

High-Throughput Quantification of Antibody-Dependent Phagocytosis using Live-Cell Analysis

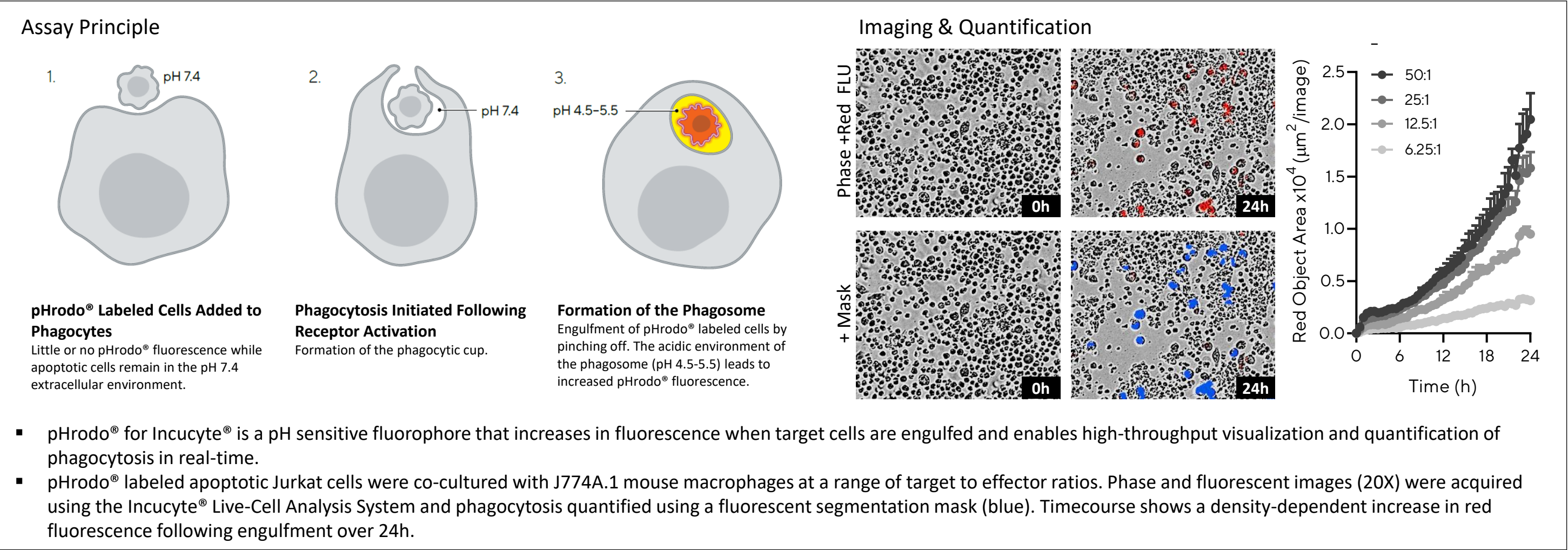
J. Trigg*, G. Lovell, N. Bevan, and T. Dale

Sartorius Corporation, Royston, Hertfordshire, UK
*Corresponding author: Jasmine.Trigg@sartorius.com

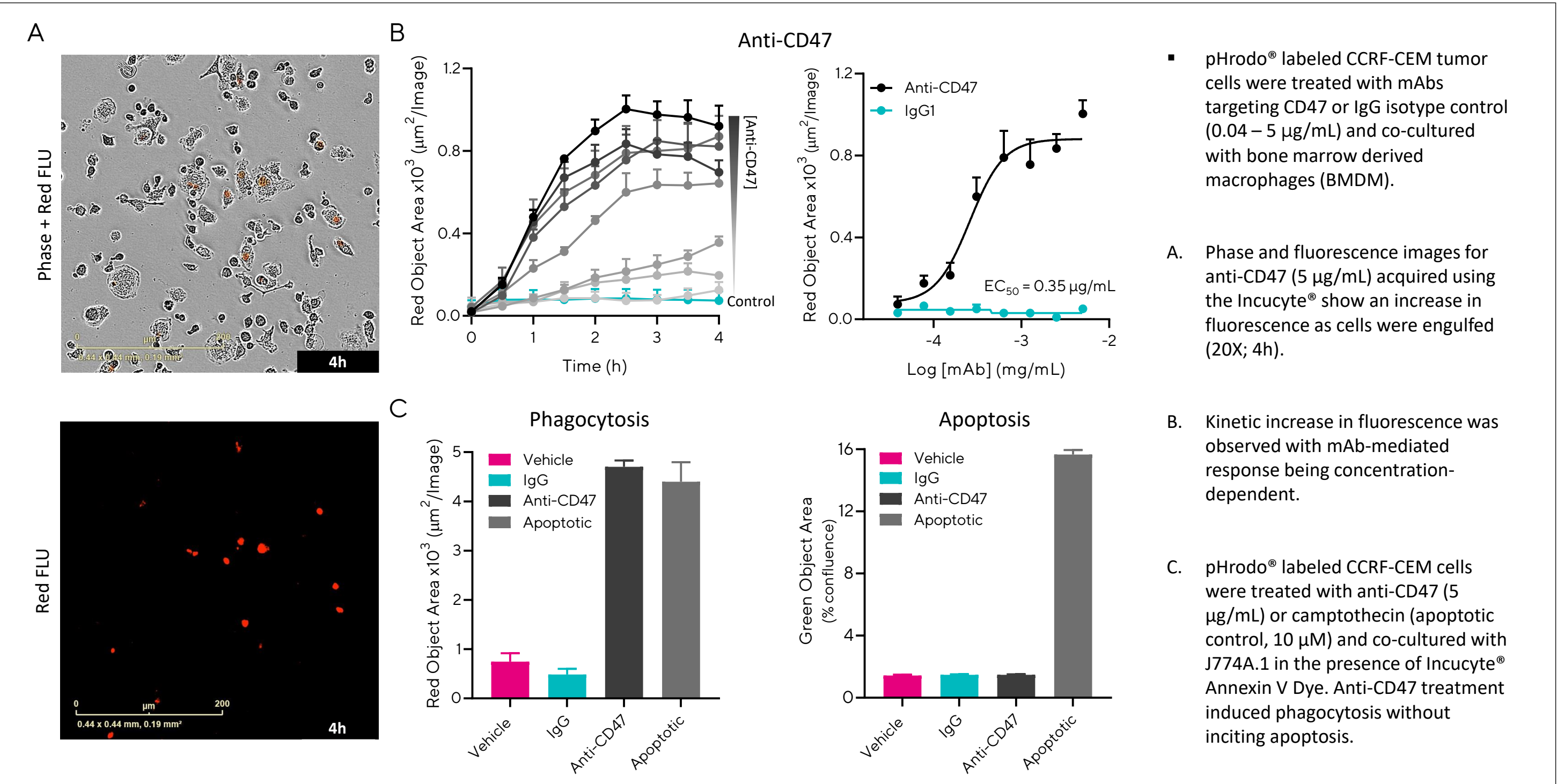
Introduction

- Phagocytosis, a specific form of endocytosis, is a critical component of innate and adaptive immune responses.
- The uptake and clearance of viable tumor cells can be promoted with monoclonal antibodies (mAbs) via antibody-dependent phagocytosis (ADCP) or through the blockade of “don’t-eat-me” signals, such as CD47. These mechanisms hold immunotherapeutic promise and are studied extensively in drug development.
- Here, we have developed and validated an *in vitro* assay for the high-throughput quantification of phagocytosis using the Incucyte® Live-Cell Analysis System.
- The Incucyte® Phagocytosis Assay combines pHrodo® for Incucyte® reagents and integrated image-based fluorescent measurements in a simple mix-and-read protocol.
- These data exemplify that live-cell analysis is a powerful tool for quantitative morphological and functional assessment of ADCP, which is amenable to screening for therapeutic candidates.

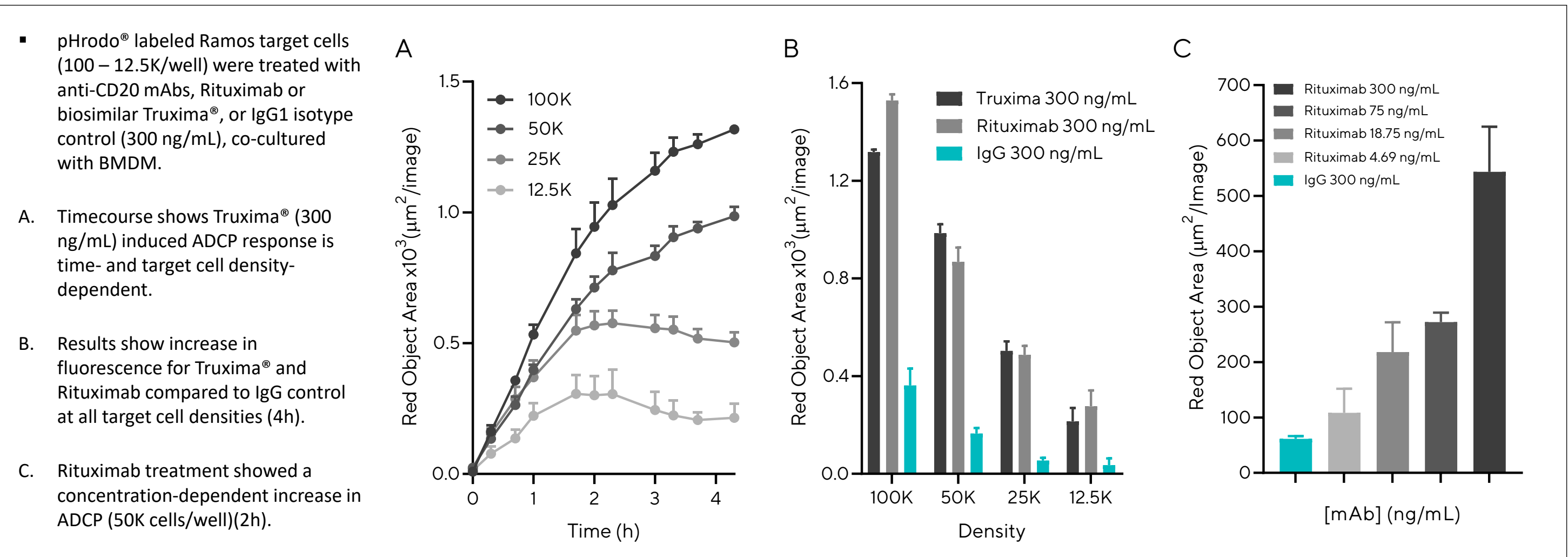
Visualize and quantify phagocytosis in real-time



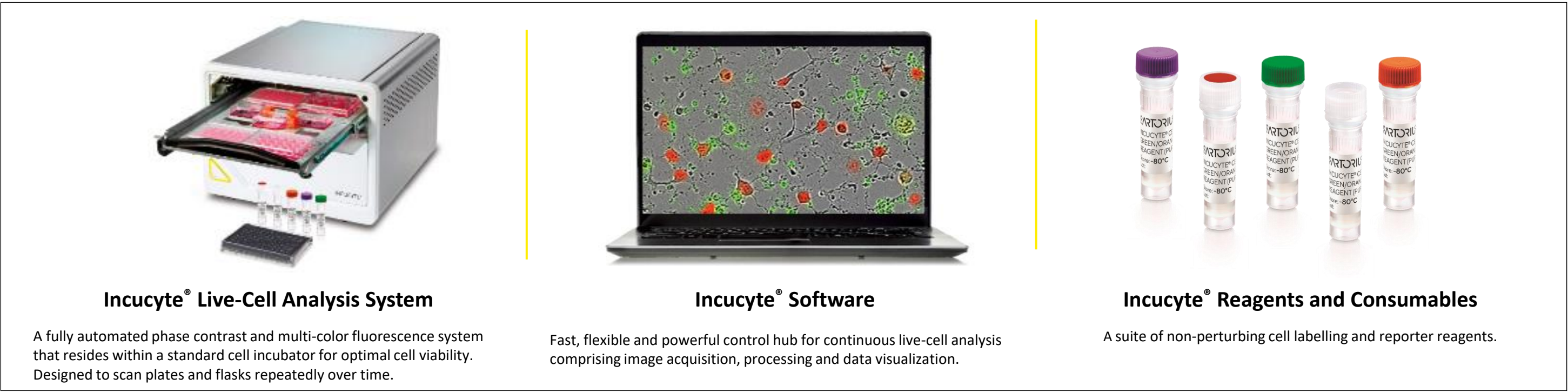
Anti-CD47 mAb stimulates macrophage-mediated phagocytosis



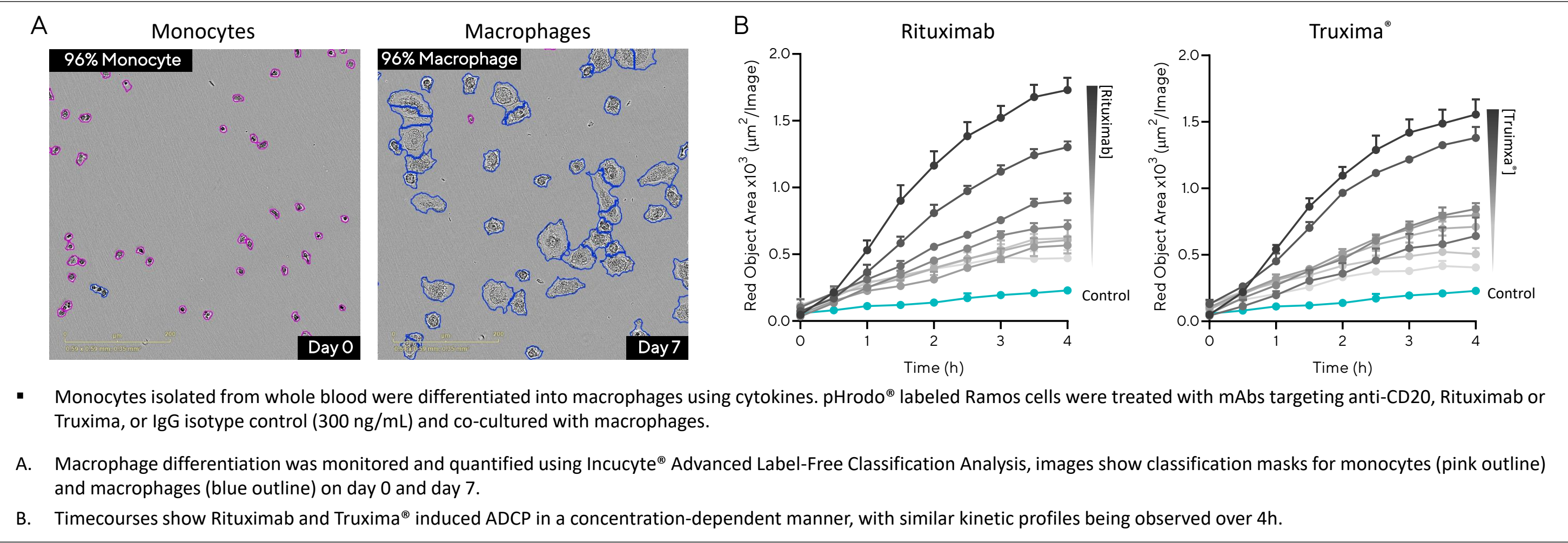
Clinical anti-CD20 mAbs Truxima® and Rituximab promote ADCP



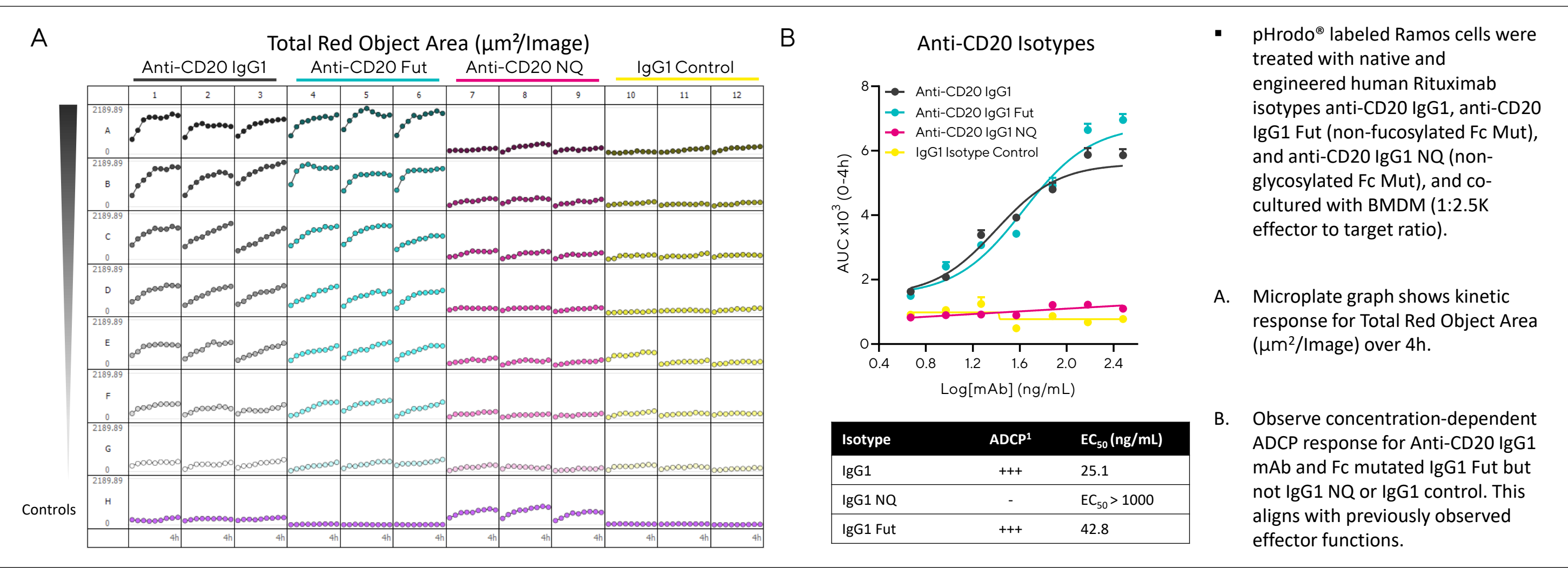
Incucyte® Live-Cell Imaging and Analysis Solutions



ADCP increases with anti-CD20 concentration in primary macrophages



High-throughput assessment of native and engineered Rituximab Fc Mutants



Target HER2 antigen expression correlates with ADCP response

