

Novel Cell-based Bioassays for Monoclonal Antibody and Bispecific Molecules in PD-1 Blockade Monotherapy and Combination Therapy

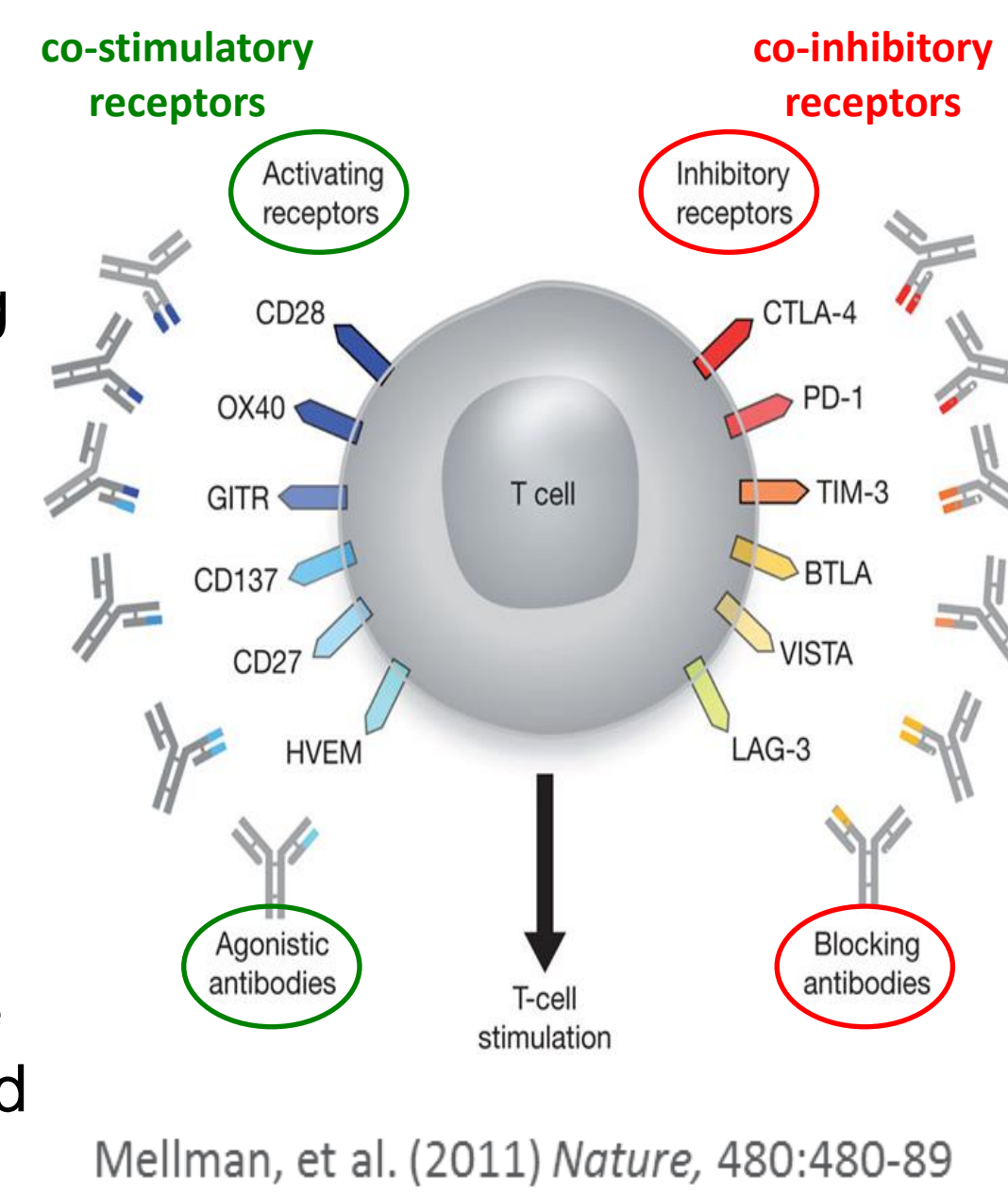
Julia Gilden, Jamison Grailer, Jun Wang, Denise Garvin, Pete Stecha, Jim Hartnett, Steven Edenson, Frank Fan, Mei Cong and Zhijie Jey Cheng

Promega Corporation, 2800 Woods Hollow Rd, Madison, WI 53711



1. Introduction

Immunotherapy targeting PD-1 has led to a paradigm shift in cancer drug discovery, due to its durable effect against a wide variety of cancers. Combining PD-1 checkpoint inhibitors with other clinically active treatments, including those targeting other immune checkpoint (IC) receptors, has also shown improved clinical results. Here, we report the development of a suite of cell-based reporter bioassays for monoclonal antibodies targeting PD-1/PD-L1, or bispecific molecules targeting PD-1/PD-L1 and a co-stimulatory receptor (e.g., 4-1BB, OX40) or an immune inhibitory receptor (e.g., CTLA-4, LAG-3, TIGIT) and show that the combination bioassays are able to measure the synergistic effect of PD-1 blockade with a second IC inhibitor receptor blockade or a costimulatory receptor activation.



Mellman, et al. (2011) *Nature*, 480:480-89

2. Assay Workflow of PD-1/PD-L1 Bioassay

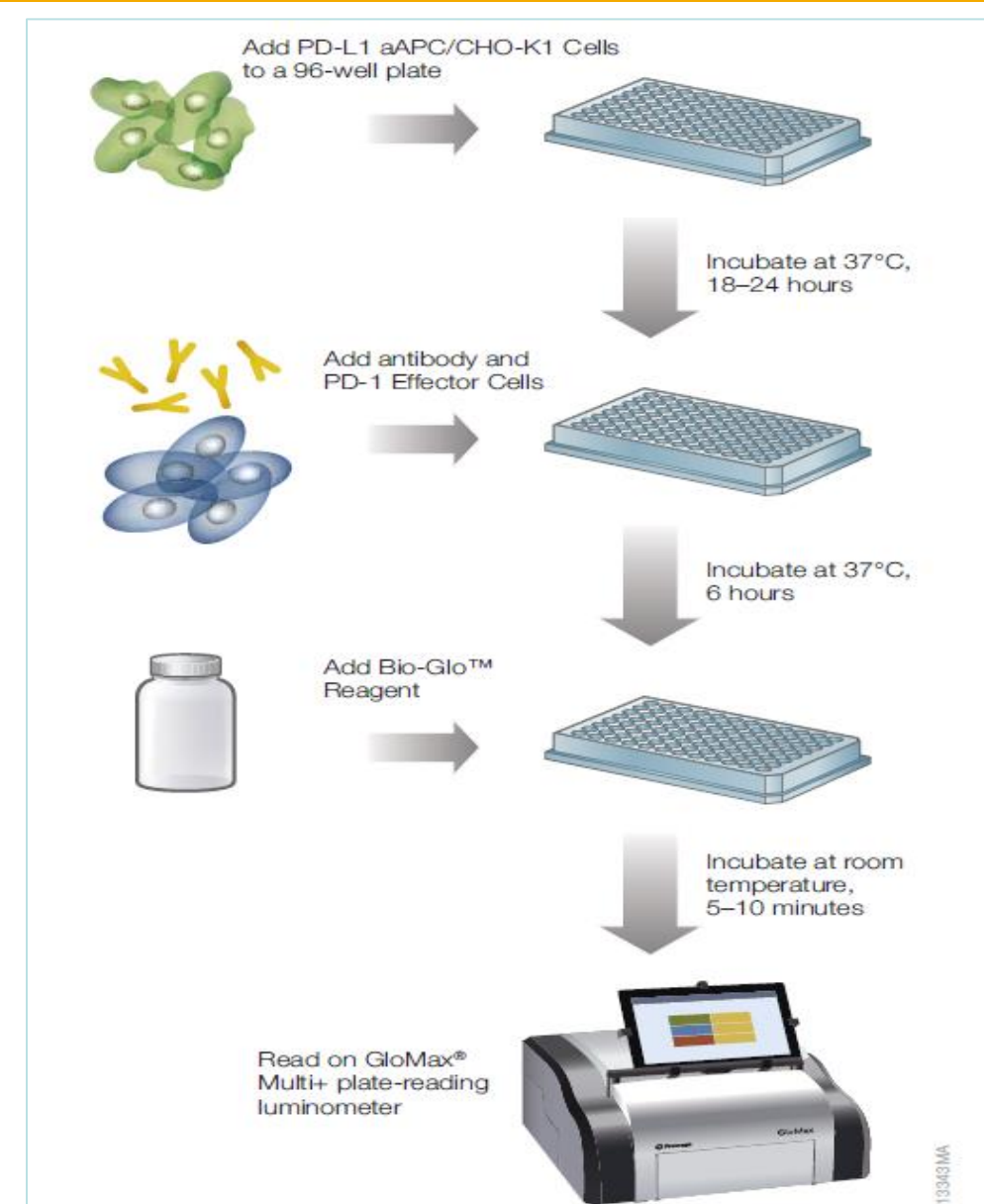
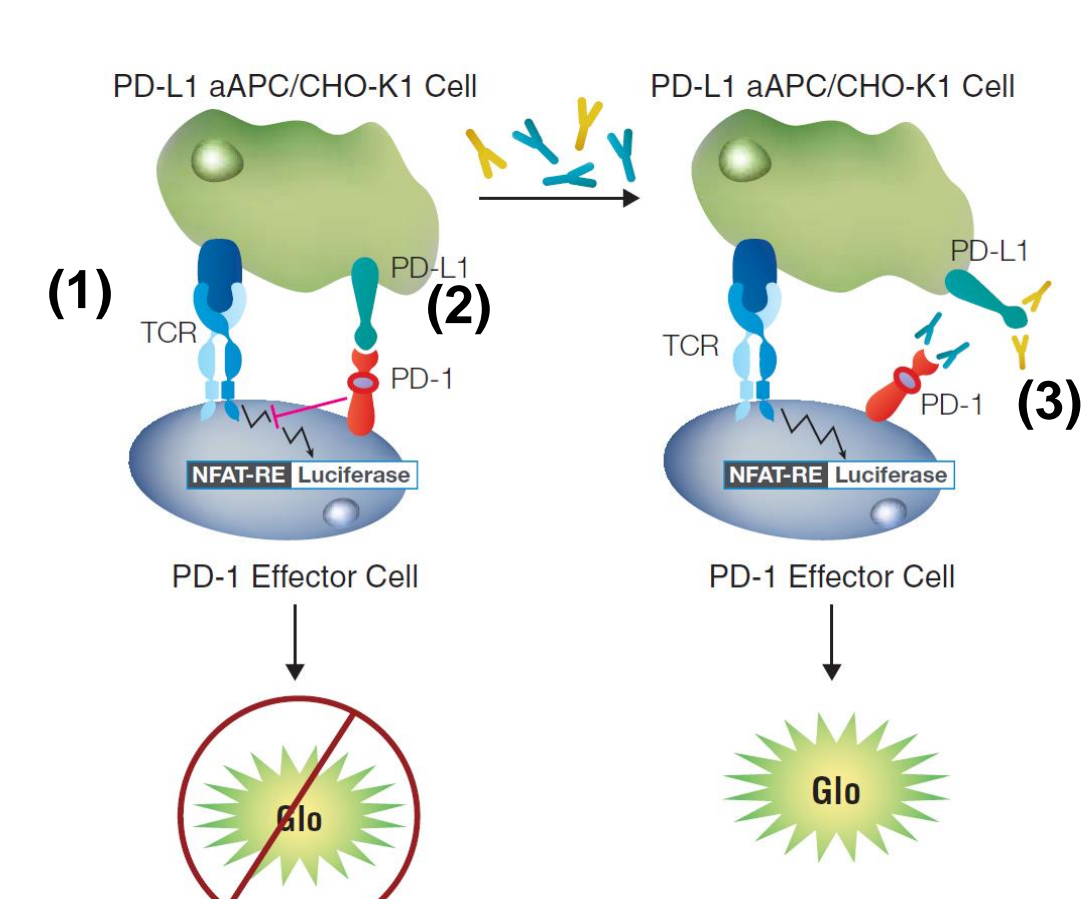
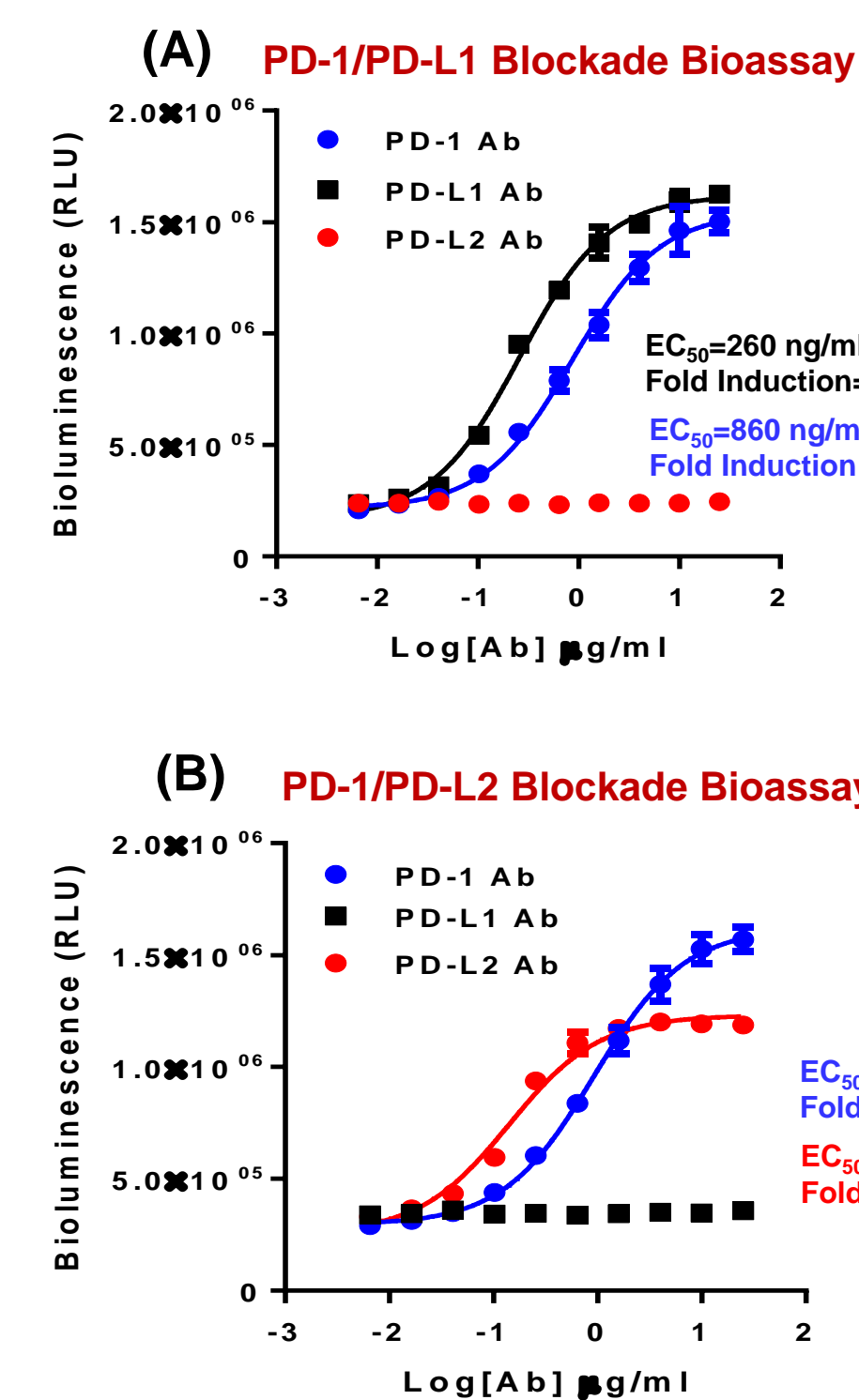


Plate PD-L1 aAPC cells
Add blocking antibody
Add PD-1 Effector cells
6-hour Induction
Add Bio-Glo™ detection reagent
Read plates

3. PD-1 Blockade Bioassays

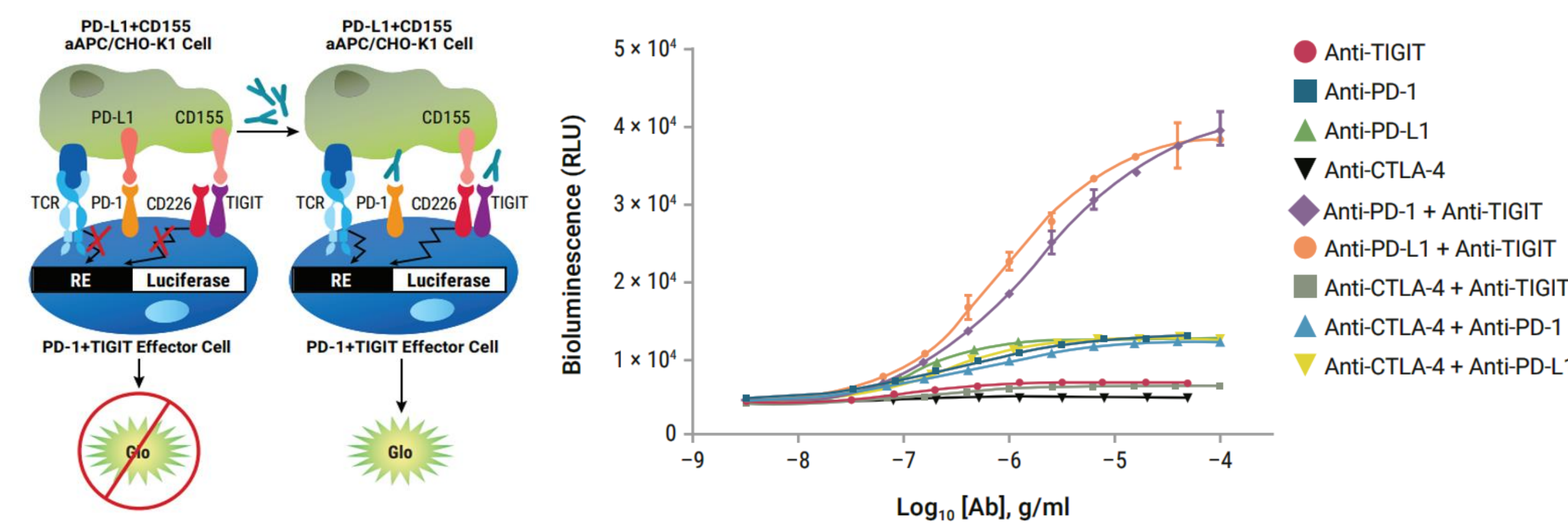


Assay Design
Luciferase activity indicating activation of T cells is:
(1) Induced by TCR engagement
(2) Inhibited by engagement of PD-1/PD-L1 or PD-1/PD-L2
(3) Restored by Ab-mediated blockade of PD-1/PD-L1 or PD-1/PD-L2



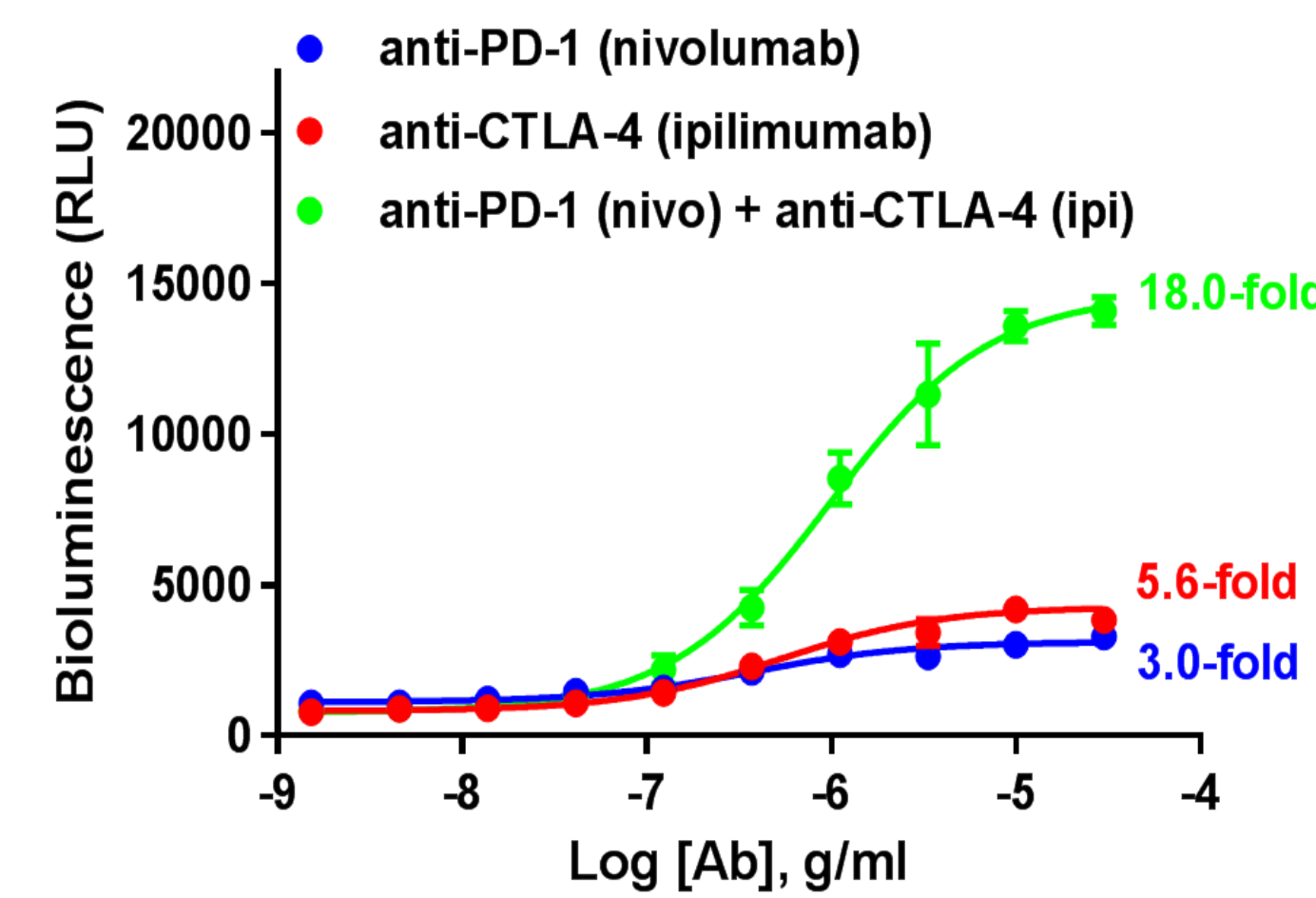
TCR-mediated luciferase activity is specially activated in PD-1/PD-L1 Bioassay (A) with anti-PD-1 or PD-L1 blocking Abs, and in PD-1/PD-L2 Bioassay (B) with anti-PD-1 or PD-L2 blocking Abs.

4. PD-1+TIGIT Combination Bioassay



Assay Design
PD-L1 and CD155 are co-expressed on aAPC cells. PD-1 and TIGIT are co-expressed on effector cells. Luciferase activity indicating activation of T cells is:
(1) Induced by engagement of TCR/CD226
(2) Inhibited by engagement of PD-1/PD-L1 and TIGIT/CD155
(3) Restored by mAb-mediated blockade of PD-1/PD-L1 and/or TIGIT/CD155.

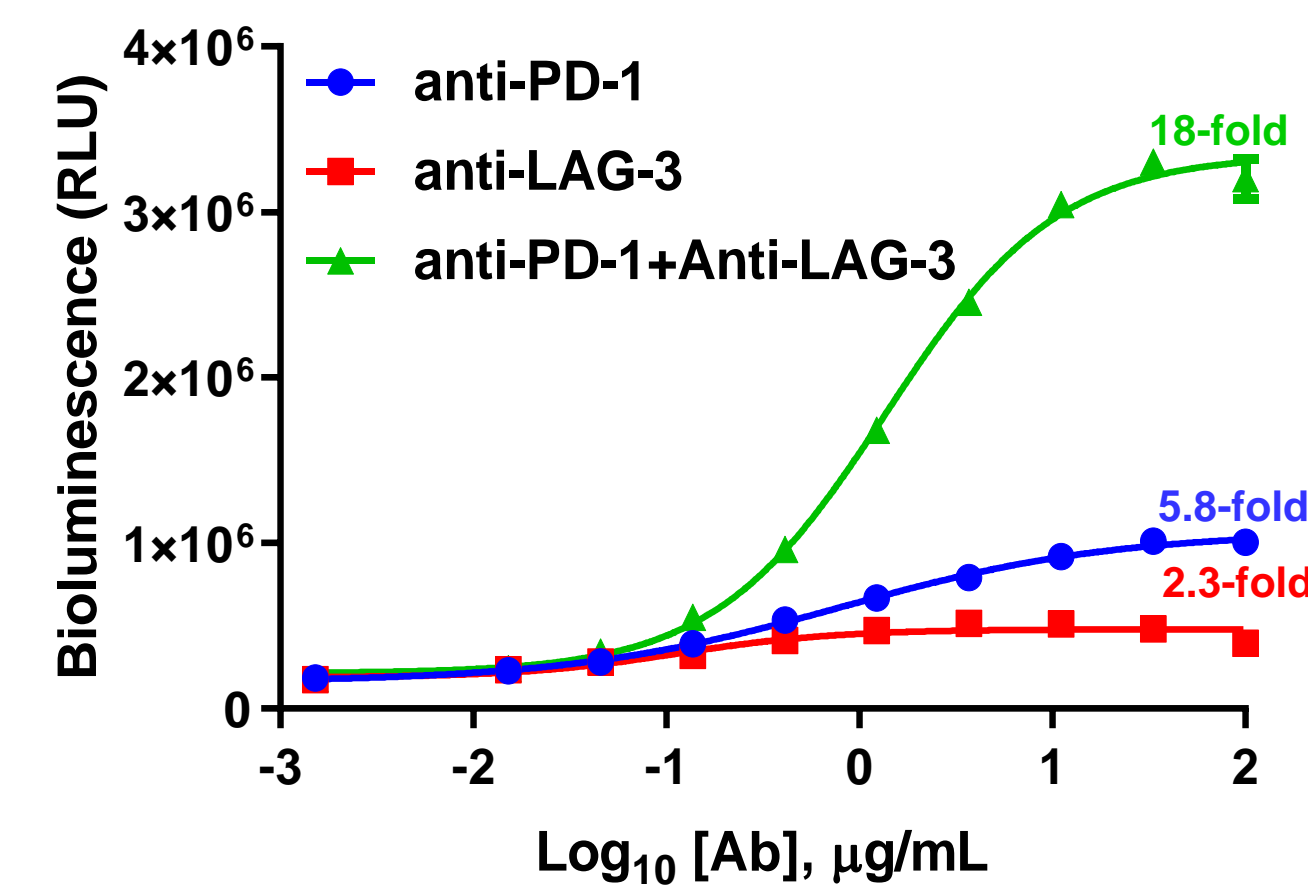
5. PD-1+CTLA-4 Combination Bioassay



Assay Design
PD-L1 and CD80 are co-expressed on aAPC cells. PD-1 and CTLA-4 are co-expressed on effector cells. Luciferase activity indicating activation of T cells is:
(1) Induced by engagement of TCR/CD28
(2) Inhibited by engagement of PD-1/PD-L1 and CTLA-4/CD80
(3) Restored by mAb-mediated blockade of PD-1/PD-L1 and/or CTLA-4/CD80

Anti-PD-1 or anti-CTLA-4 blocking Ab alone induced a 3.0- and 5.6-fold increase, while a combination of both Abs induced an 18-fold increase in luciferase activity.

6. PD-1+LAG-3 Combination Bioassay

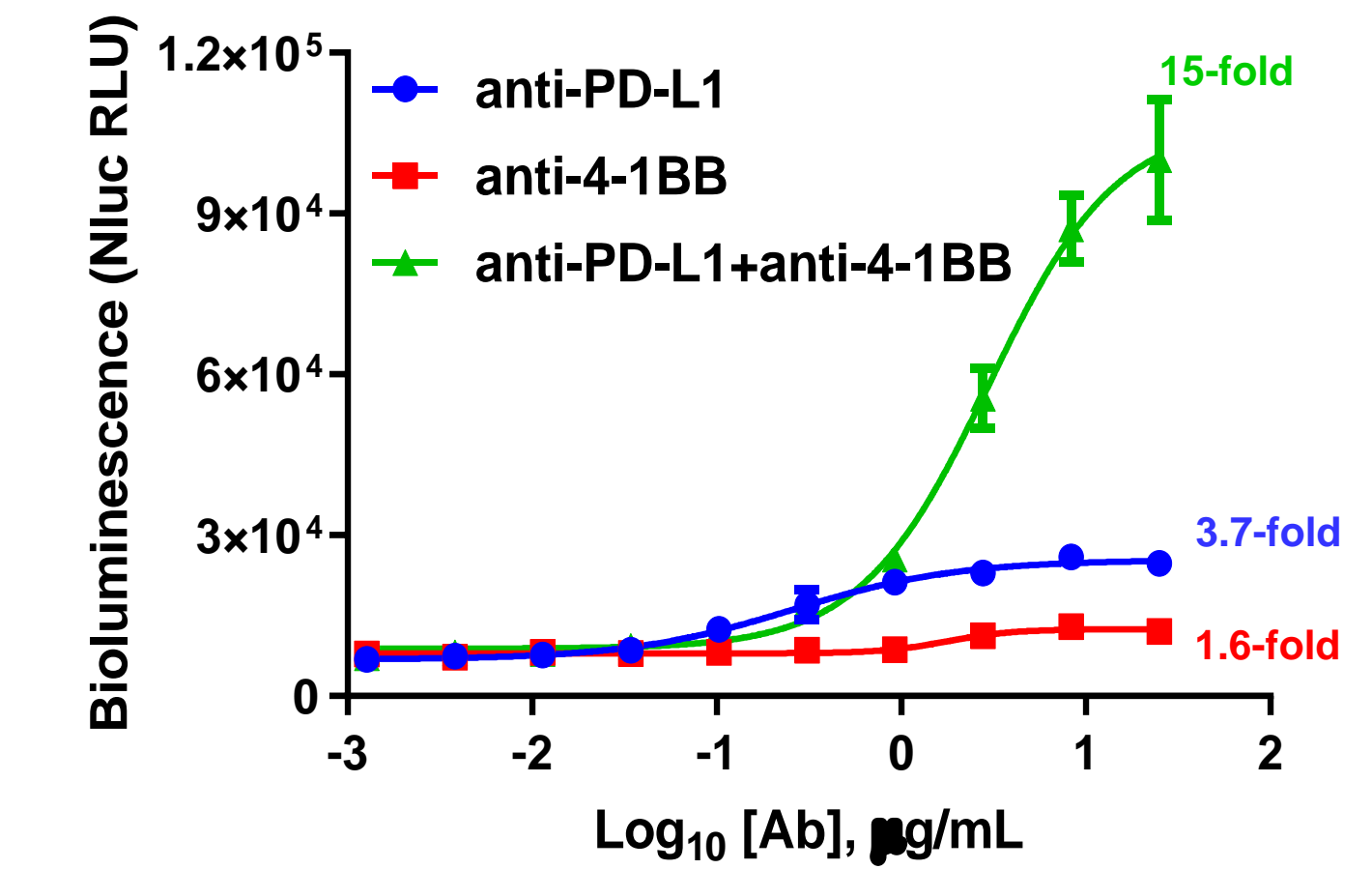


Assay Design
PD-L1 and MHC-II are co-expressed on aAPC cells. PD-1 and LAG-3 are co-expressed on effector cells. Luciferase activity indicating activation of T cells is:
(1) Induced by engagement of TCR
(2) Inhibited by engagement of PD-1/PD-L1 and LAG-3/MHC-II
(3) Restored by mAb-mediated blockade of PD-1/PD-L1 and/or LAG-3/MHC-II

Anti-PD-1 or anti-LAG-3 blocking Ab alone induced a 5.8- and 2.3-fold increase, while a combination of both Abs induced an 18-fold increase in luciferase activity.

Abstract# 1875

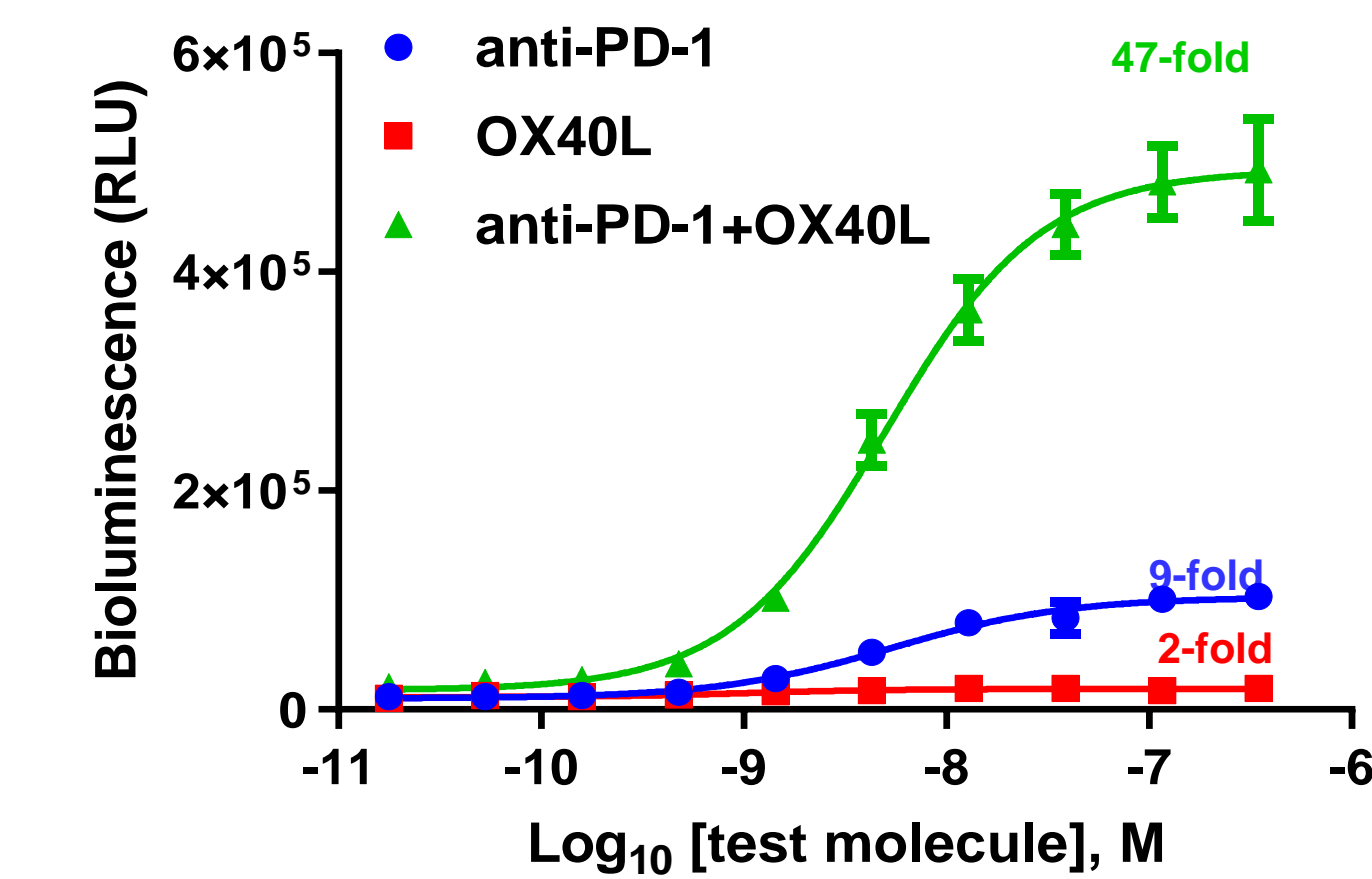
7. PD-1+4-1BB Combination Bioassay



Assay Design
PD-L1 is expressed on aAPC cells; PD-1 and 4-1BB are co-expressed on effector cells. Luciferase activity indicating activation of T cells is:
(1) Induced by activation of TCR
(2) Inhibited by engagement of PD-1/PD-L1
(3) Restored by Ab-mediated blockade of PD-1/PD-L1 and enhanced by 4-1BB agonist antibody

Anti-PD-L1 blocking Ab or anti-4-1BB agonist Ab alone induced a 3.7- and 1.6-fold increase, while a combination of both Abs induced a 15-fold increase in luciferase activity.

8. PD-1+OX40 Combination Bioassay



Assay Design
PD-L1 is expressed on aAPC cells and PD-1 and OX40 are co-expressed on effector cells. Luciferase activity indicating activation of T cells is:
(1) Induced by activation of TCR
(2) Inhibited by engagement of PD-1/PD-L1
(3) Restored by Ab-mediated blockade of PD-1/PD-L1 and enhanced by OX40 ligand

Anti-PD-1 blocking Ab or OX40L alone induced a 9- and 2-fold increase, while a combination of PD-1 Ab and OX40L induced a 47-fold increase in luciferase activity.

9. Conclusions

MOA-based reporter bioassays targeting PD-1 and a second immune checkpoint receptor overcome the limitations of primary cell-based assays and 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-mediated luciferase activity that reflects the native biology of T cell activation.
- Ability to measure the potency for immune checkpoint-targeted antibody alone or in combination.

Consistent and reliable measure of antibody activity

- Demonstrated precision, accuracy, reproducibility, robustness
- Functional performance suitable for development into potency, stability, and NAb assays

Easy-to-implement

- Rapid and convenient workflow
- All assays can be used as "Thaw-and-use" cell format, no cell culture required
- Amenable to standard 96-well and 384-well plate formats