

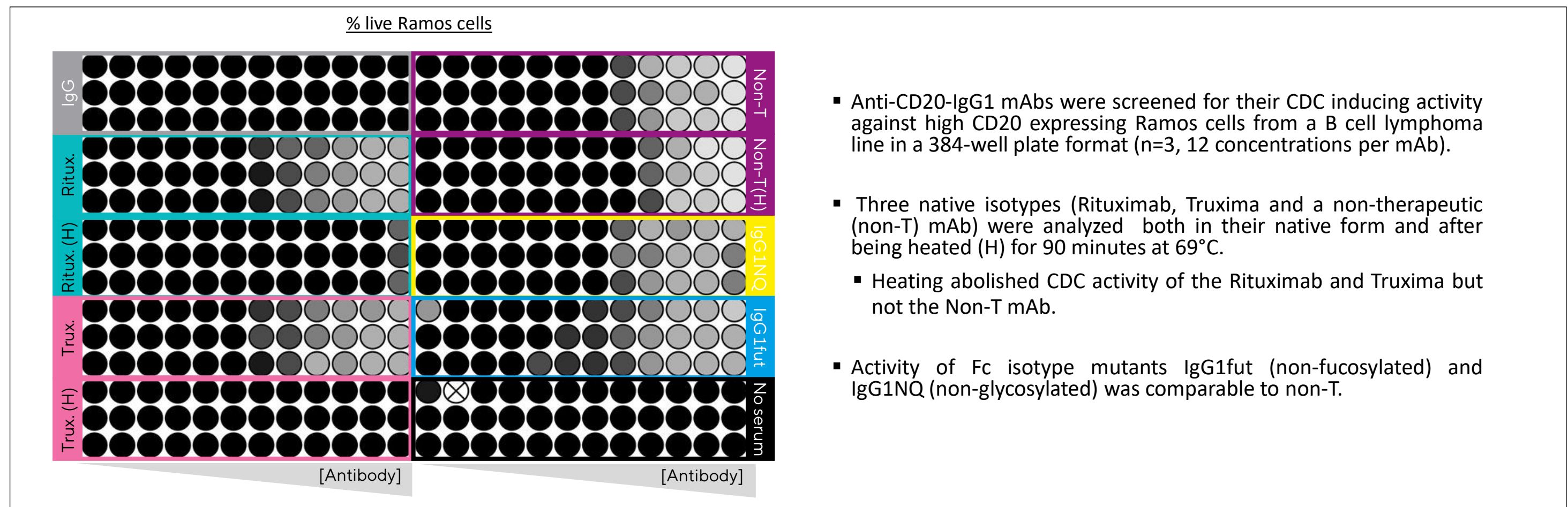
Quantifying *in vitro* complement-dependent cytotoxicity (CDC) using advanced flow cytometry

K. McBain<sup>1\*</sup>, J. Lovchik<sup>2</sup> & N. Bevan<sup>1</sup>  
Sartorius Corporation, <sup>1</sup>Royston, UK; <sup>2</sup>Albuquerque, NM  
\*Corresponding author: [Kirsty.McBain@sartorius.com](mailto:Kirsty.McBain@sartorius.com)

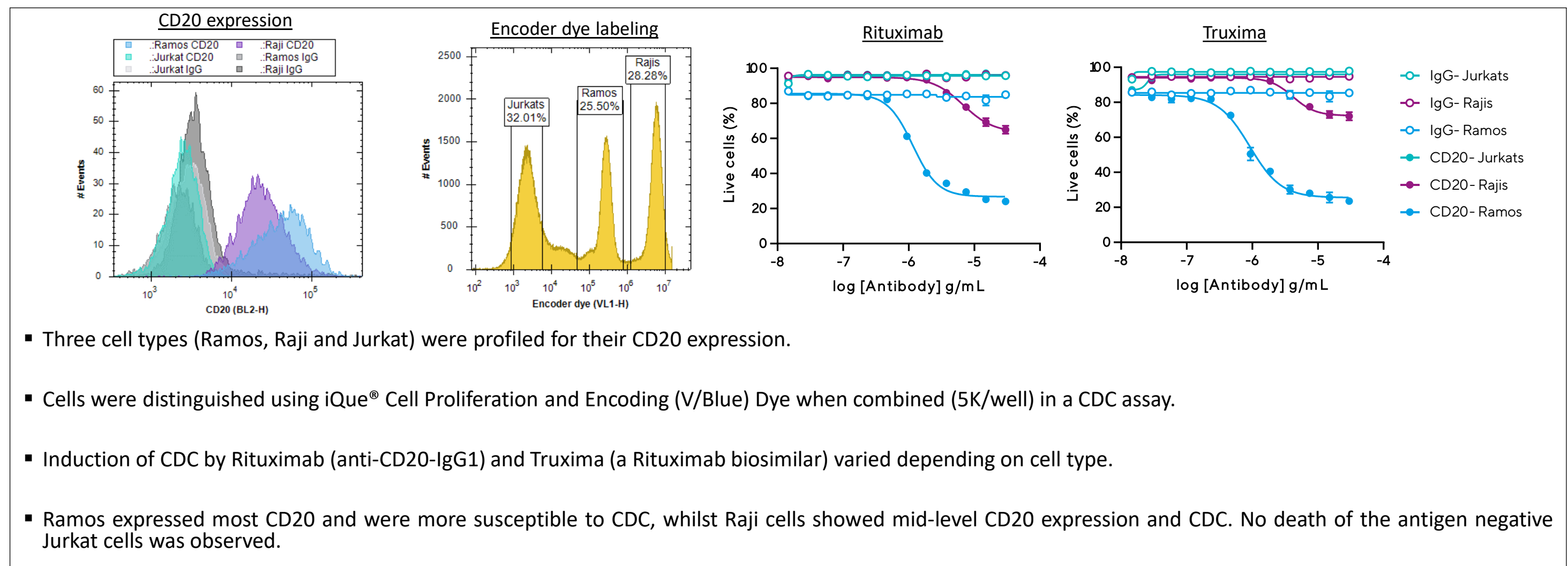
Introduction

- In recent decades, rapid progression of monoclonal antibody (mAb) development has culminated in approvals for over 100 therapeutics in the United States<sup>1</sup>.
- mAbs enhance the immune response through three key Fc-mediated functions: antibody-dependent cellular cytotoxicity (ADCC), antibody-dependent cellular phagocytosis (ADCP) and complement-dependent cytotoxicity (CDC).
- CDC is initiated when a mAb triggers activation of the complement cascade, which ultimately results in tumor cell lysis.
- We present a streamlined assay for measuring CDC using the iQue® Advanced Flow Cytometry Platform.
- Target cells are cultured with mAbs of interest and human serum, then labeled using iQue® Cell Membrane Integrity (R/Red) Dye to quantify cell death.
- Combining the high-throughput of the iQue® with rapid data analysis using the integrated iQue Forecyt® software streamlines antibody characterization and drug discovery processes.

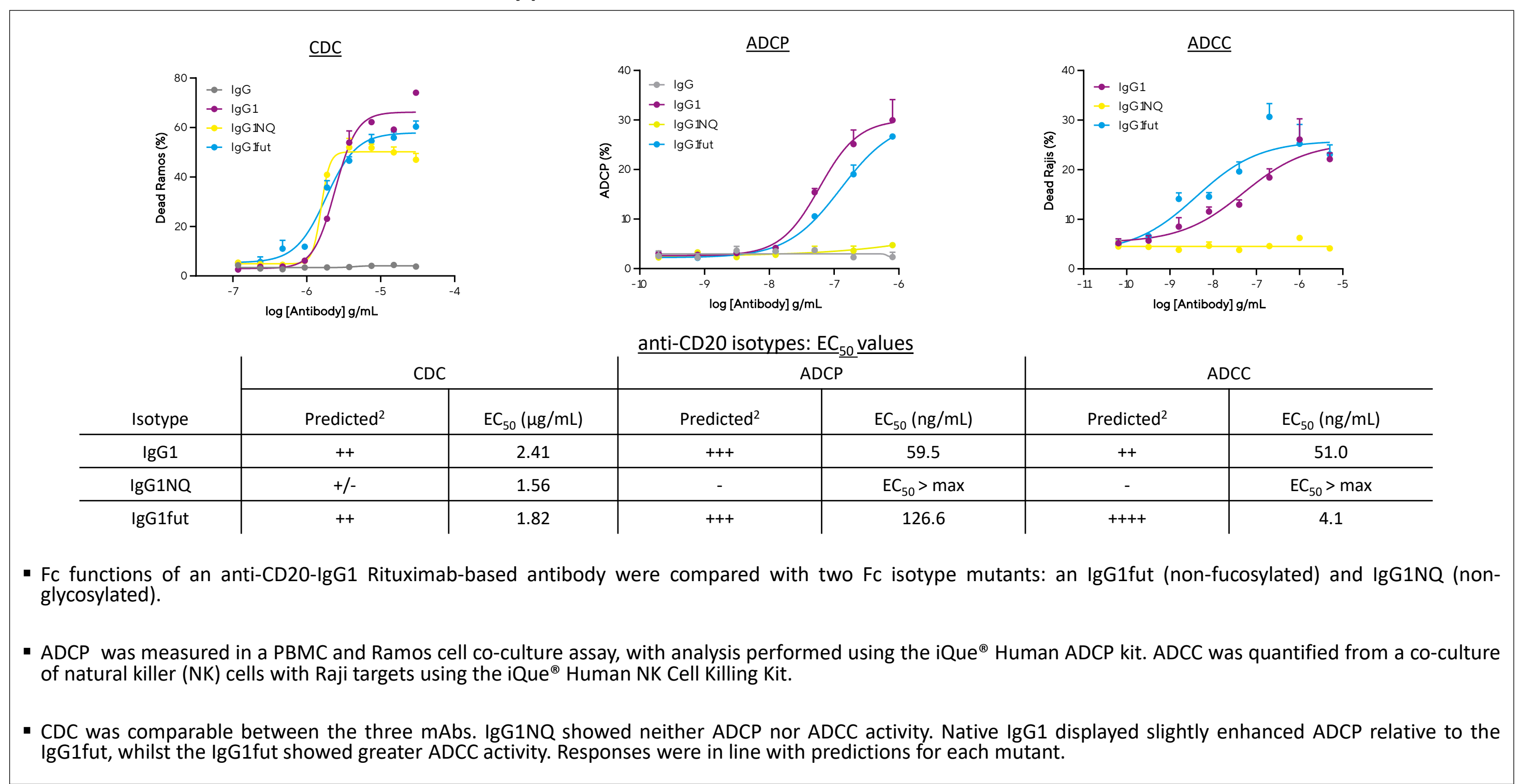
High-throughput comparison of anti-CD20 CDC activity



Induction of CDC correlates with target antigen expression



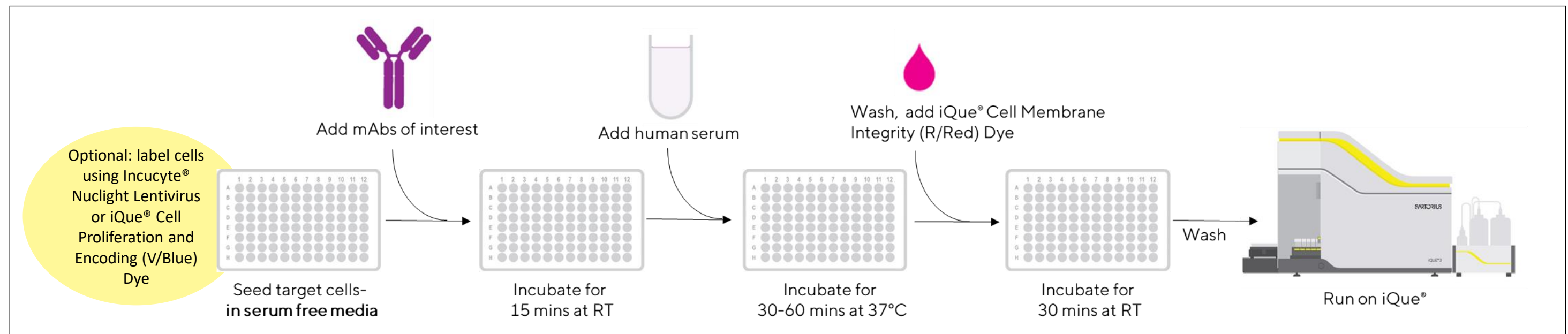
Fc function differs between isotype mutants



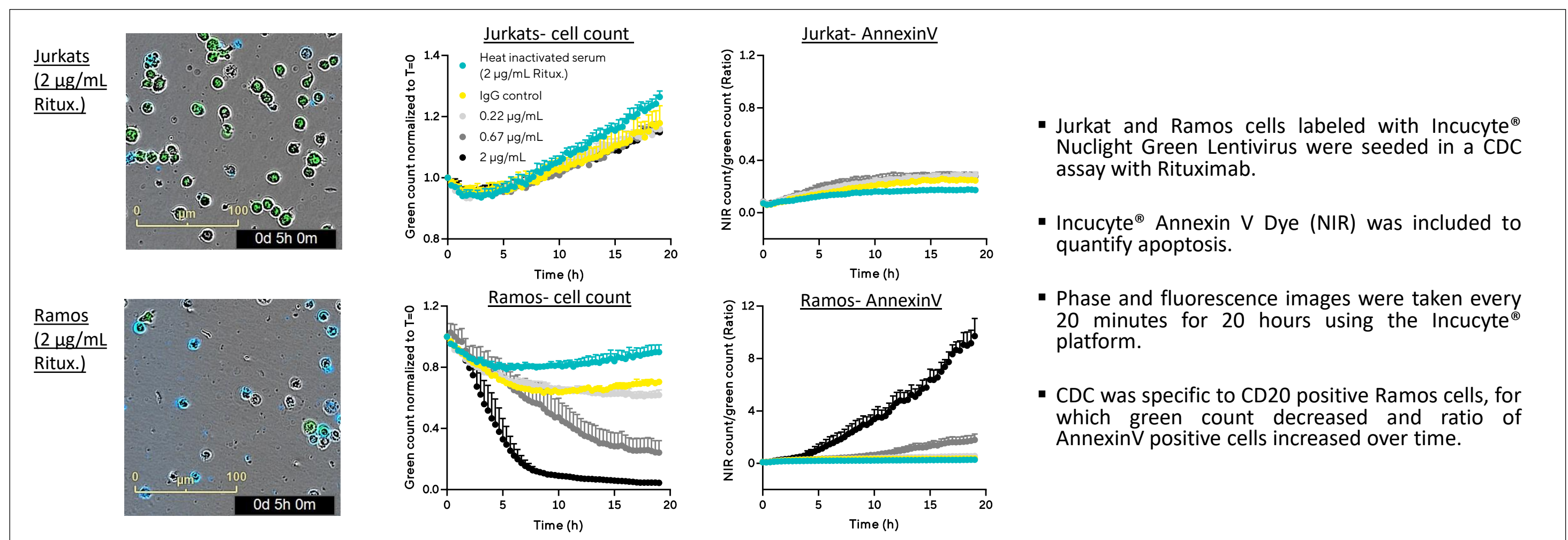
iQue® System



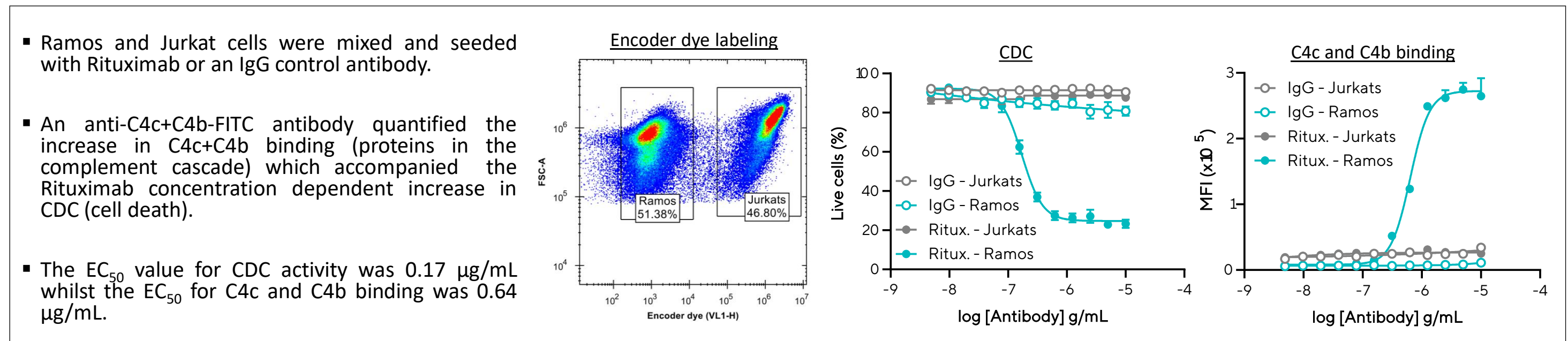
Assay Workflow



Incucyte® images quantified CDC over time



Complement protein binding during CDC



Adherent cells are resistant to CDC

