

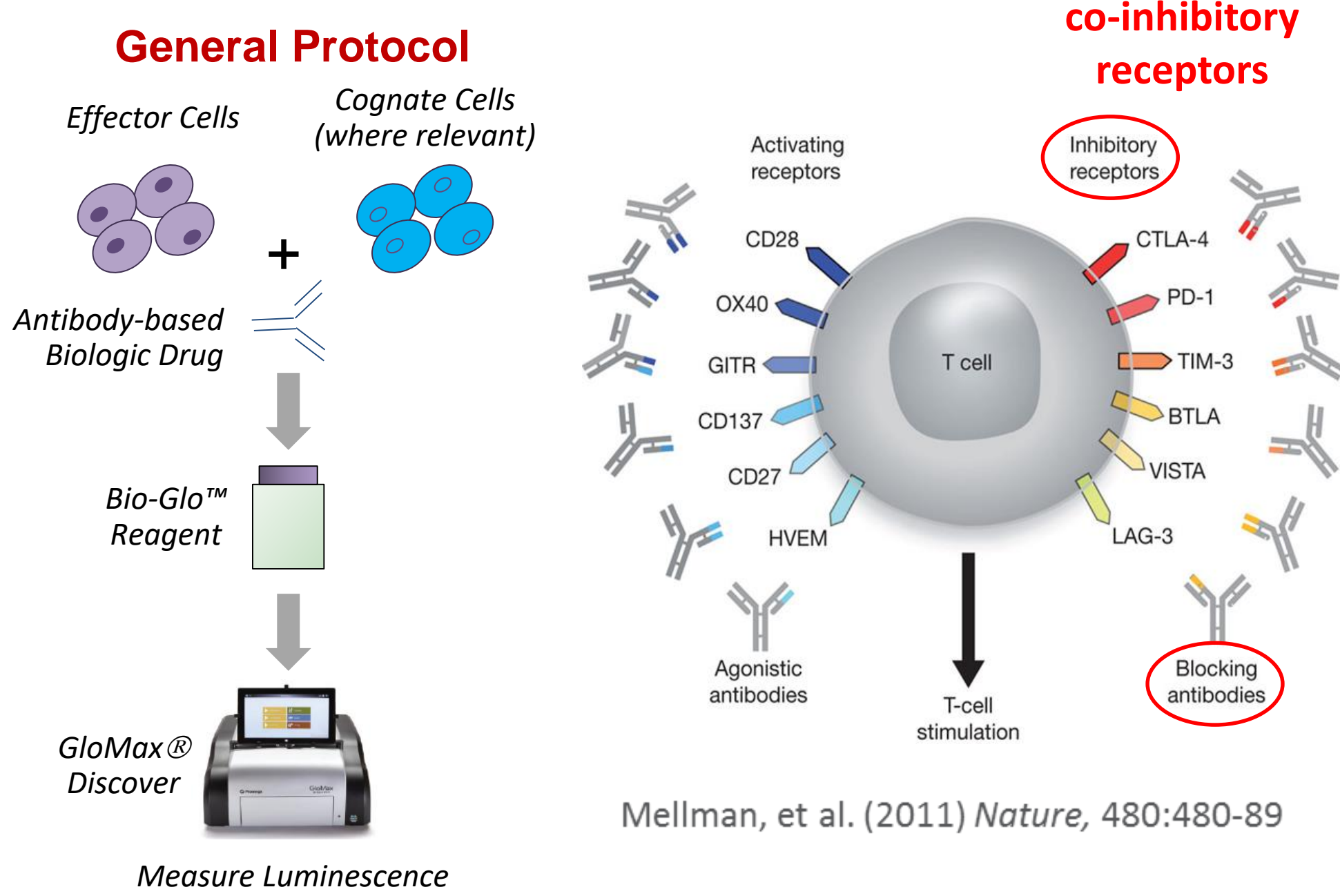
# Releasing the Brakes: Quantitative Cell-based Bioassays to Advance Individual and Combination Immune Checkpoint Immunotherapy

Jamison Grailer, Julia Gilden, Pete Stecha, Denise Garvin, Jim Hartnett, Frank Fan, Mei Cong and Zhi-jie Jey Cheng  
Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711

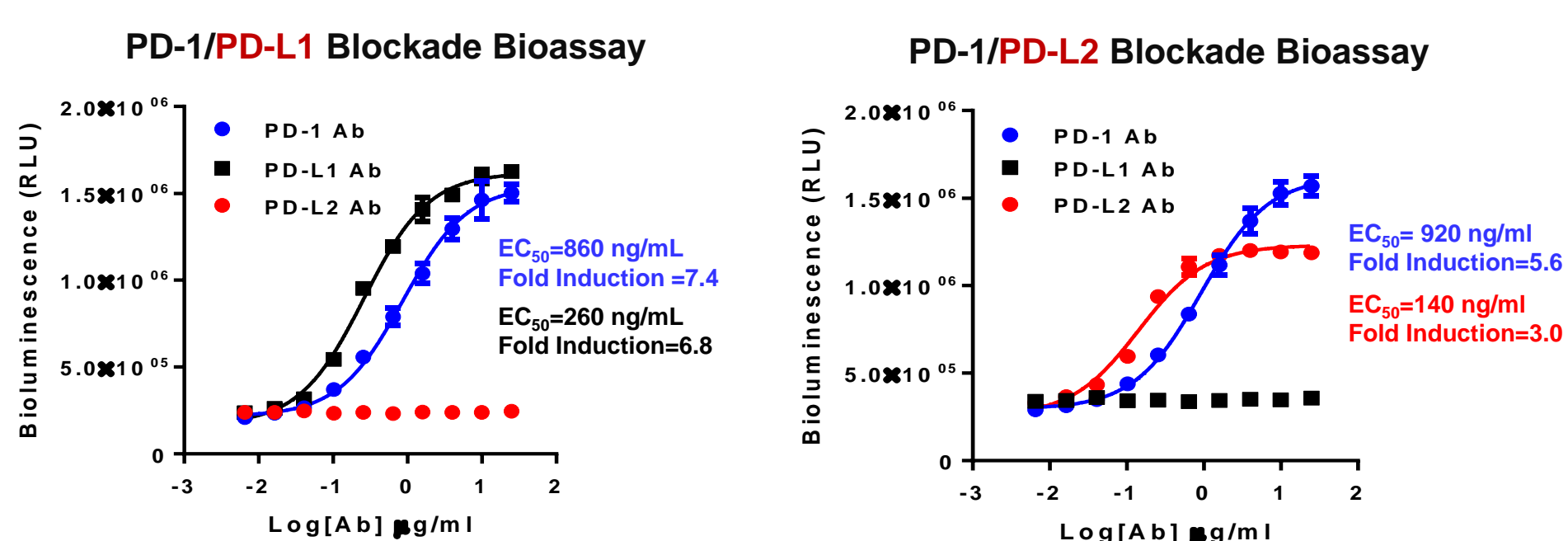
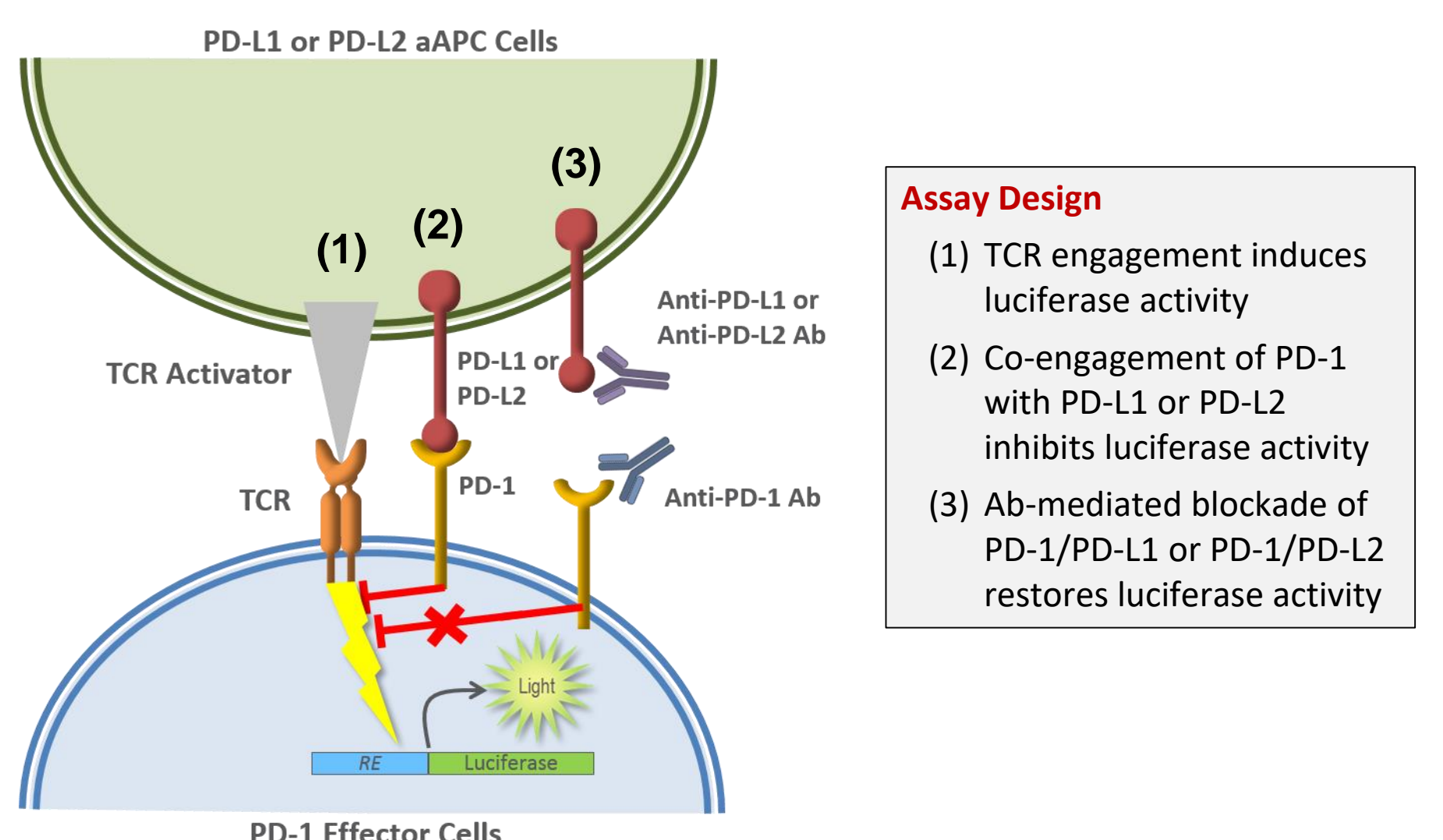


## 1. Introduction

A major challenge in the development of antibody-based biologics drugs is access to quantitative and reproducible functional bioassays. Existing methods rely on primary cells and measurement of complex functional endpoints that are cumbersome, variable, and often fail to yield data quality required for drug development in a quality-controlled environment. We have developed a portfolio of functional cell-based reporter bioassays to measure the activity of biologics drugs designed to target immune checkpoint receptors including co-inhibitory (e.g. PD-1, CTLA-4, LAG-3) and co-stimulatory (e.g. 4-1BB, GITR, OX40) receptors. These bioassays consist of stable cell lines that express luciferase under the precise control of receptor-mediated intracellular signals. Here we describe the application of MOA-based immune checkpoint co-inhibitory receptor bioassays for biologics drug discovery, development, potency and stability studies.

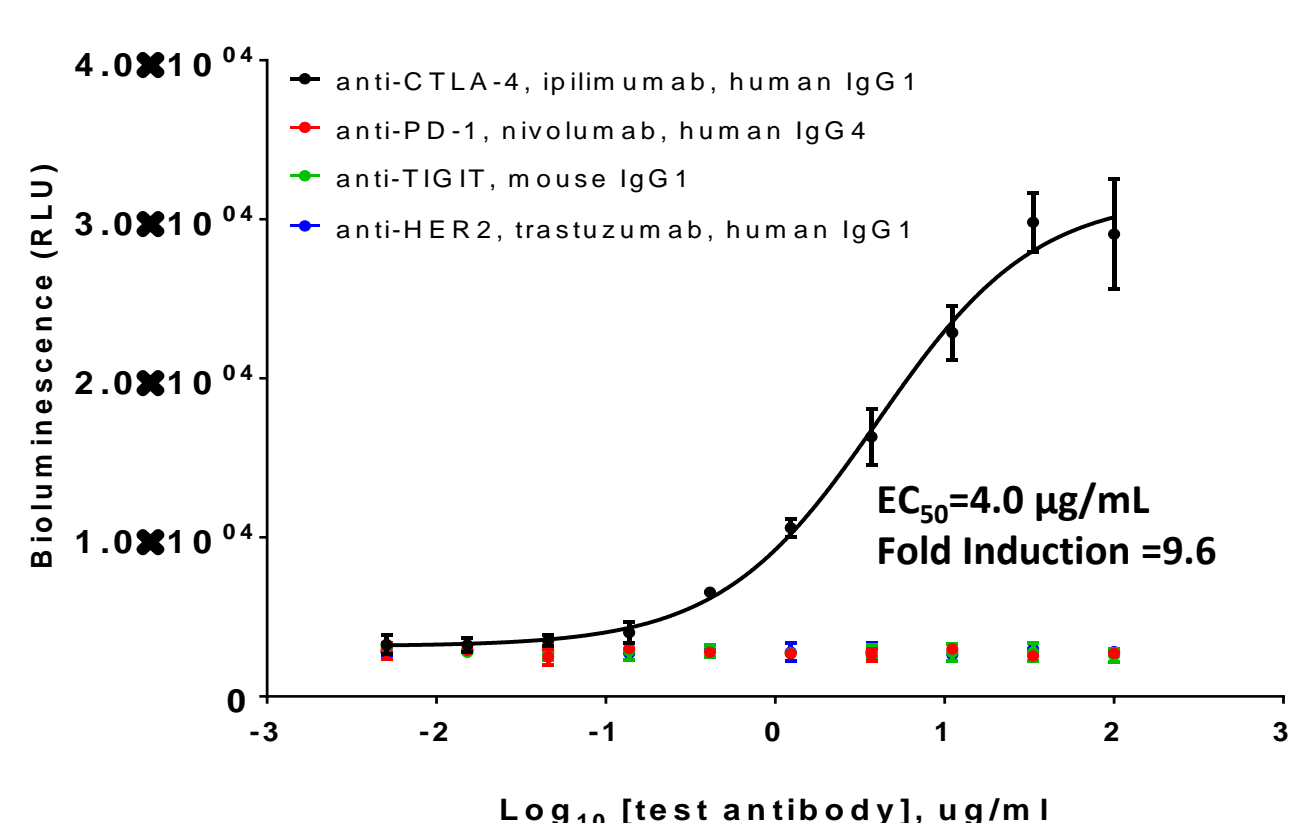
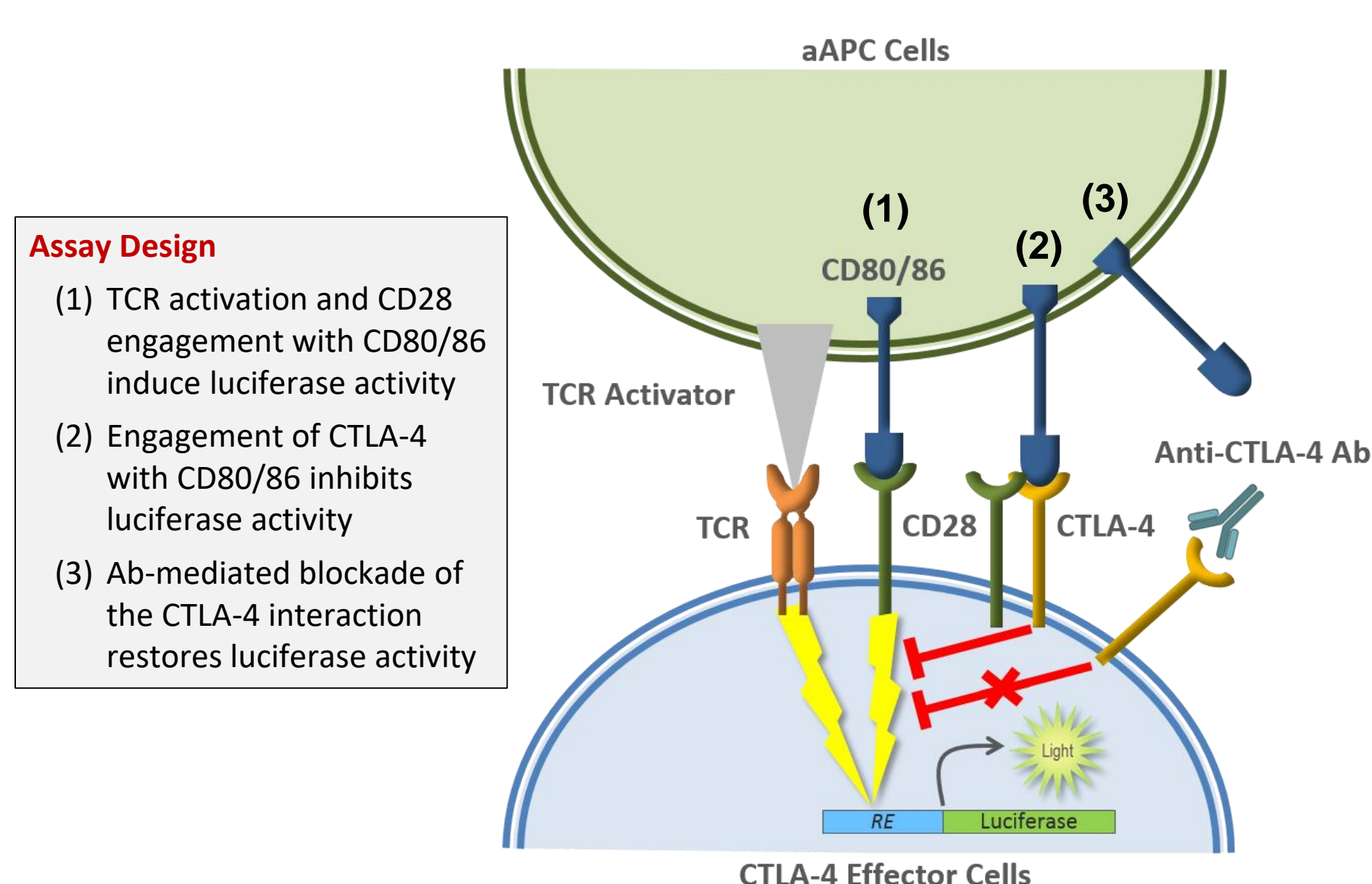


## 2. PD-1 Blockade Bioassays



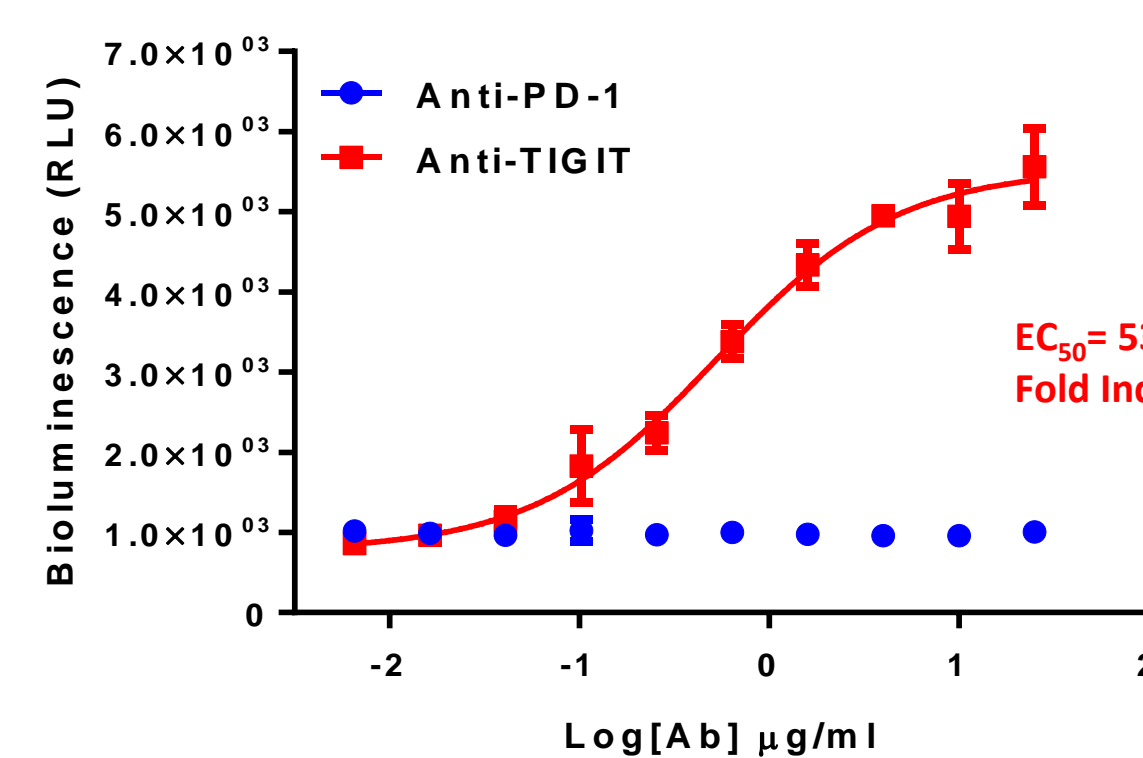
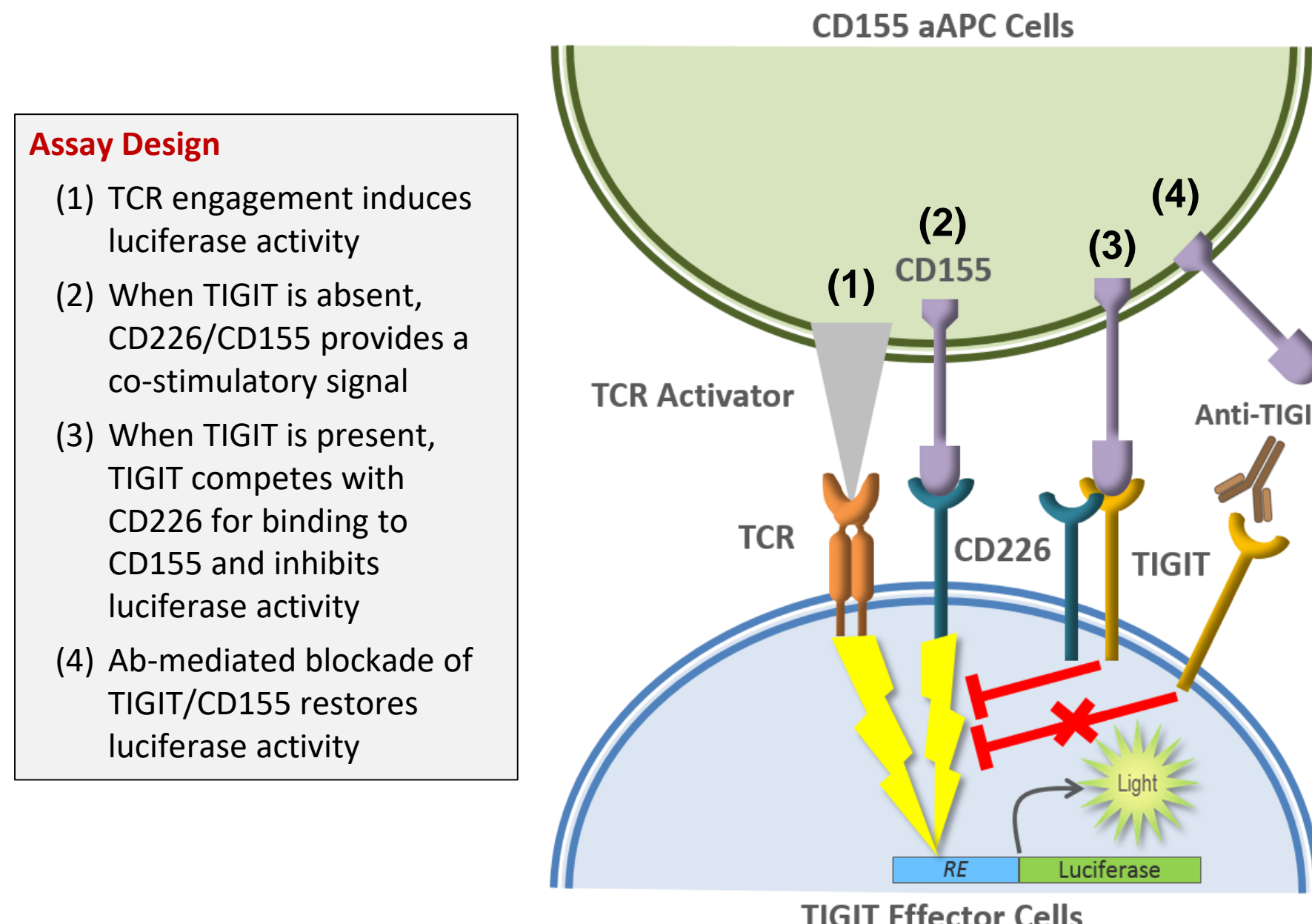
**Left:** TCR-mediated luciferase activity is recovered in the PD-1/PD-L1 bioassay with anti-PD-1 or anti-PD-L1 blocking Abs, but not with anti-PD-L2 blocking Ab.  
**Right:** TCR-mediated luciferase activity is recovered in the PD-1/PD-L2 bioassay with anti-PD-1 or anti-PD-L2 blocking Abs, but not with anti-PD-L1 blocking Abs.  
All Abs shown here are research grade.

## 3. CTLA-4 Blockade Bioassay



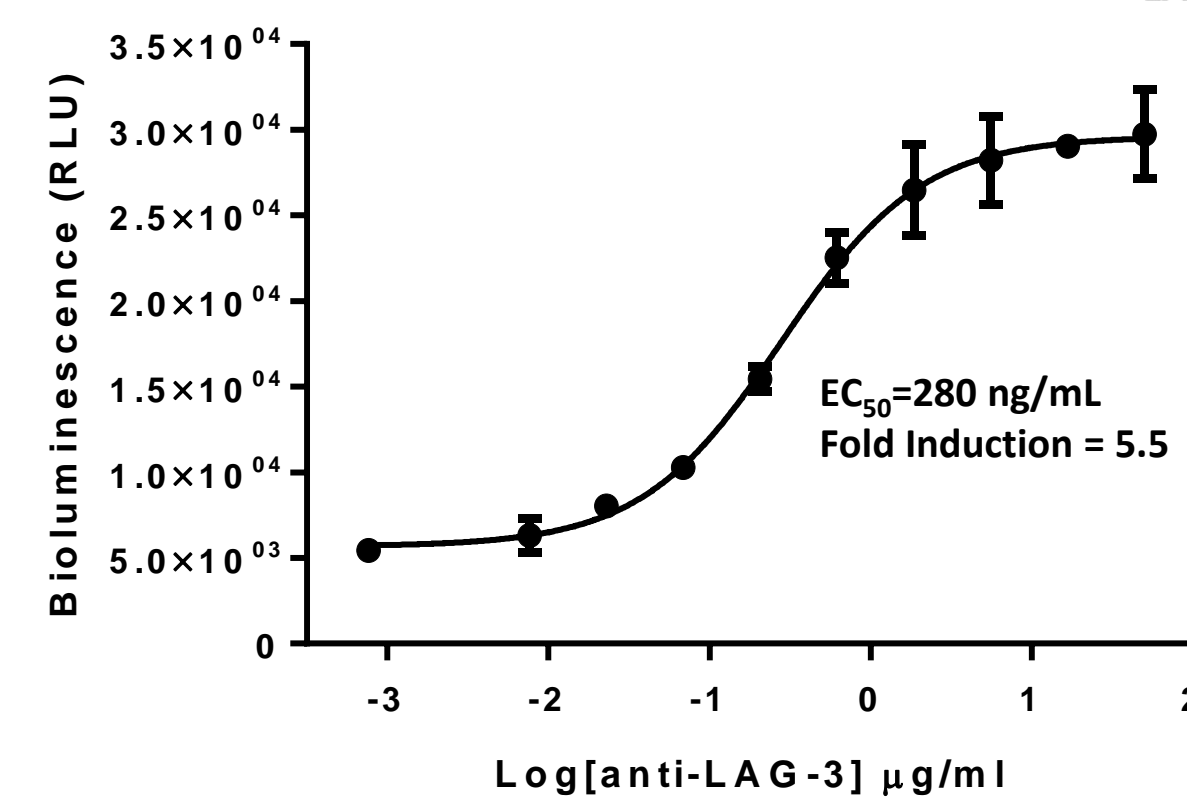
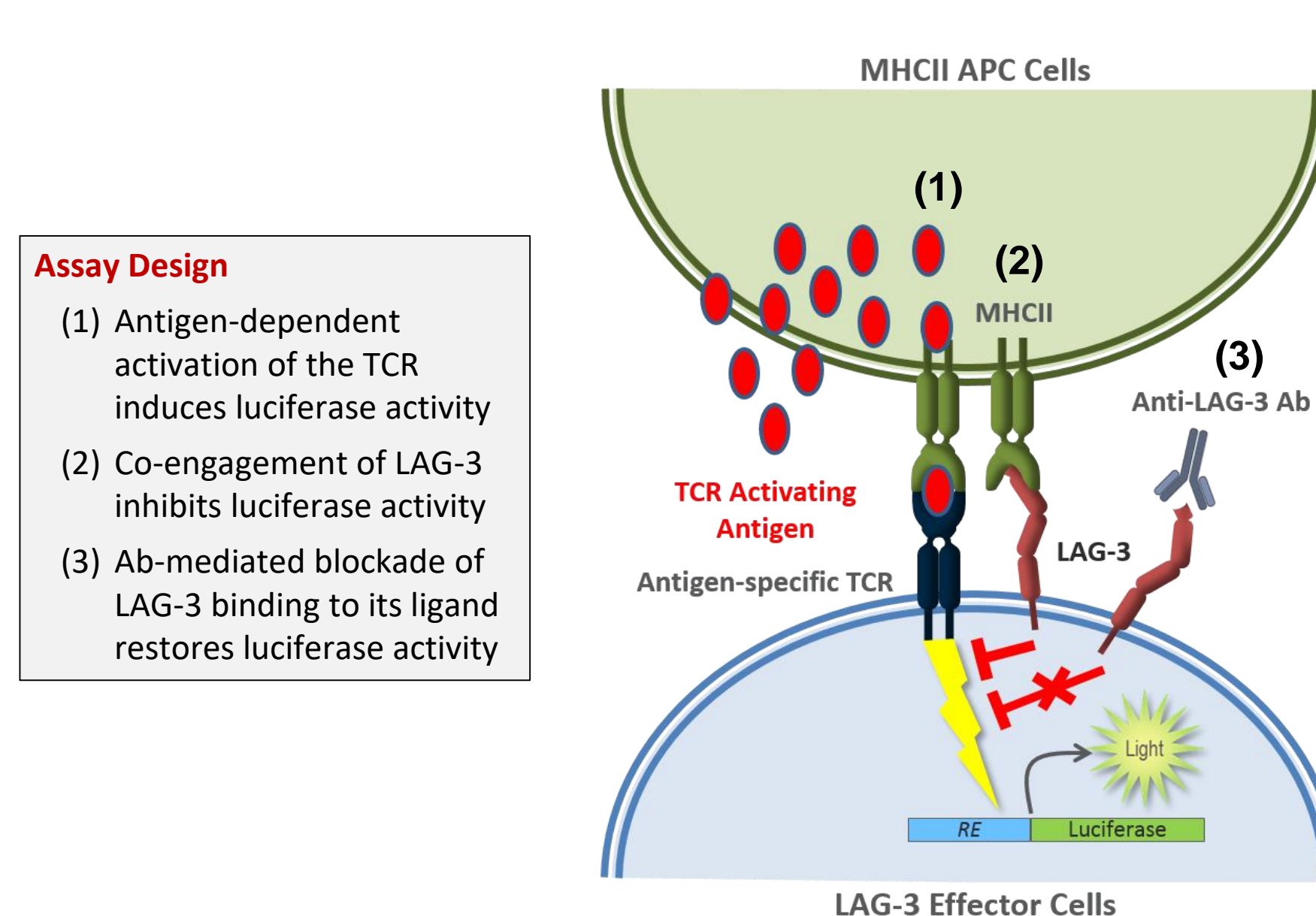
TCR and CD28-mediated luciferase activity is recovered in the CTLA-4 bioassay with an anti-CTLA-4 blocking Ab (ipilimumab), but not with anti-HER2 (trastuzumab), anti-PD-1 (nivolumab) or anti-TIGIT (clone MBSA-43) blocking Abs.

## 4. TIGIT/CD155 Blockade Bioassay



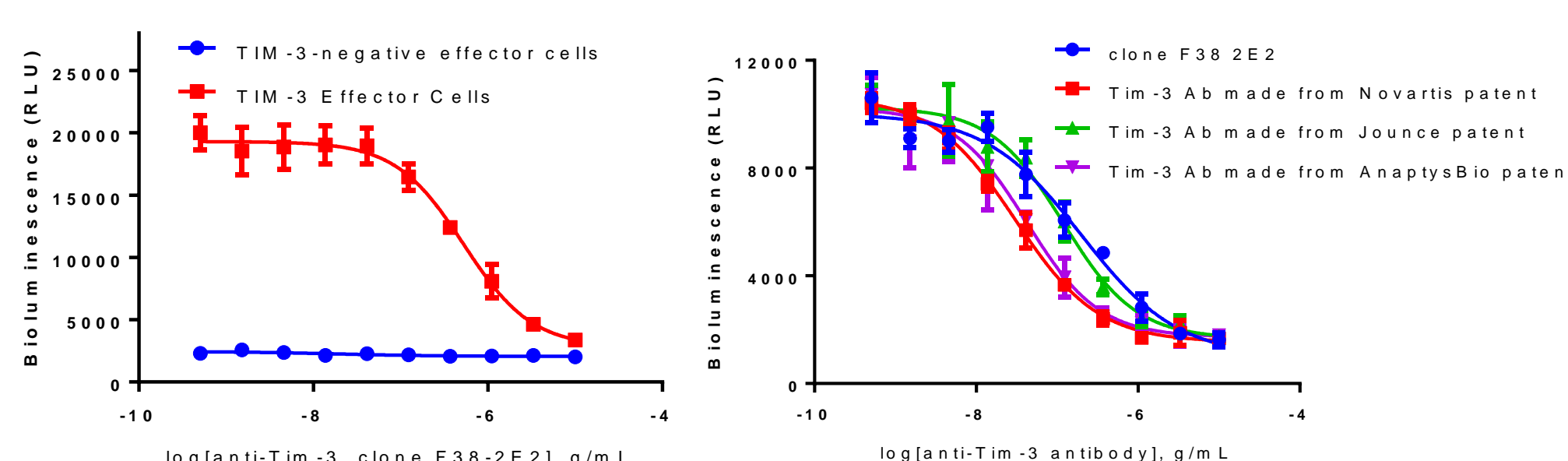
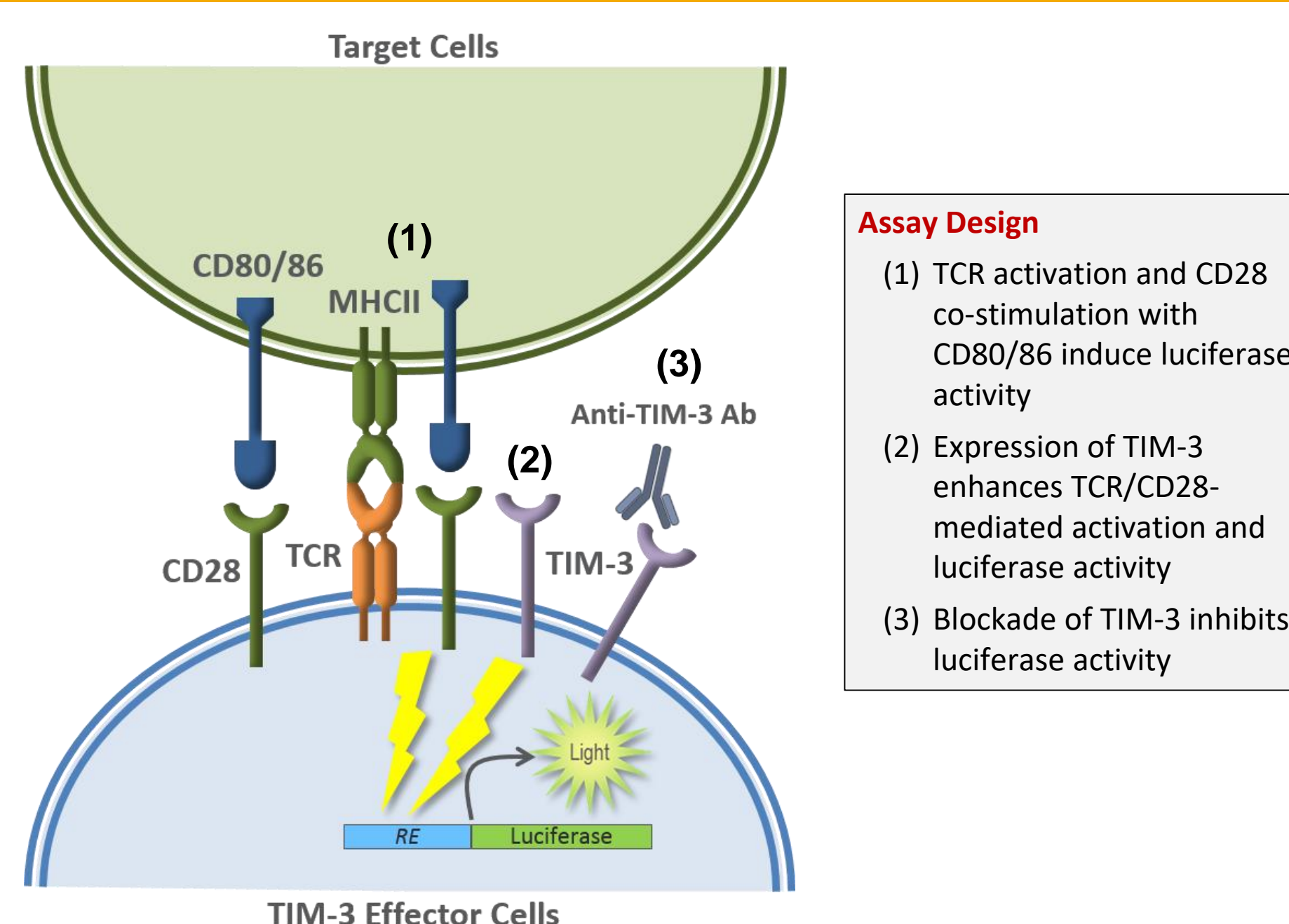
TCR- and CD226-mediated luciferase activity is recovered in the TIGIT/CD155 bioassay with an anti-TIGIT blocking Ab (clone MBSA-43), but not with an anti-PD-1 blocking Ab (nivolumab).

## 5. LAG-3/MHCII Blockade Bioassay



TCR-mediated luciferase activity is recovered in the LAG-3 bioassay with an anti-LAG-3 blocking Ab (clone 17B4).

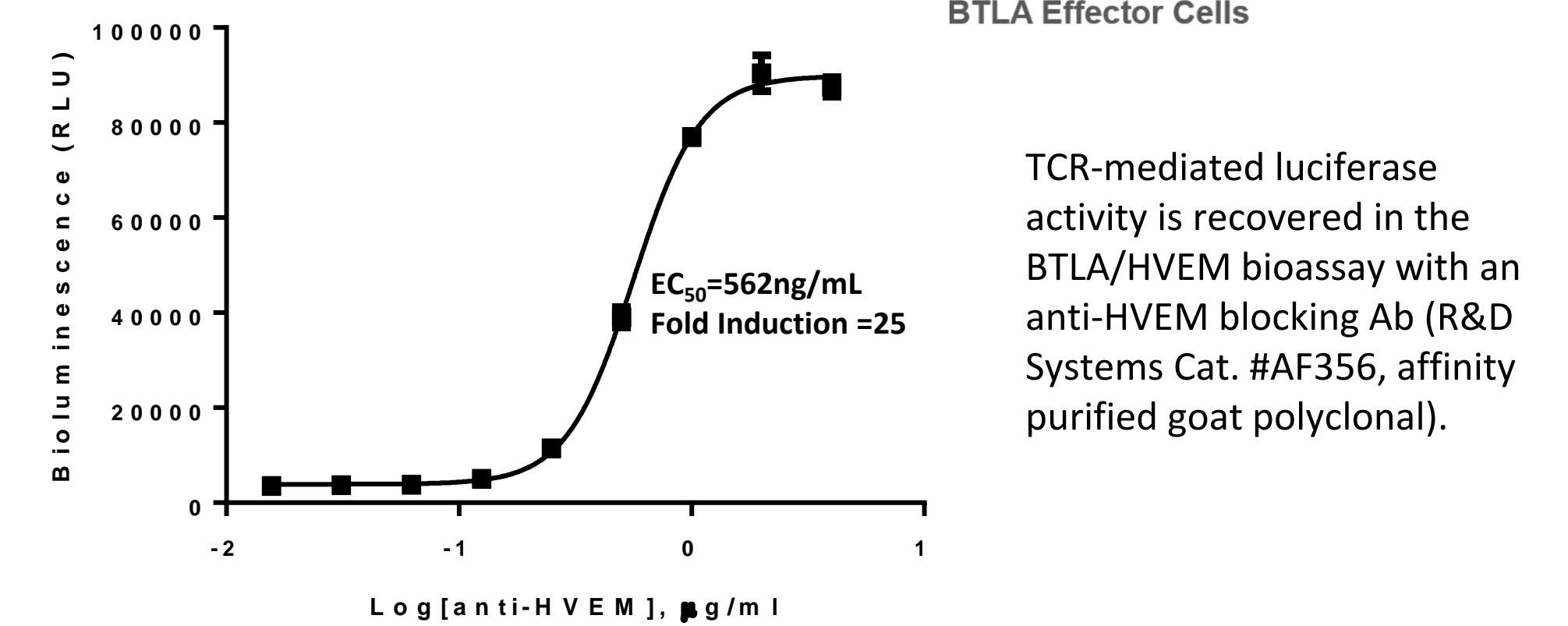
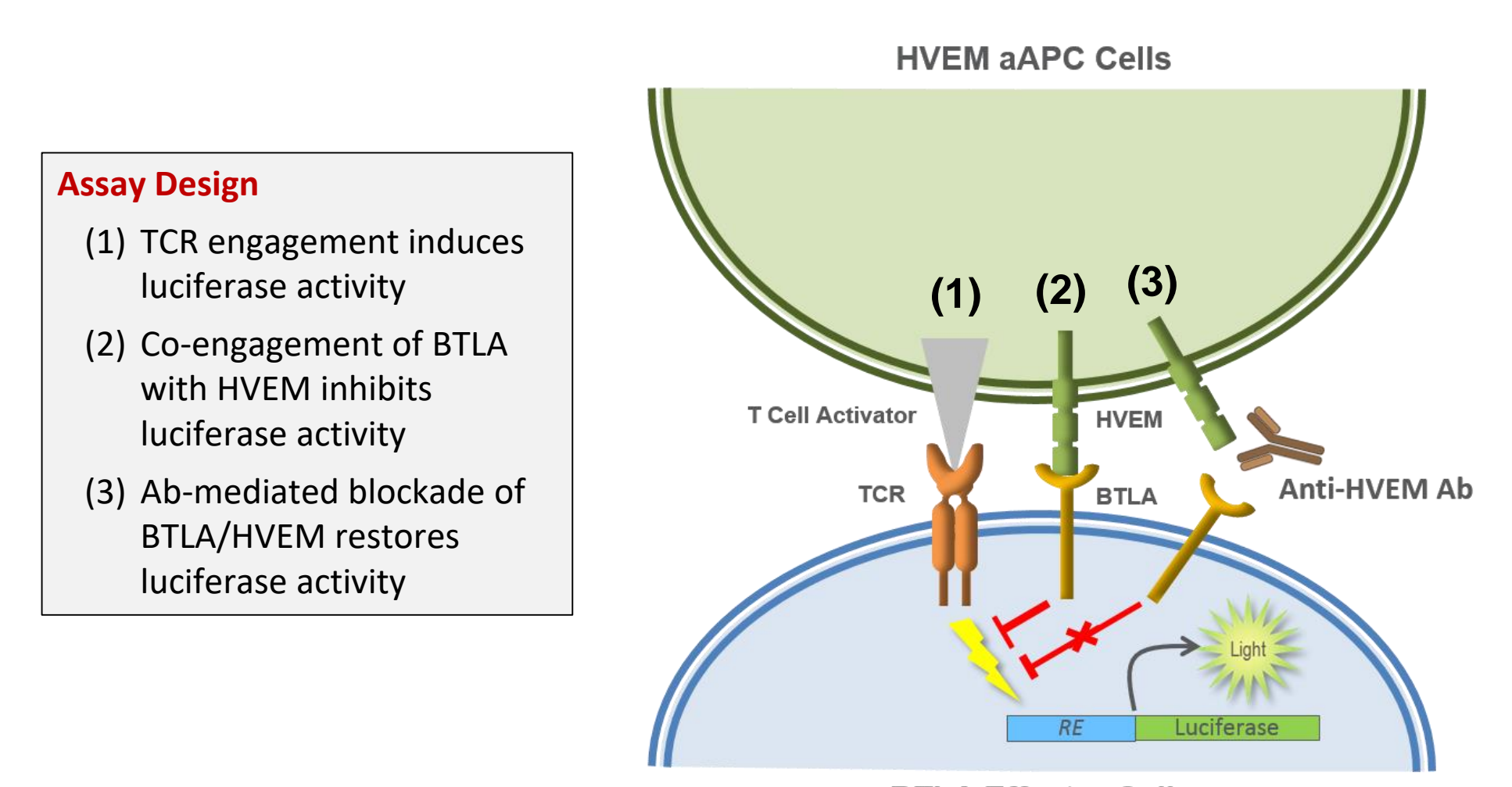
## 6. TIM-3 Bioassay



**Left:** TCR/CD28-mediated luciferase activity is enhanced with expression of TIM-3 in effector cells, which is inhibited in a dose-dependent manner with the anti-TIM-3 blocking Ab.

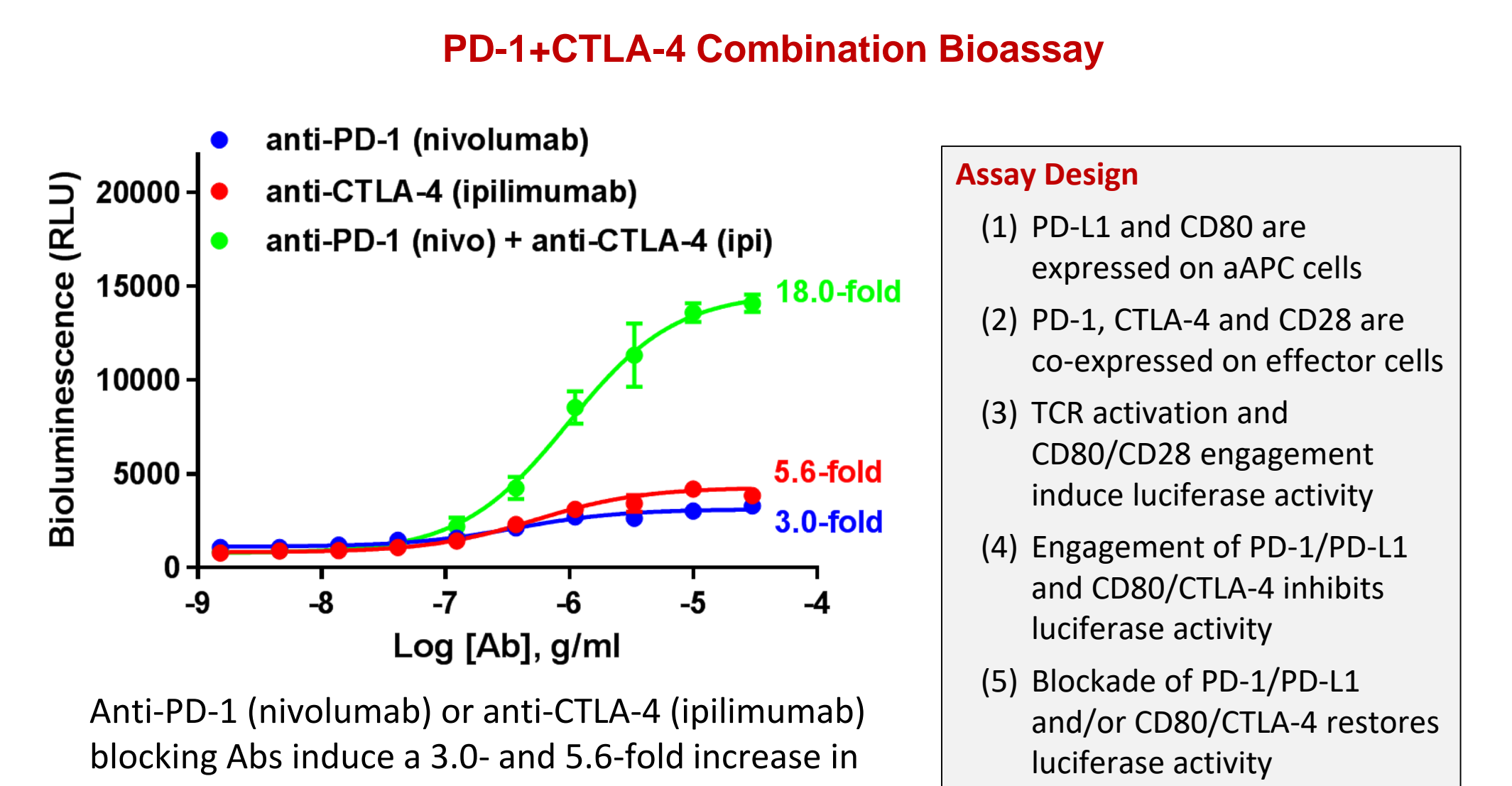
**Right:** TIM-3-mediated luciferase activity is inhibited in a dose-dependent manner with clone F38-2E2 and TIM-3 blocking antibodies made from Novartis, Jounce, and AnaptyBio.

## 7. BTLA/HVEM Blockade Bioassay

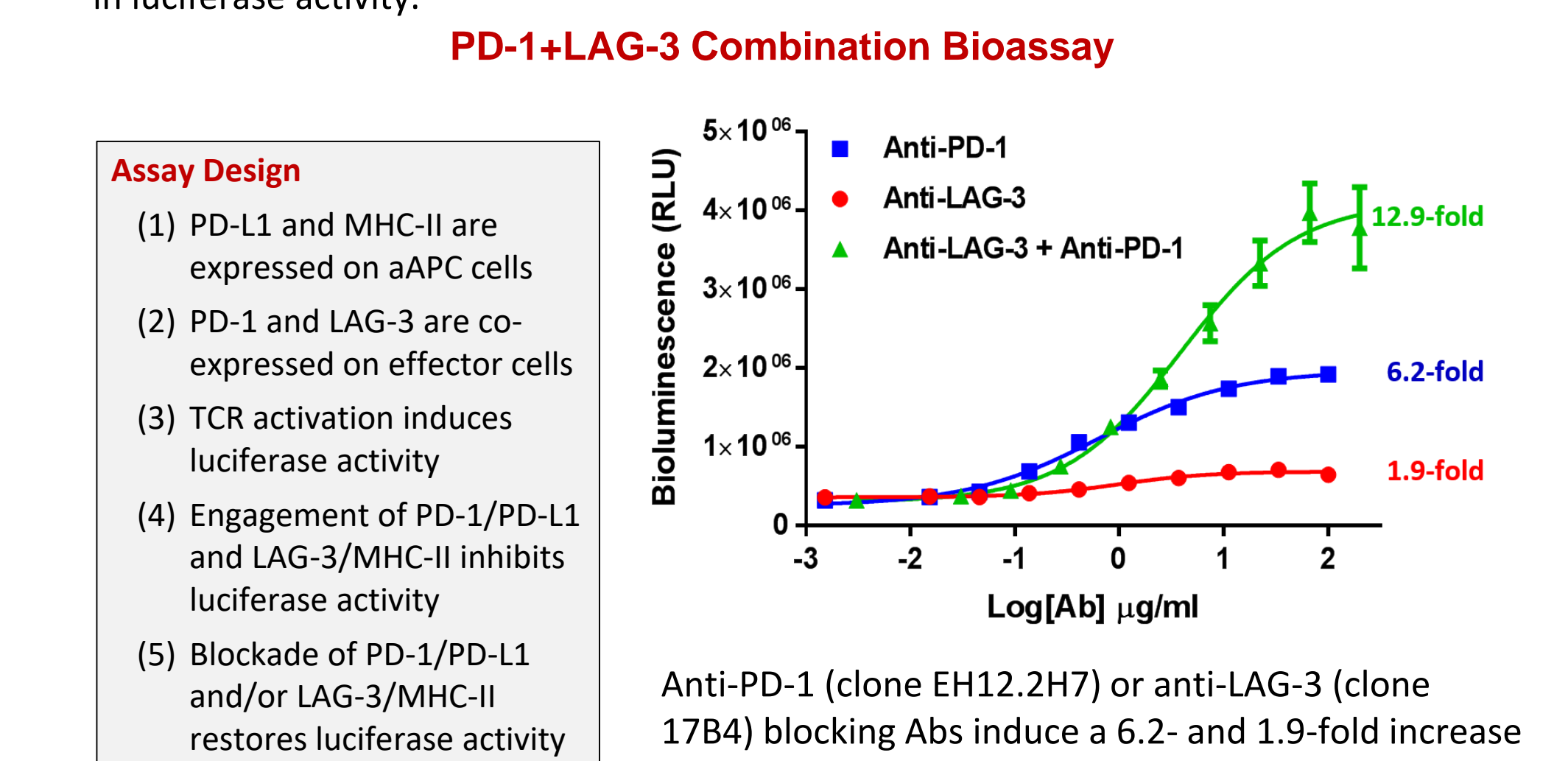


An HVEM/LIGHT co-stimulatory bioassay is also available (data not shown).

## 8. Combination Bioassays



Anti-PD-1 (nivolumab) or anti-CTLA-4 (ipilimumab) blocking Abs induce a 3.0- and 5.6-fold increase in luciferase activity when measured individually. A combination of both Abs induces an 18-fold increase in luciferase activity.



Anti-PD-1 (clone EH12.2H7) or anti-LAG-3 (clone 17B4) blocking Abs induce a 6.2- and 1.9-fold increase in luciferase activity, respectively. A combination of both Abs induces an 12.9-fold increase in activity.

Similar results were achieved using the PD-1+TIGIT Combination Bioassay (data not shown).

## 9. Conclusions

Cell-based reporter bioassays overcome the limitations of primary cell-based assays for functional characterization of antibody and other biologics drugs targeting individual or combination immune checkpoint receptors. Here we show a portfolio of MOA-based bioassays for co-inhibitory immune checkpoint receptors 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, stability, and NAb assays

### Easy-to-implement

- Rapid and convenient workflow
- Amenable to standard 96-well and 384-well plate formats