

## Visualization and quantification of phagocytosis using live-cell analysis and advanced flow cytometry

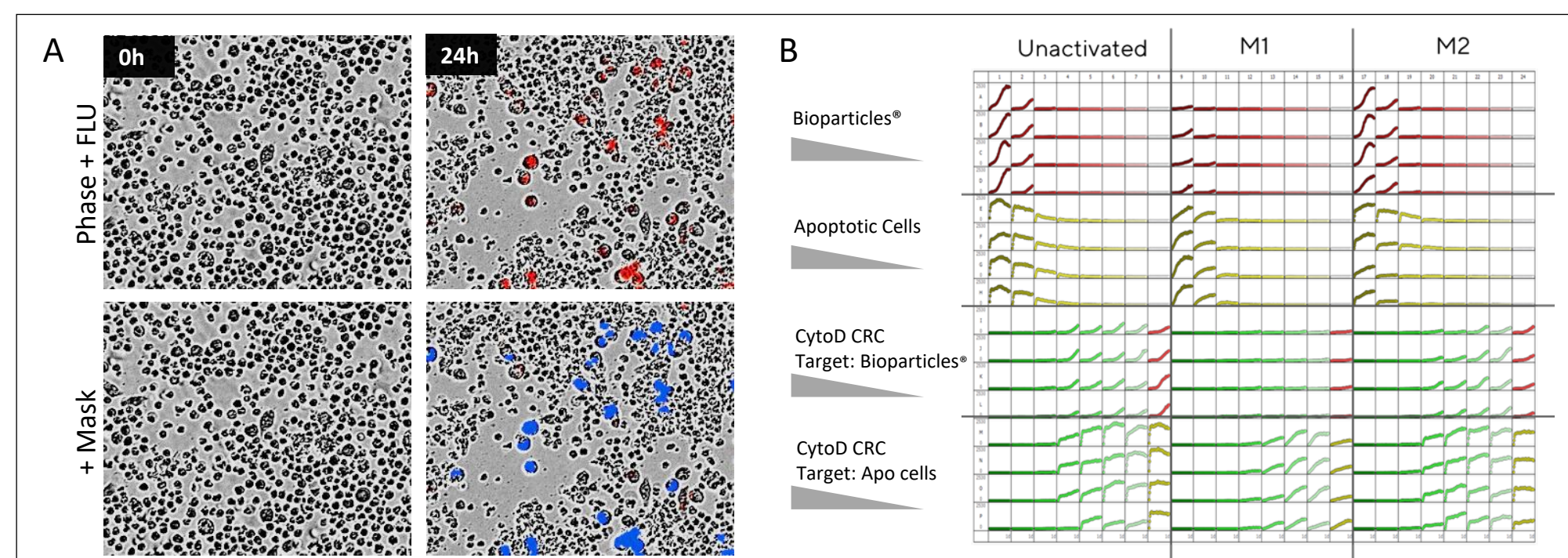
J. Trigg<sup>1</sup>, K. Mcbain<sup>1</sup>, C. Weldon<sup>2</sup>, G. Lovell<sup>1</sup>, N. Bevan<sup>1</sup>, C. Szybut<sup>1</sup>, T. Dale<sup>1</sup> and M. Carter<sup>2</sup>

<sup>1</sup>BioAnalytics, Sartorius (Essen BioScience), Royston, Hertfordshire, UK; <sup>2</sup>Sartorius, Albuquerque, NM  
\*Corresponding author: [Jasmine.Trigg@sartorius.com](mailto:Jasmine.Trigg@sartorius.com)

### Summary & Impact

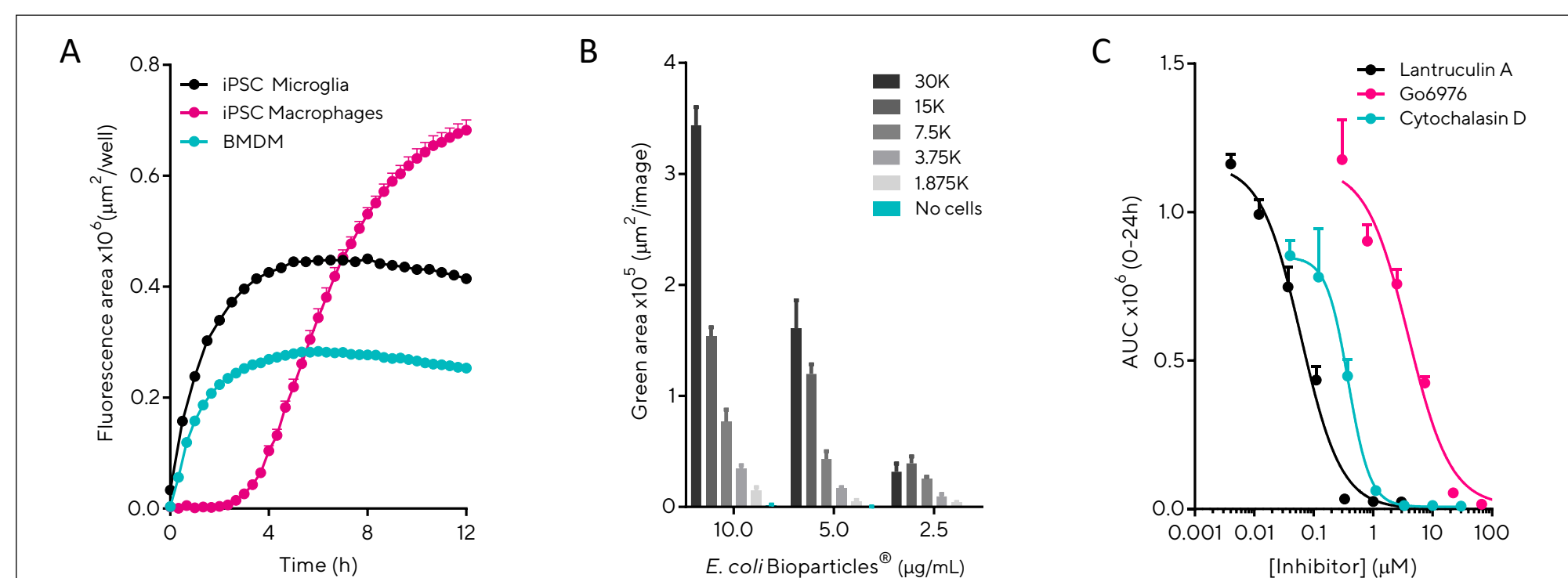
- Phagocytosis, a specific form of endocytosis, is a critical component of innate and adaptive immune responses.
  - Viable cells can be targeted with pro-phagocytic agents, such as monoclonal antibodies (mAbs), that promote engulfment and clearance through antibody-dependent cellular phagocytosis (ADCP), or by blockage of "don't-eat-me" signals (e.g., CD47). Pro-phagocytic mAbs hold great promise as therapeutics.
  - Here, we have developed and validated *in vitro* assays for quantification of phagocytosis in 96-/384-well plates using the Incucyte® Live-Cell Analysis System and/or iQue® Advanced Flow Cytometer.
- Advanced Flow Cytometer.
- The Incucyte® Phagocytosis Assay combines pHrodo® for Incucyte® reagents and integrated image-based fluorescent measurements in a simple mix-and-read protocol.
  - The iQue® Human ADCP Kit measures co-localization between live target cells and CD14+ effector cells to provide a readout for ADCP response.
  - These data exemplify that live-cell analysis, alongside advanced flow cytometry, is a powerful tool for quantitative morphological and functional assessment of phagocytosis, which is amenable to screening for therapeutic agents.

### Visualization & quantification of phagocytosis in real-time



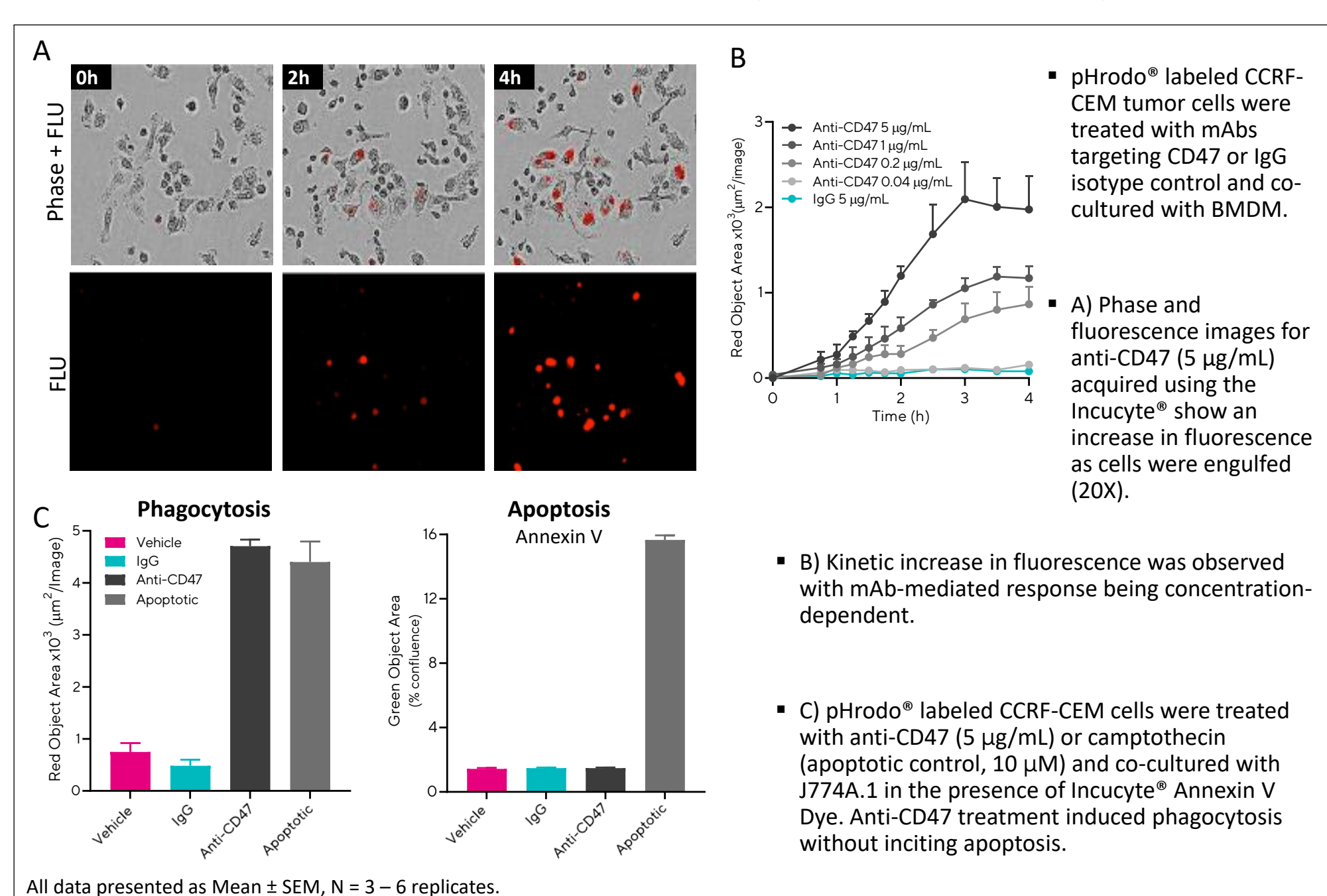
- A) pHrodo® labeled apoptotic Jurkats were co-cultured with J774A.1 macrophages. Phase and fluorescent images (20X) were acquired using the Incucyte® Live-Cell Analysis System and revealed an increase in red fluorescence following engulfment over time quantified using a fluorescent segmentation mask (blue).
- B) iPSC microglia were seeded, rested or polarized, then exposed to increasing densities of pHrodo® Bioparticles® and pHrodo® labeled apoptotic cells (Rows A - H) or increasing concentrations of Cytochalasin D and a single density of Bioparticles® or apoptotic cells (Rows I - P). Microplate view shows change in fluorescence area over 24h for all wells.

### Phagocytosis is cell type- & Bioparticle® density-dependent



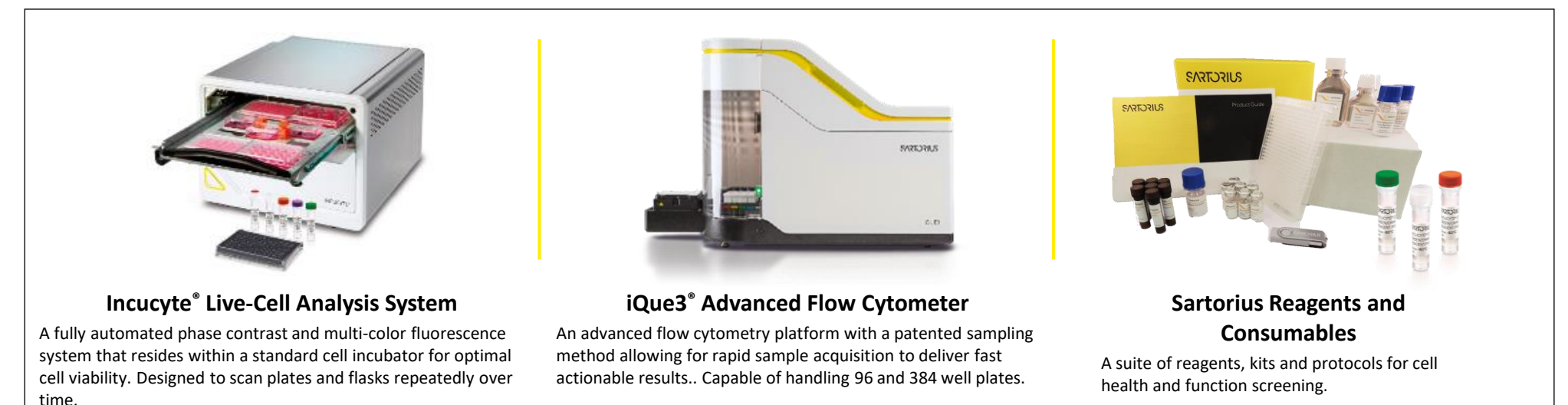
- Various phagocytes were treated with pHrodo® Green E. coli Bioparticles® and engulfment was monitored using the Incucyte®.
- A) Time-course shows rapid, cell-type dependent Bioparticle® uptake with clearance within 4h (BMDM or Microglia) or 12h (macrophages).
- B) Density range of macrophages were treated with 3 densities of Bioparticles®. Fluorescence area at 6h was compared and showed both effector cell and bioparticle density-dependent engulfment.
- C) Macrophages were treated with cytoskeletal inhibitors prior to Bioparticle® addition. Cytochalasin D (actin inhibitor), Lantrunculin A (actin stabilizer), and Go6976 (PKC inhibitor) showed concentration-dependent inhibition of phagocytosis ( $\text{IC}_{50}$  values of 0.38, 95% CI [0.30, 1.23], 0.06 95% CI [0.04, 0.08], and 3.93 95% CI [2.05, 5.80], respectively).

### Anti-CD47 mAb promotes macrophage-mediated phagocytosis

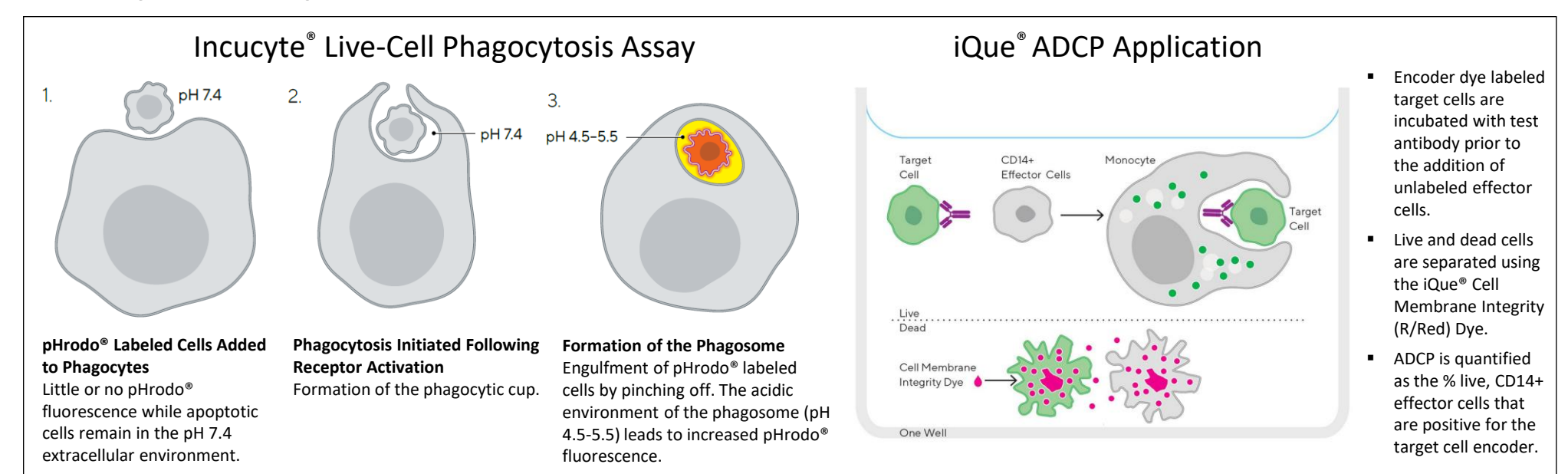


All data presented as Mean  $\pm$  SEM, N = 3 - 6 replicates.

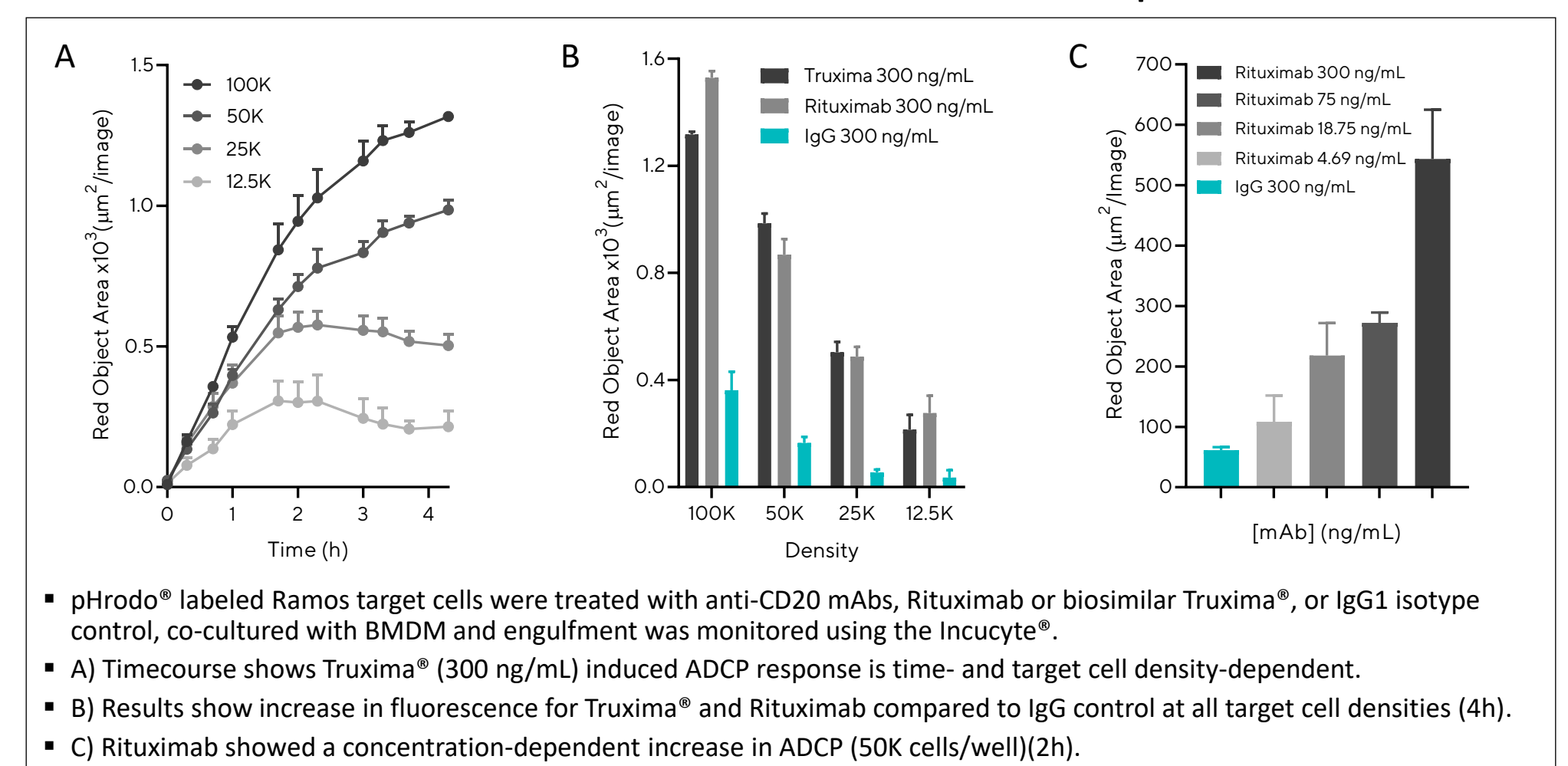
### Incucyte® & iQue3® Systems



### Assay Principles

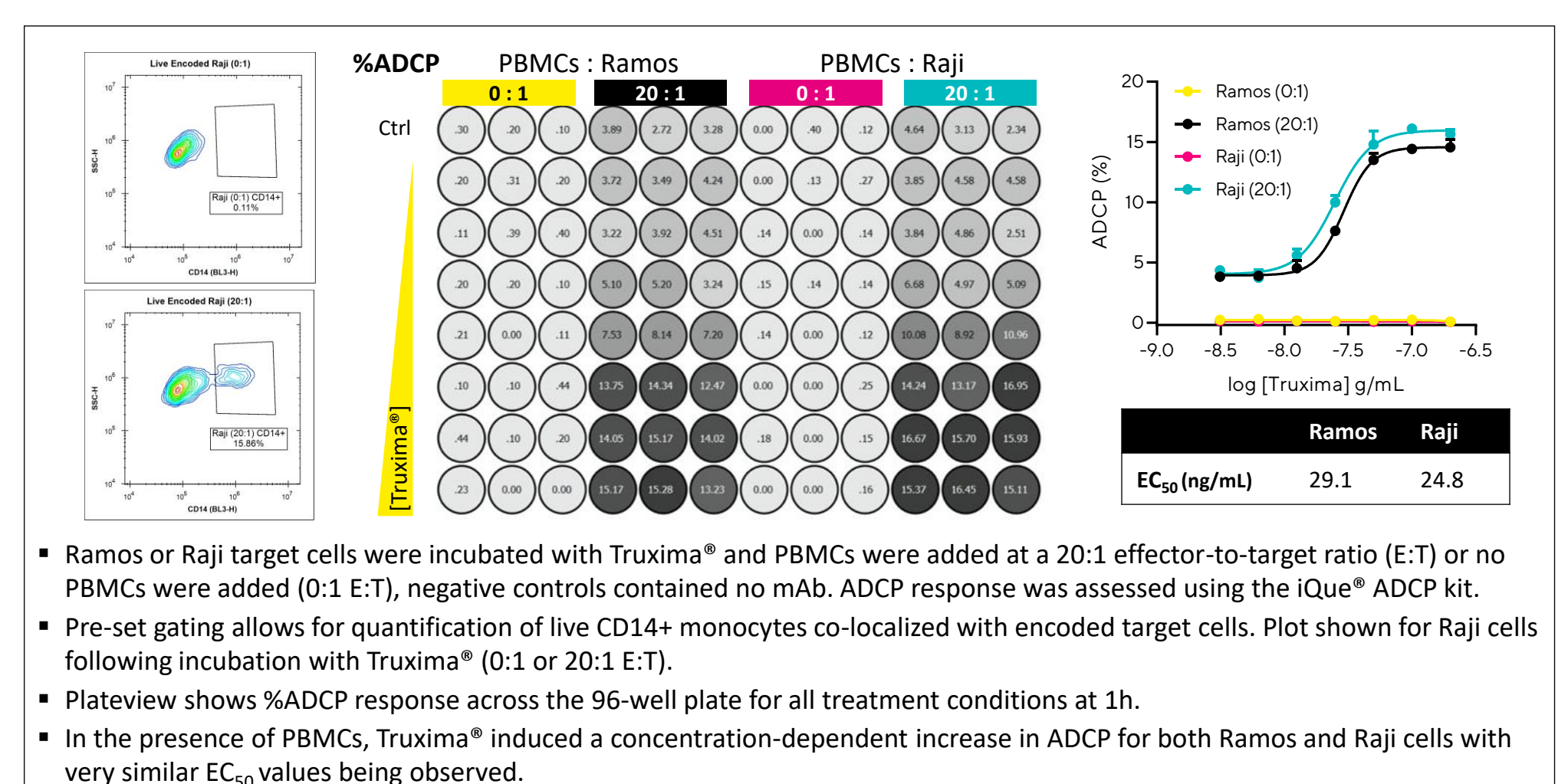


### Clinical anti-CD20 mAbs Truxima® & Rituximab promote ADCP



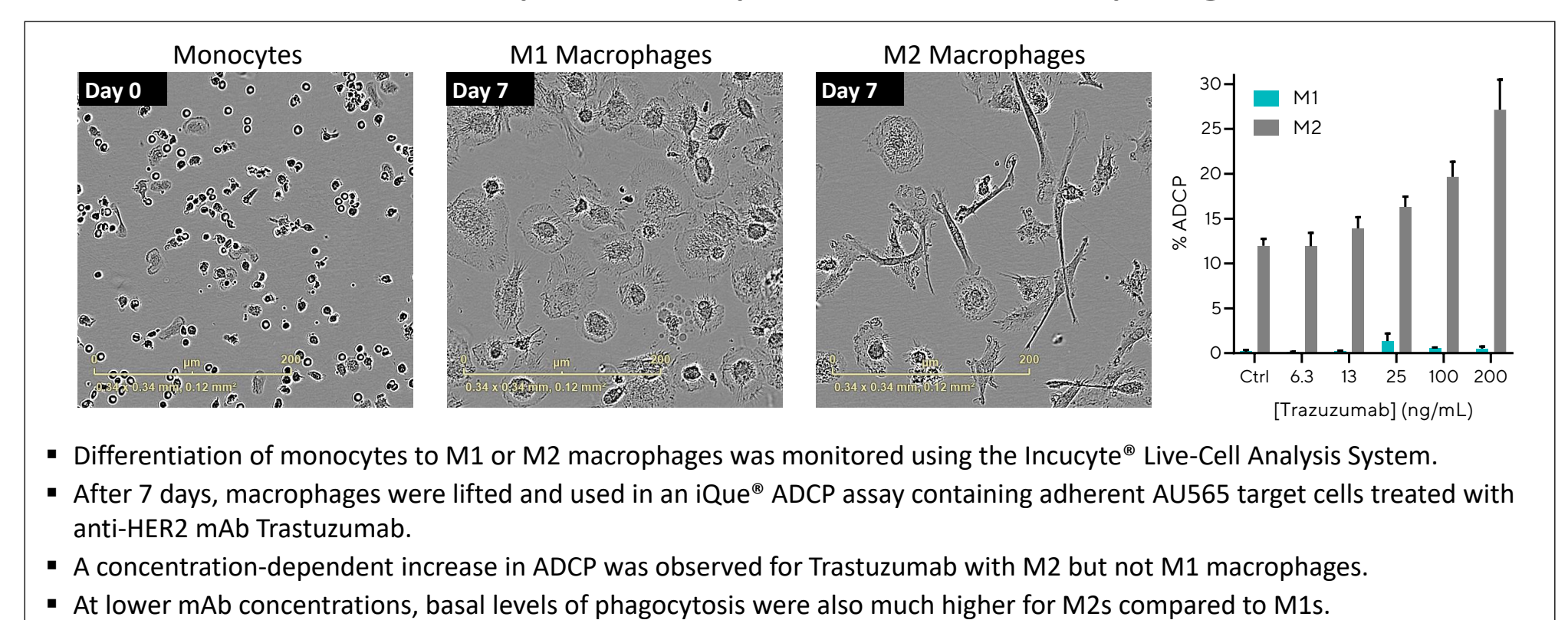
- pHrodo® labeled Ramos target cells were treated with anti-CD20 mAbs, Rituximab or biosimilar Truxima®, or IgG1 isotype control, co-cultured with BMDM and engulfment was monitored using the Incucyte®.
- A) Timecourse shows Truxima® (300 ng/mL) induced ADCP response is time- and target cell density-dependent.
- B) Results show increase in fluorescence for Truxima® and Rituximab compared to IgG control at all target cell densities (4h).
- C) Rituximab showed a concentration-dependent increase in ADCP (50K cells/well)(2h).

### ADCP increases with Truxima® concentration in B cells



- Ramos or Raji target cells were incubated with Truxima® and PBMCs were added at a 20:1 effector-to-target ratio (E:T) or no PBMCs were added (0:1 E:T), negative controls contained no mAb. ADCP response was assessed using the iQue® ADCP kit.
- Pre-set gating allows for quantification of live CD14+ monocytes co-localized with encoded target cells. Plot shown for Raji cells following incubation with Truxima® (0:1 or 20:1 E:T).
- Plateau shows %ADCP response across the 96-well plate for all treatment conditions at 1h.
- In the presence of PBMCs, Truxima® induced a concentration-dependent increase in ADCP for both Ramos and Raji cells with very similar EC<sub>50</sub> values being observed.

### Differential ADCP response of polarized macrophages



- Differentiation of monocytes to M1 or M2 macrophages was monitored using the Incucyte® Live-Cell Analysis System.
- After 7 days, macrophages were lifted and used in an iQue® ADCP assay containing adherent AU565 target cells treated with anti-HER2 mAb Trastuzumab.
- A concentration-dependent increase in ADCP was observed for Trastuzumab with M2 but not M1 macrophages.
- At lower mAb concentrations, basal levels of phagocytosis were also much higher for M2s compared to M1s.