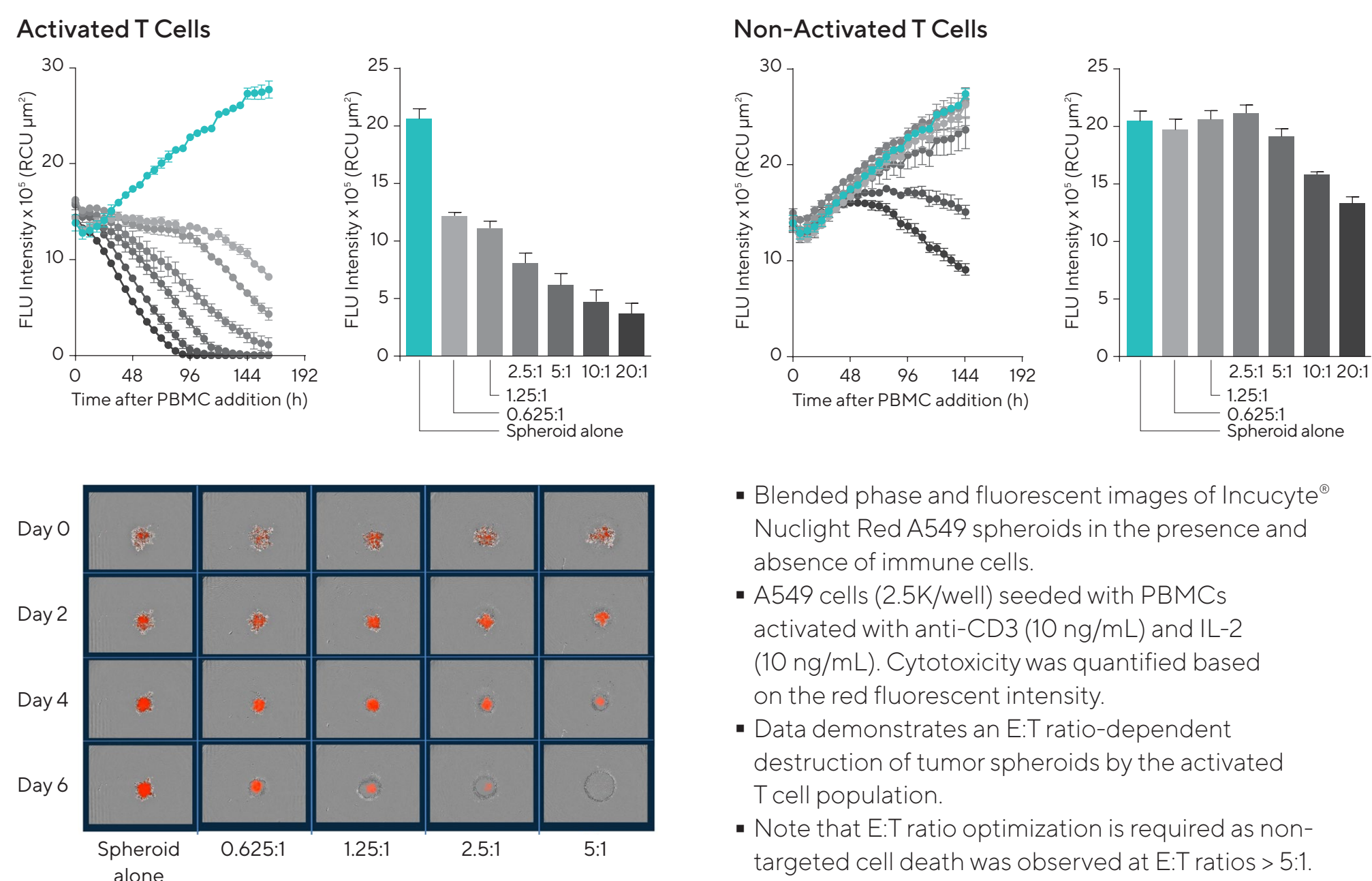
M. Oliver¹, K. Patel¹, N. Holtz², E. Endsley², T. Dale¹ and D. Trezise¹

1. Essen BioScience Ltd., a Sartorius Company, Units 2 & 3 The Quadrant, Newark Close, Royston Hertfordshire SG8 5HL UK
2. Essen BioScience Inc., a Sartorius Company, 5700 Pasadena Avenue NE, Albuquerque, NM 87113 USA

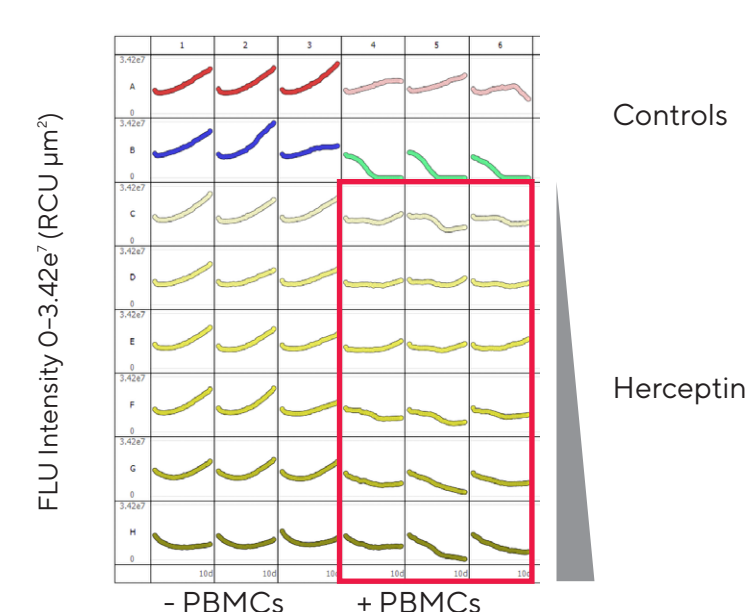
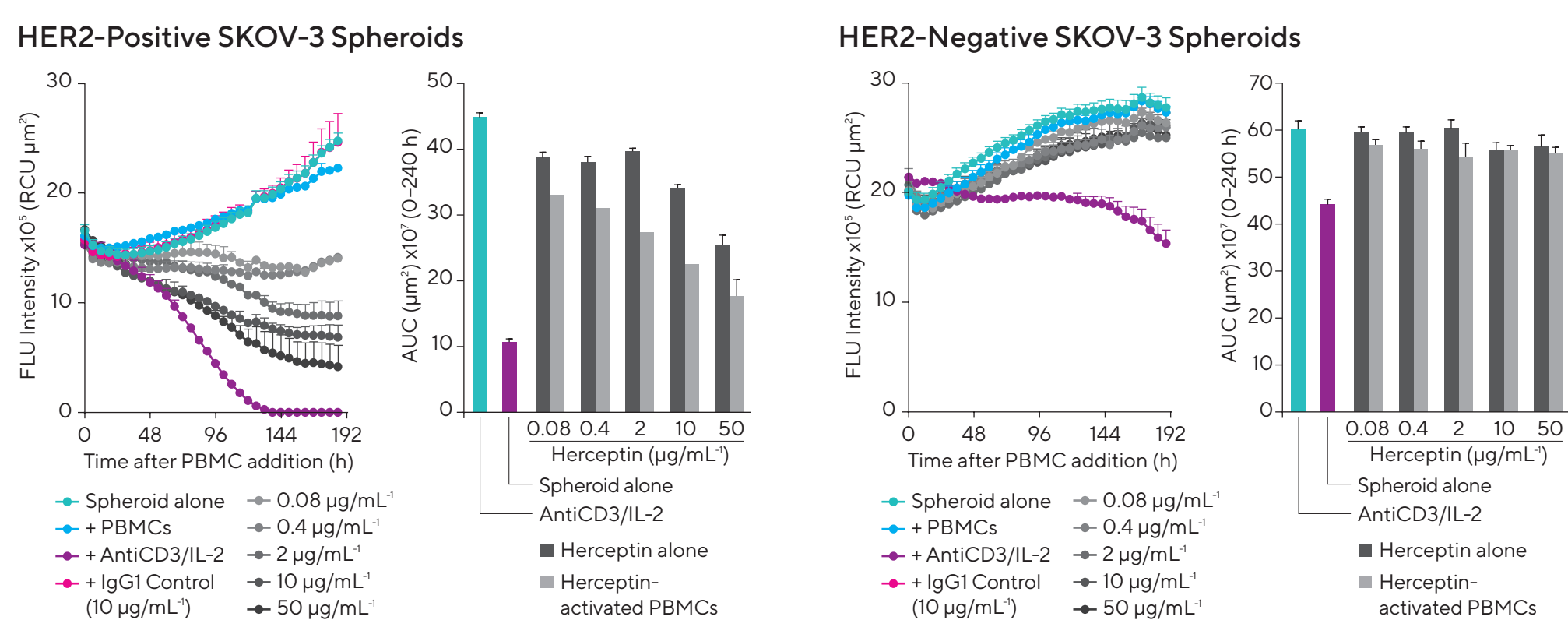
Summary and Impact

Immunotherapies such as checkpoint inhibitors, CAR-Ts and immune-targeting Abs have great promise for cancer treatment. Translational cell-based assays are required to optimize these approaches. Here we describe image-based, immune cell-killing assays of 3D tumor spheroids, geared for assessing the efficacy of novel immune-modulators. Human tumor 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 Imaging and Analysis. This method is exemplified using a range of immune cell types (PBMCs, T cells, NK-cells) and activators, including anti-CD3 and IL-2. In an ADCC format, Herceptin induced a concentration-dependent specific killing of HER2 expressing tumors. 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



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 tumor 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 center 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® 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.

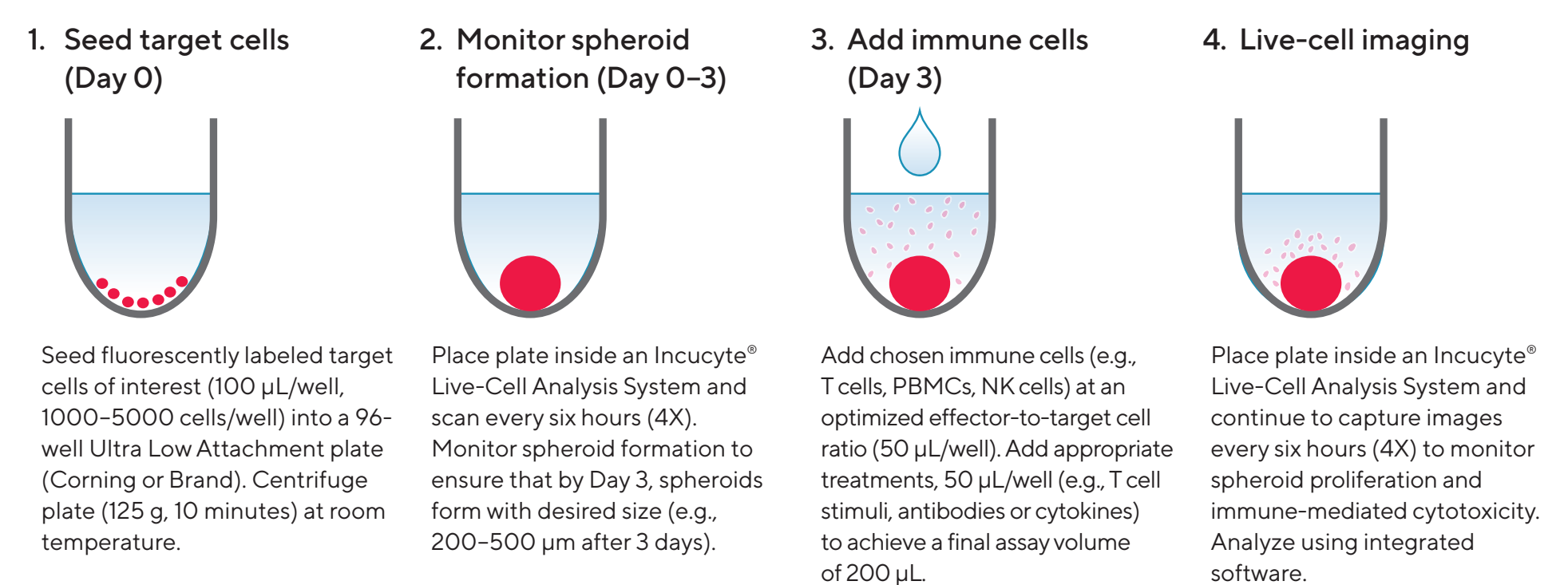
Incucyte® Software

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

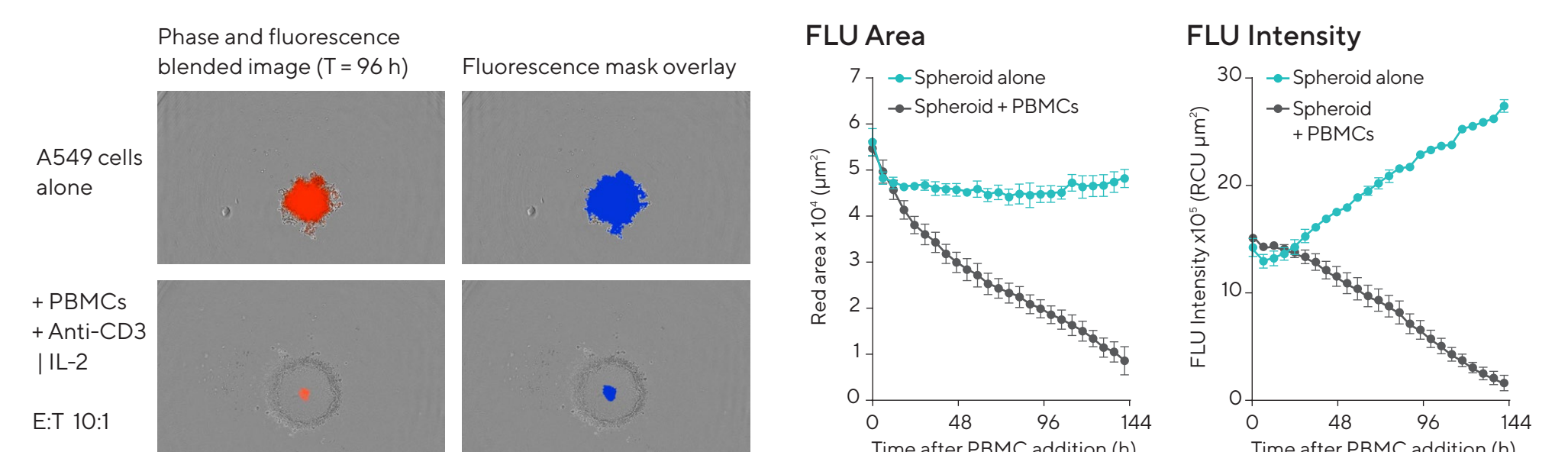
Incucyte® Reagents and Consumables

A suite of non-perturbing cell labeling and reporter reagents. Includes nuclear-targeted fluorescent proteins for cell counting, no-wash reagents for evaluating cell health, morphology, and function.

96-Well 3D Immune Cell Killing Assay Workflow

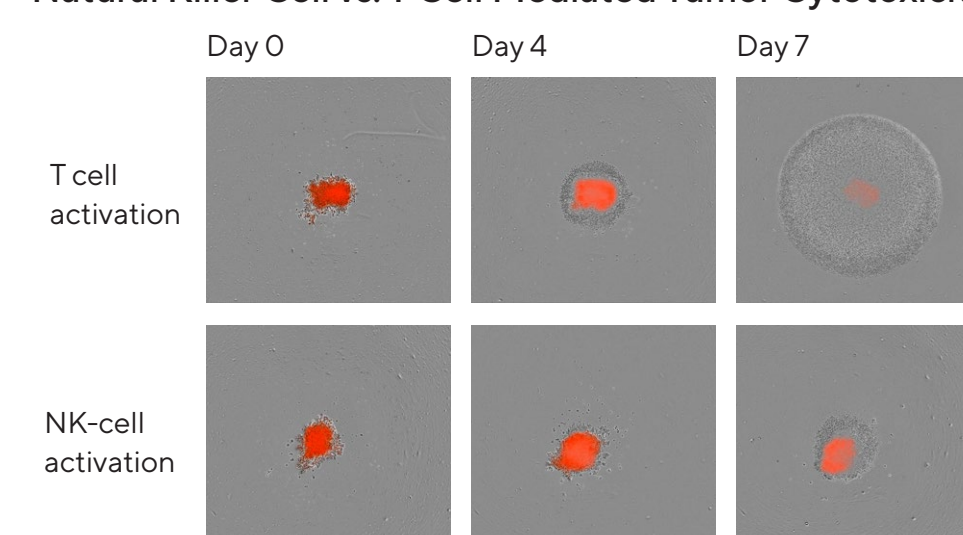


Fluorescence as a Measure of Spheroid Cytotoxicity



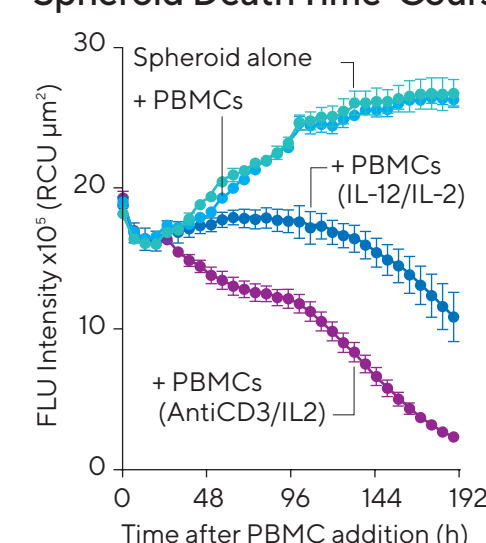
Activator-Dependent Tumor Cytotoxicity

Natural Killer Cell vs. T Cell Mediated Tumor Cytotoxicity

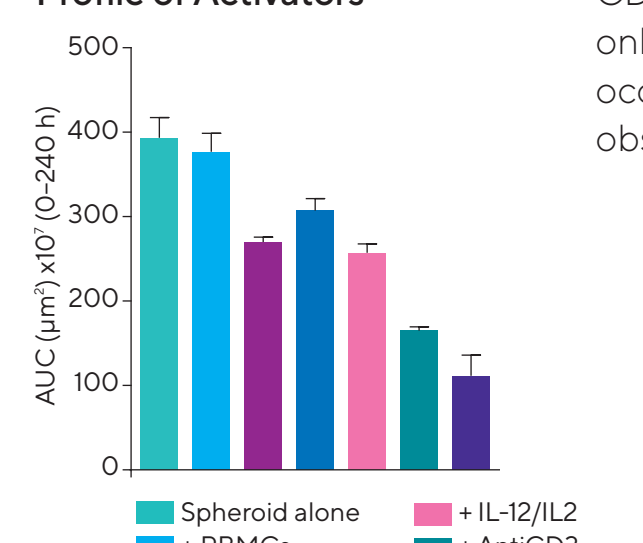


- Incucyte® Nuclight Red A549 cells were treated with isolated PBMCs activated with anti-CD3 | IL-2 (10 ng mL⁻¹) or IL-12 | IL-2 (10 ng mL⁻¹) targeting T cell or NK-cell populations, respectively.
- Activating different populations of PBMCs, resulted in differential effects on tumor spheroid destruction.
- In contrast to activated NK-cells, T cell activated populations exhibit increased proliferation.
- T cell mediated spheroid destruction occurred more rapidly compared to NK-cell mediated death.
- Data illustrates that T cell mediated spheroid destruction is significantly driven by the presence of anti-CD3 antibody.
- The frequencies of cell types within PBMC populations vary from donor to donor. Typically, CD3+ T cells account for 45-70% of PBMCs, while only 5-20% are NK cells. This variability in occurrence could explain the differential effects observed in tumor spheroid cytotoxicity.

Spheroid Death Time-Course



Profile of Activators



Incucyte® Nuclight Red A549 cells (2.5K/well). Optimized E:T ratios of 2.5:1