

## Validation of Novel Continuous Live-Cell Assays for Immune Cell Activation and Killing of Blood Cell Cancers

Nicola Bevan<sup>1</sup>, Hinnah Campwala<sup>1</sup>, Clare Szybut<sup>1</sup>, Kalpana Patel<sup>1</sup>, Dan Appledorn<sup>2</sup>, Tim Dale<sup>1</sup>, Derek Trezise<sup>1</sup>

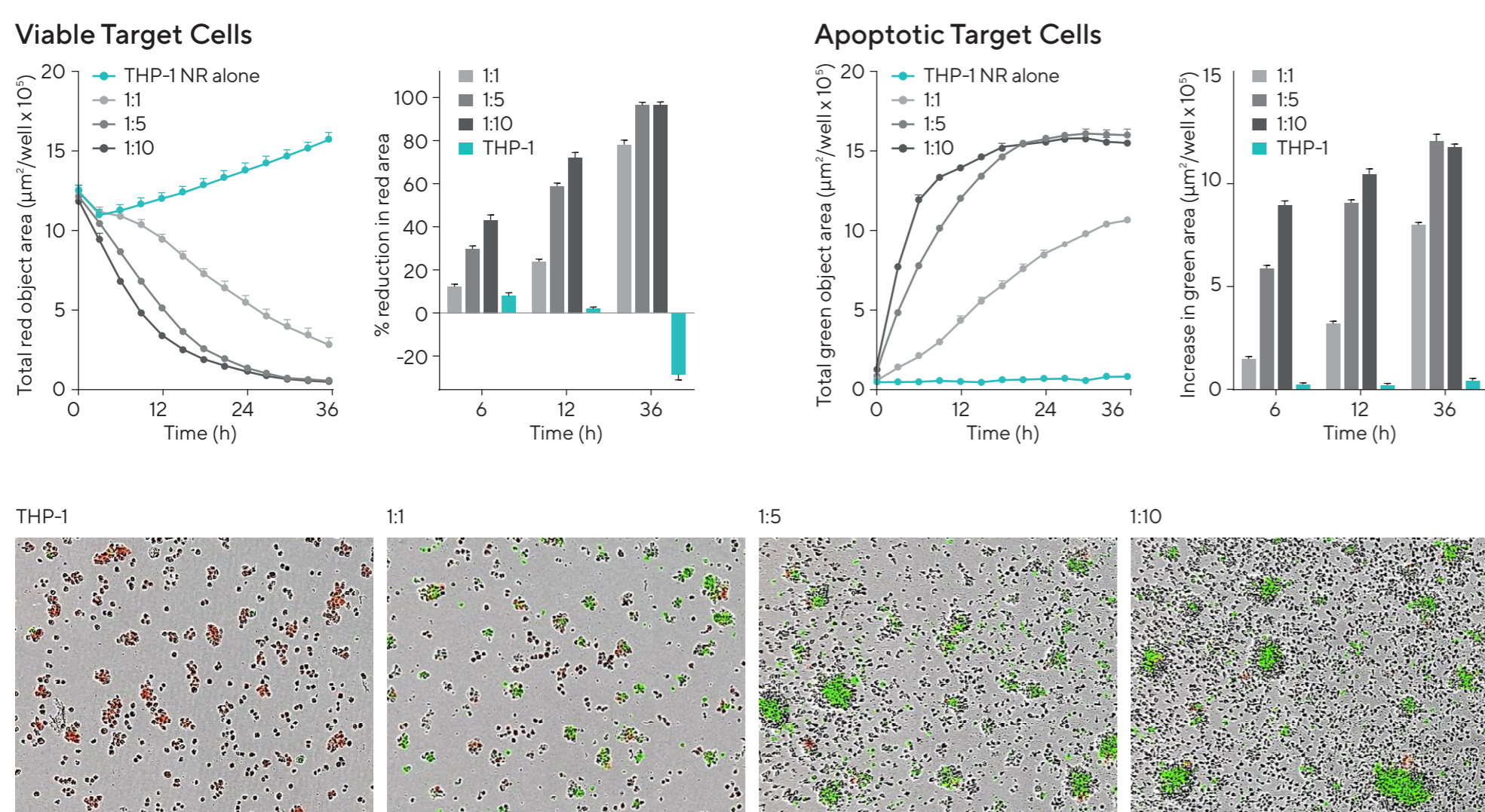
1. Essen BioScience Ltd., a Sartorius Company, Units 2 & 3 The Quadrant, Newark Close, Royston Hertfordshire SG8 5HL UK

2. Essen BioScience Inc., a Sartorius Company, 300 West Morgan Road, Ann Arbor, MI 48108 USA

### Overview

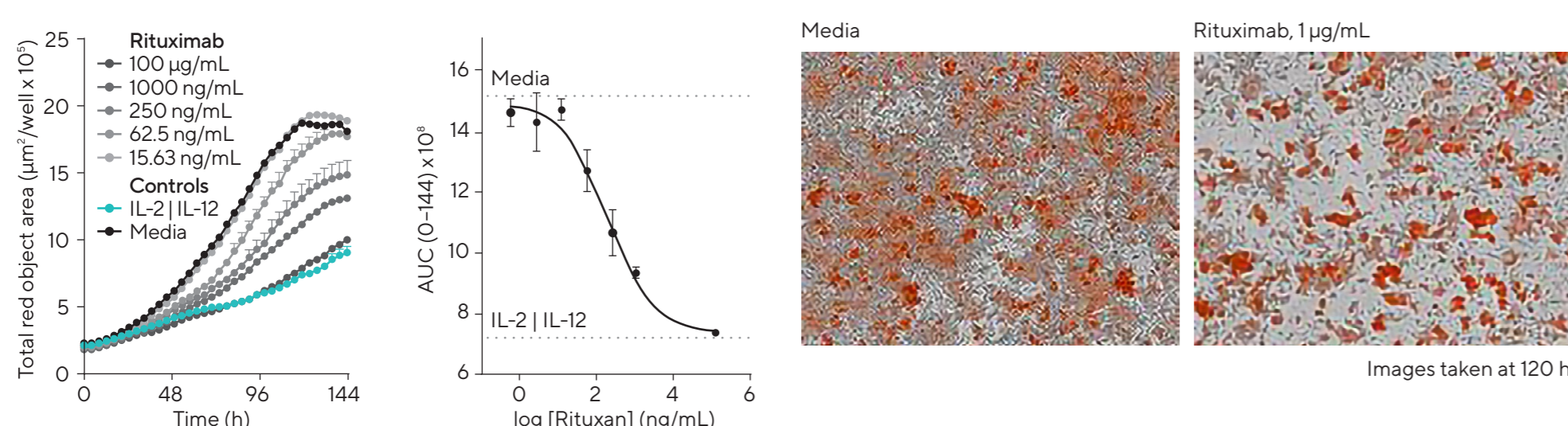
The blood cancers leukemia, lymphoma and myeloma are expected to cause the deaths of an estimated 56,840 people in the US in 2020. New immunological approaches afford great promise for improved therapies. Here, we describe novel high throughput live-cell image-based assays for immune cell activation and killing of blood cancer cells that are geared towards screening for new treatments for these malignancies. The data examples illustrate how continuous live-cell analysis can be used with suspension cells, to provide both additional biological insight (full time-courses, clustering information, morphology) and enhanced productivity (automation, miniaturization) compared to other techniques. We believe this approach will be a valuable addition to the technology toolbox for academic and industrial immuno-oncology researchers.

### Immune Cell Killing of Leukemic Monocytes



- THP-1 Nuclight® Red labeled monocytes (10 K/well) were seeded on 96-well PLO-coated flat-bottom plates in the absence or presence of pre-activated PBMCs (IL-2/α-CD-3 10 ng/mL, 4 days) and Incucyte® Annexin V green reagent
- Immune cell killing assay—note the time-dependent decrease in the number of red target cells and an increase in the Annexin V signal indicating cell death with increasing effector cell number
- Images show the aggregation of effector and target cells associated with cell death (increase in green fluorescence)

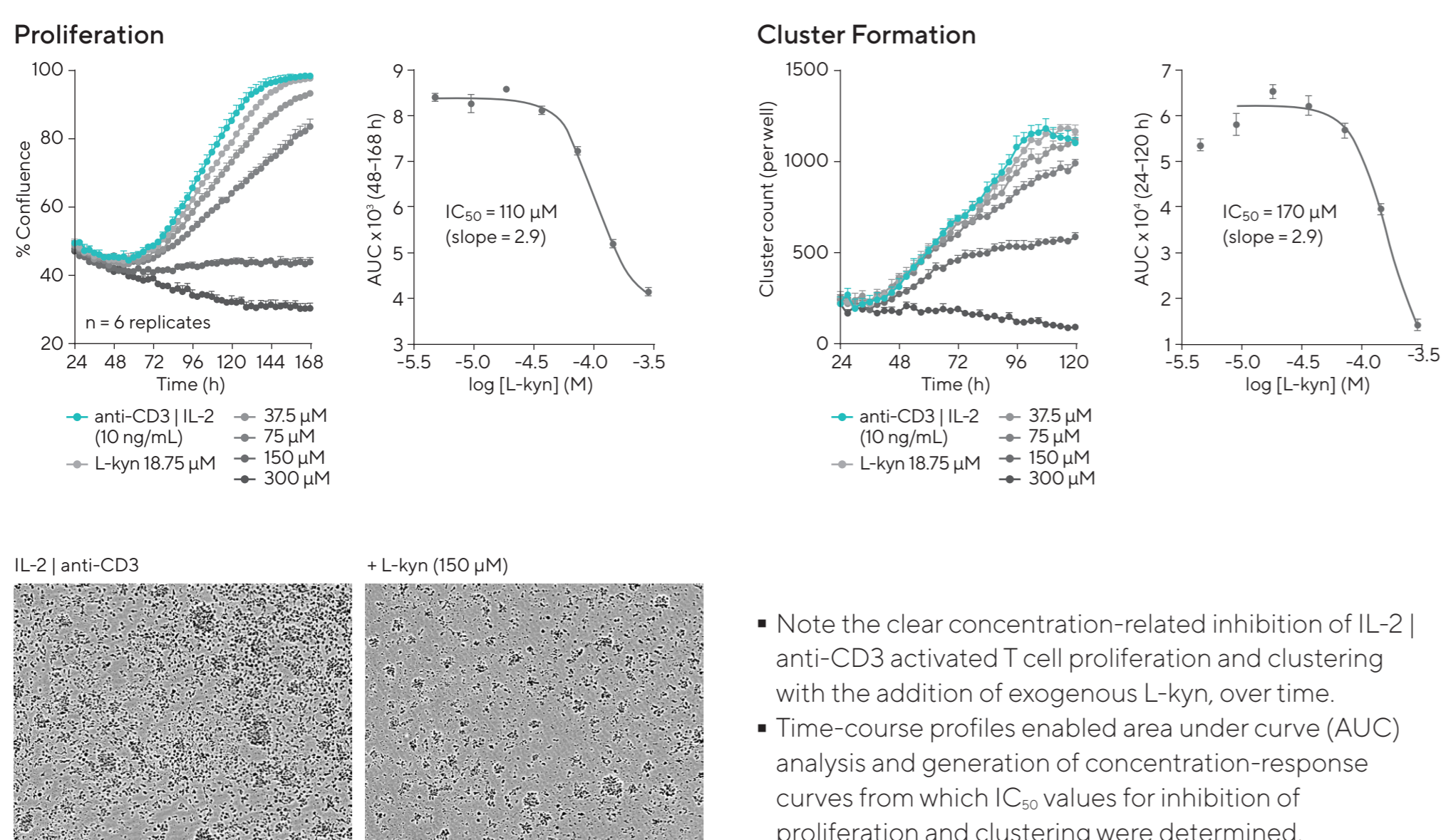
### Antibody-Dependent Cell-Mediated Cytotoxicity of B Cells



- CD20-positive Ramos B lymphocytes (10K/well), expressing Nuclight® Red, were seeded on 96-well PLO-coated flat bottom plates in combination with PBMCs (1:10 Target:Effector ratio)
- Rituximab induced a concentration-dependent ADCC (reduction in red fluorescence); IC<sub>50</sub> 0.22 µg/mL
- IL-2 | IL-12 activation of the NK cell population was used as a positive control

### L-Kynurenine Inhibits T Cell Proliferation and Clustering

- L-kynurenine (L-kyn) is a metabolite formed from the catabolism of L-tryptophan by the enzymes IDO and TDO
- Some cancers increase L-kyn production in a bid to block antigen-driven T cell proliferation and induce T cell death, thus allowing cancer cells to escape immune surveillance
- Inhibitors of IDO and/or TDO are, therefore, promising therapeutic targets for the treatment of cancer



- Note the clear concentration-related inhibition of IL-2 | anti-CD3 activated T cell proliferation and clustering with the addition of exogenous L-kyn, over time.
- Time-course profiles enabled area under curve (AUC) analysis and generation of concentration-response curves from which IC<sub>50</sub> values for inhibition of proliferation and clustering were determined.

### 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.

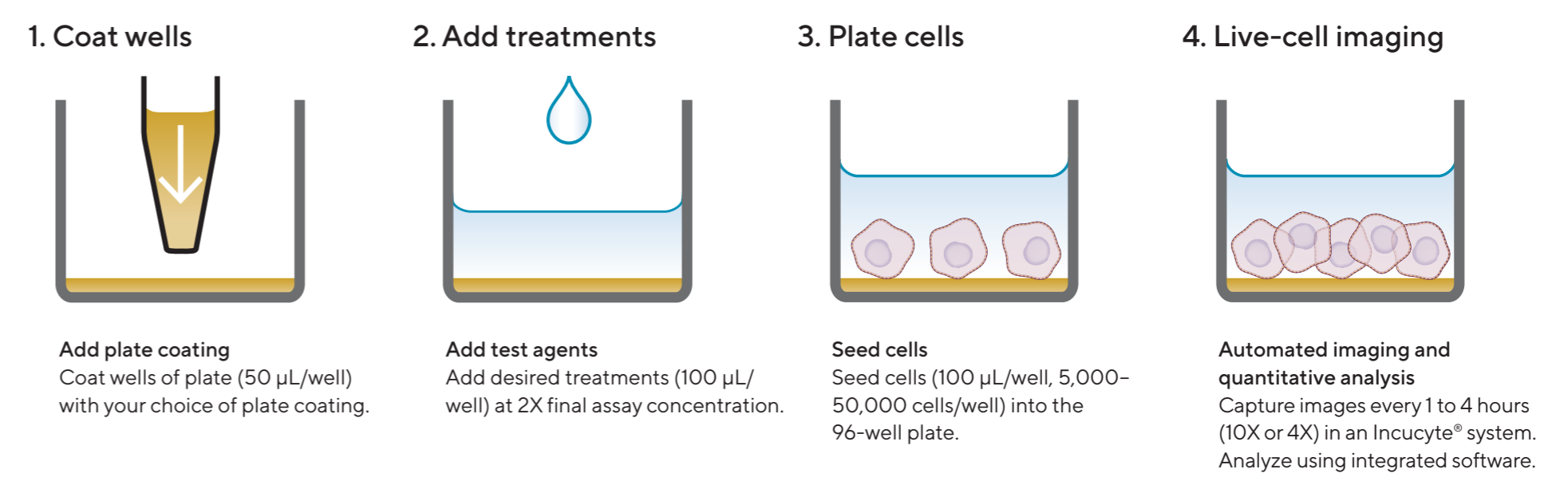
#### Incucyte® Software

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

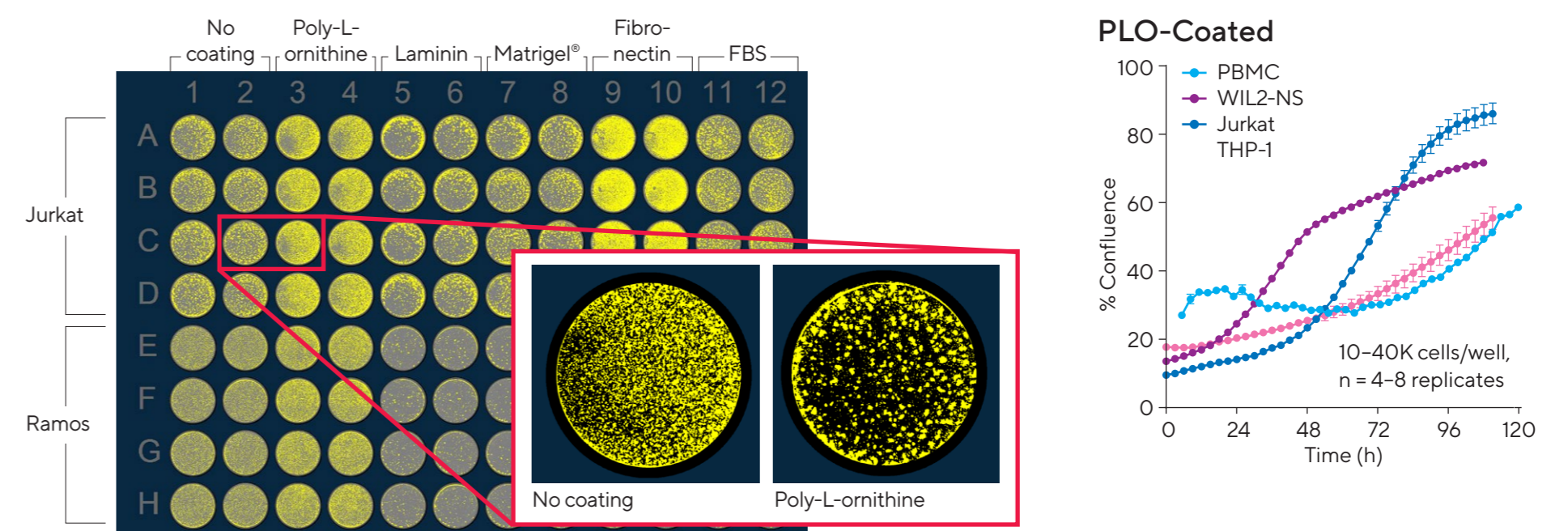
#### Incucyte® Reagents and Consumables

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.

### Non-Adherent Cell Assay Methodology

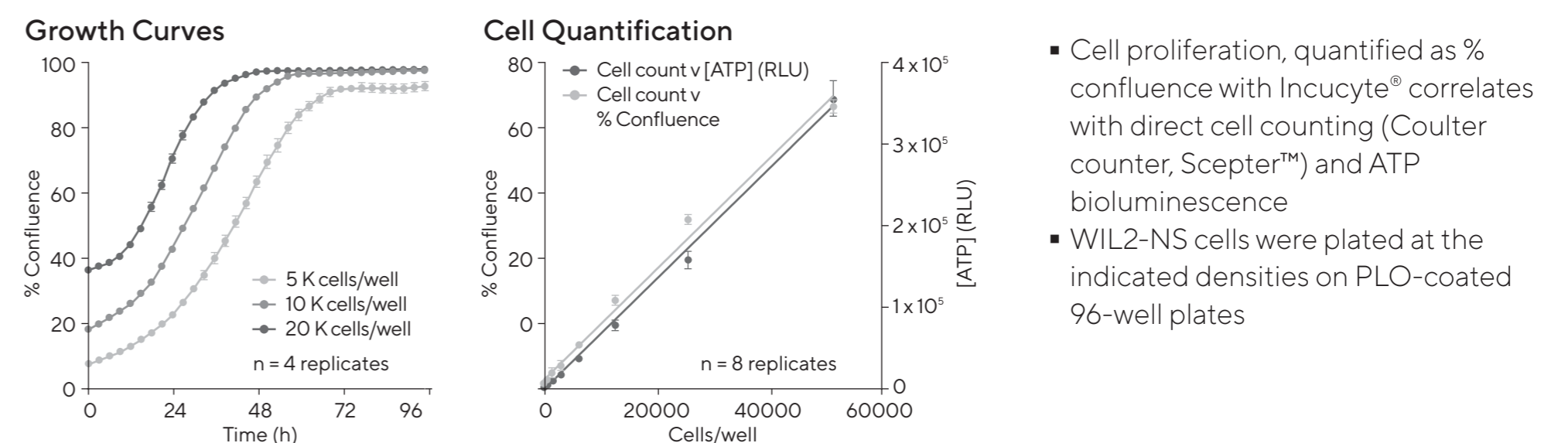


### Plate-Coatings: Impact on Distribution of Cells



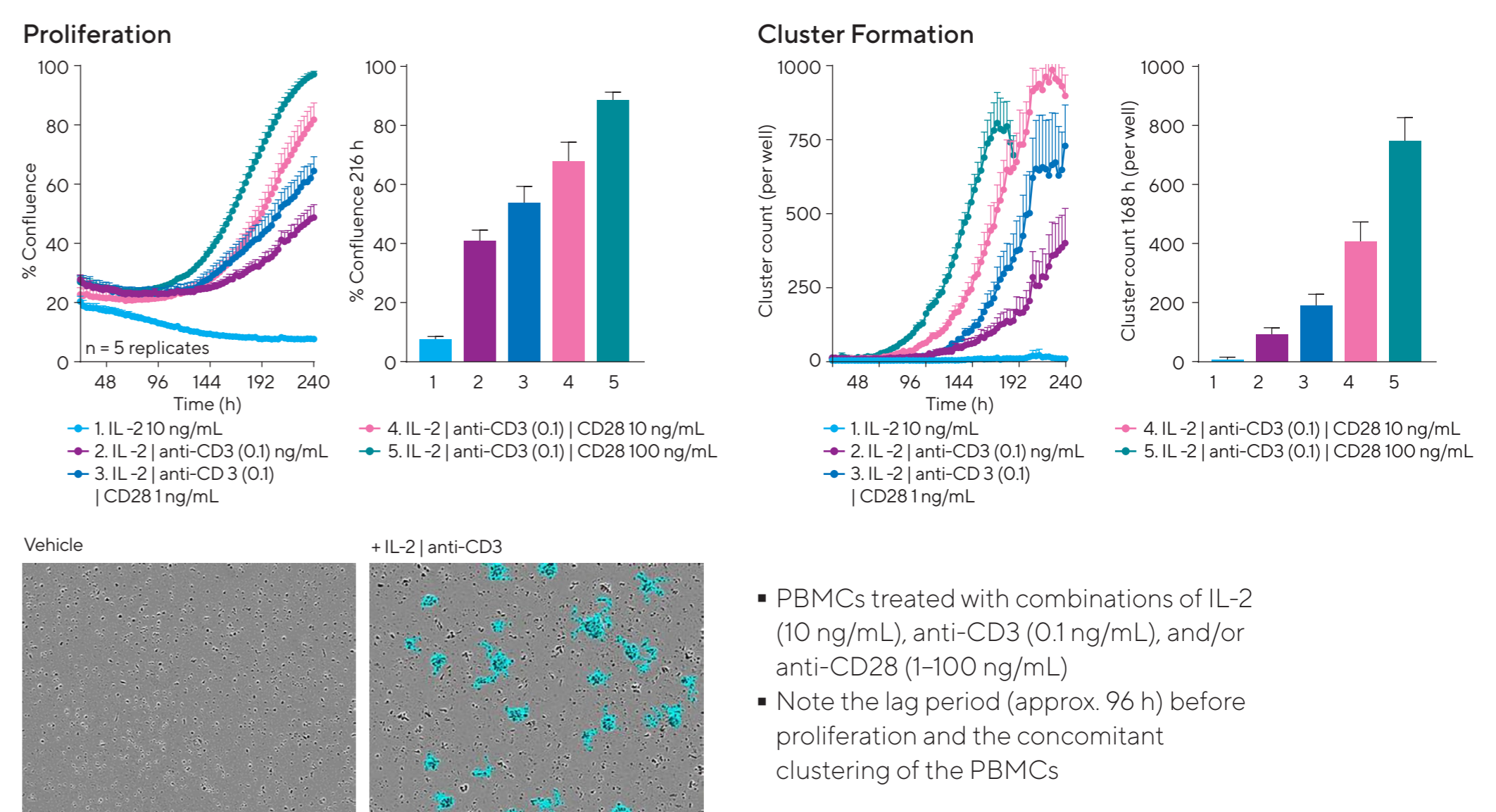
- Coating of plates with Poly-L-ornithine (PLO) facilitates uniform distribution of cells within each well
- Fibronectin is also suitable for most cell types but may enhance cell proliferation per se
- Yellow = phase confluence mask

### % Confluence as a Measure of Cell Proliferation



- Cell proliferation, quantified as % confluence with Incucyte® correlates with direct cell counting (Coulter counter, Scepter™) and ATP bioluminescence
- WIL2-NS cells were plated at the indicated densities on PLO-coated 96-well plates

### T Cell Activation is Stimulus and Concentration-Dependent



- PBMCs treated with combinations of IL-2 (10 ng/mL), anti-CD3 (0.1 ng/mL), and/or anti-CD28 (1-100 ng/mL)
- Note the lag period (approx. 96 h) before proliferation and the concomitant clustering of the PBMCs