

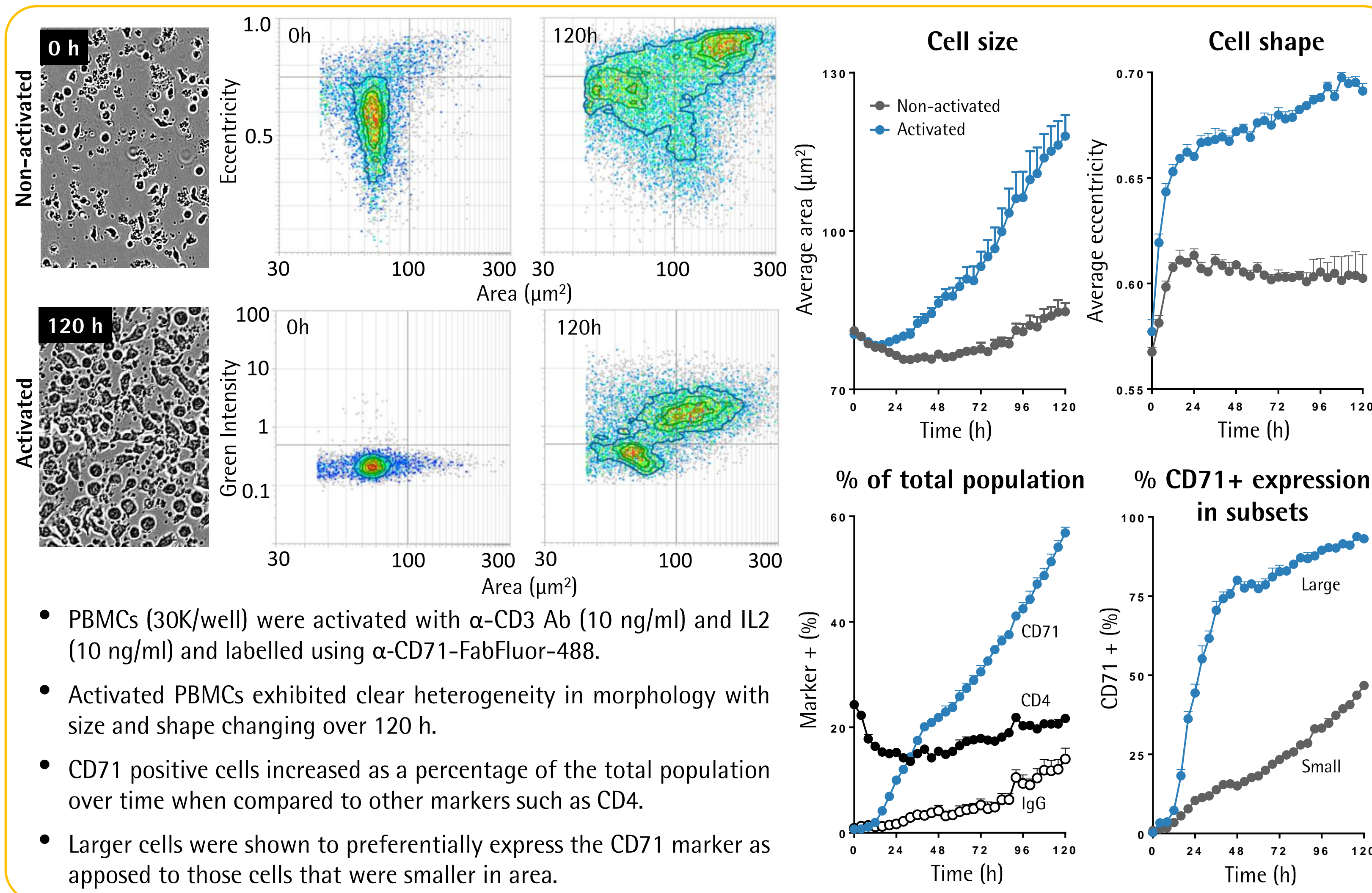
Quantifying immune cell subsets in living cultures over time using IncuCyte® live-cell analysis

C. Szybut, N. Bevan, H. Campwala, L. Kelsey, N. Dana, T. Jackson, N. Holtz, E. Endsley, T. Dale, D. Trezise
Essen BioScience, Welwyn Garden City, AL7 3AX UK & Ann Arbor, MI, 48108, USA

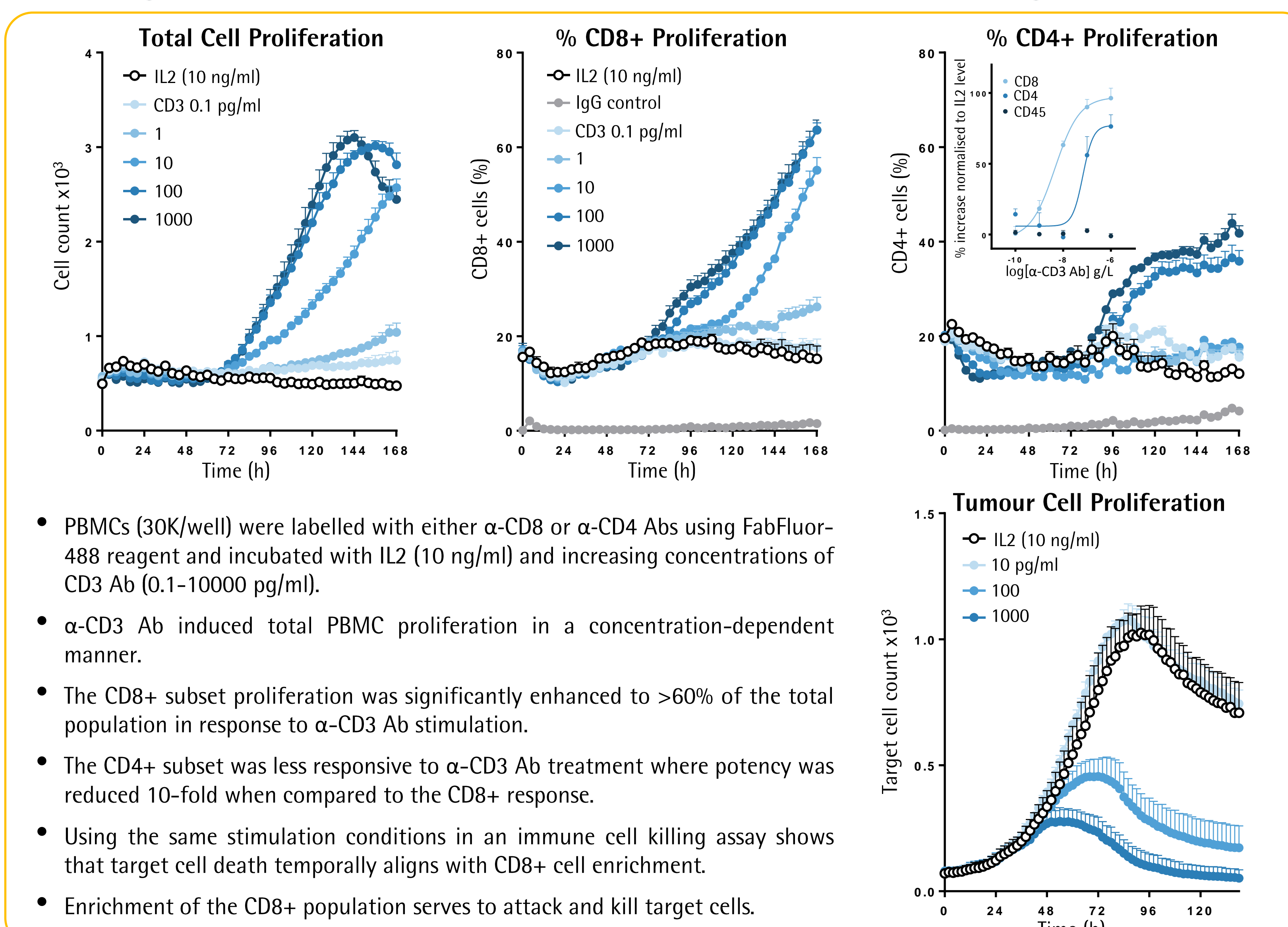
Summary & Impact

- Heterogeneity exists in all cellular populations, ranging from the cell types present to differences at the genetic level or stage of cell cycle. This heterogeneity plays an important role in how populations react in response to therapeutics and biological stimuli.
- To date IncuCyte® analysis has been solely based on population-averaged measures whereby object (cell) data is consolidated into an aggregate metric.
- However, effects on subpopulations can sometimes be masked by larger numbers of 'non-responsive' cells or similar sized populations may produce opposite responses that result in a net zero result.
- Analysis at the cell-by-cell level promises valuable and additional biological insight beyond which whole population measures may deliver.
- The IncuCyte cell-by-cell software module and FabFluor-488 labelling protocol provides automated image capture and analysis in real time in order to provide an integrated solution for monitoring at the cell-by-cell level and increase biological insight.
- Here, we present data validating the cell-by-cell analysis segmentation and classification and use examples of PBMC activation to show parameter changes over time. Importantly, these changes can be related to function.

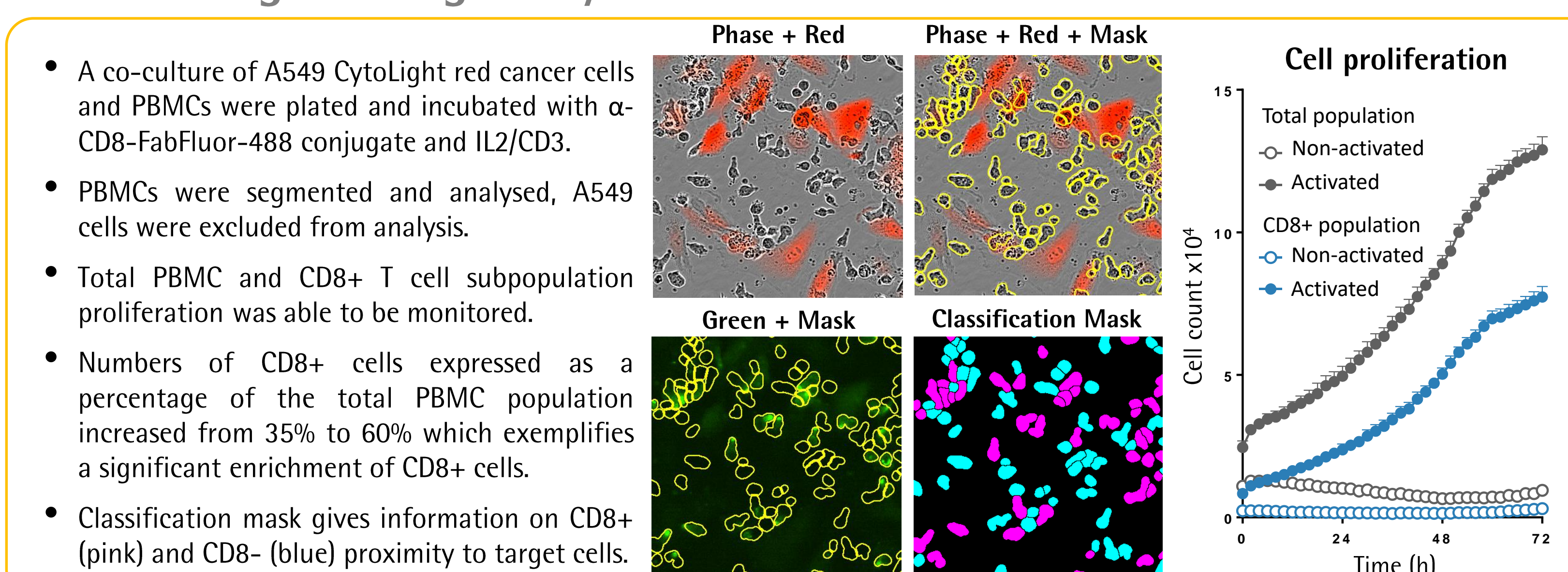
Monitoring PBMC activation: morphology & protein expression



Linking PBMC subset proliferation to immune cell killing



Monitoring heterogeneity in the tumour microenvironment

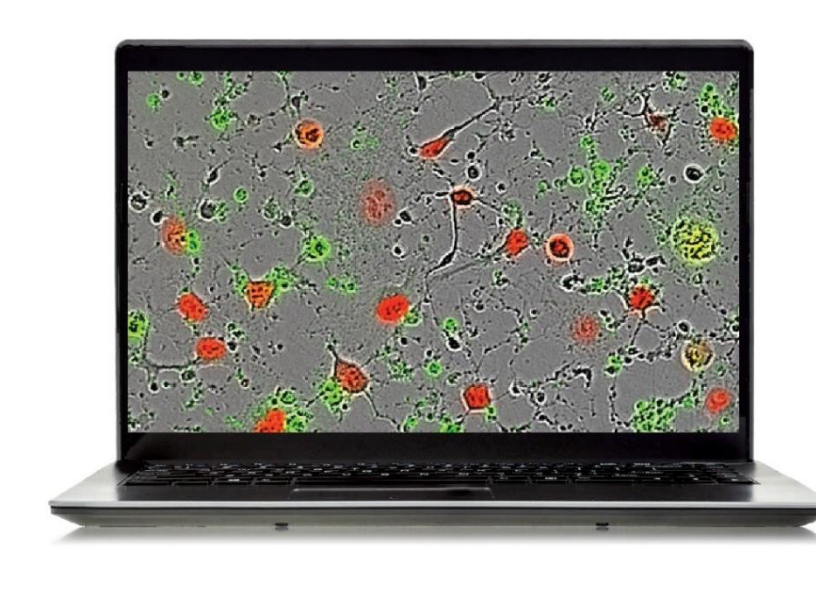


IncuCyte® System for Continuous Live-cell Analysis: Methodology



IncuCyte® S3 Live-Cell Analysis System

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



IncuCyte® Software

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

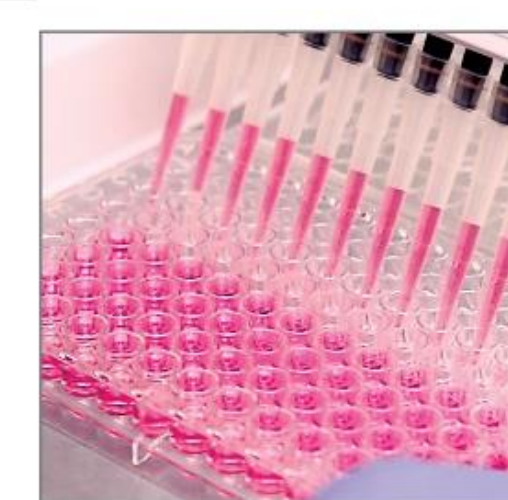
NEW: IncuCyte® S3 Cell-by-Cell Software Module



IncuCyte® Reagents and Consumables

A suite of non-perturbing cell labeling and reporter reagents. Includes nuclear-targeted GFP and RFPs for cell counting plus no-wash cell health reagents for apoptosis and cytotoxicity.

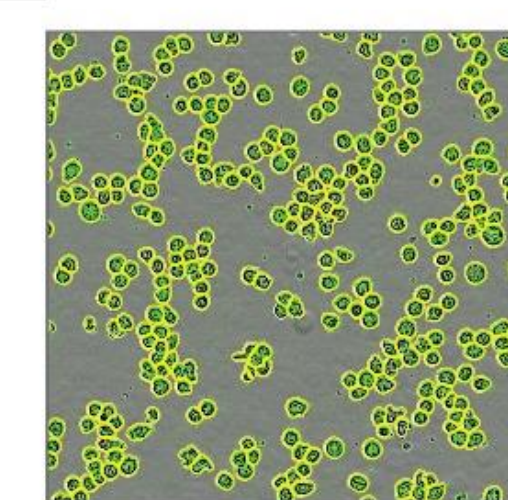
1 PREPARE ASSAY PLATES 2 ACQUIRE IMAGES 3 IDENTIFY OBJECTS 4 MEASURE and CLASSIFY 5 VISUALIZE DATA



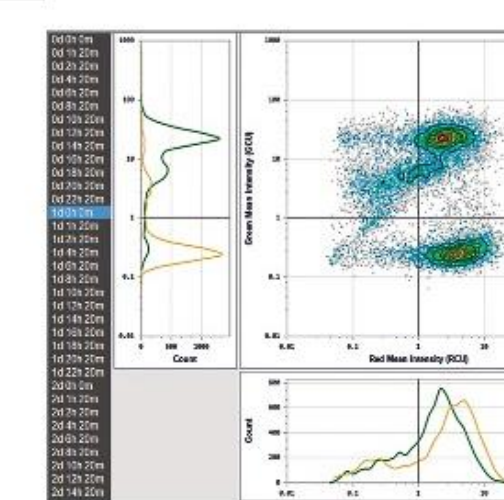
Seed cells, add treatments and optional non-perturbing reagents.



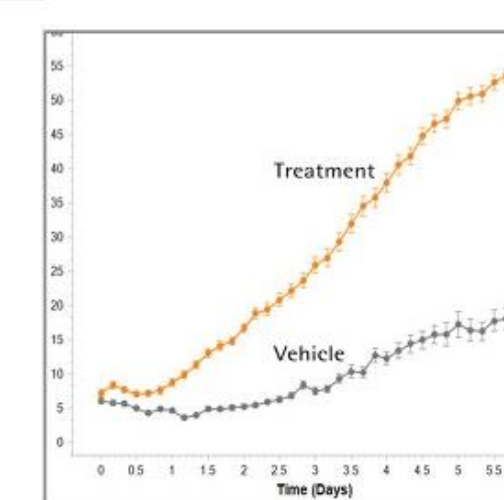
Acquire images with 20x objective, at appropriate time interval and length for assay.



Segment image and create a mask that delineates cell boundaries.

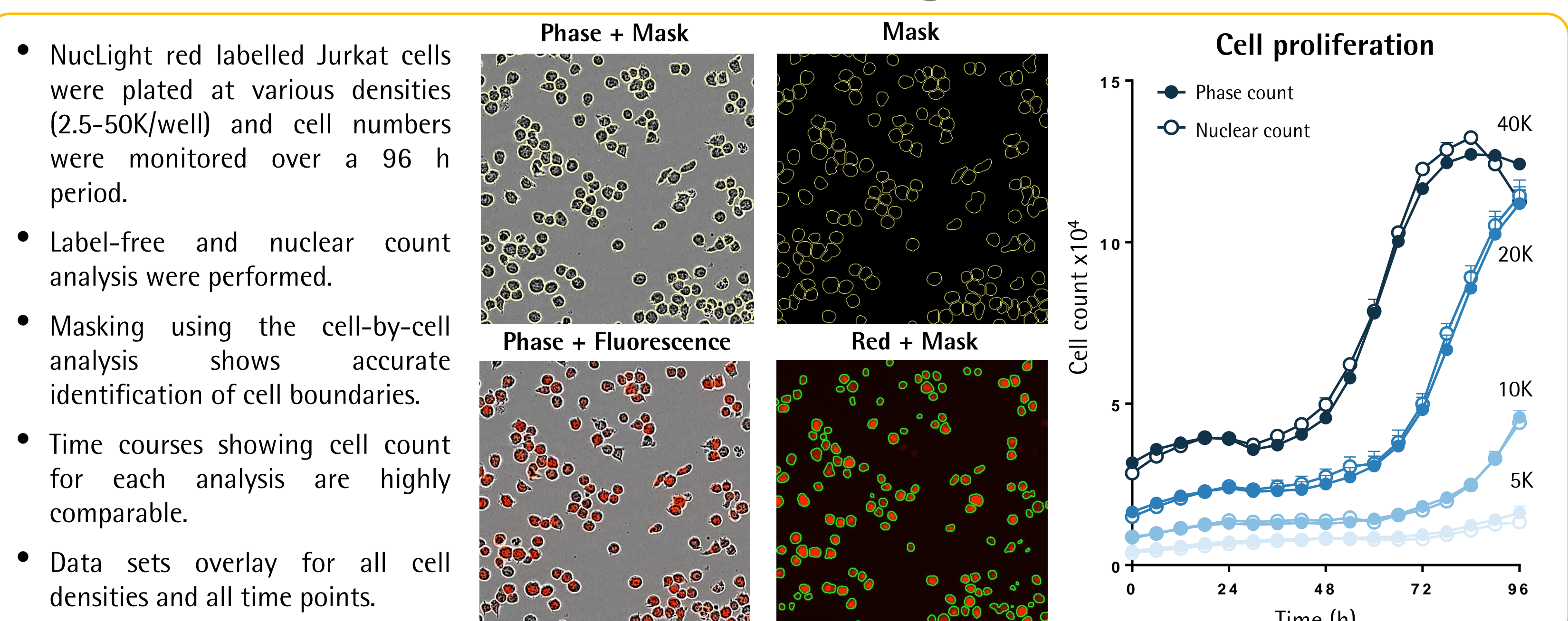


Generate metrics for each cell in the mask. Classify cells into subsets based on shape, size or fluorescence properties.

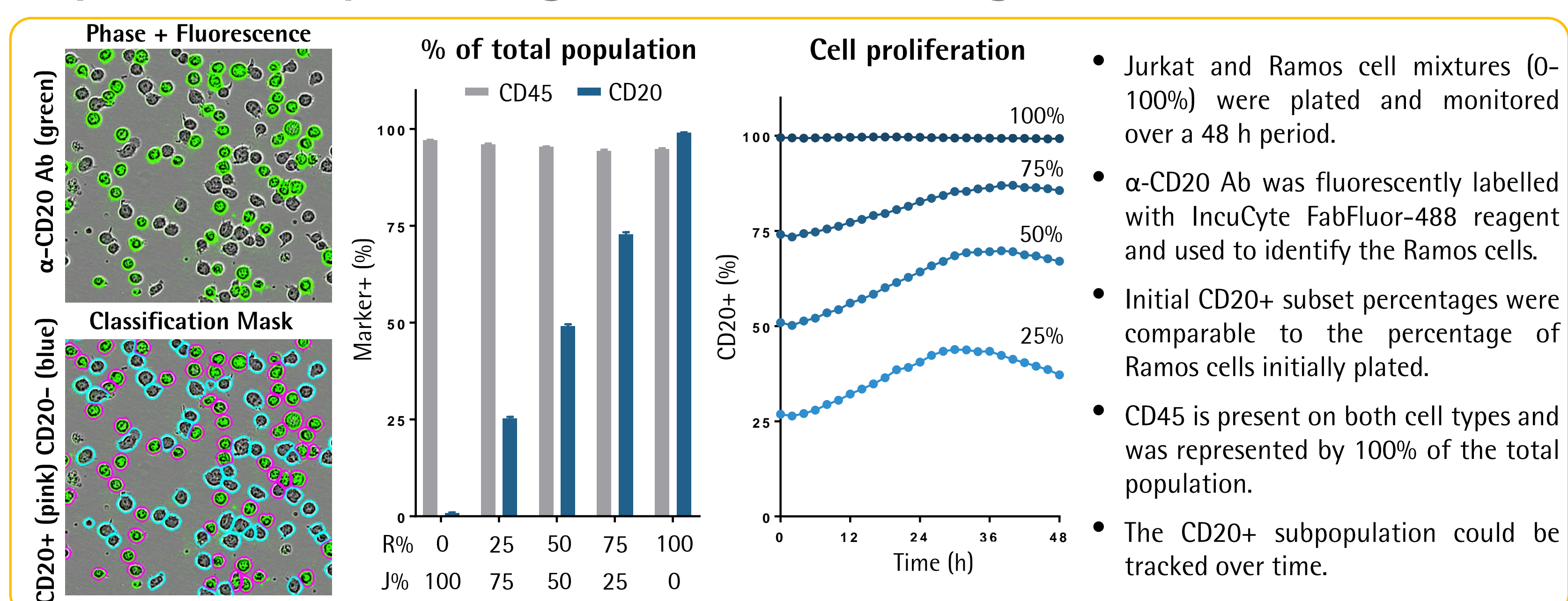


Graph changes in subsets over time.

Label-free, accurate non-adherent cell segmentation



Specific, non-perturbing fluorescent labelling



Track cell health of identified subsets

