

Measurement of Fc-mediated ADCC and CDC of anti-TNF α and anti-VEGF Therapeutic Antibodies using Reporter-based Bioassays and Engineered TNF α ⁺ and VEGF⁺ Target Cells

Denise Garvin, Rich Moravec, Jamison Grailer, Jim Hartnett, Chris Heid, Frank Fan, Mei Cong and Zhi-jie Jey Cheng
Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711



1. Introduction

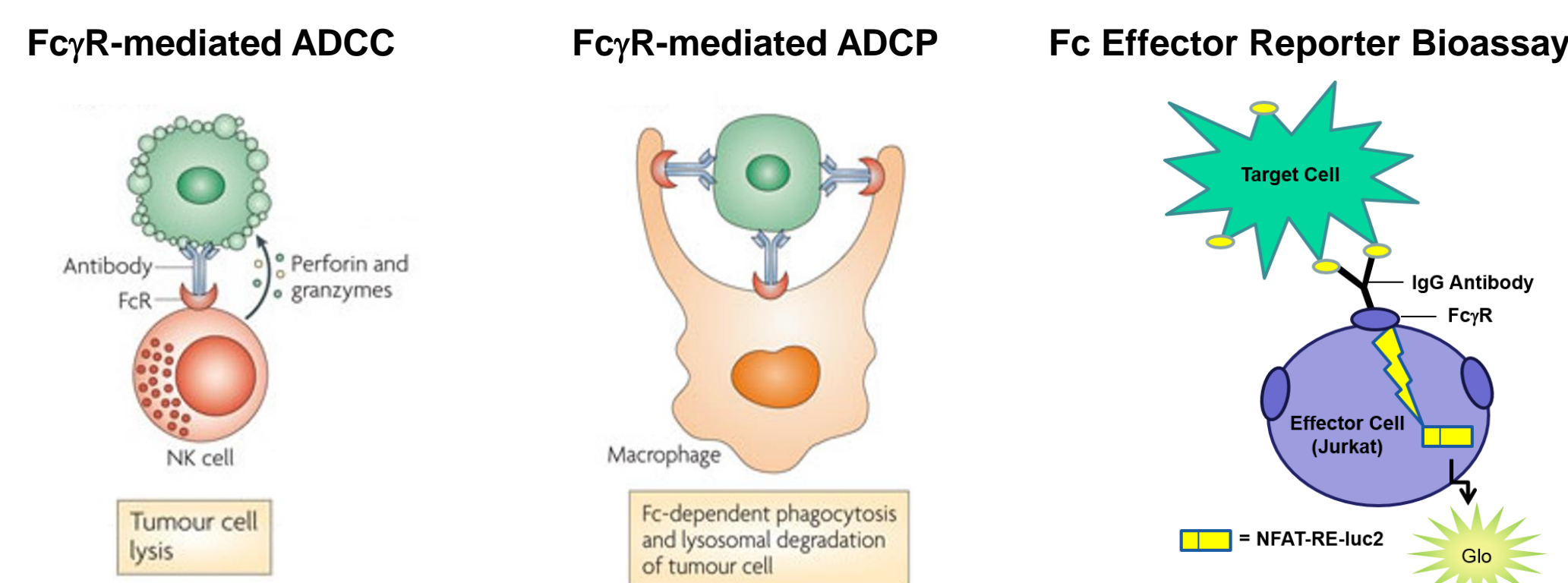
Fc-mediated effector functions are critical to the efficacy and safety of therapeutic antibodies. Measurement of Fc-mediated antibody-mediated cellular cytotoxicity (ADCC) and complement-dependent cytotoxicity (CDC) during antibody drug discovery and development is not only important for antibodies that harness ADCC and/or CDC as their primary mechanism of action (e.g. rituximab, trastuzumab), but also for antibodies designed to target and block soluble ligands such as TNF α and VEGF.

We previously reported the development of a cell-based reporter bioassay platform which has been used to measure ADCC and ADCP mediated through Fc γ RI, Fc γ RIIa and Fc γ RIIIa. These reporter bioassays exhibit the specificity, accuracy, precision, and robustness necessary for qualification according to ICH guidelines and have been used extensively to characterize and measure the potency of antibody-based biologics drugs that target cell surface immune receptors. In the current study, we sought to evaluate Fc-mediated ADCC and CDC activities of therapeutic antibodies designed to target and block soluble ligands including TNF α and VEGF.

To measure ADCC activity of anti-TNF α and anti-VEGF blocking antibodies, we developed engineered target cells that express either membrane-bound TNF α or VEGF. When used as target cells with reporter-based effector cells expressing a relevant Fc γ R, ADCC activity of adalimumab (anti-TNF α) and bevacizumab (anti-VEGF) was detected in a specific and dose-dependent manner. Similarly, when used in a luminescence-based CDC assay, the engineered target cells elicited an appropriate Fc γ R-mediated response. The assay signals demonstrated IgG isotype specificity as IgG4 variants showed minimal activity in both ADCC and CDC assays. Our results demonstrate that the combined use of cell-based reporter bioassays with target cells engineered to express membrane-bound soluble ligands can provide a simple, specific, and quantitative platform to measure Fc-mediated effector functions of therapeutic antibodies targeting soluble ligands.

2. MOA-based ADCC & ADCP Reporter Bioassays

Surrogate Measure of In Vivo Biology

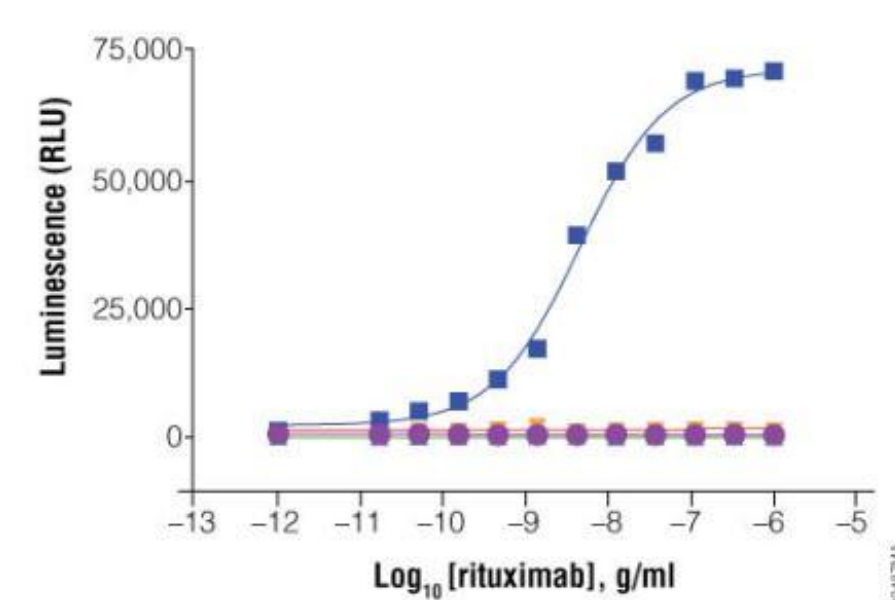
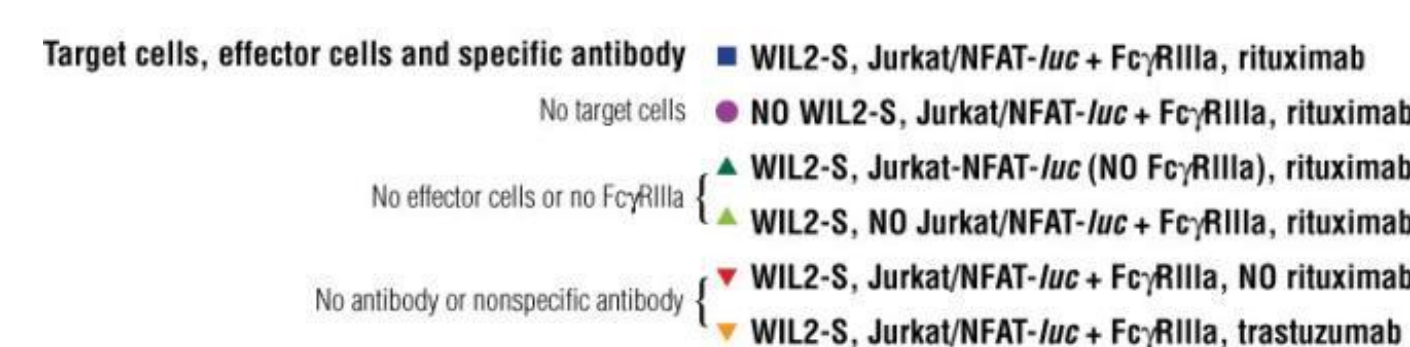


Fc γ R Reporter Bioassays are Specific

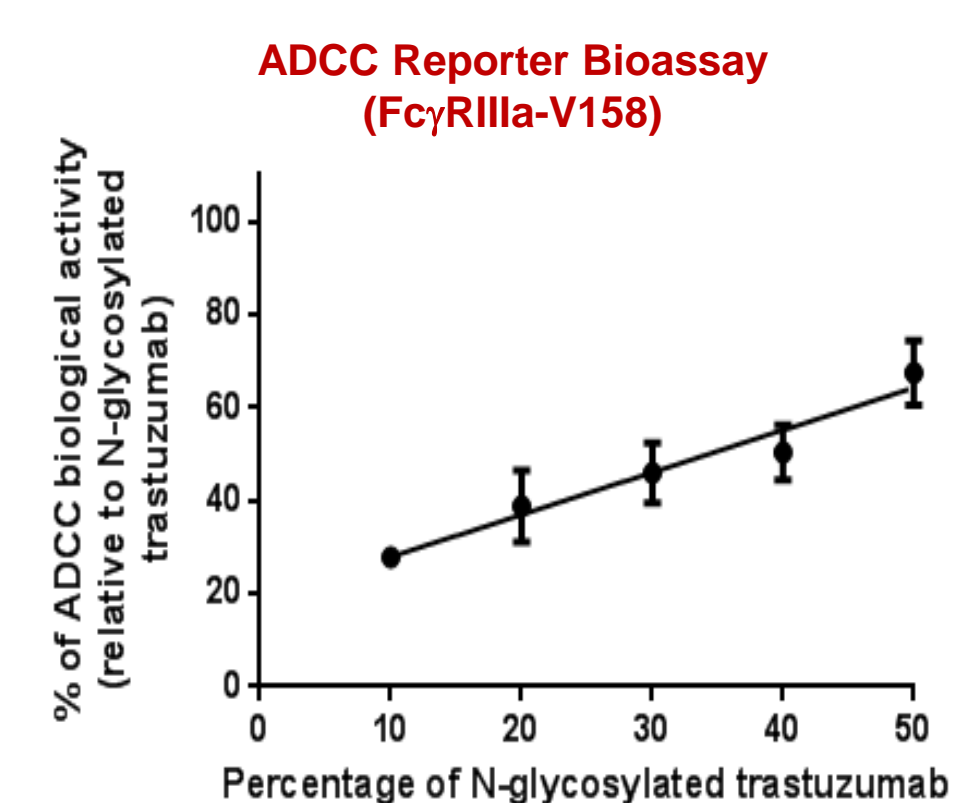
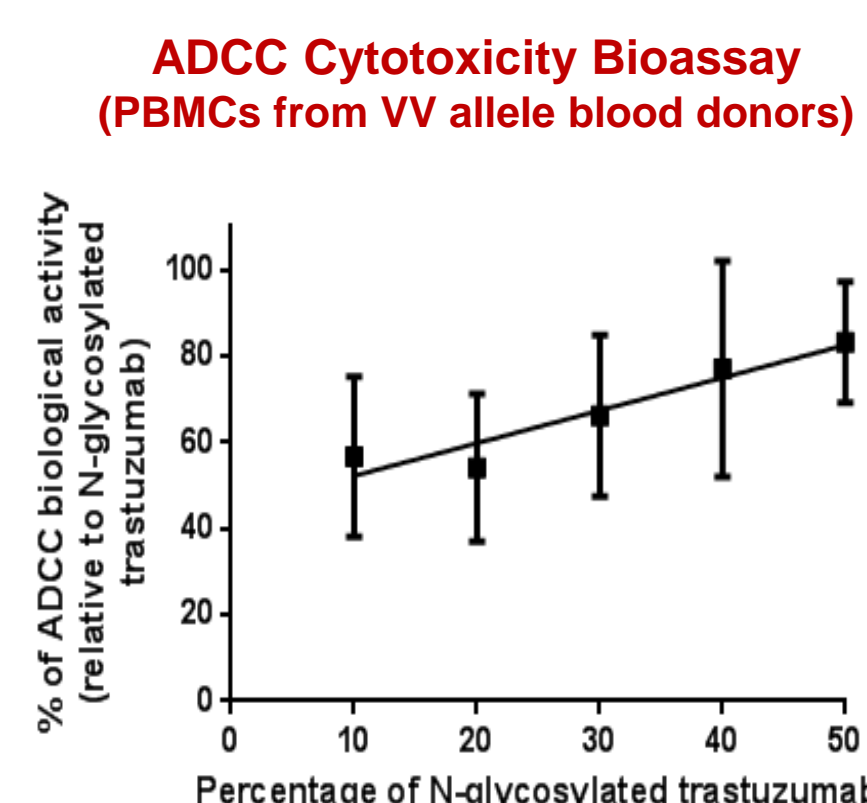
Variations of the ADCC Reporter Bioassay (Fc γ RIIIa-V158) were used to measure the Fc effector function of either Rituximab (anti-CD20) or Trastuzumab (anti-Her2) to demonstrate assay specificity.

Assay signal is dependent on:

- Target cells expressing Ab-targeted antigen
- Effector cells expressing Fc γ R
- Antigen-specific IgG antibody
- Similar results were obtained using human Fc γ RIIIa, Fc γ RIIa, Fc γ RI and mouse Fc γ RIII and Fc γ RIV bioassays.



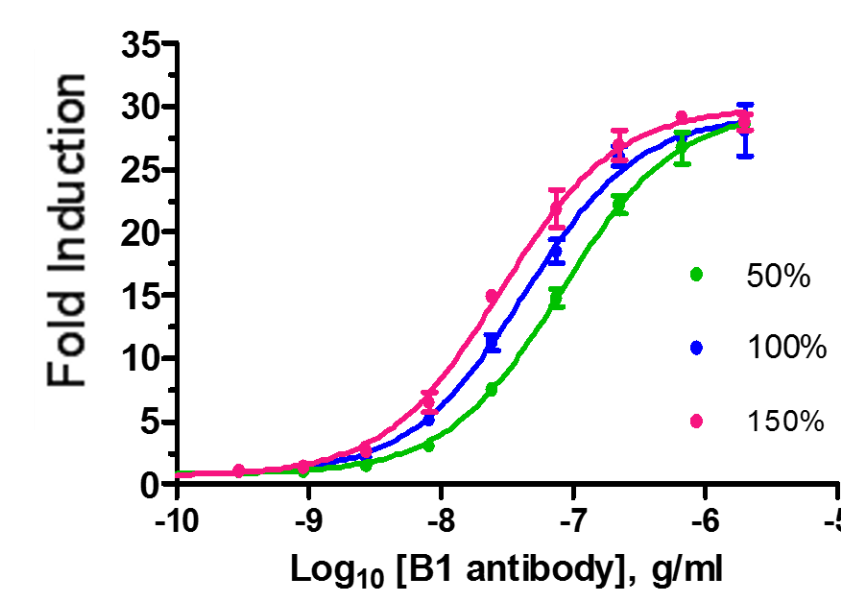
3. ADCC Reporter Bioassay Correlates with Primary Cell-based ADCC Cytotoxicity Assay



Antibody (trastuzumab) potency was measured using a primary cell-based ADCC assay (left panel) and the ADCC Reporter Bioassay (Fc γ RIIIa-V158, right panel). The surrogate ADCC Reporter Bioassay shows similar potency ranking but much less variability (as indicated by the smaller error bars).

4. Fc γ R Reporter Bioassays Measure Antibody Relative Potency and Stability

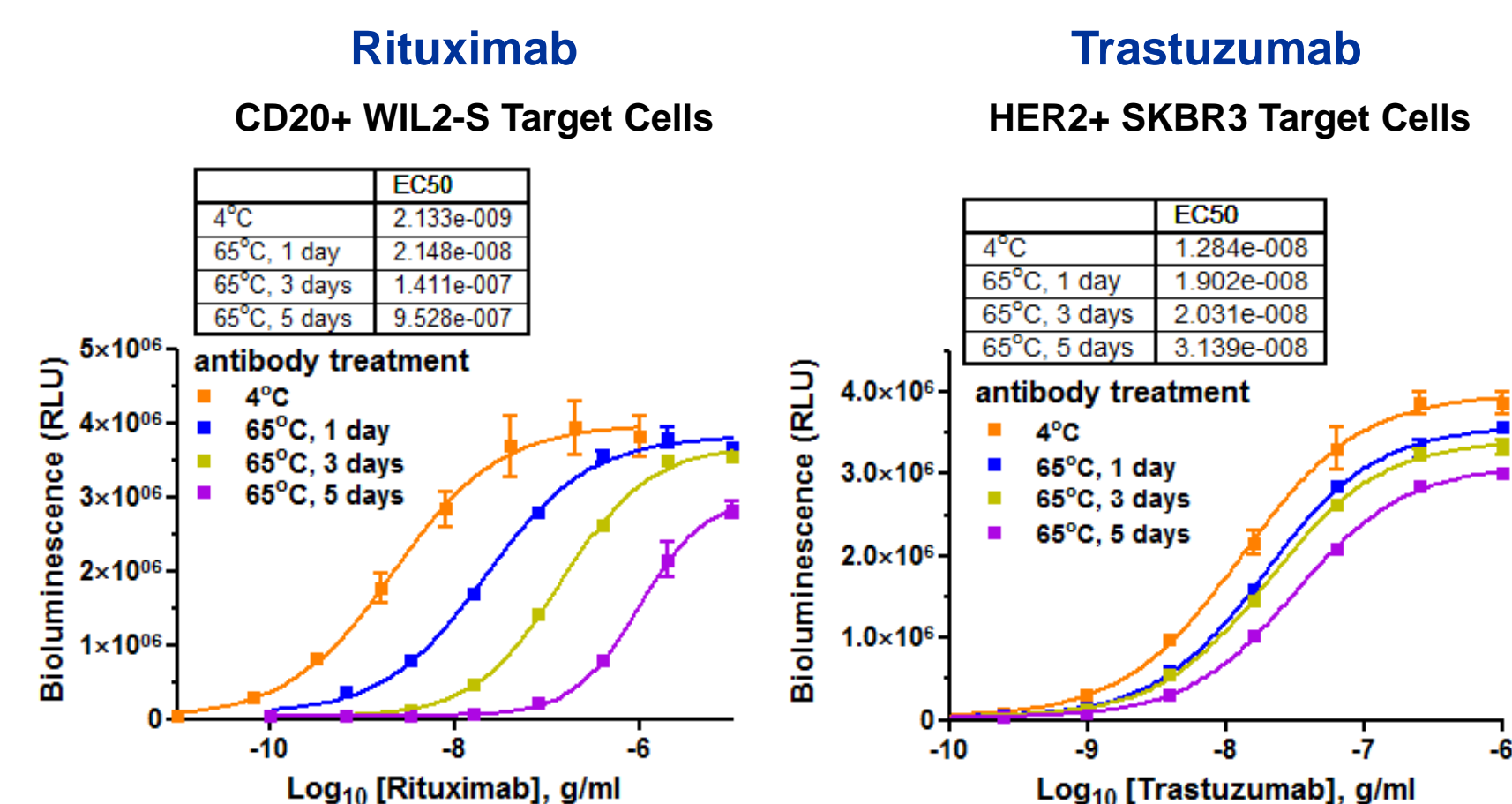
Potency Determination



The relative potency of anti-CD20 Ab preparations were measured using the ADCC Reporter Bioassay (Fc γ RIIIa-V158).

Excellent accuracy, precision, and linearity in the range of 50-150% relative potency (per ICH guidelines) was demonstrated (shown here and data not shown).

Stability Indication

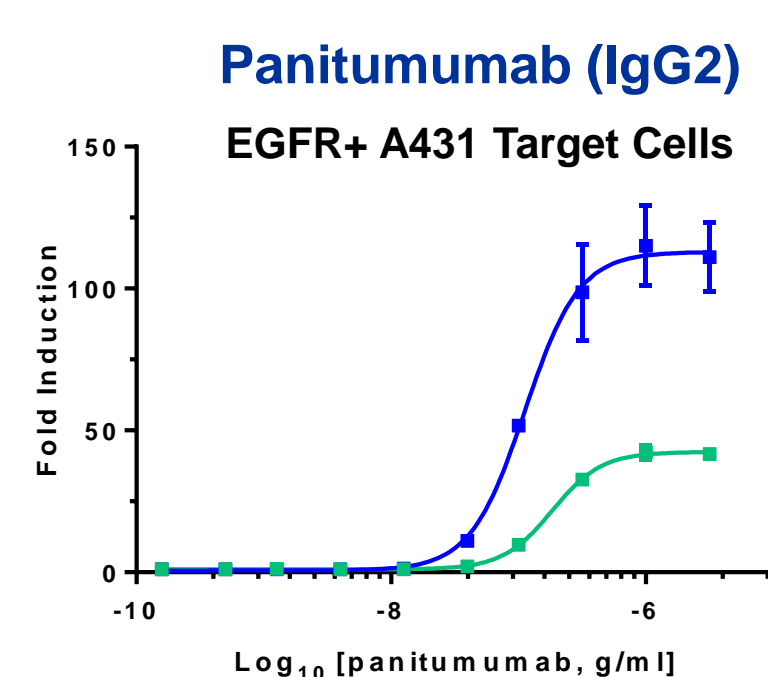


The ADCC Reporter Bioassay (Fc γ RIIIa-V158) was used in a stability study of Rituximab and Trastuzumab following heat denaturation at 65°C for the indicated number of days.

Similar results were obtained using human Fc γ RIIIa, Fc γ RIIa, Fc γ RI and mouse Fc γ RIII and Fc γ RIV bioassays.

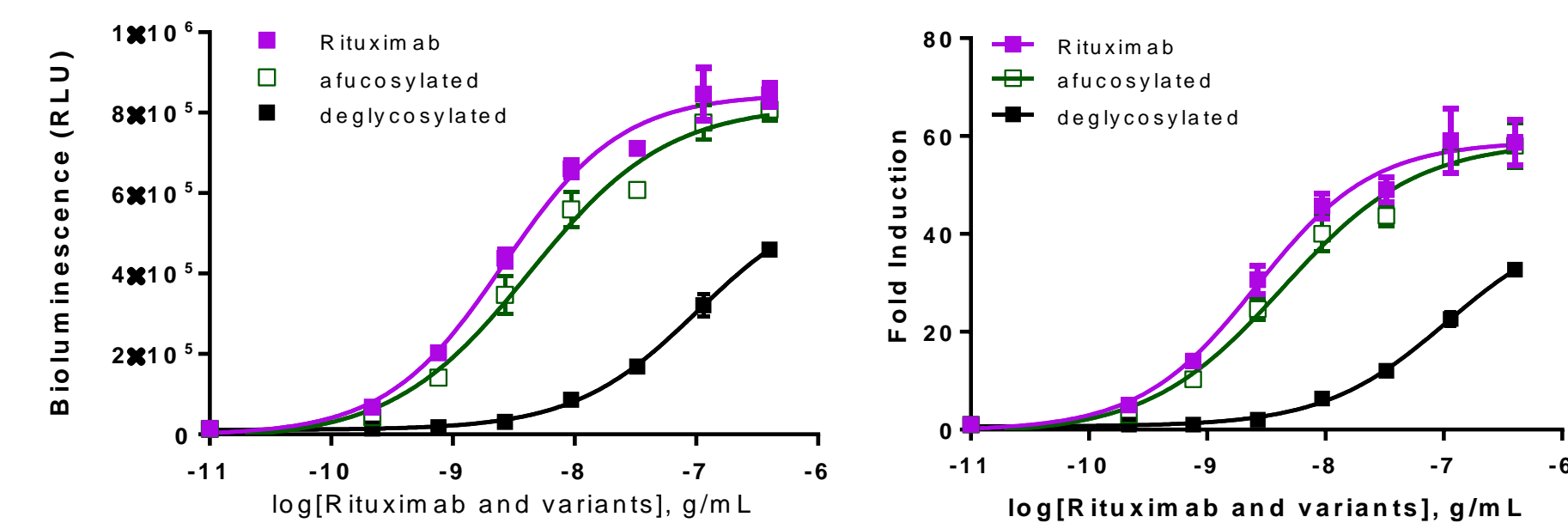
5. Application of Fc γ RIIIa & Fc γ RI ADCP Bioassays to Characterize Antibody MOA and Glycosylation

Fc γ RIIIa-mediated MOA



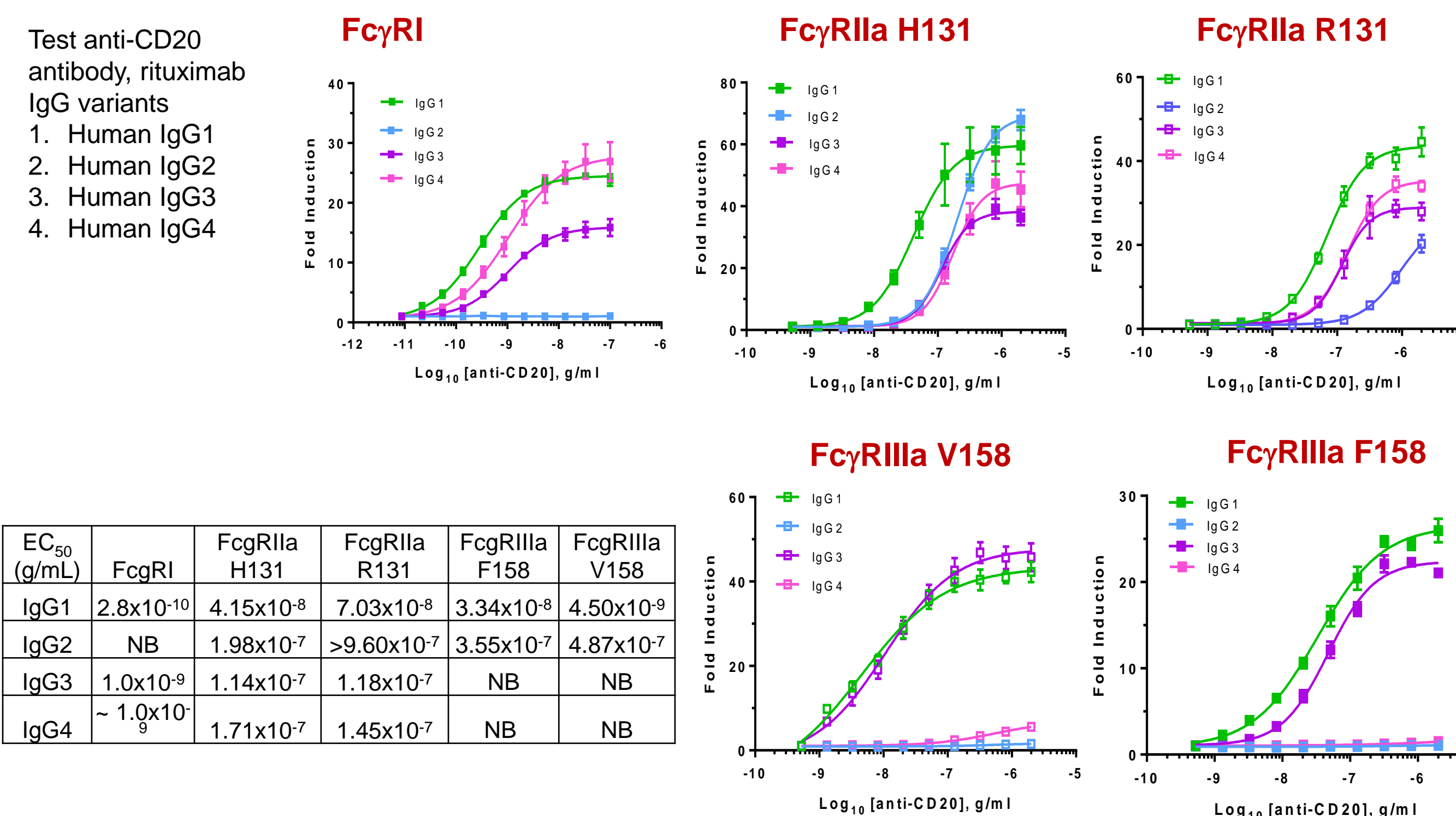
Legend: Fc γ RIIIa-H ADCP Reporter Bioassay (blue line), Fc γ RIIIa-R ADCP Reporter Bioassay (green line).

Antibody Deglycosylation Measured by Fc γ RI



CD20+ Raji Target Cells were assayed with Rituximab preparations that were either untreated, afucosylated or deglycosylated, as indicated. Deglycosylated antibody showed a significant decrease in bioluminescence (left panel) and fold induction (right panel), as expected.

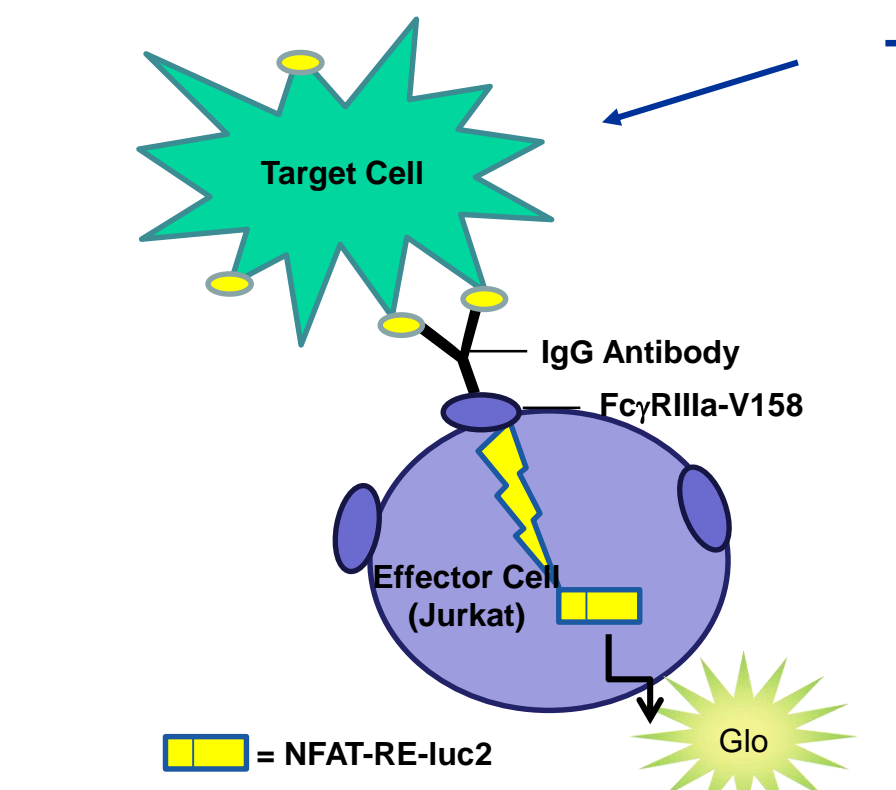
6. Fc γ R bioassays show appropriate IgG isotype Specificity



EC ₅₀ (g/mL)	Fc γ RI	Fc γ RIIIa H131	Fc γ RIIIa R131	Fc γ RIIIa F158	Fc γ RIIIa V158
IgG1	2.8x10 ⁻¹⁰	4.15x10 ⁻⁸	7.03x10 ⁻⁸	3.34x10 ⁻⁸	4.50x10 ⁻⁹
IgG2	NB	1.98x10 ⁻⁷	>9.60x10 ⁻⁷	3.55x10 ⁻⁷	4.87x10 ⁻⁷
IgG3	1.0x10 ⁻⁹	1.14x10 ⁻⁷	1.18x10 ⁻⁷	NB	NB
IgG4	~1.0x10 ⁻⁹	1.71x10 ⁻⁷	1.45x10 ⁻⁷	NB	NB

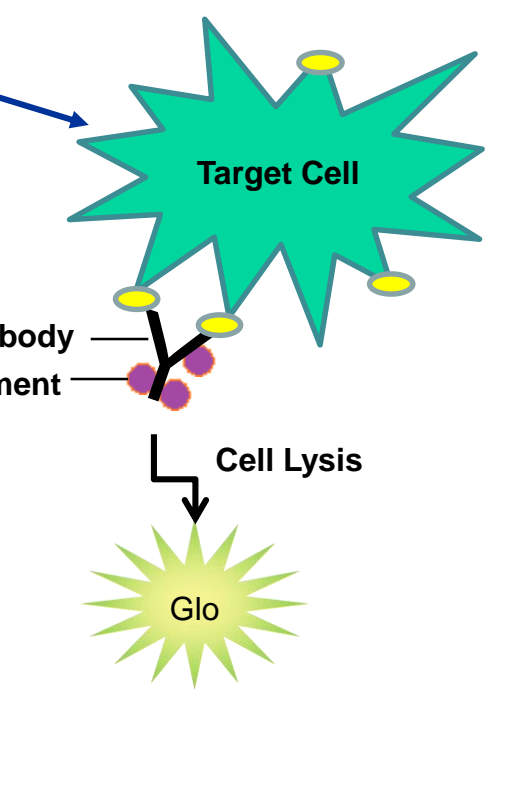
7. Anti-TNF α & Anti-VEGF ADCC & CDC Assays using Engineered Target Cells

ADCC Reporter Bioassay (Fc γ RIIIa-V158)



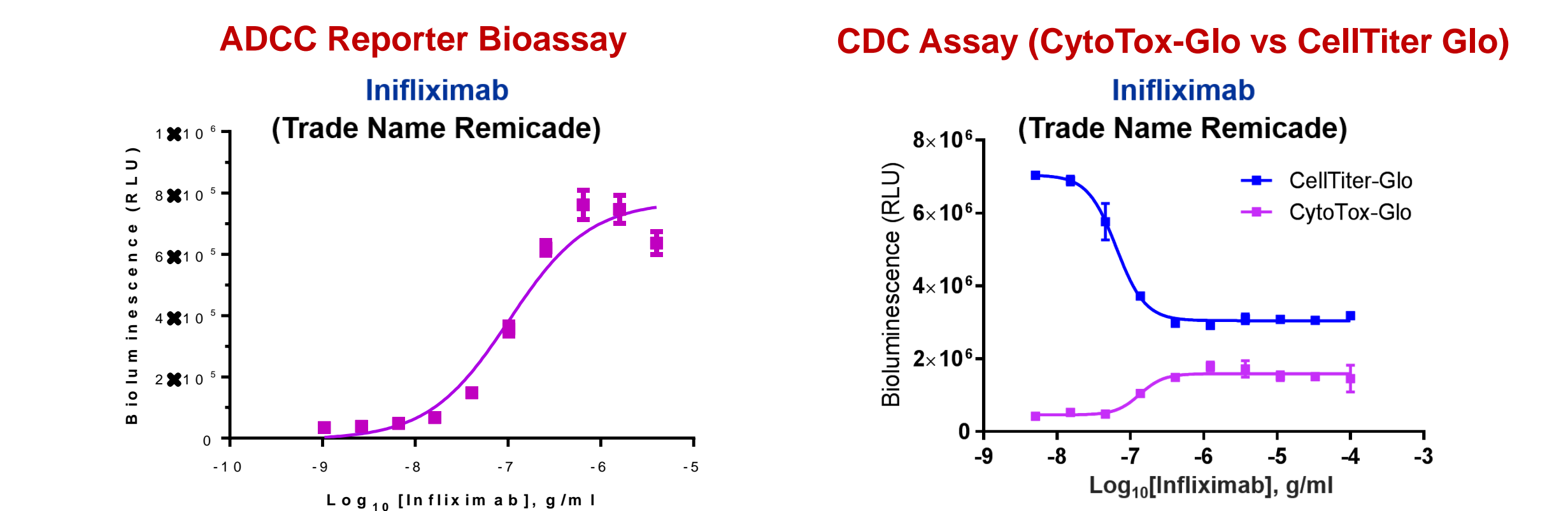
- Target cell-bound antibody binds to Fc γ RIIIa-V158
- Fc γ RIIIa-V158 signals through NFAT pathway in ADCC Effector Cells
- Luciferase activity measured using Bio-Glo reagent

CDC Assay

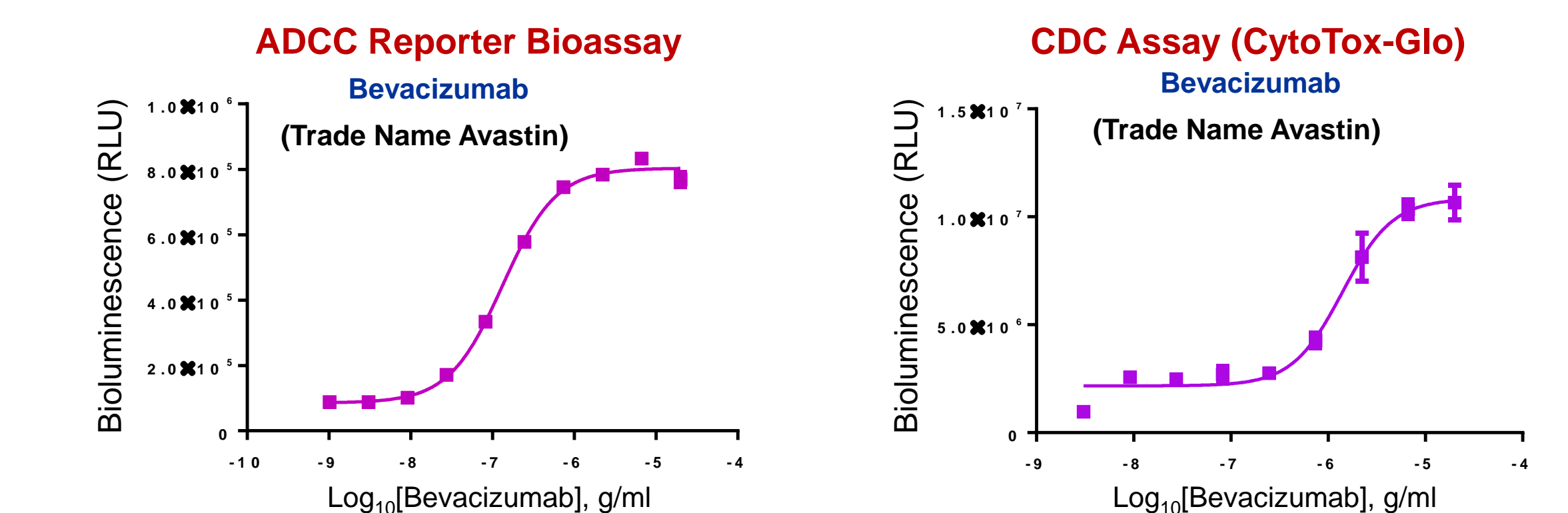


- Complement binds to target cell-bound antibody
- Activation of the complement cascade results in cell lysis
- Cell viability or cytotoxicity of target cells is measured using CellTiter-Glo (loss-of-signal) or CytoTox-Glo (gain-of-signal)

Measure anti-TNF α mAb-induced ADCC & CDC Activity



Measure anti-VEGF mAb-induced ADCC & CDC Activity



8. Conclusions

Functional cell-based assays for the measurement of ADCC and CDC are critical in the development of therapeutic antibodies. Here we show application of Fc γ R reporter bioassays, engineered target cells, cell viability and cytotoxicity assays to elucidate and characterize the Mechanism of Action (MOA) of antibody biologics for their Fc Effector Function.

Biologically relevant measurement of antibody Fc effector function

- ADCC reporter bioassays with specific Fc γ R-induced expression of NFAT-RE driven luciferase
- Specific Ab-mediated target cell killing via CDC
- Appropriate IgG isotype specificity in ADCC and CDC assays

Consistent and reliable measure of antibody activity

- Demonstrated precision, accuracy, reproducibility, robustness

Easy-to-implement

- Multiple product formats meet diverse workflows; commercial kits include necessary reagents
- Rapid protocols for standard 96-well plate format