

## Characterisation of Bi-specific T-cell Engager (BiTE) Antibody-mediated Cell Killing Using a Combined Live-cell and Flow Cytometry Workflow

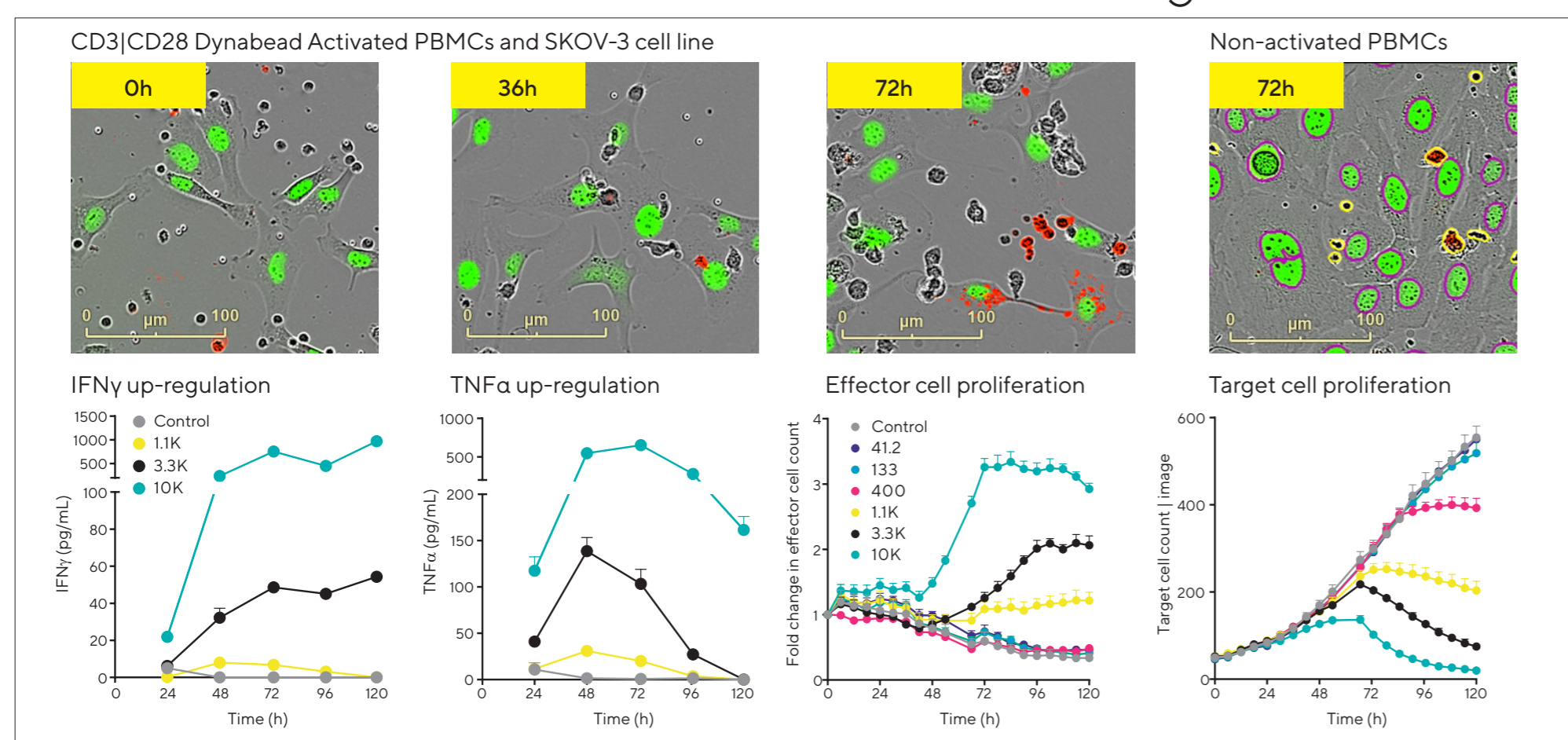
C. Szybut, L. Kelsey, K. McBain & T. Dale

Sartorius, Welwyn Garden City, AL7 3AX UK

### Summary & Impact

- The development of novel immunotherapies depends on the ability to evaluate immune cell effects on tumor cells.
- Here we describe a workflow combining live-cell analysis and flow cytometry enabling complementary parameters to be quantified.
- To demonstrate the potential insight gained by combining platforms, we have characterised immune cell killing induced by a  $\alpha$ -hCD3  $\times$  CD19 BiTE antibody.
- Phase and fluorescent images were captured with Incucyte<sup>®</sup> and segmented objects were quantified.
- Cytokine secretion and T-cell activation markers were evaluated using the iQue<sup>®</sup> platform.
- This combined approach has the potential to facilitate a greater understanding of the complicated biology involved in immune cell activation and killing of tumor cells.

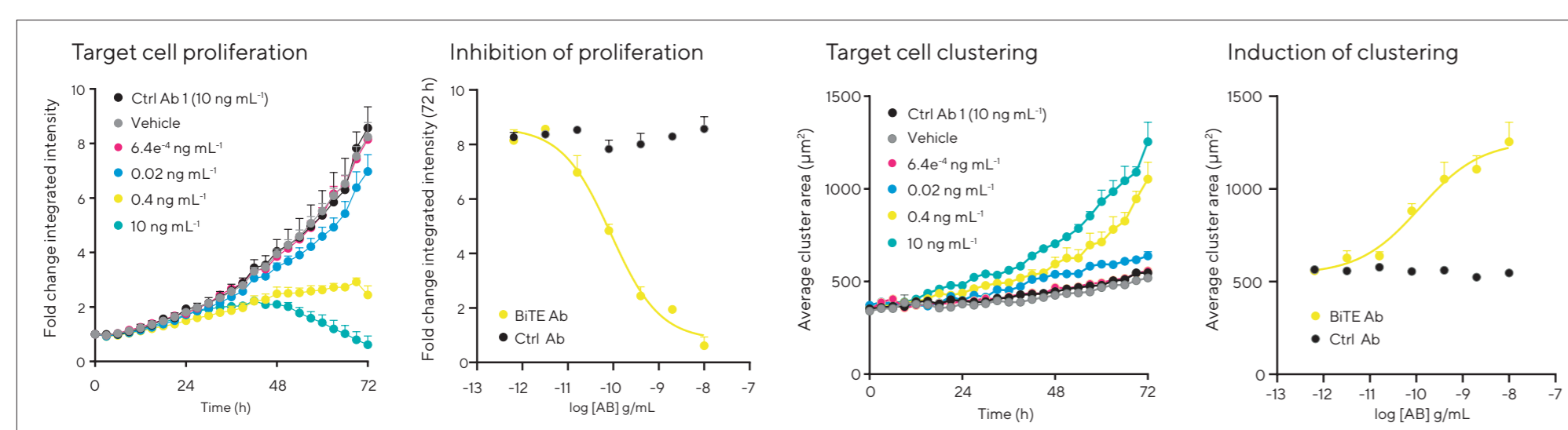
### Timeline of Immune Cell Activation and Killing



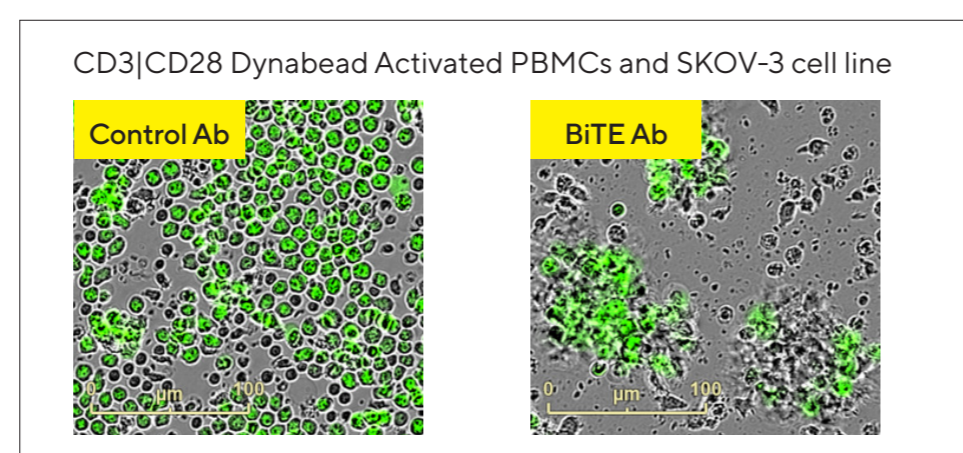
- Adherent SKOV-3 NucLight Green target cells (nuclear segmentation, pink mask) were seeded (4K | well) with non-labelled PBMCs (Cell-by-Cell segmentation, yellow mask) at a 1:5 target to effector cell ratio.
- Cell health was monitored using Annexin V Red reagent and PBMCs were activated with CD3 | CD28 Dynabeads<sup>®</sup> (41-10,000 beads) for 5 days, images were captured and analysed using Incucyte<sup>®</sup>.
- Supernatant samples were taken daily (10  $\mu$ L) and secreted cytokines IFN $\gamma$  and TNF $\alpha$  were quantified using the QBeads component from the T-Cell Activation Cell and Cytokine Profiling Kit.
- Cytokine levels were shown to increase in both a time- and concentration-dependent manner.
- Effector cell population increased over time with the top three bead densities eliciting the greatest response 48h post bead addition.
- CD3 | CD28 activation yielded a time- and concentration-dependent decrease in SKOV-3 cell numbers with the maximal amount of beads inducing near complete abolition (>95% at 120h).
- This timeline of events details how activation of the T-cell subpopulation ultimately leads to immune cell-mediated cytotoxicity.

### BiTE Induced Spatial and Proliferative Changes in Target Cells

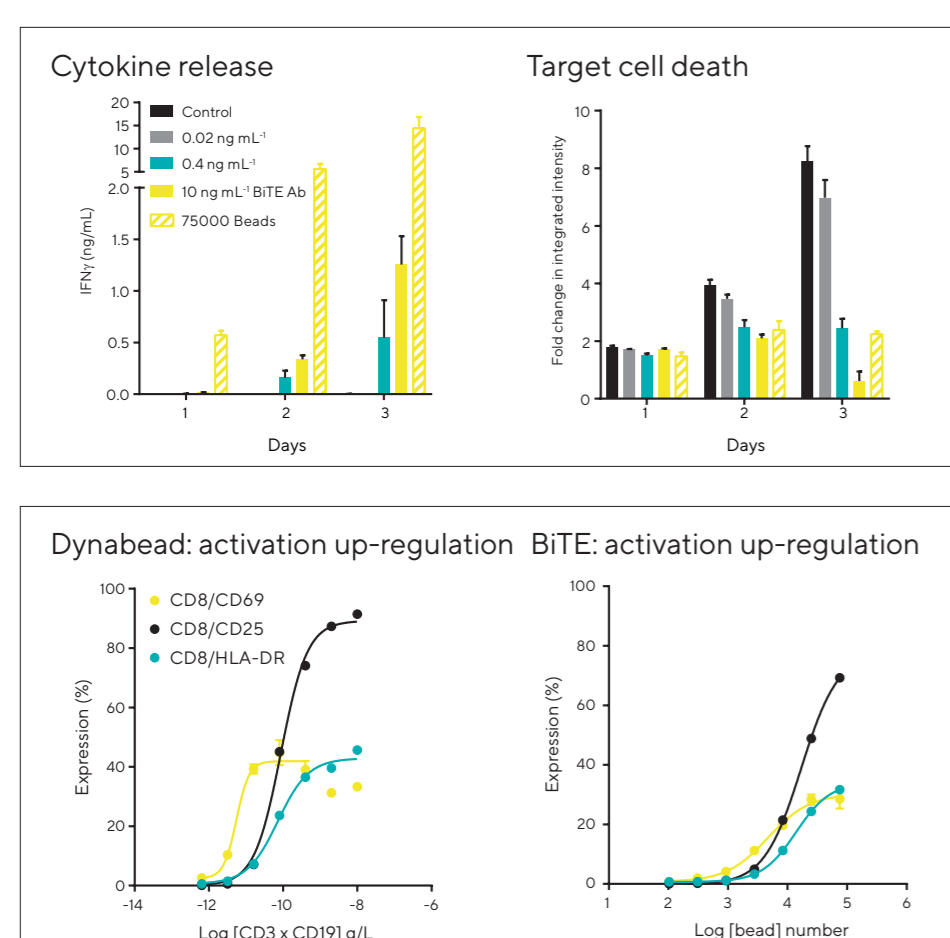
N2A cells were pre-treated with staurosporine (24 hrs), labeled with the Incucyte<sup>®</sup> pHrodo<sup>®</sup> Orange Cell Labeling Kit, and added to pre-plated primary rat microglia (Brain Bits, 20,000 K/well). N2A cells alone have minimal fluorescence (left image). Engulfment of labeled apoptotic N2A cells by microglia causes an increase in orange fluorescence (fluorescent image) that is segmented using automated Incucyte<sup>®</sup> Live-Cell Analysis Software (fluorescent image with mask). A kinetic N2A density-dependent response is quantified over time (right).



- Ramos NucLight Green target cells were seeded (15K | well) with non-labelled PBMCs (1:5 target to effector cell ratio) and activated with BITE antibody ( $\alpha$ -hCD3  $\times$  CD19, 0.6 pg-10 ng mL<sup>-1</sup>) or control antibody ( $\alpha$ -hCD3  $\times$   $\beta$ Gal, 0.6 pg-10 ng mL<sup>-1</sup>).
- Representative images at 72h in the presence of control or BITE Ab.
- Phase and fluorescence images acquired using Incucyte<sup>®</sup> were quantified, taking into account both the brightness and size of green objects (GCU  $\times$   $\mu$ m<sup>2</sup>) and expressed as fold change in integrated intensity.
- Inclusion of BITE antibody induced a concentration-dependent increase in target cell cytotoxicity as well as target cell clustering between 24 and 72h.



### Differential Profiles Between BiTE- and CD3 | CD28-induced Activation



- Cytokine samples taken daily (10  $\mu$ L) from the BITE ( $\alpha$ -hCD3  $\times$  CD19) co-culture were compared with CD3 | CD28 Dynabead activation.
- BITE Ab elicited a time- and concentration-dependent increase in IFN $\gamma$  levels, reaching a maximal concentration of 1.3  $\pm$  0.3 ng mL<sup>-1</sup>.
- Although CD3 | CD28 activation yielded a significantly higher IFN $\gamma$  level (14.4  $\pm$  2.5 ng mL<sup>-1</sup>), BITE stimulation resulted in greater target cell death (93% compared to 72% following Dynabeads).

- After 3 days cells were lifted and labelled using the TCA-kit antibody cocktail and membrane integrity dye prior to analysis using the iQue<sup>®</sup> platform.
- Inclusion of BITE antibody induced a clear left shift in CD69 expression pattern when compared to bead activation, with low concentrations of BITE (20 pg mL<sup>-1</sup>) capable of producing almost exclusive expression of the CD69 early activation marker (EC<sub>50</sub> value 5.5 pg mL<sup>-1</sup>).

### Incucyte<sup>®</sup> System and iQue<sup>®</sup>3 Combined



#### Incucyte<sup>®</sup> S3 Live-Cell Analysis System

A fully automated phase contrast and two-color fluorescence system that resides within a standard cell incubator for optimal cell viability. Designed to scan plates and flasks repeatedly over time.



#### iQue<sup>®</sup>3 Flow Cytometer

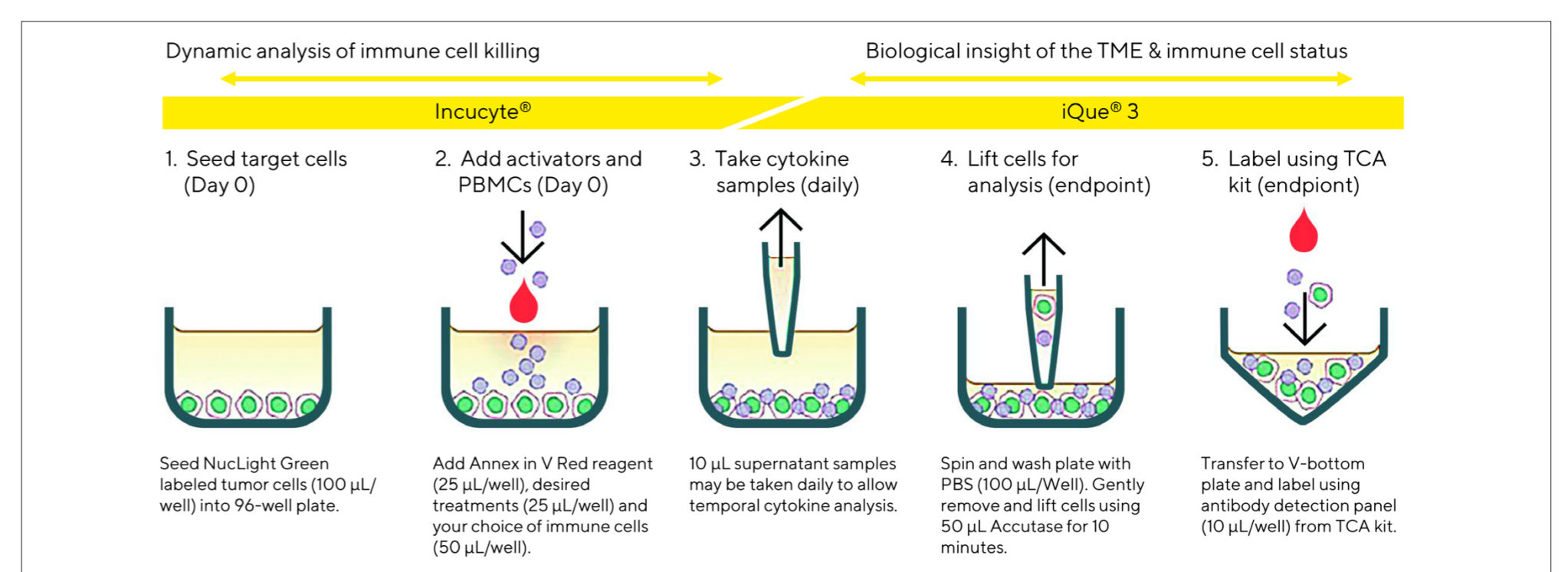
An advanced flow cytometry platform with a patented sampling method allowing for rapid sample acquisition which delivers fast actionable results. Capable of handling 96 and 384 well plates.



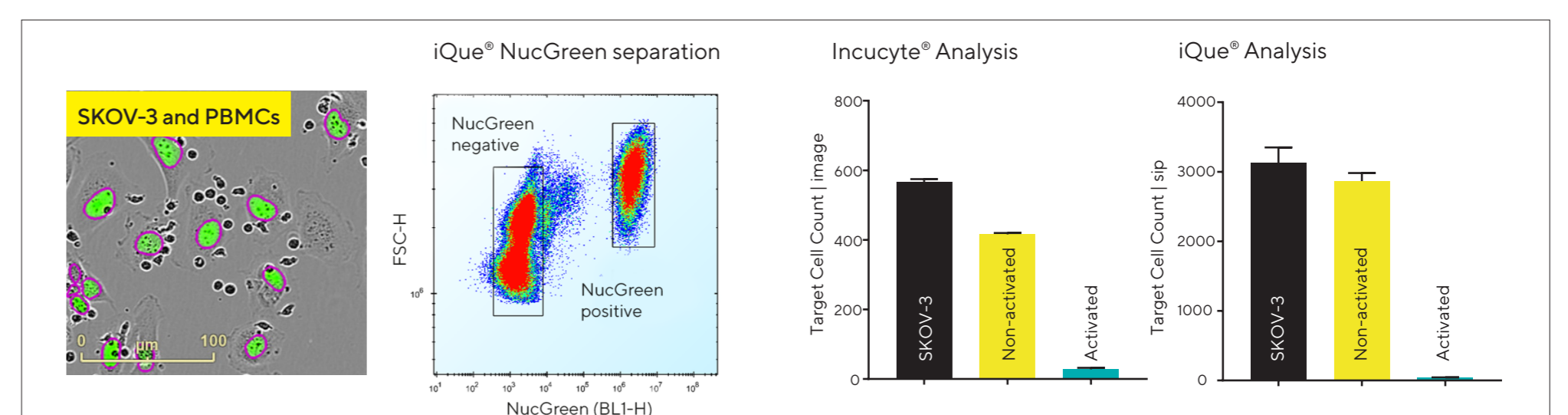
#### Sartorius Reagents

A suite of reagents, kits and protocols for cell health and function screening.

### Live-cell Analysis and Flow Cytometry Workflow: Methodology

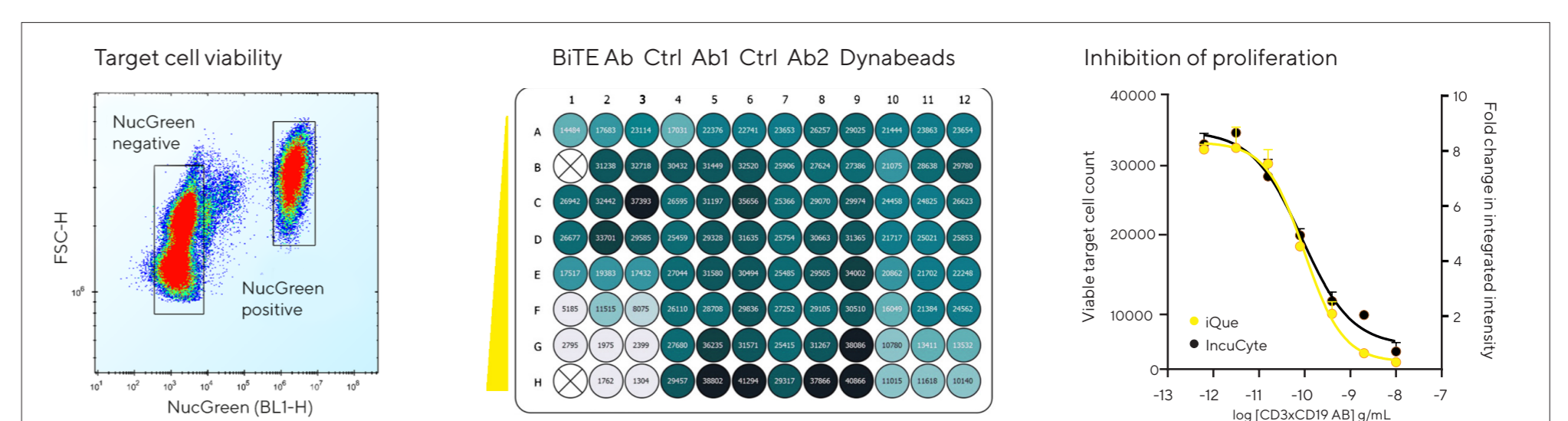


### Reagents and Kits are Compatible Across Both Platforms



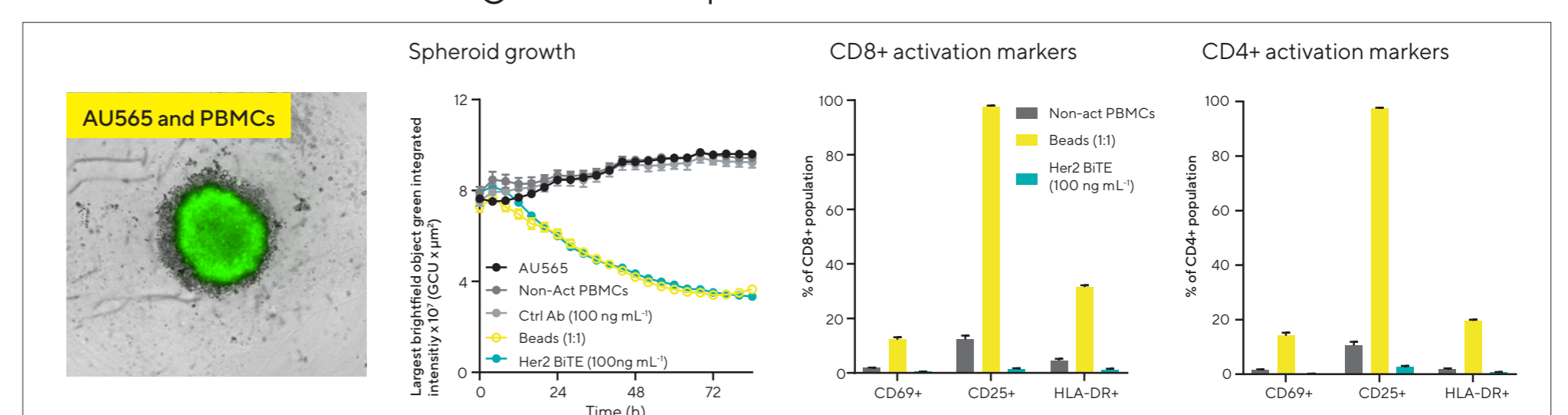
- NucLight Green cell lines were seeded with PBMCs (label free) in the presence of an activator e.g. CD3 | CD28 Dynabeads, BITE Ab.
- Incucyte<sup>®</sup> target cell quantification of adherent SKOV-3 cells (nuclear segmentation, pink mask) describes the effect of activated PBMCs, where target cell numbers were significantly reduced in comparison to the target cells alone.
- Cells were then lifted from the assay plate and analysed using the iQue<sup>®</sup>. Gating using Forecy software enabled separation of the target and effector cell types based on positive or negative green fluorescent populations.
- Cell count analysis from iQue<sup>®</sup> analysis showed activated PBMCs induced high levels of cytotoxicity, this trend was highly comparable across both platforms.

### Quantitative Pharmacology Demonstrates Comparable Results



- Ramos NucLight Green target cells were seeded (15K | well) with non-labelled PBMCs (1:5 target to effector cell ratio) and activated with BITE ( $\alpha$ -hCD3  $\times$  CD19, 0.6 pg-10 ng mL<sup>-1</sup>) or control Abs ( $\alpha$ -hCD3  $\times$   $\beta$ Gal and  $\alpha$ -hCD19  $\times$   $\beta$ Gal, 0.6 pg-10 ng mL<sup>-1</sup>).
- Viable and non-viable target cells were identified using a cell membrane integrity dye (Sartorius) and gated populations were quantified.
- Live events were displayed using a heat map where increasing shades of purple equated to increased number of target cells.
- Activation with BITE and CD3 | CD28 Ab induced concentration-dependent decrease in target cell numbers.
- Importantly proliferation potency concentrations were comparable between the iQue<sup>®</sup> and Incucyte<sup>®</sup> (IC50 values 99 pg ml<sup>-1</sup>; 88 pg ml<sup>-1</sup> respectively).

### Immune Cell Killing of 3D Spheroids



- AU565 NucLight Green labelled breast cancer target cells were seeded in a round bottom ULA plate (4K | well) and allowed to form spheroids over 3 days.
- Spheroids were subsequently treated with PBMCs (1:5 target to effector cell ratio) which were either non-activated or in well-activated with an  $\alpha$ -hCD3  $\times$  Her-2 BITE Ab or CD3 | CD28 Dynabeads.
- Spheroid growth in the presence of BITE-activated (100 ng ml<sup>-1</sup>) PBMCs was inhibited by 64  $\pm$  0.02% (comparable to Dynabead activation).
- Interestingly, the BITE Ab did not elicit a classic T-cell activation response suggesting an alternative mechanism of action.