

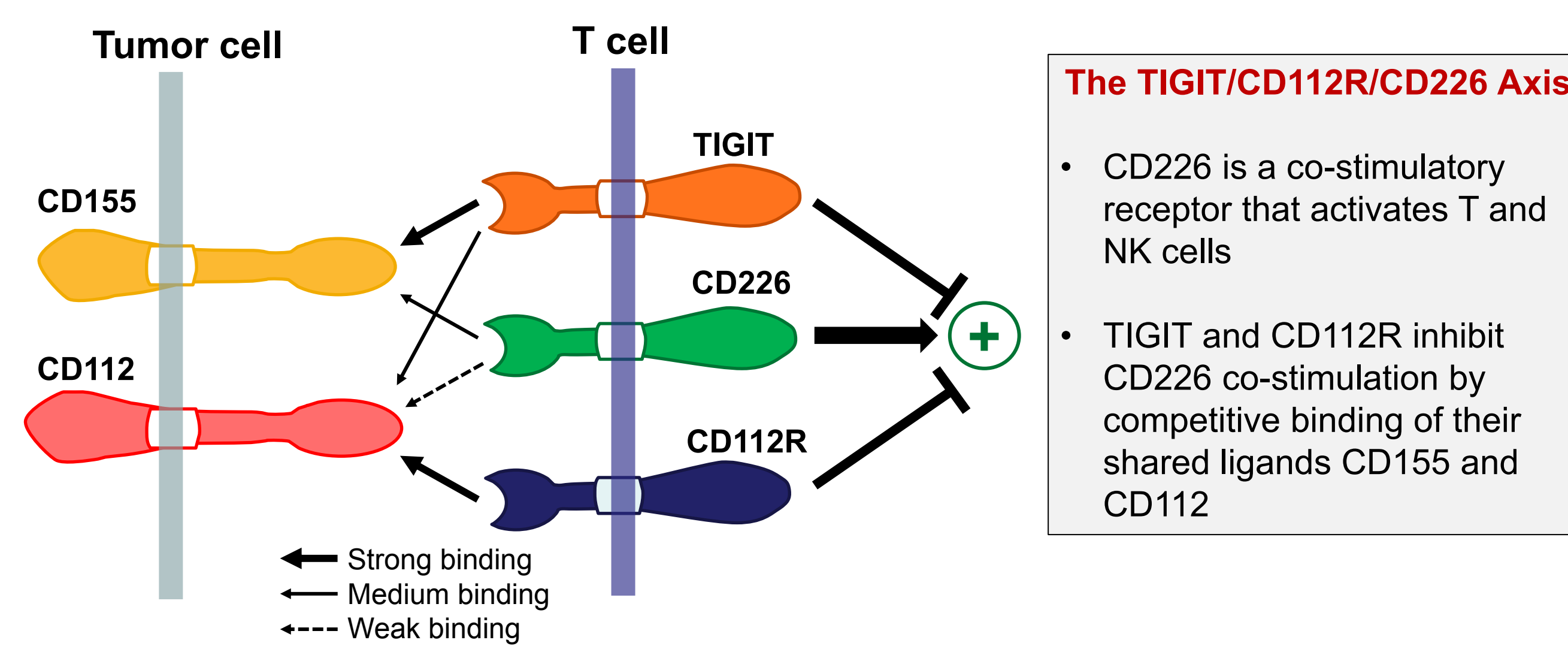
MoA-Based Potency Bioassays for Immunotherapy Programs Targeting the TIGIT/CD112R/CD226 Axis



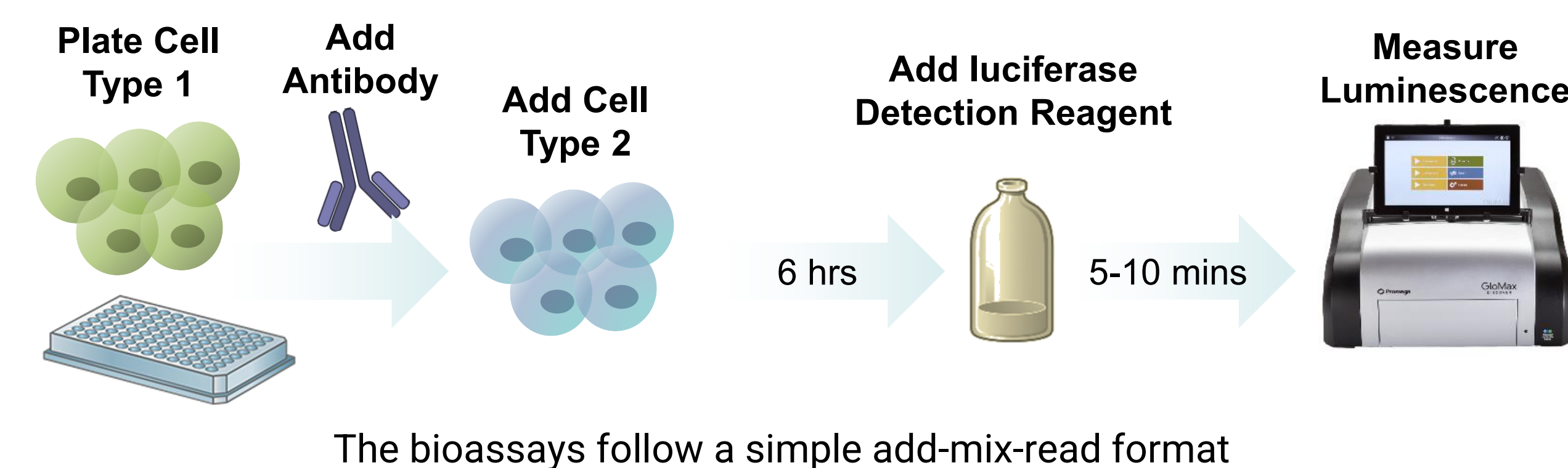
Jamison Grailer, Denise Garvin, Jim Hartnett, Kai Hillman, Frank Fan, Mei Cong, and Zhi-Jie Jey Cheng
Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711

1. Introduction

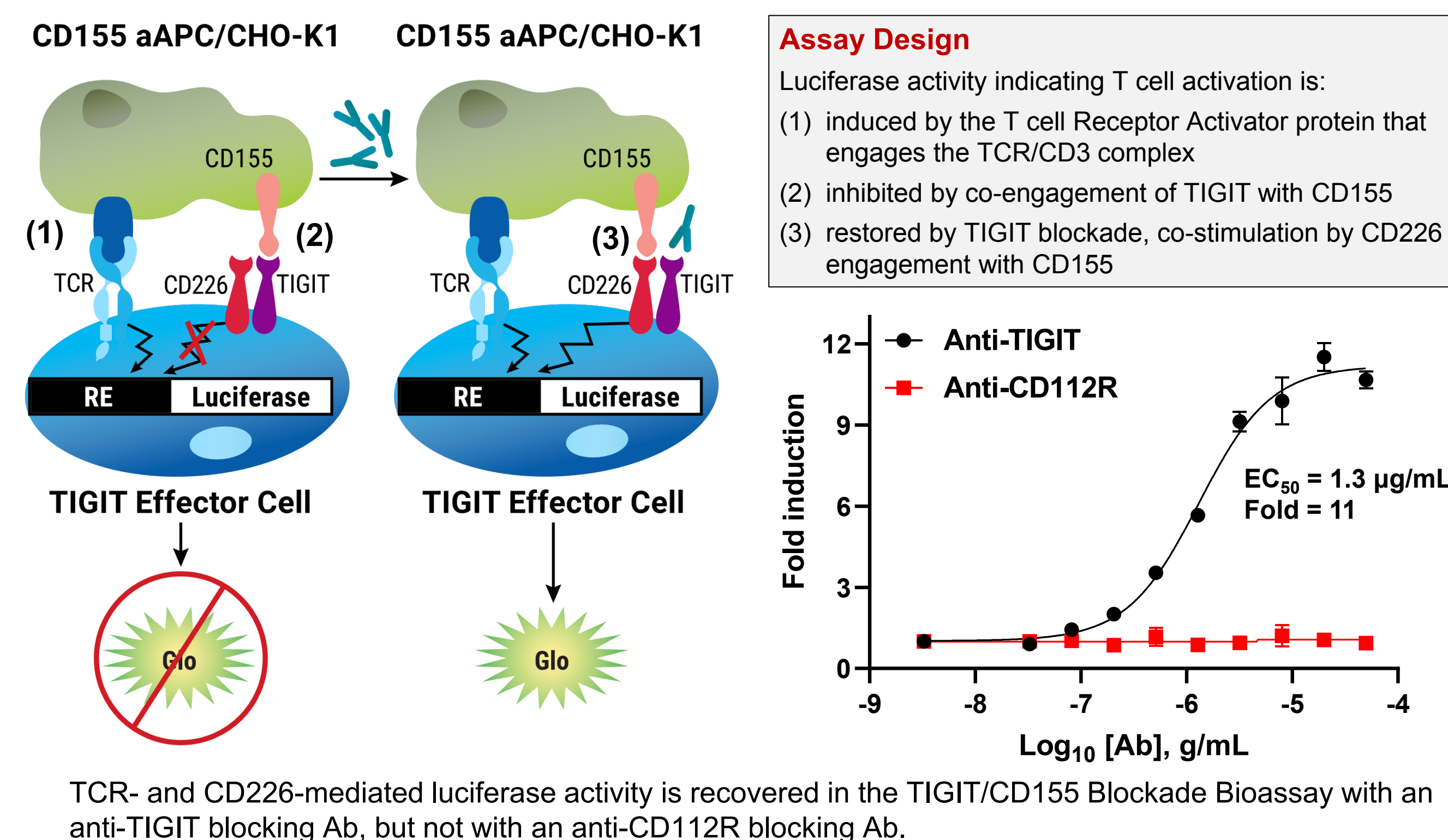
- TIGIT, CD112R and CD226 are immune checkpoints that have gained attention for cancer immunotherapy due to promising preclinical results
- We developed a portfolio of cell-based reporter bioassays to measure the potency of biologics designed to target the TIGIT/CD112R/CD226 axis



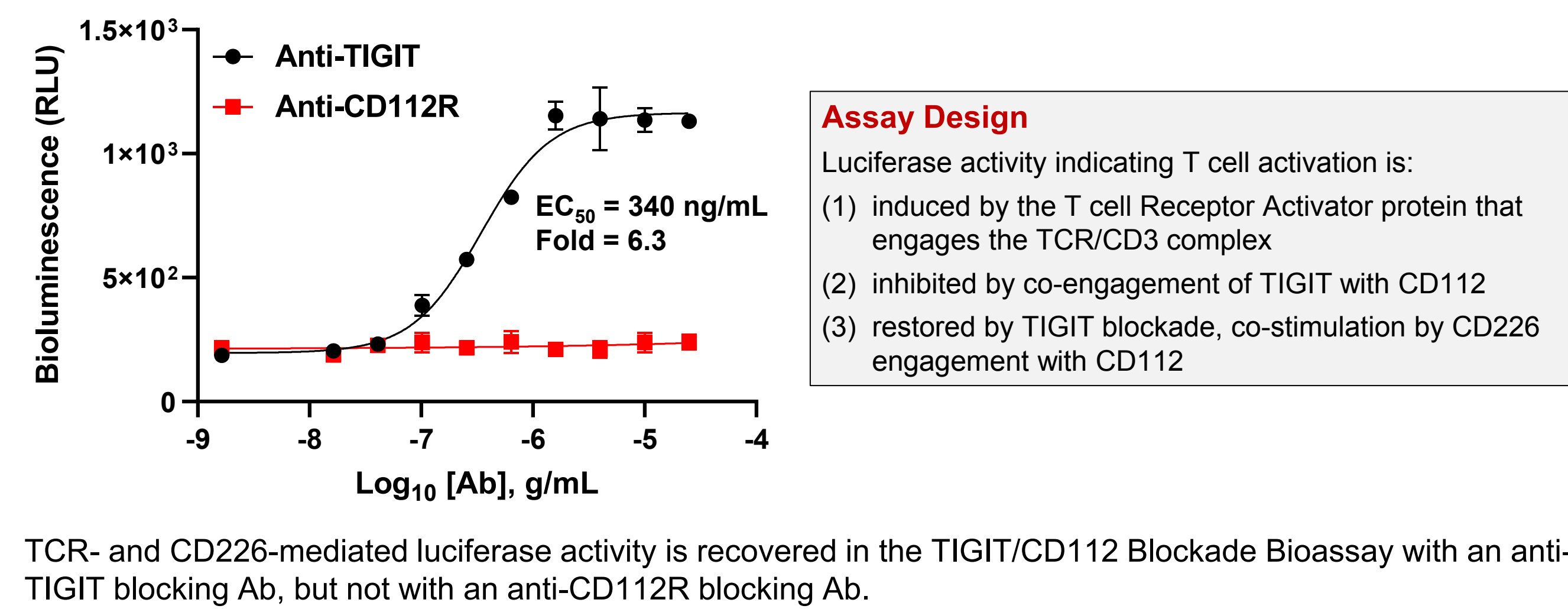
2. General Bioassay Workflow



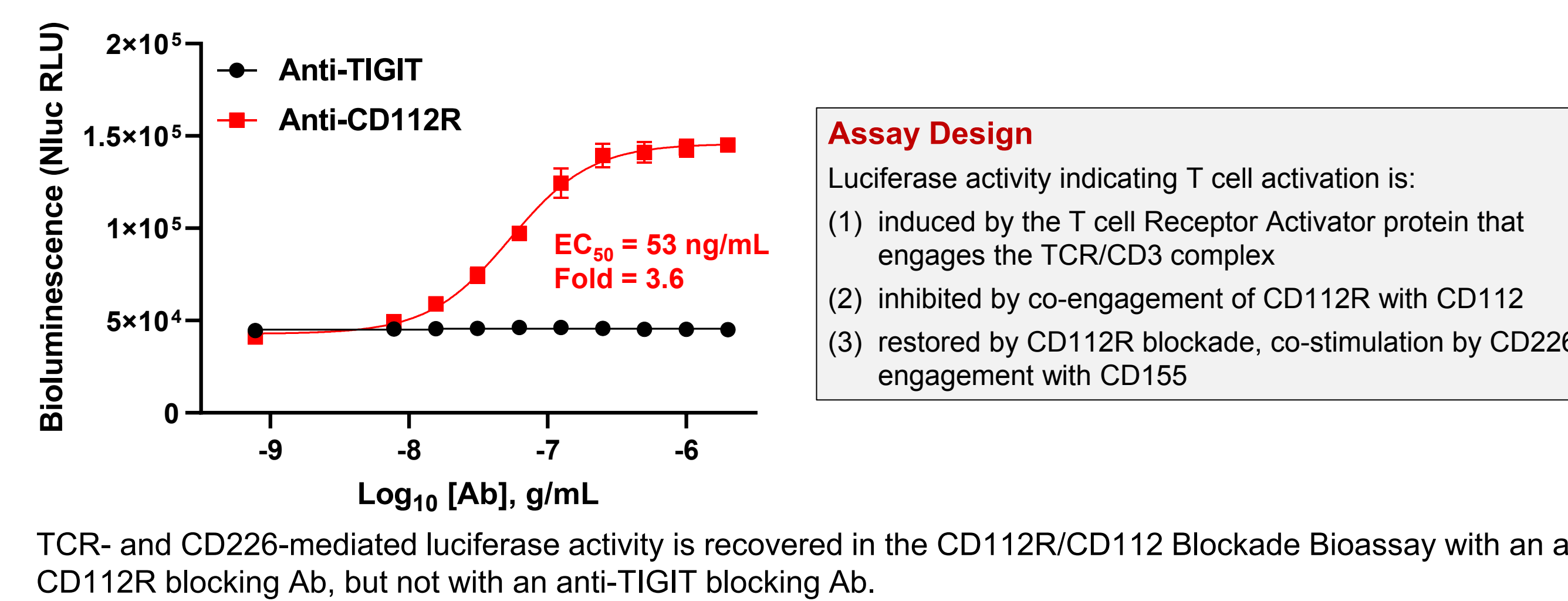
3. The TIGIT/CD155 Blockade Bioassay Measures the Potency of TIGIT/CD155 Blocking Antibodies



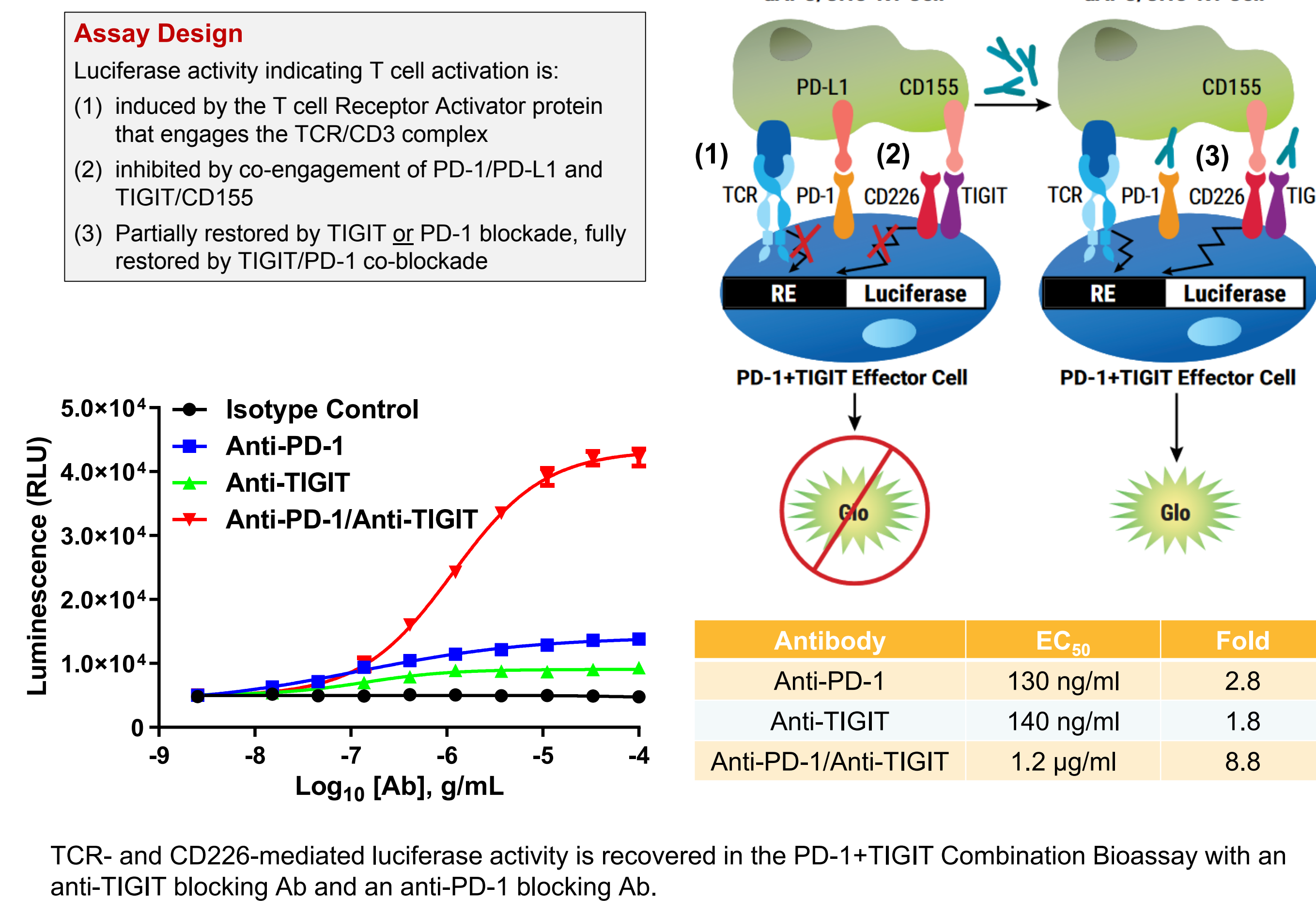
4. The TIGIT/CD112 Blockade Bioassay Measures the Potency of TIGIT/CD112 Blocking Antibodies



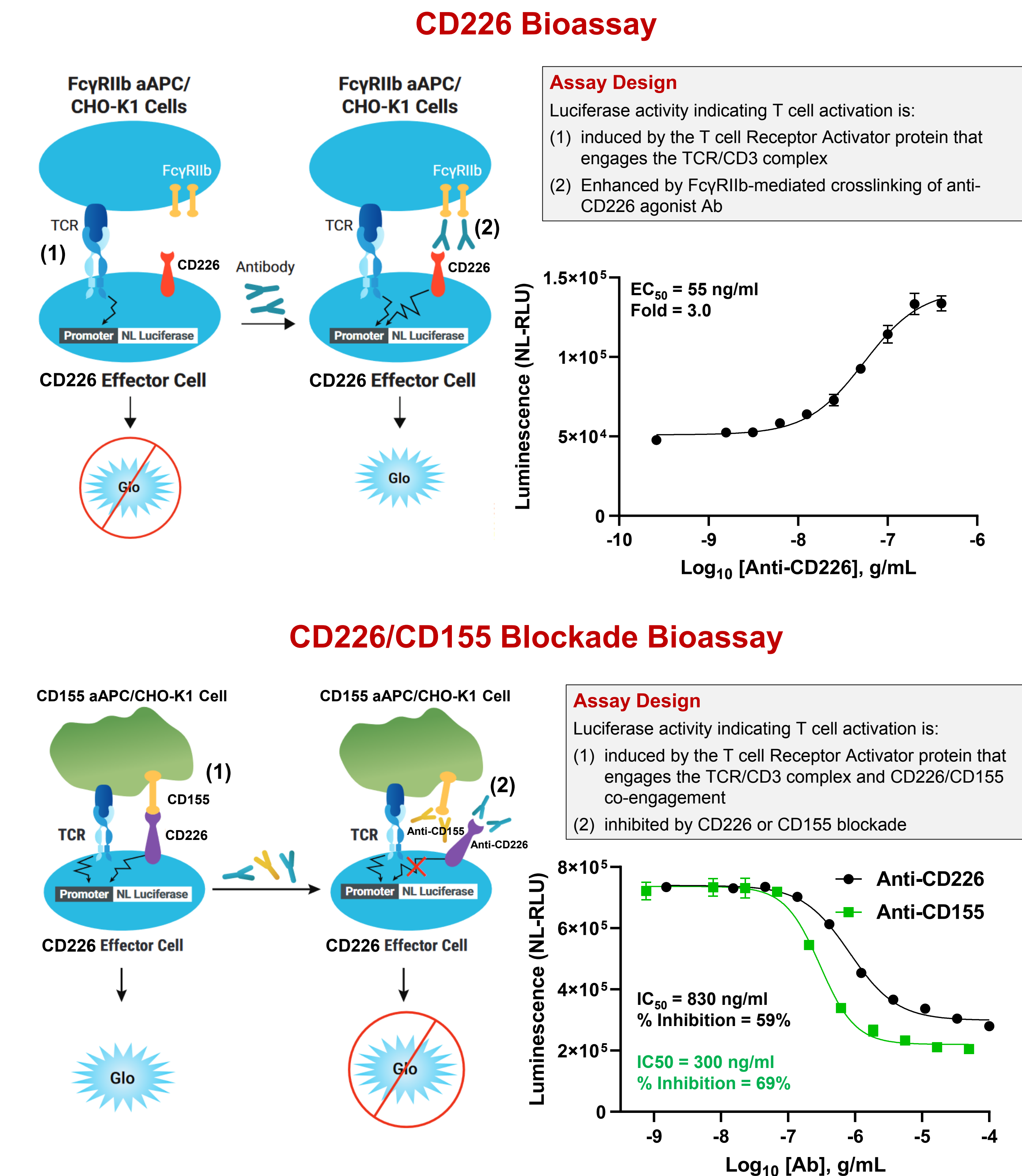
5. The CD112R/CD112 Blockade Bioassay Measures the Potency of CD112R Blocking Antibodies



6. The PD-1+TIGIT Combination Bioassay Measures Synergetic Effects of Blocking Antibodies for PD-1 and TIGIT



7. The CD226 Bioassays Measure the Potency of CD226 Agonist or Antagonist Antibodies



8. Conclusions

Here we show a portfolio of MOA-based bioassays for the TIGIT/CD112R/CD226 axis that can be used for antibody screening, characterization, potency and stability studies. These bioassays provide the following:

Biologically relevant measurement of antibody MOA

- Specific immune checkpoint-regulated expression of luciferase that reflects the native biology of T cell activation
- Demonstrated ability to measure the potencies of immune checkpoint-targeted antibodies

Consistent and reliable measure of antibody activity

- Demonstrated precision, accuracy, reproducibility, robustness
- All assays can be used as "Thaw-and-use" cell format, no cell culture required
- Functional performance suitable for development into potency and stability assays

Easy-to-implement

- Rapid and convenient workflow, amenable to 96- and 384-well plates