

A human FcγRIIa (H131 variant)-specific reporter assay for evaluating antibody activity in antibody-dependent phagocytosis (ADCP) pathway activation

Teresa Surowy, Rich Moravec, Denise Garvin, Aileen Paguio and Frank Fan

R&D, Promega Corporation, 2800 Woods Hollow Road, Madison, WI 53711, USA. Email: teresa.surowy@promega.com



1. Introduction

Antibody-dependent cellular phagocytosis (ADCP) is an important mechanism of action of therapeutic antibodies that is effective at killing tumor cells in addition to antibody-dependent cell-mediated cytotoxicity (ADCC). Antibody drug development is beginning to address and realize the potential of this mechanism of action more attentively. Additionally, viral vaccine development is realizing the potential of an ADCP response for vaccine effectiveness.

However, even more so than classic ADCC bioassay, classic ADCP assays are tedious and highly variable. They are FACS-based rather than plate-based, and rely on isolation of primary cells, typically monocytes, that then need to be differentiated into phagocytic macrophages in culture for many days, and the need for multiple labeling reactions of effector and target cells. We sought to develop a plate-based, less tedious, faster and less variable ADCP bioassay. With this, understanding, manipulating and quantifying ADCP in both therapeutic antibody and vaccine development could be improved.

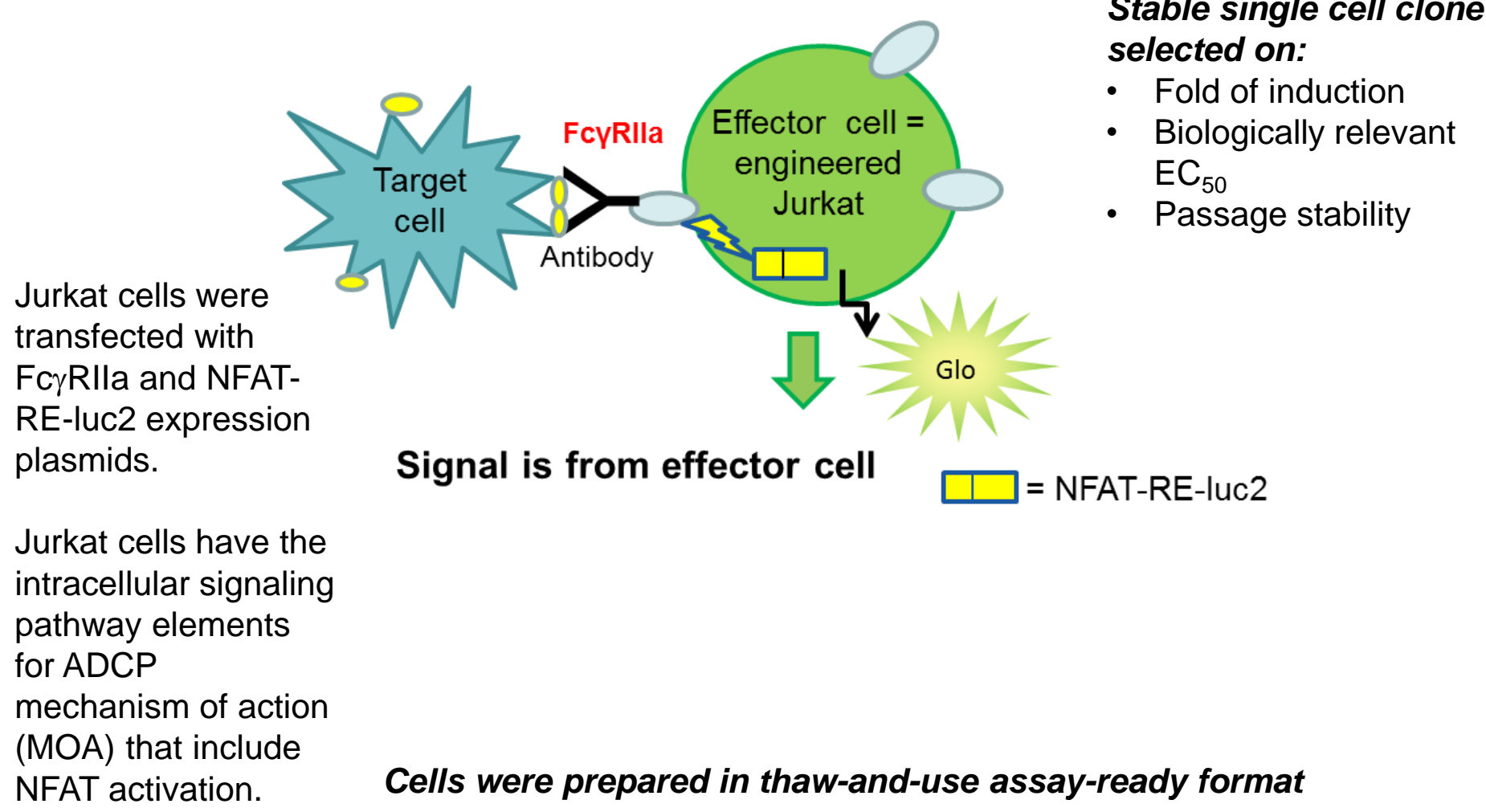
Macrophage effector cells and other myeloid-derived phagocytes mediating ADCP are more dependent on Fcγ activating receptors beyond the NK-cell expressing and ADCC-predominant FcγRIIIa. The FcγRIIa receptor has been reported to have the dominant role in ADCP by macrophages.

We describe here our development and evaluation of a human FcγRIIa (H131 variant)-specific reporter bioassay for quantifying antibody bioactivity in ADCP pathway activation. The H131 variant is able to bind all human IgG isotypes. The assay uses dual-engineered Jurkat effector cells that express the receptor and mediate FcγRIIa-dependent NFAT activation of reporter luciferase expression only when receptors are bound by target cell-bound antibodies.

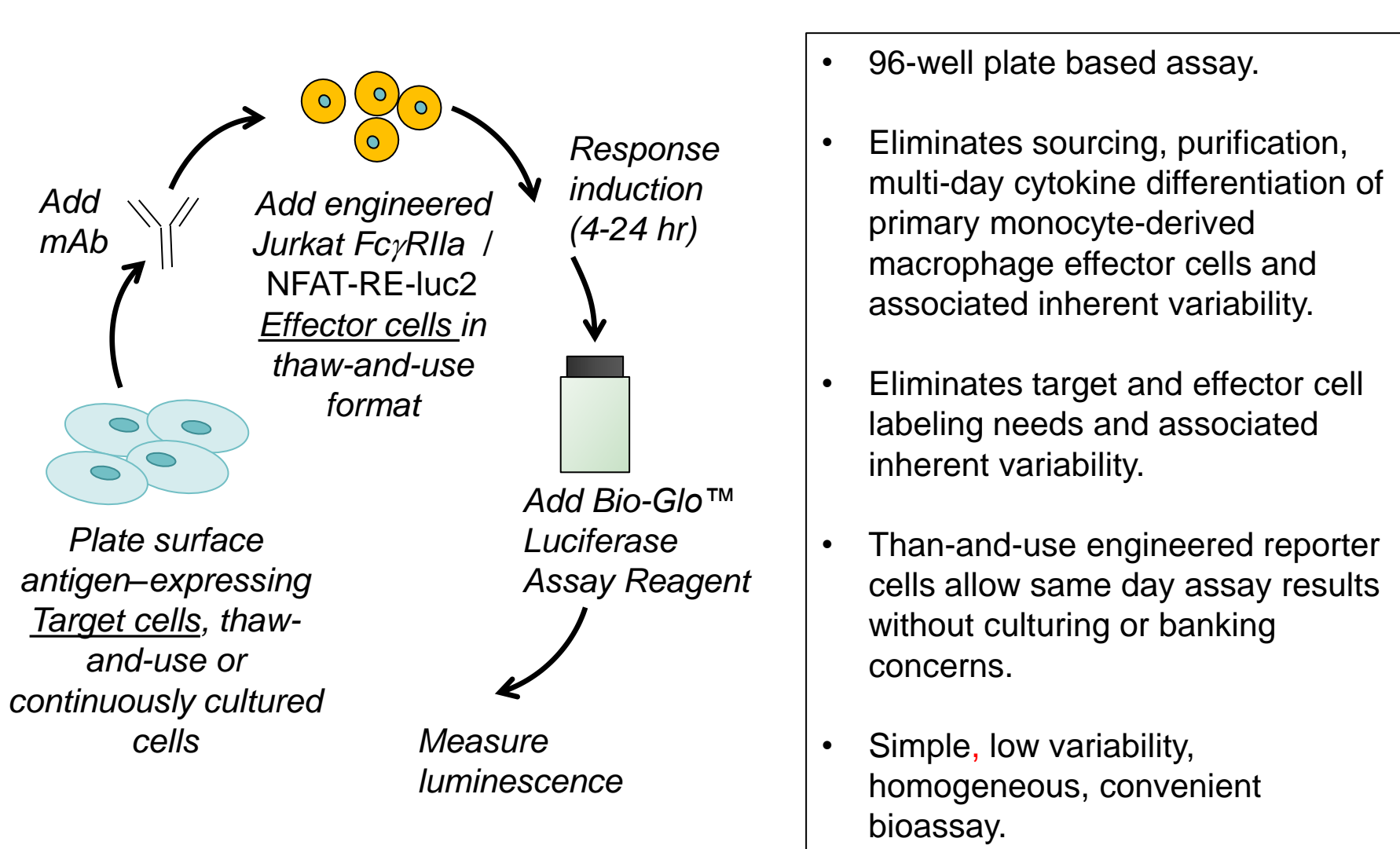
We confirmed surface expression of the FcγRIIa (H131 variant) in our effector cells via FACS and passage stability of the cell line response for cell bank development. We developed and optimized the ADCP reporter bioassay using the effector cells in thaw-and-use format (immediate use, no culture), for ease of use and assay consistency. Rituximab, a chimeric IgG1 antibody drug with known ADCP activity through FcγRIIa, was used in much of the bioassay development. Using a series of rituximab isotype variants, we demonstrate that all four human IgG isotypes are active in the assay with expected relative activities. The assay is also specific, induction being dependent on the presence of FcγRIIa (H131 variant) and target-cell bound antibody. We further demonstrate that the assay can quantify the biological activity of other antibodies such as anti-EGFR panitumumab (an IgG2), anti-Her2 trastuzumab and an anti-GD2 antibody via this receptor. In a bioassay qualification study quantifying relative potency, the assay was accurate, precise and demonstrated good linearity and range. We also demonstrated the assay was stability-indicating using heat stressed rituximab and that it was robust.

2. Generation of engineered FcγRIIa/NFAT-RE-luc2 effector cells

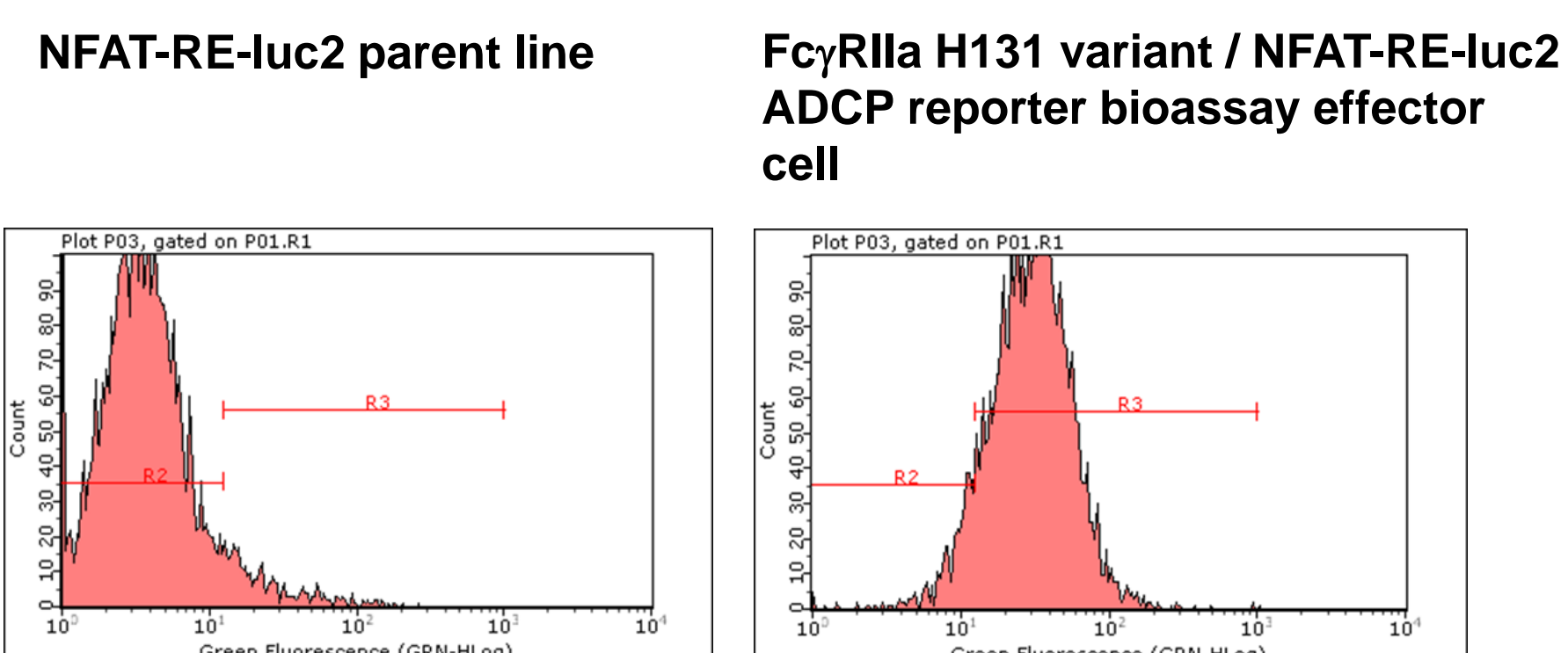
Reporter-based FcγRIIa ADCP bioassay



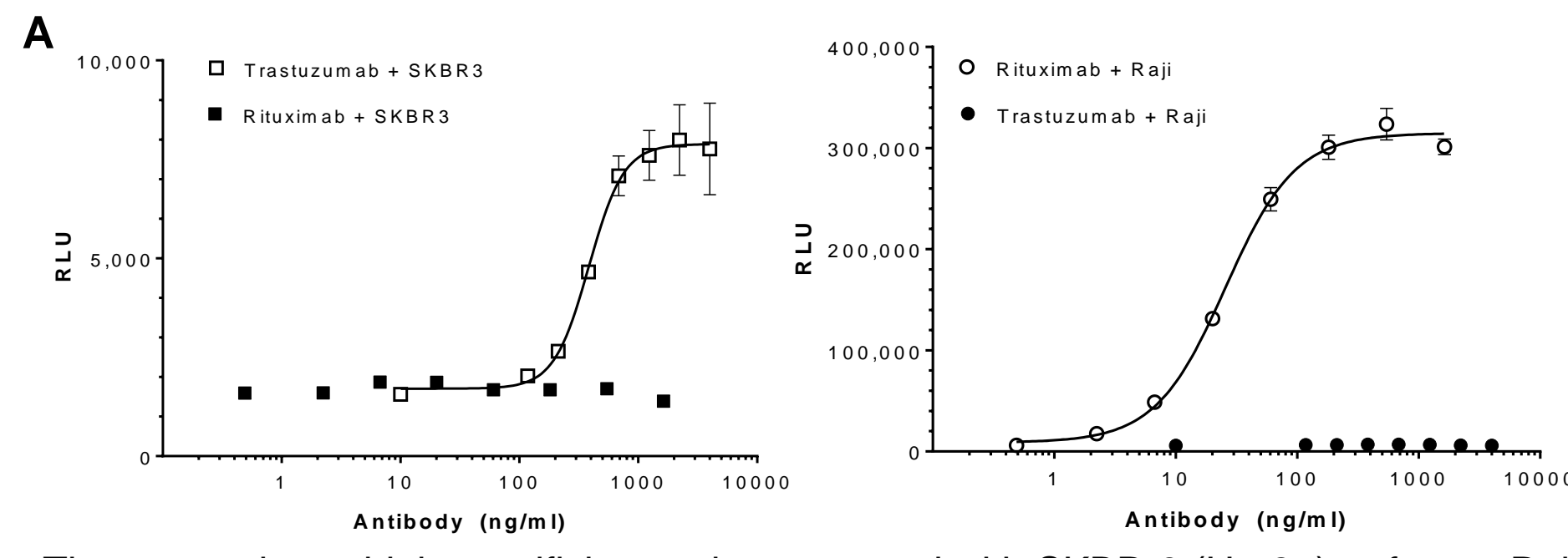
3. Protocol and advantages of the bioassay



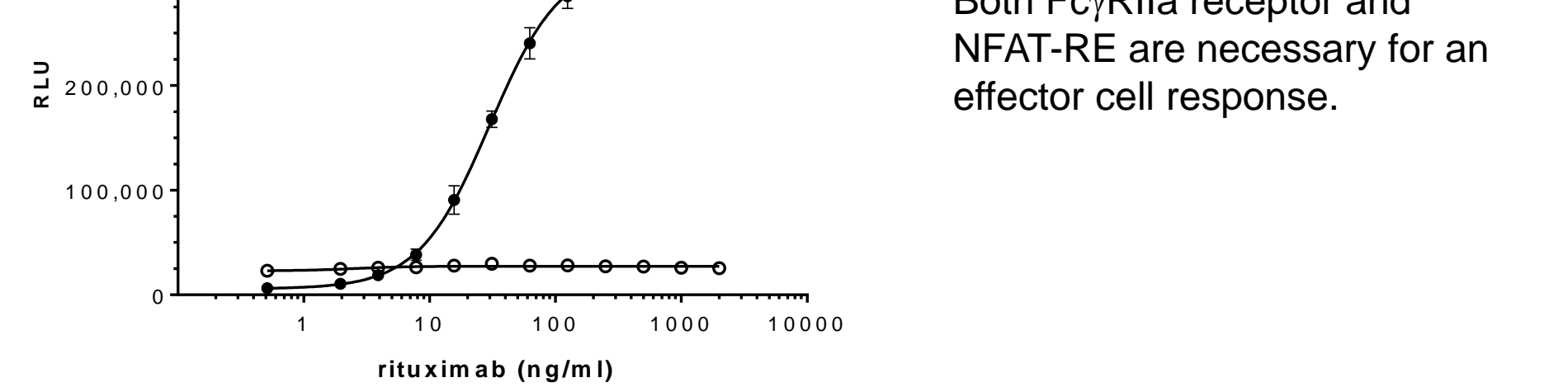
4. Confirmation of cell surface FcγRIIa expression by FACS analysis



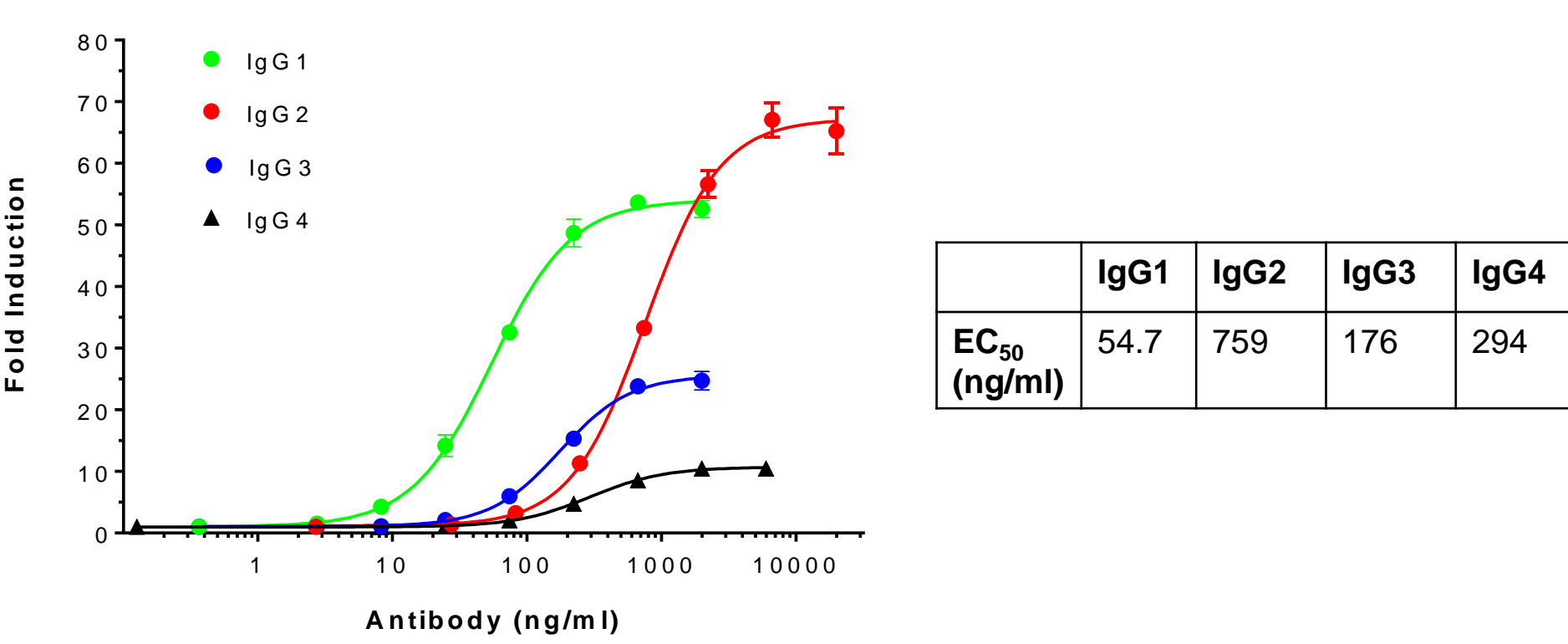
5. The bioassay is specific



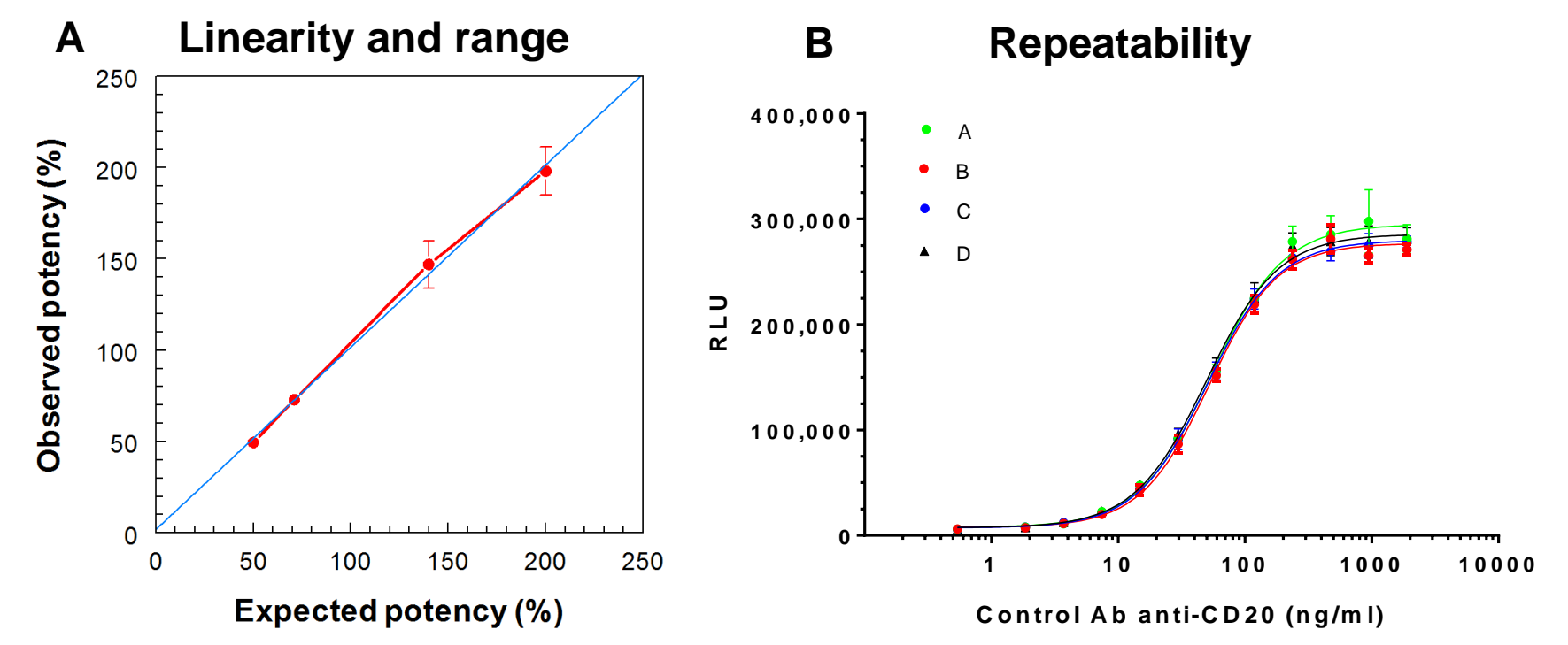
6. FcγRIIa (H131) bioassay response shows appropriate IgG isotype bioactivities



7. Bioassay can quantify bioactivity in ADCP pathway activation of a variety of drug and research antibodies



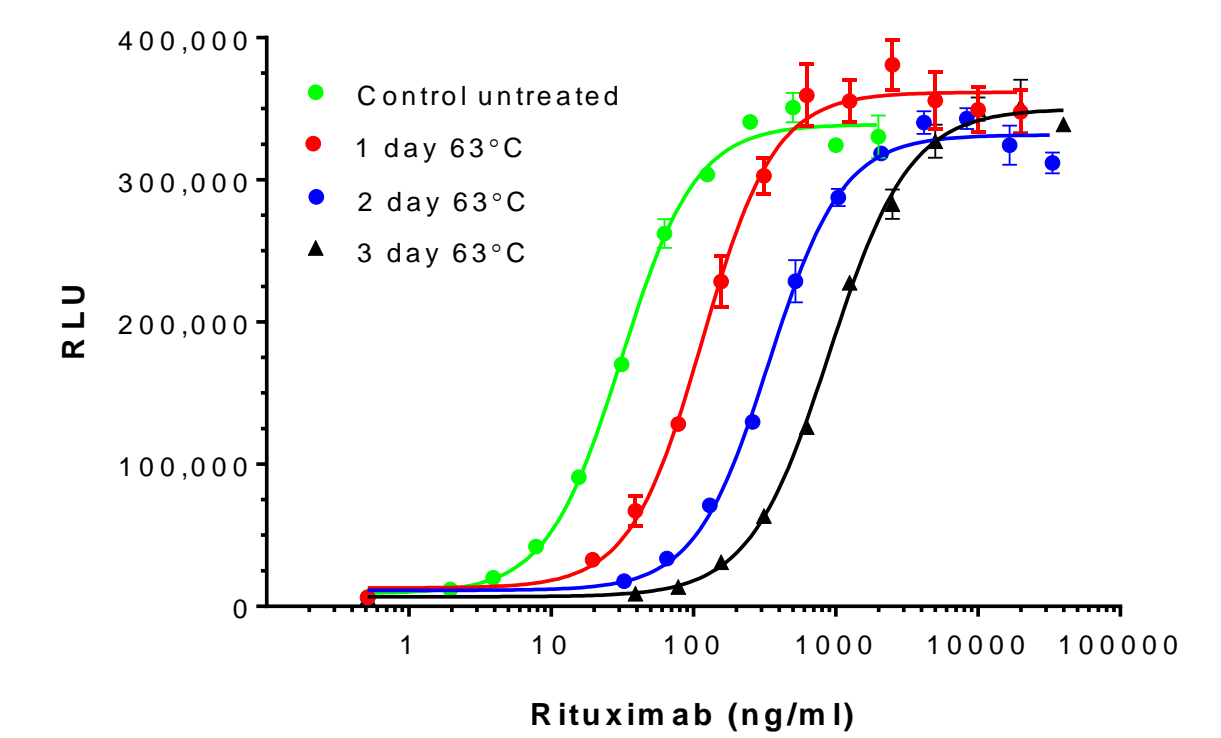
8. Potency qualification using ADCP Control Ab, Anti-CD20



Potency qualification (3-day) with ADCP Control Ab, anti-CD20 (IgG1) using frozen thaw-and-use FcγRIIa (H131) effector and Raji target cells. Assay was performed on three days using 35K effector cells + 15K Raji target cells, with data analyzed using JMP 11 software.

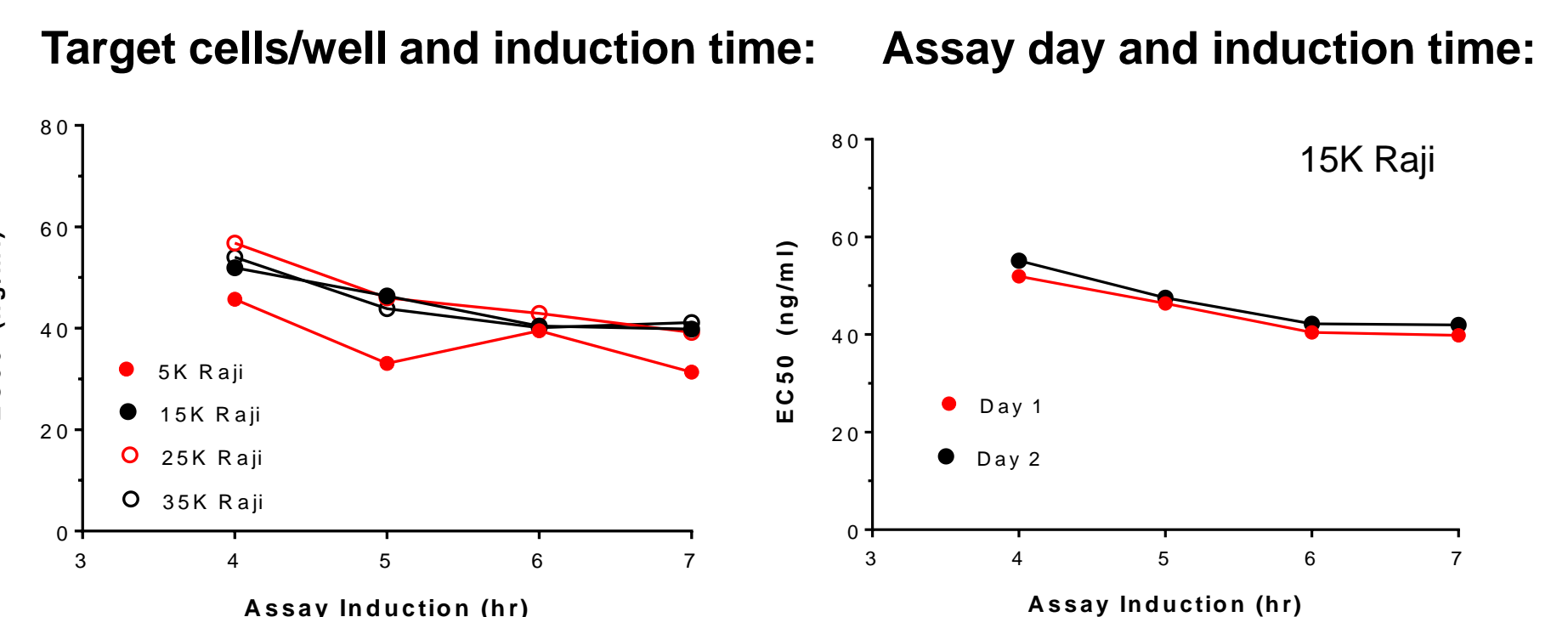
A. Linearity and range. **B.** Example within-one-day repeatability (100% reference sample EC₅₀ values on 4 plates: 43.6, 46.4, 46.7, and 47.1 ng/ml).

9. Bioassay demonstrates stability indicating properties using heat stressed rituximab



Rituximab was heat stressed at 63°C for 1-3 days and compared to control unheated sample. Results obtained following a 6 hour induction using frozen thaw-and-use FcγRIIa (H131) effector and Raji target cells.

10. ADCP reporter bioassay is robust



Summary

- We have developed an FcγRIIa-specific ADCP reporter assay that quantifies ADCP MOA pathway activation for the predominant Fcγ receptor mediating macrophage ADCP.
- The luminescent 96-well plate assay is simple, homogeneous, convenient, and fast. It should provide a much less variable and tedious method of assessing ADCP pathway MOA than FACS analysis using monocyte-derived macrophages differentiated in culture.
- It is able to quantify ADCP pathway activation for a variety of research and drug antibodies and demonstrates appropriate human IgG isotype relative activities.
- The bioassay is specific, robust and stability-indicating. In a qualification study it demonstrated excellent linearity, recovery and intra- and intermediate assay precision, indicating potential application for plate-based lot release testing.