

# Development of a bioluminescent cell-based bioassay to measure Fc receptor functionality in antibody-dependent cell-mediated cytotoxicity

Zhi-jie J. Cheng, Shan Chung\*, Denise Garvin, Rich Moravec, Aileen Paguio, Valerie Quarmbly\*, Frank Fan and Teresa K. Surowy

Department of Research, Promega Corporation, Madison, WI 53711. email: [zey.cheng@promega.com](mailto:zey.cheng@promega.com);

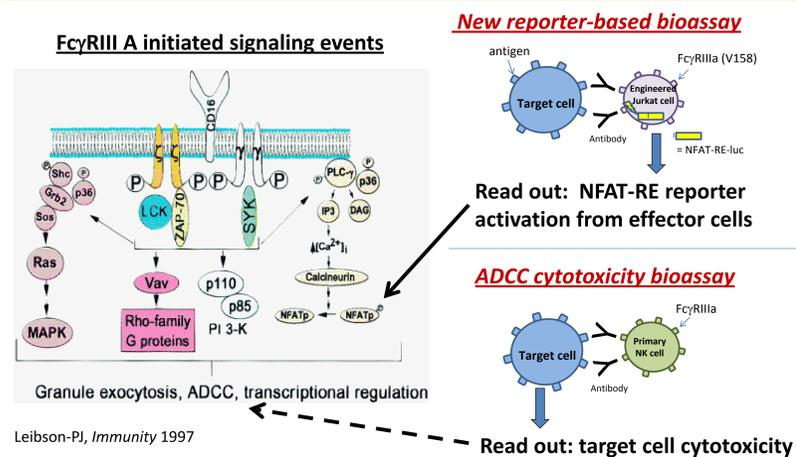
\* BioAnalytical Sciences, Genentech, Inc., South San Francisco, CA 94080



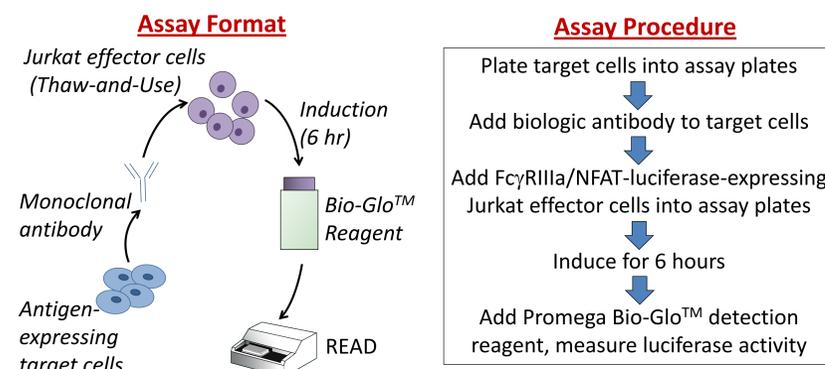
## 1. Abstract

Primary peripheral blood mononuclear cells (PBMCs) are routinely used in traditional bioassays to quantify antibody drug potency in antibody-dependent cellular cytotoxicity (ADCC). These bioassays are labor intensive and have high inherent assay variability. Here, we report the development of a bioluminescent cell-based bioassay which measures activation of effector cells via cross-linking of FcγRIIIA with target cell-bound antibodies. For this, Jurkat T-cell stable cell lines that stably express NFAT-luciferase reporter and human FcγRIIIA were generated to replace primary PBMCs as effector cells in ADCC bioassay. Effector cells were also developed in a frozen, thaw-and-use format to minimize assay variability due to cell culture and handling. The resultant bioassay using this format demonstrated good assay precision and accuracy in bioassay qualification. Bioassay using the engineered effector cells is robust, specific and is able to quantify the potencies of rituximab and trastuzumab, two monoclonal antibody drugs for cancer. When used to measure effects of Fc glycosylation on effector functions of therapeutic antibodies, a linear correlation was observed between relative antibody activity and the extent of Fc glycosylation. Thus, the bioluminescent cell-based reporter bioassay provides a simple and robust approach to measure potency of therapeutic antibodies in ADCC with high precision.

## 2. Introduction

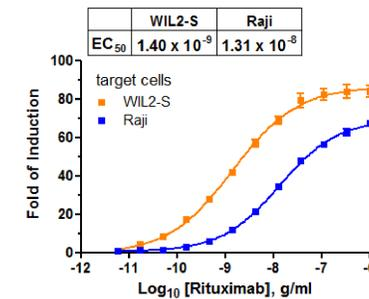


## 3. Reporter bioassay method and assay format

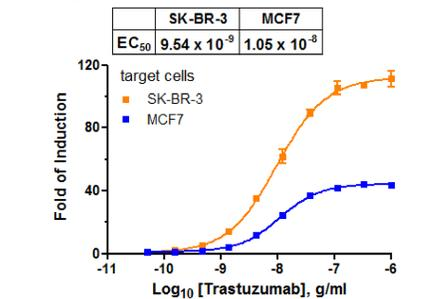


## 4. Measuring Potency for Biologic Drugs Rituximab (anti-CD20) and Trastuzumab (anti-Her2)

### A. CD20<sup>+</sup> B cell lines (suspension) as target cells



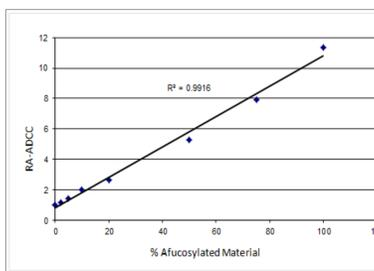
### B. Her2<sup>+</sup> breast cancer cell lines (adherent) as target cells



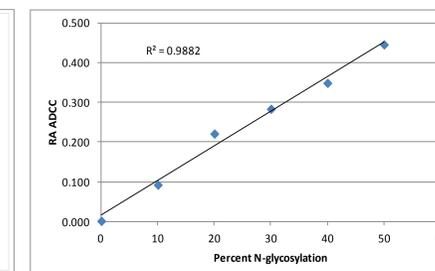
Quantifying the potencies of Rituximab (trade name: Rituxan, anti-CD20 chimeric monoclonal antibody drug) or Trastuzumab (trade name: Herceptin, anti-Her2 humanized monoclonal antibody drug) with appropriate target cell systems.

## 5. Fc effector functionality in new ADCC reporter bioassay correlates with antibody afucosylation and N-glycosylation

### A. Relative ADCC reporter activity vs percentage of antibody afucosylation



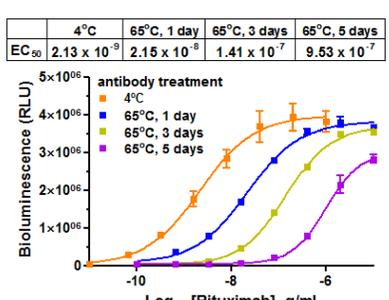
### B. Relative ADCC reporter activity vs percentage of Rituximab N-glycosylation



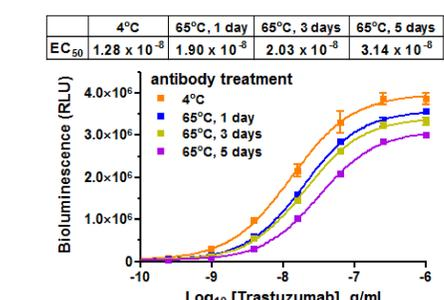
- Linear correlation shown between percentage of afucosylated antibody in blended antibody samples and relative luciferase reporter activity in ADCC reporter assay
- Linear correlation shown between percentage of N-glycosylated antibody in blended Rituximab samples and relative reporter activity in ADCC reporter assay

## 6. Stability indicating properties of new reporter bioassay for monoclonal antibody biologic drugs

### A. Potencies of heat-treated Rituximab

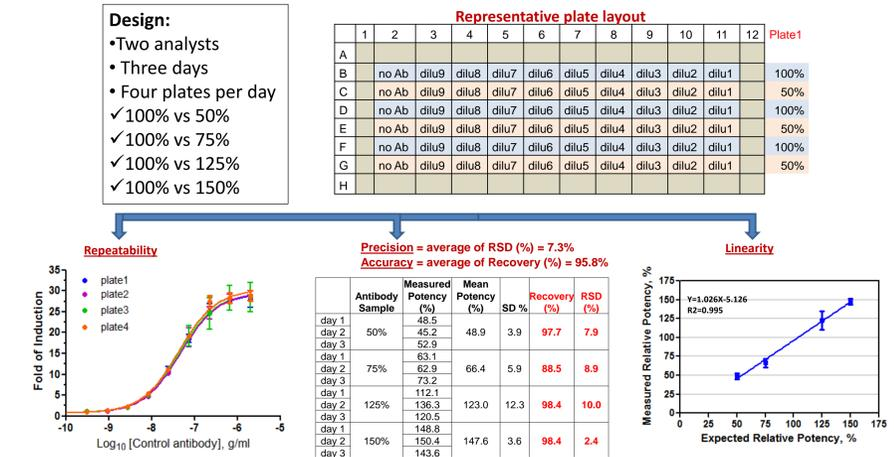


### B. Potencies of heat-treated Trastuzumab



Quantifying the potencies of heat-treated Rituximab (A) or Trastuzumab (B) samples using ADCC reporter bioassay.

## 7. Bioassay Qualification: Repeatability, Precision, Accuracy and Linearity



## 8. ADCC Bioassay kit configurations

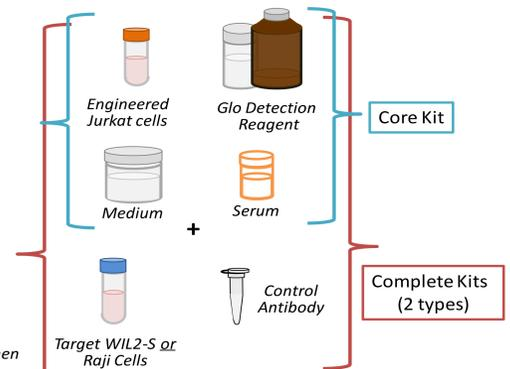
### Core Kit

- Engineered Jurkat Effector cells (NFAT-RE-luc2 / FcγRIIIA); frozen, thaw-and-use
  - Bio-Glo™ Luciferase Assay System
  - RPMI Medium
  - Low IgG Serum
- Use with your own Abs and target cells

### Complete Kits (2 types)

- Above PLUS:
- Target WIL2-S cells; frozen, thaw-and-use or Target Raji cells; frozen thaw-and-use
- Control Ab (CD20)

Use as an assay run control, or when WIL2-S or Raji target cells are suitable or your target is CD20



## 9. Summary

- We developed a bioluminescent cell-based reporter bioassay which can be used to quantify biologic potency of therapeutic monoclonal antibodies based on Fc effector functionality as a result of target cell-bound antibody activation of FcγRIIIA receptor on engineered effector cells.
- Replacement of primary NK cells with the engineered Jurkat effector cells eliminates the high bioassay variability typically seen in assays using NK cells from blood donors.
- The bioassay was used to assess the impact of glycan afucosylation and N-glycosylation of antibodies on their Fc effector functionality. Linear correlations obtained between the extent of antibody afucosylation or N-glycosylation and relative antibody bioactivity show that this assay, like traditional ADCC bioassays, is able to appropriately quantify Fc effector functionality of monoclonal antibodies.
- Use of thaw-and-use effector and target cells improves assay precision due to tight control over consistency of cell preparations for the assay.
- The assay was qualified as a bioassay for therapeutic monoclonal antibodies.