

TECHNICAL MANUAL

# 4-1BB Bioassay

Instructions for Use of Products  
**JA2351 and JA2355**

**Note:** This Technical Manual includes a protocol for FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255) for use as needed.

# 4-1BB Bioassay

All technical literature is available at: [www.promega.com/protocols/](http://www.promega.com/protocols/)  
 Visit the web site to verify that you are using the most current version of this Technical Manual.  
 E-mail Promega Technical Services if you have questions on use of this system: [techserv@promega.com](mailto:techserv@promega.com)

1. Description .....	2
2. Product Components and Storage Conditions .....	8
3. Before You Begin .....	10
3.A. Materials to Be Supplied by the User .....	10
4. Assay Protocol for FcγRIIb-Dependent Antibodies.....	11
4.A. Preparing Assay Buffer, Bio-Glo™ Reagent and Antibody Samples .....	11
4.B. Plate Layout Design.....	13
4.C. Preparing and Plating FcγRIIb CHO-K1 Cells.....	13
4.D. Preparing Antibody Serial Dilutions.....	14
4.E. Preparing 4-1BB Effector Cells and Setting Up Assay.....	16
4.F. Adding Bio-Glo™ Reagent .....	16
4.G. Data Analysis .....	16
5. Assay Protocol for Ligand or FcγRIIb-Independent Antibodies.....	17
5.A. Preparing Assay Buffer, Bio-Glo™ Reagent and Antibody Samples .....	17
5.B. Plate Layout Design.....	19
5.C. Preparing Ligand or Antibody Serial Dilutions .....	19
5.D. Preparing 4-1BB Effector Cells and Setting Up Assay.....	21
5.E. Adding Bio-Glo™ Reagent .....	21
5.F. Data Analysis .....	22
6. Troubleshooting .....	22
7. References .....	23
8. Appendix.....	24
8.A. Representative Assay Results with FcγRIIb-Dependent Antibody.....	24
8.B. Representative Assay Results with Crosslinked Ligand or FcγRIIb-Independent Antibody .....	25
8.C. Related Products.....	26
9. Summary of Changes .....	29

## 1. Description

The human immune system is regulated by a complex network of inhibitory and stimulatory receptors that facilitate the elimination of pathogens, while maintaining tolerance to self-antigens. Stimulatory immune checkpoint receptors appear to have a significant role in cancer progression and autoimmune disease. Several co-stimulatory immune checkpoint receptors such as Glucocorticoid-induced TNF receptor family-related protein (GITR), 4-1BB, OX40, CD40 and Inducible T-cell Costimulator (ICOS) have been identified. Activating these receptors with ligands or agonist antibodies has emerged to be the next generation of immunotherapeutic strategy to enhance anti-tumor immune responses and promote immune-mediated tumor rejection (1–3).

4-1BB (CD137/TNFRSF9), a member of the tumor necrosis factor receptor superfamily, is an inducible co-stimulatory receptor expressed on T cells, natural killer (NK) cells and innate immune cell populations (4). When present on the cell surface, 4-1BB interacts with 4-1BB ligand (4-1BBL) and induces subsequent cell proliferation and production of interferon gamma (IFN $\gamma$ ) and IL-2, particularly in T and NK cells (5,6).

Current methods used to measure the activity of biologic drugs targeting 4-1BB rely on primary human T cells and measurement of functional endpoints such as cell proliferation, cell surface marker expression and IFN $\gamma$  and interleukin-2 (IL-2) production. These assays are laborious and highly variable due to their reliance on donor primary cells, complex assay protocols and nonqualified assay reagents. As a result, these assays are difficult to establish in a quality-controlled, drug-development setting.

The 4-1BB Bioassay<sup>(a-e)</sup> (Cat.# JA2351, JA2355), is a bioluminescent cell-based assay that overcomes the limitations of existing assays and can be used to measure the potency and stability of ligands or agonist antibodies that can bind and activate 4-1BB (6,7). The assay consists of a genetically engineered Jurkat T cell line that expresses human 4-1BB and a luciferase reporter driven by a response element that can respond to 4-1BB ligand/agonist antibody stimulation. The 4-1BB Effector Cells<sup>(a,c,d)</sup> are provided in thaw-and-use format as cryopreserved cells that can be thawed, plated and used in an assay without the need for cell culture and propagation.

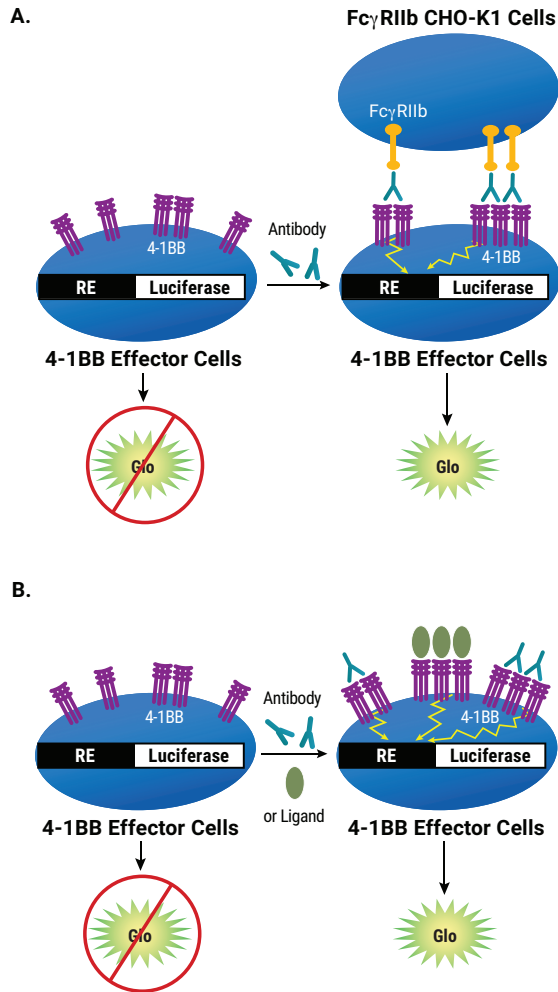
The antibodies used to stimulate T cell activation via interaction with co-stimulatory receptors on the effector cells, can be crosslinked via binding to Fc $\gamma$ RIIb receptors. This crosslinking increases avidity, stabilizes the interaction of the antibody to the receptor, enhances receptor clustering and downstream signaling (8,9).

The 4-1BB Bioassay can be conducted with FcγRIIb CHO-K1 Cells<sup>(a,c,d)</sup> (Cat.# JA2251, JA2255) to test whether agonist antibodies activate 4-1BB in an FcγRIIb-dependent manner. FcγRIIb CHO-K1 Cells may be required to crosslink agonist antibodies but are not required for testing ligands. It is recommended that, when screening for agonist antibodies of co-stimulatory immune checkpoints, you perform the assay both with and without FcγRIIb CHO-K1 Cells to ascertain the need for these cells in enhancing the effect of the agonist antibodies raised against the co-stimulatory immune checkpoint targets.

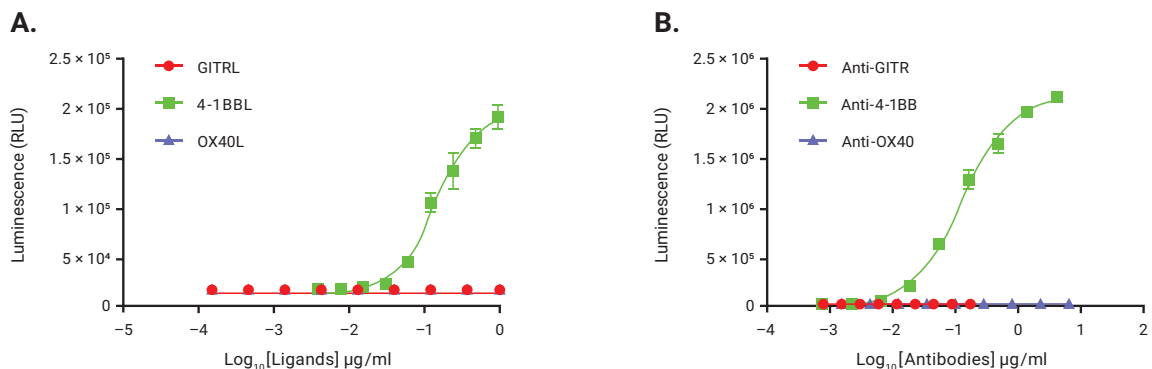
Induction of the 4-1BB Effector Cells with a 4-1BB ligand or agonist antibody results in response element-mediated luminescence (Figure 1). The bioluminescent signal is quantified using the Bio-Glo™ Luciferase Assay System and a standard luminometer such as the GloMax® Discover System (see Section 8.C, Related Products).

The 4-1BB Bioassay reflects the mechanism of action (MOA) of biologics designed to activate 4-1BB. Specifically, 4-1BB-mediated luminescence is detected following the incubation of 4-1BB agonist antibodies or 4-1BB ligand (Figure 2). The bioassay is prequalified according to ICH guidelines and shows the precision, accuracy and linearity required for routine use in potency and stability studies (Table 1 and Figure 3). The assay can be performed in a one-day or two-day time frame depending on antibody properties. The bioassay workflow is simple and robust, and is compatible with both 96-well and 384-well plate formats used for antibody screening in early drug discovery (Figure 4). In addition, the bioassay can be used with up to 100% human serum (in ligand or antibody samples) with minimal impact on fold induction (Figure 5), indicating potential for further development into a neutralizing antibody bioassay.

1. Description (continued)



**Figure 1. Representation of the 4-1BB Bioassay. Panel A.** Assay with Fc $\gamma$ RIIb-dependent agonist antibody. The bioassay consists of two engineered cell lines, 4-1BB Effector Cells and Fc $\gamma$ RIIb CHO-K1 Cells. In the presence of Fc $\gamma$ RIIb CHO-K1 Cells, the anti-4-1BB antibody can be crosslinked, thereby inducing 4-1BB pathway-activated luminescence. **Panel B.** Assay with Fc $\gamma$ RIIb-independent agonist antibody or ligand. The bioassay consists of one engineered cell line, 4-1BB Effector Cells. In the absence of agonist antibody or 4-1BB ligand, the 4-1BB receptor is not activated and luminescence signal is low. The addition of agonist antibody or 4-1BB ligand induces the 4-1BB pathway-activated luminescence, which can be detected in a dose-dependent manner.

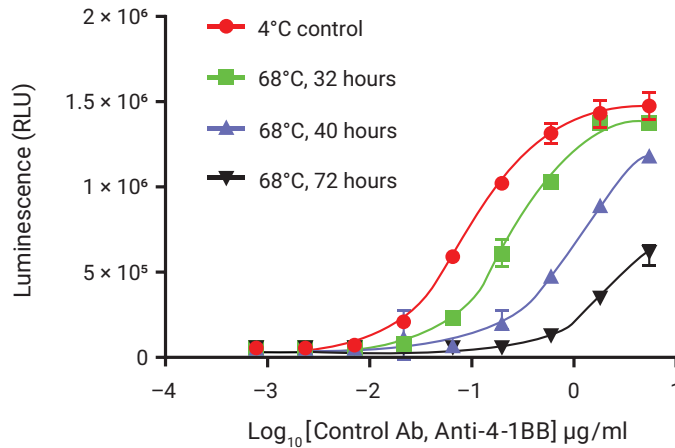


**Figure 2. The 4-1BB Bioassay reflects the mechanism of action (MOA) and shows specificity of biologics designed to activate 4-1BB.** **Panel A.** 4-1BB Effector Cells were induced, respectively, with a serial titration of ligands: GITRL, 4-1BBL and OX40L as indicated. **Panel B.** 4-1BB Effector Cells were induced with a serial titration of anti-GITR antibody, Control Ab, Anti-4-1BB (Cat. # K1161) and anti-OX40 antibody, as indicated, in the presence of FcγRIIb CHO-K1 Cells (Cat. # JA2251, JA2255). After a 6-hour incubation, Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were generated using thaw-and-use cells and fitted to a 4-parameter logistic curve using GraphPad Prism® software.

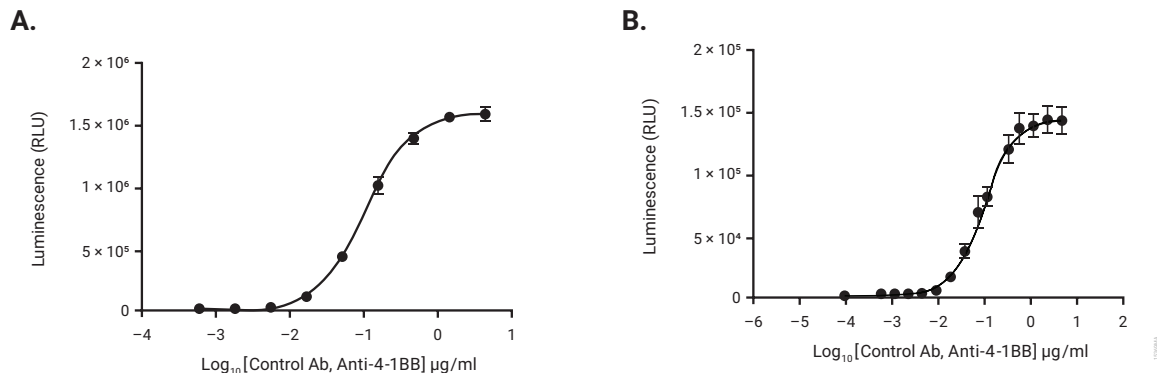
**Table 1. The 4-1BB Bioassay Shows Precision, Accuracy and Linearity.**

Parameter	Results	
	% Expected Relative Potency	% Recovery
Accuracy	50	106.0
	70	96.2
	140	97.3
	200	103.2
Repeatability (% CV)	100% (Reference)	3.5
Intermediate Precision (% CV)		6.6
Linearity (r <sup>2</sup> )		0.996
Linearity (y = mx + b)		y = 1.025x - 1.835
<p>A 50–200% theoretical potency series of recombinant human IL-15 was analyzed in triplicate in three independent experiments performed on three days by two analysts (for a total of six independent experiments). Bio-Glo™ Reagent was added and luminescence quantified using the GloMax® Discover System. Data were analyzed and relative potencies calculated after parallelism determination using JMP® software. Data were generated using thaw-and-use cells.</p>		

1. Description (continued)



**Figure 3. The 4-1BB Bioassay is stability-indicating.** Samples of Control Ab, Anti-4-1BB (Cat.# K1161), were maintained at 4°C (control) or heat-treated at 68°C for indicated times, and then analyzed using the 4-1BB Bioassay with FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255). After a 6-hour incubation, Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were generated using thaw-and-use cells and fitted to a 4-parameter logistic curve using GraphPad Prism® software.



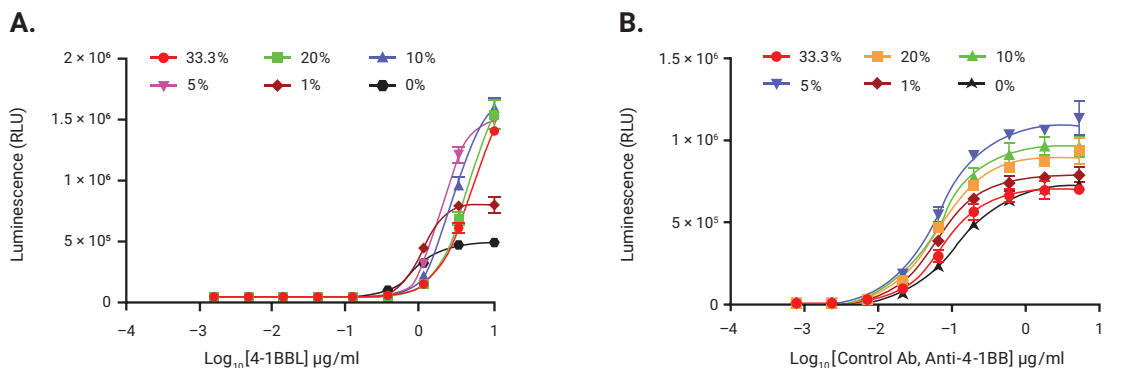
**Figure 4. The 4-1BB Bioassay is amenable to 384-well plate format and compatible with laboratory automation. Panel A.**

The 4-1BB Bioassay was performed in 96-well plates as described in this technical manual using Control Ab, Anti-4-1BB.

**Panel B.** The 4-1BB Bioassay was performed in 384-well format using a Mantis<sup>®</sup> liquid handler to dispense the cells and Echo<sup>®</sup> Acoustic liquid handler for antibody handling. On the day before assay, FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255) were plated at 8 × 10<sup>3</sup> cells/10µl/well. On the day of the assay, Control Ab, Anti-4-1BB (Cat.# K11161), was serially diluted and added to the plate at 0.2µl/well. Finally, 4-1BB effector cells were added at 1.0 × 10<sup>4</sup> cells/10µl/well. After a 6-hour incubation, 20µl of Bio-Glo<sup>™</sup> Reagent was added, and luminescence was quantified using the GloMax<sup>®</sup> Discover System. Data were fitted to a 4-parameter logistic curve using GraphPad Prism<sup>®</sup> software. The EC<sub>50</sub> values were 0.13µg/ml and 0.095µg/ml, and the fold inductions were 103 and 58 for 96-well and 384-well format, respectively. Data were generated using thaw-and-use cells.



## 1. Description (continued)



**Figure 5. The 4-1BB Bioassay is tolerant to human serum. Panel A.** 4-1BB ligand (4-1BBL, R&D Systems Cat. #2295-4L-025/CF) was analyzed in the presence of increasing concentrations of pooled normal human serum (0–100% in the ligand sample), resulting in final assay concentration of human serum (0–33.3%). **Panel B.** Control Ab, Anti-4-1BB was analyzed in the presence of Fc $\gamma$ RIIb CHO-K1 Cells (Cat. # JA2251, JA2255) and increasing concentrations of pooled normal human serum (0–100% in the antibody sample), resulting in final assay concentration of human serum (0–33.3%). After a 6-hour incubation, Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were fitted to a 4-parameter logistic curve using GraphPad Prism® software. Data were generated using thaw-and-use cells. The 4-1BB Bioassay is tolerant to serum with this human serum pool. A different human serum pool showed similar effects on the assay (data not shown).

## 2. Product Components and Storage Conditions

PRODUCT	SIZE	CAT. #
4-1BB Bioassay	1 each	JA2351

Not for Medical Diagnostic Use.

Each kit contains sufficient reagents for 120 assays using the inner 60 wells of two 96-well plates. Includes:

- 1 vial 4-1BB Effector Cells (0.5ml)
- 36ml RPMI 1640 Medium
- 4ml Fetal Bovine Serum
- 1 vial Bio-Glo™ Luciferase Assay System (lyophilized)
- 10ml Bio-Glo™ Luciferase Assay Buffer

## 2. Product Components and Storage Conditions (continued)

PRODUCT	SIZE	CAT.#
<b>4-1BB Bioassay 5X</b>	<b>5 each</b>	<b>JA2355</b>

Not for Medical Diagnostic Use.

Each kit contains sufficient reagents for 600 assays using the inner 60 wells of two 96-well plates. Includes:

- 5 vials 4-1BB Effector Cells (0.5ml per vial)
- 5 × 36ml RPMI 1640 Medium
- 5 × 4ml Fetal Bovine Serum
- 5 vials Bio-Glo™ Luciferase Assay System (lyophilized)
- 5 × 10ml Bio-Glo™ Luciferase Assay Buffer

### Items Available Separately

PRODUCT	SIZE	CAT.#
<b>Fc<math>\gamma</math>R1Ib CHO-K1 Cells</b>	<b>1 each</b>	<b>JA2251</b>

Not for Medical Diagnostic Use.

Each kit contains sufficient reagents for 120 assays using the inner 60 wells of two 96-well plates. Includes:

- 1 vial Fc $\gamma$ R1Ib Cells (0.5ml)
- 36ml RPMI 1640 Medium

PRODUCT	SIZE	CAT.#
<b>Fc<math>\gamma</math>R1Ib CHO-K1 Cells</b>	<b>5 each</b>	<b>JA2255</b>

Not for Medical Diagnostic Use.

Each kit contains sufficient reagents for 600 assays using the inner 60 wells of two 96-well plates. Includes:

- 5 vials Fc $\gamma$ R1Ib Cells (0.5ml per vial)
- 5 × 36ml RPMI 1640 Medium

**Note:** The 4-1BB Bioassay components are shipped separately because of different temperature requirements. The 4-1BB Effector Cells are shipped on dry ice. The Bio-Glo™ Luciferase Assay Substrate and Buffer and Fetal Bovine Serum are shipped on dry ice, separately from the cells. The RPMI 1640 Medium is shipped at ambient temperature.

**Storage Conditions:** Upon arrival, immediately transfer the cell vials to below  $-140^{\circ}\text{C}$  (freezer or liquid nitrogen vapor phase) for long-term storage. Do not store cell vials submerged in liquid nitrogen. **Do not** store cell vials at  $-80^{\circ}\text{C}$  because this will decrease cell viability and cell performance.

Store Bio-Glo™ Luciferase Assay Substrate, Bio-Glo™ Luciferase Assay Buffer and Fetal Bovine Serum at  $-30^{\circ}\text{C}$  to  $-10^{\circ}\text{C}$ . Avoid multiple freeze-thaw cycles of the serum.

For optimal performance, use reconstituted Bio-Glo™ Reagent on the day of preparation. However, once reconstituted, Bio-Glo™ Reagent can be stored at  $-30^{\circ}\text{C}$  to  $-10^{\circ}\text{C}$  for up to 6 weeks.

Store RPMI 1640 Medium at  $+2^{\circ}\text{C}$  to  $+10^{\circ}\text{C}$  protected from fluorescent light.

### 3. Before You Begin

**Please read through the entire protocol to become familiar with the components and the assay procedure before beginning.**

Remove the product label from the box containing vials with cells or note the catalog number and kit lot number from the label. This information can be used to download documents for the specified product from the website such as Certificate of Analysis.

The 4-1BB Bioassay is intended to be used with user-provided antibodies or ligands designed to activate 4-1BB. Control Ab, Anti-4-1BB (Cat.# K1161) and FcγRIIb CHO-K1 Cells Cat.# JA2251, JA2255) are available separately for use in assay optimization and routine quality control. We strongly recommend including 4-1BB Ligand or Control Ab, Anti-4-1BB, along with FcγRIIb CHO-K1 Cells as a positive control in the first few assays to gain familiarity with the assay. Data generated using these reagents are shown in Figures 2–5 and Sections 8.A and B, Representative Assay Results.

The 4-1BB Effector Cells are provided in frozen, thaw-and-use format and are ready to be used without any additional cell culture or propagation. When thawed and diluted as instructed, the cells will be at the appropriated concentration for the assay. The cells are sensitive, and care should be taken to follow cell thawing and plating procedures as described in Section 4.E.

The 4-1BB Bioassay produces a bioluminescent signal and should work with all major luminometers or luminescence plate readers for the detection of luminescence. Bioassay development and performance data included in this Technical Manual were generated using the GloMax<sup>®</sup> Discover System (see Section 8.C, Related Products). An integration time of 0.5 second/well was used for all readings.

#### 3.A. Materials to Be Supplied by the User

- user-defined anti-4-1BB antibodies or other biologics samples
- solid-white, flat-bottom 96-well assay plates (e.g., Corning<sup>®</sup> Cat.# 3917) for plating and reading luminescence
- sterile clear V-bottom 96-well plate with lid (e.g., Corning<sup>®</sup> Cat.# 3896 or Linbro Cat.# 76-223-05) for preparing antibody dilutions
- pipettes (single-channel and 12-channel; for best results use both manual and electronic pipettes as needed)
- sterile 15ml and 50ml conical tubes
- sterile reagent reservoirs (e.g., Corning<sup>®</sup> Cat.# 4870)
- 37°C, 5% CO<sub>2</sub> incubator
- 37°C water bath
- plate reader with glow luminescence measuring capability or luminometer (e.g., GloMax<sup>®</sup> Discover System or equivalent system)
- **optional:** FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255; If using this assay for the first time and/or testing an Ab that might be dependent on crosslinking FcγRIIb, use this cell line.)
- **optional:** Control Ab, Anti-4-1BB (Cat.# K1161)
- **optional:** control 4-1BB ligand/TNFSF9 with His tag (R&D Systems Cat.# 2295-4L-025/CF)
- **optional:** anti-His tag Ab for crosslinking control 4-1BB ligand (Biolegend Cat.# 652502)

#### 4. Assay Protocol for FcγRIIb-Dependent Antibodies

This assay protocol requires two engineered cell lines: 4-1BB Effector Cells and FcγRIIb CHO-K1 Cells. The FcγRIIb CHO-K1 Cells are provided in Thaw-and-Use format (Cat.# JA2251, JA2255) and CPM format (Cat.# JA2252; see *FcγRIIb CHO-K1 Cells Propagation Model Technical Manual*, #TM569, for details).

The procedure below illustrates the use of the 4-1BB Bioassay to test two FcγRIIb-dependent antibody samples against a reference sample in a single assay run using the FcγRIIb CHO-K1 Cells thaw-and-use format. Each test and reference antibody is run in triplicate, in a 10-point dilution series, in a single 96-well assay plate using the inner 60 wells. Other experimental and plate layouts are possible but may require further optimization.

**Note:** When preparing test and reference antibodies, choose an appropriate starting concentration and dilution scheme to achieve a complete dose-response curve with proper upper and lower asymptotes and sufficient points on the slope. For reference, we use 10µg/ml as a starting concentration (1X) and threefold serial dilution when testing Control Ab, Anti-4-1BB.

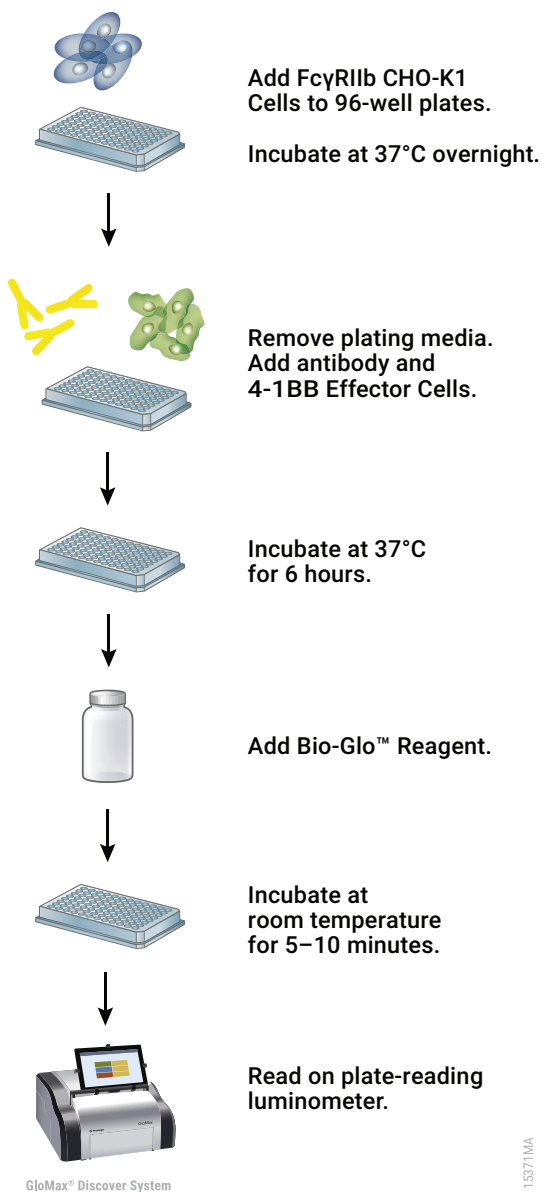
##### 4.A. Preparing Assay Buffer, Bio-Glo™ Reagent and Antibody Samples

1. **FcγRIIb CHO-K1 Cell Recovery Medium:** On the day before the assay, prepare 30ml of cell recovery medium (95% RPMI 1640/5% FBS) in a 50ml conical tube. Thaw the FBS overnight at 4°C or in a 37°C water bath on the day of use. Add 1.5ml of FBS to 28.5ml of RPMI 1640 medium. Mix well and warm to 37°C before use.
2. **Assay Buffer:** On the day of the assay, prepare an appropriate amount of assay buffer (99% RPMI 1640/1% FBS). Thaw the FBS overnight at 4°C or in a 37°C water bath on the day of use. Mix well and warm to 37°C before use. For reference, 30ml of assay buffer is typically sufficient for 120 wells in a 96-well assay format using the inner 60 wells.

**Note:** The recommended assay buffer contains 1% FBS. This concentration of FBS works well for the Control Ab, Anti-4-1BB, that we tested.

3. **Bio-Glo™ Reagent:** For reference, 10ml of Bio-Glo™ Reagent is sufficient to assay 120 wells in a 96-well assay format. Thaw the Bio-Glo™ Luciferase Assay Buffer at 4°C overnight or in a room temperature water bath on the day of assay. Equilibrate the Bio-Glo™ Luciferase Assay Buffer to ambient temperature, protected from light. Transfer all of the Bio-Glo™ Luciferase Assay Buffer into the amber bottle containing the Bio-Glo™ Luciferase Assay Substrate and mix by inversion until the Substrate is thoroughly dissolved. Equilibrate and store the reconstituted Bio-Glo™ Reagent at ambient temperature (22–25°C) protected from light before adding to assay plates. The Bio-Glo™ Reagent can be stored at room temperature after reconstitution with only ~18% loss in luminescence after 24 hours.
4. **Test and Reference Samples:** Using assay buffer as the diluent, prepare starting dilutions (dilu1, 3X final concentration) of two test antibodies (180µl each) and one reference antibody (400µl) in 1.5ml tubes. Store the tubes containing antibody starting dilutions appropriately before making antibody serial dilutions.

**Note:** If you are using Control Ab, Anti-4-1BB, as a reference antibody in your assay, prepare a 400µl starting dilution with 30µg/ml of anti-4-1BB antibody (dilu1, 3X final concentration) by adding 12µl of anti-4-1BB stock (1,000µg/ml) to 388µl of assay buffer. Store the antibody starting dilution on ice until ready to use in the assay.



**Figure 6. Schematic protocol for 4-1BB Bioassay with Fc $\gamma$ RIIb-dependent antibody.**

#### 4.B. Plate Layout Design

For the protocol described here, use the plate layout illustrated in Figure 7 as a guide. The protocol describes serial replicate dilutions (n = 3) of test and reference antibody to generate two 10-point dose-response curves for each plate.

Recommended Plate Layout Design													
	1	2	3	4	5	6	7	8	9	10	11	12	
A	B	B	B	B	B	B	B	B	B	B	B	B	Assay Buffer (B)
B	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Ab
C	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Ab
D	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Ab
E	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Ab
F	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Ab
G	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Ab
H	B	B	B	B	B	B	B	B	B	B	B	B	Assay Buffer (B)

**Figure 7. Example plate layout showing nonclustered sample locations of test antibody and reference antibody dilution series and wells containing assay buffer (denoted by “B”) alone.**

#### 4.C. Preparing and Plating FcγRIIb CHO-K1 Cells

The thaw-and-use FcγRIIb CHO-K1 cells (not included in this kit) are sensitive, and care should be taken to follow the cell thawing and plating procedures **exactly** as described. Do not overmix or overwarm the cell reagents. No additional cell culture or cell manipulation is required or recommended. We recommend that you thaw and dilute a maximum of two vials of thaw-and-use cells at one time.



Follow institutional guidelines for handling, including use of personal protective equipment (PPE), and waste disposal for biohazardous material.

#### 4.C. Preparing and Plating FcγRIIb CHO-K1 Cells (continued)

**Note:** Perform the following steps in a sterile cell culture hood.

1. On the day before performing the assay, prepare 30ml of FcγRIIb CHO-K1 cell recovery medium (RPMI 1640/5% FBS) as described in Section 4.A
2. Warm the recovery medium in a 37°C water bath for 15 minutes and transfer 14.5ml of recovery medium to a 15ml conical tube.
3. Remove one vial of thaw-and-use FcγRIIb CHO-K1 Cells from -140°C storage and transfer to the lab bench on dry ice. Thaw the vial in a 37°C water bath until cells are just thawed (about 2 minutes). While thawing, gently agitate and visually inspect.
4. Gently mix the cell suspension in the vial by pipetting, and then transfer 0.5ml of cells to the tube containing 14.5ml of recovery medium. Mix well by gently inverting 1–2 times.
5. Transfer the cell suspension to a sterile reagent reservoir. Immediately, using a multichannel pipette, dispense 100μl of cell suspension to each well of the inner 60 wells of two 96-well, white, flat-bottom assay plates.
6. Add 100μl of assay medium per well to outside wells of the assay plates.
7. Cover the assay plates with a lid, and incubate the cells overnight in a 37°C, 5% CO<sub>2</sub> incubator.

#### 4.D. Preparing Antibody Serial Dilutions

The instructions described here are for preparing a single stock of threefold serial dilutions of a single antibody for analysis in triplicate (120μl of each antibody dilution provides a sufficient volume for analysis in triplicate). Alternatively, you can prepare three independent stocks of serial dilutions to generate triplicate samples. To prepare threefold serial dilutions, you will need 400μl of reference antibody at 3X the highest antibody concentration in your dose-response curve. You will need 180μl of each test antibody at 3X the highest antibody concentration for test antibody dose-response curves. For other dilution schemes, adjust the volumes accordingly.

**Note:** If you are using Control Ab, Anti-4-1BB, as a control in the assay, follow the instructions below to prepare threefold serial dilutions. A threefold serial dilution for test antibodies is listed as an example below as well.

1. On the day of assay, prepare an appropriate amount of assay buffer as described in Section 4.A.
2. Warm the assay buffer in a 37°C water bath for 15 minutes and transfer 9.5ml of assay buffer to a 15ml conical tube. Set tube aside to be used in Section 4.E.
3. To a sterile, clear V-bottom 96-well plate, add 180μl of reference antibody starting dilution (dilu1, 3X final concentration) to wells A11 and B11 (see Figure 8).

Recommended Plate Layout for Antibody Dilutions Prepared from a Single Antibody Stock												
	1	2	3	4	5	6	7	8	9	10	11	12
A		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	Reference Ab
B		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	Reference Ab
C												
D												
E		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	Test Ab 1
F												
G		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	Test Ab 2
H												

**Figure 8. Example plate layout showing antibody serial dilutions.**

- Add 180µl of test antibodies 1 and 2 starting dilution (dilu1, 3X final concentration) to wells E11 and G11, respectively (see Figure 8).
  - Add 120µl of assay buffer to other wells in these four rows, from column 10 to column 2.
  - Transfer 60µl of the antibody starting dilutions from column 11 into column 10. Mix well by pipetting. Avoid creating bubbles.
  - Repeat equivalent threefold serial dilutions across the columns from right to left through column 3. Do not dilute into column 2.
- Note:** Wells A2, B2, E2 and G2 contain 120µl of assay buffer without antibody as a negative control.
- Take the 96-well assay plates containing FcγRIIb CHO-K1 Cells out of the incubator and remove 95µl of medium per well from the inner 60 wells using manual multichannel pipette.
  - Using an electronic multichannel pipette, immediately add 25µl of the appropriate antibody dilution (see Figure 8) to the preplated FcγRIIb CHO-K1 Cells according to the plate layout in Figure 7.
  - Cover the assay plate with a lid and keep at ambient temperature (22–25°C) while preparing the 4-1BB Effector Cells.



#### 4.E. Preparing 4-1BB Effector Cells and Setting Up Assay

The thaw-and-use 4-1BB Effector cells included in this kit are sensitive, and care should be taken to follow the cell thawing and plating procedures **exactly** as described. Do not overmix or overwarm the cell reagents. No additional cell culture or cell manipulation is required or recommended. We recommend that you thaw and dilute a maximum of two vials of thaw-and-use cells at one time.

1. Remove one vial of thaw-and-use 4-1BB Effector Cells from  $-140^{\circ}\text{C}$  storage and transfer to the bench on dry ice. Thaw the vial in a  $37^{\circ}\text{C}$  water bath until cells are just thawed (about 2 minutes). While thawing, gently agitate and visually inspect.
2. Gently mix cell suspension in the vial by pipetting, and then transfer 0.5ml of cells to the 15ml conical tube containing 9.5ml of assay buffer (see Section 4.D). Mix well by gently inverting the tube 1–2 times.
3. Transfer the 4-1BB cell suspension to a sterile reagent reservoir. Immediately, using a multichannel pipette, dispense  $50\mu\text{l}$  of cell suspension to each well of the inner 60-wells of the plates containing Fc $\gamma$ RIIb CHO-K1 Cells and anti-4-1BB antibody.
4. Add  $75\mu\text{l}$  of assay buffer to the outside wells of the 96-well assay plates.
5. Place a lid on the plates and incubate the plates in a  $\text{CO}_2$  incubator at  $37^{\circ}\text{C}$  for 6 hours.

#### 4.F. Adding Bio-Glo™ Reagent

**Note:** Bio-Glo™ Reagent should be at ambient temperature ( $22\text{--}25^{\circ}\text{C}$ ) when added to assay plates.

1. Remove the assay plates from the incubator and equilibrate to ambient temperature for 10–15 minutes.
2. Using a manual multichannel pipette, add  $75\mu\text{l}$  of Bio-Glo™ Reagent to the inner 60 wells of the assay plates, taking care not to create bubbles.
3. Add  $75\mu\text{l}$  of Bio-Glo™ Reagent to wells B1, C1 and D1 of each assay plate to measure the background signal.
4. Incubate at ambient temperature for 5–15 minutes.  
**Note:** Varying the incubation time will affect the raw relative light unit (RLU) values but should not significantly change the  $\text{EC}_{50}$  value and fold induction.
5. Measure luminescence using a luminometer or luminescence plate reader.

#### 4.G. Data Analysis

1. Determine the plate background by calculating the average RLU from wells B1, C1 and D1.
2. Calculate fold induction = 
$$\frac{\text{RLU (induced-background)}}{\text{RLU (no antibody control-background)}}$$
3. Graph data as RLU versus  $\text{Log}_{10}$  [antibody] and fold induction versus  $\text{Log}_{10}$  [antibody]. Fit curves and determine the  $\text{EC}_{50}$  value of antibody response using appropriate curve fitting software (such as GraphPad Prism® software).

## 5. Assay Protocol for Ligand or FcγRIIb-Independent Antibodies

This assay protocol illustrates the use of the 4-1BB Bioassay to test two antibody samples against a reference sample in a single assay run. Each test and reference antibody is run in triplicate, in a 10-point dilution series, in a single 96-well assay plate using the inner 60 wells. Other experimental and plate layouts are possible but may require further optimization.

**Note:** When preparing test antibodies or ligand, choose an appropriate starting concentration and dilution scheme to achieve a full dose-response curve with proper upper and lower asymptotes and sufficient points on the slope. For reference, we use 1 µg/ml as a starting concentration (1X) and 2-fold dilution when testing 4-1BB ligand.

### 5.A. Preparing Assay Buffer, Bio-Glo™ Reagent and Antibody Samples

1. **Assay Buffer:** On the day of the assay, prepare an appropriate amount of assay buffer (99% RPMI 1640/1% FBS). Thaw the FBS overnight at 4°C or in a 37°C water bath on the day of use. Mix well and warm to 37°C before use. For reference, 30ml of assay buffer is typically sufficient for 120 wells in a 96-well assay format using the inner 60 wells.

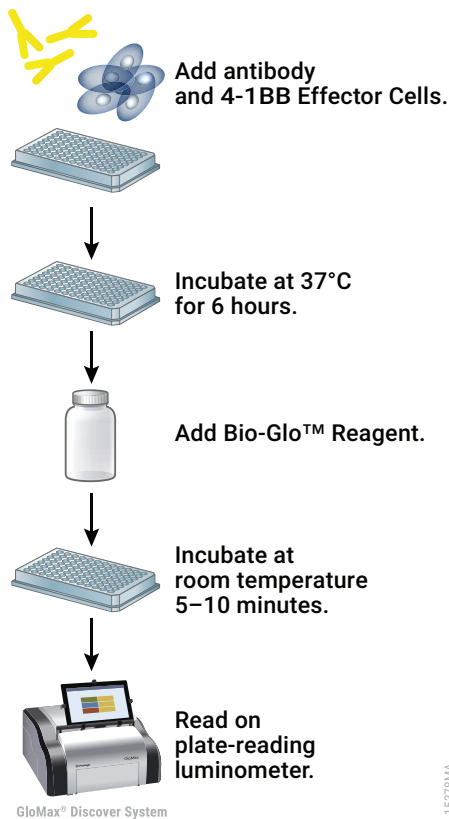
**Note:** The recommended assay buffer contains 1% FBS. This concentration of FBS works well with the crosslinked control ligand, 4-1BBL, that we tested. If you experience assay performance issues when using this assay buffer, we recommend testing different serum concentrations in the range of 0.5–10%.

2. **Bio-Glo™ Reagent:** For reference, 10ml of Bio-Glo™ Reagent is sufficient to assay 120 wells in a 96-well assay format. Thaw the Bio-Glo™ Luciferase Assay Buffer at 4°C overnight or in a room temperature water bath on the day of assay. Equilibrate the Bio-Glo™ Luciferase Assay Buffer to ambient temperature, protected from light. Transfer all of the Bio-Glo™ Luciferase Assay Buffer into the amber bottle containing the Bio-Glo™ Luciferase Assay Substrate and mix by inversion until the substrate is thoroughly dissolved. Equilibrate and store the reconstituted Bio-Glo™ Reagent at ambient temperature (22–25°C) protected from light before adding to assay plates. The Bio-Glo™ Reagent can be stored at room temperature after reconstitution with only ~18% loss in luminescence after 24 hours.
3. **Test and Reference Samples:** Using assay buffer as the diluent, prepare starting dilutions (dilu1, 3X final concentration) of two test antibodies (180 µl each) and one reference crosslinked ligand (500 µl) in 1.5ml tubes. Store the tubes containing ligand or antibody starting dilutions appropriately before making antibody serial dilutions.

#### Notes:

If you are using 4-1BBL as a reference ligand, prepare a 500 µl starting dilution with 3 µg/ml 4-1BBL (His-tagged) and 15 µg/ml crosslinking antibody, anti-His (dilu1, 3X final concentration) by adding 15 µl of 4-1BBL, stock (100 µg/ml) and 15 µl of crosslinking antibody, anti-His stock (500 µg/ml) to 470 µl of assay buffer. The final (1X) starting concentration is 1 µg/ml of 4-1BBL and 5 µg/ml of anti-His Ab. Store the antibody starting dilution on ice until ready to use in the assay.

To streamline assay setup, prepare antibody or ligand serial dilutions prior to harvesting and plating cells.



**Figure 9. Schematic protocol for 4-1BB Bioassay with 4-1BB Ligand or FcγRIIb-independent antibody.**

## 5.B. Plate Layout Design

For the protocol described here, use the plate layout illustrated in Figure 10 as a guide. The protocol describes serial replicate dilutions ( $n = 3$ ) of test antibody and reference ligand to generate two 10-point dose-response curves for each plate.

Recommended Plate Layout Design													
	1	2	3	4	5	6	7	8	9	10	11	12	
A	B	B	B	B	B	B	B	B	B	B	B	B	Assay Buffer (B)
B	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Ligand
C	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Ab
D	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Ligand
E	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Ab
F	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Ligand
G	B	no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Ab
H	B	B	B	B	B	B	B	B	B	B	B	B	Assay Buffer (B)

**Figure 10. Example plate layout showing nonclustered sample locations of test antibody and reference ligand dilution series and wells containing assay buffer (denoted by “B”) alone.**

## 5.C. Preparing Ligand or Antibody Serial Dilutions

The instructions described here are for preparation of a single stock of twofold serial dilutions of a crosslinked ligand for analysis in triplicate (120 $\mu$ l of each crosslinked ligand dilution provides a sufficient volume for analysis in triplicate). Alternatively, you can prepare three independent stocks of serial dilutions to generate triplicate samples. To prepare twofold serial dilutions for crosslinked ligand, you will need 500 $\mu$ l of reference crosslinked ligand at 3X the highest concentration in your dose response curve. To prepare threefold serial dilutions for test antibodies, you will need 180 $\mu$ l of each test antibody at 3X the highest antibody concentration in each of the test antibody dose-response curves. For other dilution schemes, adjust the volumes accordingly.

**Note:** If you are using 4-1BBL as a control in the assay, use the following instructions to prepare twofold serial dilutions. A threefold serial dilution for test antibodies is shown as an example on the next page.

**5.C. Preparing Ligand or Antibody Serial Dilutions (continued)**

1. On the day of assay, prepare an appropriate amount of assay buffer as described in Section 5.A.
2. Warm the assay buffer in a 37°C water bath for 15 minutes and transfer 9.5ml of assay buffer to a 15ml conical tube. Set the tube aside to be used in Section 5.D.
3. To a sterile, clear V-bottom 96-well plate, add 240µl of reference crosslinked ligand starting dilution (dilu1, 3X final concentration) to wells A11 and B11 (see Figure 11).
4. Add 180µl of test antibodies 1 and 2 starting dilution (dilu1, 3X final concentration) to wells E11 and G11, respectively (Figure 11).
5. Add 120µl of assay buffer to other wells in these four rows, from column 10 to column 2.
6. Transfer 120µl of the crosslinked ligand or 60µl test antibody starting dilutions from column 11 into column 10. Mix well by pipetting. Avoid creating bubbles.
7. Repeat equivalent twofold or threefold serial dilutions across the columns from right to left through column 3. Do not dilute into column 2.
8. Cover the assay plate with a lid and keep at ambient temperature (22–25°C) while preparing the 4-1BB Effector Cells.

**Note:** Wells A2, B2, E2 and G2 contain 120µl of assay buffer without antibody as a negative control.

Recommended Plate Layout for Antibody Dilutions Prepared from a Single Antibody Stock													
	1	2	3	4	5	6	7	8	9	10	11	12	
A		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		Reference Ligand
B		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		Reference Ligand
C													
D													
E		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		Test Ab 1
F													
G		no Ab	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		Test Ab 2
H													

**Figure 11. Example plate layout showing antibody serial dilutions.**

#### 5.D. Preparing 4-1BB Effector Cells and Setting Up Assay

The thaw-and-use 4-1BB Effector Cells included in this kit are sensitive, and care should be taken to follow the cell thawing and plating procedures exactly as described. Do not overmix or overwarm the cell reagents. No additional cell culture or cell manipulation is required or recommended. We recommend that you thaw and dilute a maximum of two vials of thaw-and-use cells at one time.

1. Remove one vial of thaw-and-use 4-1BB Effector cells from  $-140^{\circ}\text{C}$  storage and transfer to the bench on dry ice. Thaw the vial in a  $37^{\circ}\text{C}$  water bath until cells are just thawed (about 2 minutes). While thawing, gently agitate and visually inspect.
2. Gently mix the cell suspension in the vial by pipetting, and then transfer 0.5ml of cells to the 15ml conical tube containing 9.5ml of assay buffer (from Section 5.C). Mix well by gently inverting the tube 1–2 times.
3. Transfer the 4-1BB cell suspension to a sterile reagent reservoir. Immediately, using a multichannel pipette, dispense  $50\mu\text{l}$  of cell suspension to each well of the inner 60 wells of two 96-well, white, flat-bottom assay plates.
4. Using an electronic multichannel pipette, add  $25\mu\text{l}$  of the appropriate antibody or ligand titrations prepared in Section 5.C. to the assay plates according to the plate layout in Figure 10.
5. Add  $75\mu\text{l}$  of assay buffer to the outside wells of the 96-well assay plates.
6. Cover each assay plate with a lid and incubate in a  $37^{\circ}\text{C}$ , 5%  $\text{CO}_2$  incubator for 6 hours.

#### 5.E. Adding Bio-Glo™ Reagent

**Note:** Bio-Glo™ Reagent should be at ambient temperature ( $22\text{--}25^{\circ}\text{C}$ ) when added to assay plates.

1. Remove the assay plates from the incubator and equilibrate to ambient temperature for 10–15 minutes.
2. Using a manual multichannel pipette, add  $75\mu\text{l}$  of Bio-Glo™ Reagent to the inner 60 wells of the assay plates, taking care not to create bubbles.
3. Add  $75\mu\text{l}$  of Bio-Glo™ Reagent to wells B1, C1 and D1 of each assay plate to measure the background signal.
4. Incubate at ambient temperature for 5–15 minutes.

**Note:** Varying the incubation time will affect the raw relative light unit (RLU) values but should not significantly change the  $\text{EC}_{50}$  value and fold induction.

5. Measure luminescence using a luminometer or luminescence plate reader.

## 5.F. Data Analysis

1. Determine the plate background by calculating the average RLU from wells B1, C1 and D1.
2. Calculate fold induction = 
$$\frac{\text{RLU (induced-background)}}{\text{RLU (no antibody control-background)}}$$
3. Graph data as RLU versus  $\text{Log}_{10}$  [antibody] and fold induction versus  $\text{Log}_{10}$  [antibody]. Fit curves and determine the  $\text{EC}_{50}$  value of antibody response using appropriate curve fitting software (such as GraphPad Prism® software).

## 6. Troubleshooting

For questions not addressed here, please contact your local Promega Branch Office or Distributor. Contact information available at: [www.promega.com](http://www.promega.com). Email: [techserv@promega.com](mailto:techserv@promega.com)

### Symptoms

### Possible Causes and Comments

Low luminescence measurements (RLU readout)

Choose an instrument designed for plate-reading luminescence detection. Instruments designed primarily for fluorescence detection are not recommended. Luminometers measure and report luminescence as relative values, and actual RLU numbers will vary between instruments.

Insufficient cells per well can lead to low RLU. Handle and plate cells according to the instructions to ensure a sufficient number of viable cells per well.

If performing the assay for the first time, we recommend that you try using the Fc $\gamma$ RIIb Cells since your Ab of your interest may be dependent on crosslinking by Fc $\gamma$ RIIb. In the case of ligands, crosslinking by an antibody may be necessary.

Weak assay response (low fold induction)

Low activity of Bio-Glo™ Reagent leads to low RLU. Store and handle the Bio-Glo™ Reagent according to instructions.

Optimize the concentration range of your test sample(s) to achieve a full dose response with complete upper and lower asymptotes. The  $\text{EC}_{50}$  value obtained in the 4-1BB Bioassay may vary from the  $\text{EC}_{50}$  value obtained using other methods such as primary T cell-based assays.

Determine if the antibody used is dependent on crosslinking for performance by testing in the presence of Fc $\gamma$ RIIb Cells.

If untreated control RLU is less than 100-fold above plate reader background RLU, subtract plate background RLU from all samples before calculating fold induction.

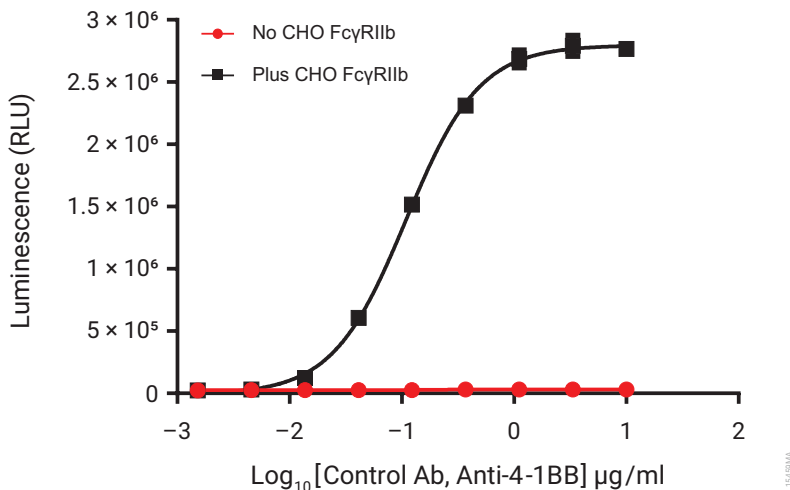
## 7. References

1. Mahoney, K.M., Rennert, P.D. and Freeman, G.J. (2015) Combination cancer immunotherapy and new immunomodulatory targets. *Nat. Rev. Drug Discov*, **14**, 561–84.
2. Melero, I. *et al.* (2015) Evolving synergistic combinations of targeted immunotherapies to combat cancer. *Nat. Rev. Cancer* **15**, 457–72.
3. Chester, C. *et al.* (2018) Immunotherapy targeting 4-1BB: Mechanistic rationale, clinical results, and future strategies. *Blood* **131**, 49–57.
4. Chester, C., Ambulkar, S. and Kohrt, H.E. (2016) 4-1BB agonism: Adding the accelerator to cancer immunotherapy. *Cancer Immunol. Immunother.* **65**, 1234–8.
5. Lee, H.W. *et al.* (2002) 4-1BB promotes the survival of CD8+ T lymphocytes by increasing expression of Bcl-xL and Bfl-1. *J. Immunol.* **169**, 4882–8.
6. Snell, L.M. *et al.* (2011) T-cell intrinsic effects of GITR and 4-1BB during viral infection and cancer immunotherapy. *Immunol. Rev.* **244**, 197–217.
7. Wang, J. *et al.* (2018) Cell-based reporter bioassays to evaluate the FcγR-dependent agonist activity of therapeutic antibodies against co-stimulatory receptors. *American Association of Cancer Research (AACR) Annual meeting*, Poster abstract #2732.
8. White, A.L. *et al.* (2011) Interaction with FcγRIIB is critical for the agonistic activity of anti-CD40 monoclonal antibody. *J. Immunol.* **187**, 1754–63.
9. Wilson, N.S. *et al.* (2011) An Fcγ receptor-dependent mechanism drives antibody-mediated target-receptor signaling in cancer cells. *Cancer Cell* **19**, 101–13.



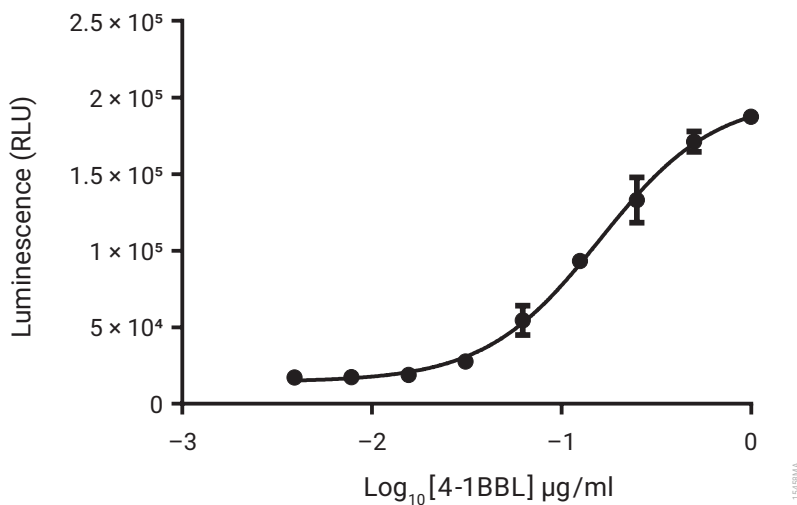
## 8. Appendix

### 8.A. Representative Assay Results with FcγRIIb-Dependent Antibody



**Figure 12. The 4-1BB Bioassay measures the activity of Control Antibody, Anti-4-1BB.** Thaw-and-use FcγRIIb CHO-K1 Cells were plated overnight. On the following day, a titration of Control Ab, Anti-4-1BB, (Section 4.C) was added, followed by addition of thaw-and-use 4-1BB Effector Cells. After 6 hours of induction at 37°C, Bio-Glo™ Luciferase Assay Reagent was added, and luminescence was determined using a GloMax® Discover System. Four-parameter logistic curve analysis was performed with GraphPad Prism® software. The EC<sub>50</sub> response determined was 0.11 µg/ml, and the fold induction was ~110.

### 8.B. Representative Assay Results with Crosslinked Ligand or FcγRIIb-Independent Antibody



**Figure 13. The 4-1BB Bioassay measures the activity of crosslinked 4-1BB Ligand.** On the assay day, thaw-and-use 4-1BB Effector Cells were plated in a 96-well plate at 50µl/well. Cells were incubated with various concentrations of crosslinked 4-1BBL (Section 5.C). After 6 hours of induction at 37°C, Bio-Glo™ Luciferase Assay Reagent was added, and luminescence was determined using a GloMax® Discover System. Four-parameter logistic curve analysis was performed with GraphPad Prism® software. The EC<sub>50</sub> response determined was ~0.16µg/ml, and the fold induction was 12.

## 8.C. Related Products

### Fc Effector Bioassays

Product	Size	Cat.#
ADCC Reporter Bioassay, Complete Kit (Raji)*	1 each	G7015
ADCC Reporter Bioassay, Core Kit*	1 each	G7010
ADCC Reporter Bioassay, F Variant, Core Kit**	1 each	G9790
ADCC Reporter Bioassay, Target Kit (Raji)*	1 each	G7016
FcγRIIIa-H ADCP Reporter Bioassay, Complete Kit**	1 each	G9901
FcγRIIIa-H ADCP Reporter Bioassay, Core Kit**	1 each	G9991
Mouse FcγRIV ADCC Bioassay, Complete Kit	1 each	M1201
Mouse FcγRIV ADCC Bioassay, Core Kit	1 each	M1211
Membrane TNFα Target Cells**	1 each	J3331
Membrane RANKL Target Cells**	1 each	J3381

\*For Research Use Only. Not for use in diagnostic procedures.

\*\*Not for Medical Diagnostic Use. Additional kit formats are available.

### Fc Effector Immunoassay

Product	Size	Cat.#
Lumit™ FcRn Binding Immunoassay	100 assays	W1151

Not for Medical Diagnostic Use. Additional kit formats are available.

### Immune Checkpoint Bioassays

Product	Size	Cat.#
CD28 Bioassay	1 each	JA6701
CD28 Blockade Bioassay	1 each	JA6101
CD40 Bioassay	1 each	JA2151
CTLA-4 Blockade Bioassay	1 each	JA3001
GITR Bioassay	1 each	JA2291
ICOS Bioassay	1 each	JA6801
ICOS Blockade Bioassay	1 each	JA6001
LAG-3/MHCII Blockade Bioassay	1 each	JA1111
OX40 Bioassay	1 each	JA2191
PD-1/PD-L1 Blockade Bioassay	1 each	J1250

### Immune Checkpoint Bioassays (continued)

Product	Size	Cat.#
PD-1+TIGIT Combination Bioassay	1 each	J2211
PD-L1 Negative Cells	1 each	J1191
TIGIT/CD155 Blockade Bioassay	1 each	J2201

Not for Medical Diagnostic Use. Additional kit formats are available.

### Macrophage-Derived Bioassays

Product	Size	Cat.#
TLR Bioassay	1 each	JA9011
SIRPα/CD47 Blockade Bioassay	1 each	JA6011
SIRPα/CD47 Blockade Bioassay, Fc-Dependent	1 each	JA4801

Not for Medical Diagnostic Use. Additional kit formats are available.

### T Cell Activation Bioassays

Product	Size	Cat.#
T Cell Activation Bioassay (IL-2)	1 each	J1651
T Cell Activation Bioassay (NFAT)	1 each	J1621
T Cell Activation Bioassay (TCRαβ-KO, CD4+)	1 each	GA1172
T Cell Activation Bioassay (TCRαβ-KO, CD8+)	1 each	GA1162
T Cell Activation Bioassay (TCRαβ-KO, CD4+, CD8+)	1 each	GA1182

Not for Medical Diagnostic Use. Additional kit formats are available.

### Cytokine and Growth Factor Bioassays

Product	Size	Cat.#
IL-2 Bioassay	1 each	JA2201
IL-6 Bioassay	1 each	JA2501
IL-12 Bioassay	1 each	JA2601
IL-15 Bioassay	1 each	JA2011
IL-23 Bioassay	1 each	JA2511
RANKL Bioassay	1 each	JA2701
VEGF Bioassay	1 each	GA2001

Not for Medical Diagnostic Use. Additional kit formats are available.

### 8.C. Related Products (continued)

#### Control Antibodies and Proteins

Product	Size	Cat.#
Control Ab, Anti-4-1BB	50µg	K1161
Control Ab, Anti-CD20	5µg	GA1130
Control Ab, Anti-CD40	50µg	K1181
Control Ab, Anti-CTLA-4	100µg	JA1020
Control Ab, Anti-LAG-3	100µg	K1150
Control Ab, Anti-OX40	50µg	K1191
Control Ab, Anti-PD-1	100µg	J1201
Control Ab, Anti-SIRPα	50µg	K1251
Control Ab, Anti-TIGIT	100µg	J2051
Control Ab, Anti-TIM-3	100µg	K1210
Recombinant VEGF ligand	10µg	J2371

#### Detection Reagents

Product	Size	Cat.#
Bio-Glo™ Luciferase Assay System	10ml	G7941
	100ml	G7940
Bio-Glo-NL™ Luciferase Assay System	10ml	J3081
	100ml	J3082
	1,000ml	J3083

Not for Medical Diagnostic Use.

#### Luminometers

Product	Size	Cat.#
GloMax® Navigator System	1 each	GM2000
GloMax® Discover System	1 each	GM3000
GloMax® Explorer System	1 each	GM3500

For Research Use Only. Not for use in diagnostic procedures.

**Note:** Additional Fc Effector, Immune Checkpoint, T Cell Activation, Cytokine, Macrophage, Primary Cell and Target Cell Killing Bioassays are available. To view and order Promega Bioassay products visit: [www.promega.com/products/reporter-bioassays/](http://www.promega.com/products/reporter-bioassays/) or e-mail: [EarlyAccess@promega.com](mailto:EarlyAccess@promega.com). For information on custom bioassay development and services visit the Promega Tailored R&D Solutions web site: [www.promega.com/custom-solutions/tailored-solutions/](http://www.promega.com/custom-solutions/tailored-solutions/).

## 9. Summary of Changes

The following changes were made to the 1/24 revision of this document:

1. Corrected the title of Table 1 to “The 4-1BB Bioassay Shows Precision, Accuracy and Linearity”.
2. Changed font and cover image.
3. Updated Section 8 and patent statements.
4. Made minor text edits.

<sup>(a)</sup>NOT FOR MEDICAL DIAGNOSTIC USE. FOR IN VITRO USE ONLY. BY USE OF THIS PRODUCT, RECIPIENT AGREES TO BE BOUND BY THE TERMS OF THIS LIMITED USE STATEMENT. If the recipient is not willing to accept the conditions of this limited use statement, and the product is unused, Promega will accept return of the unused product and provide the recipient with a full refund.

This product may not be further sold or transferred by the recipient and may be used only by the recipient, and then only for (1) research use, (2) discovery, development and monitoring of biologic drugs and vaccines, (3) quality assurance testing of biologic drugs and vaccines, and (4) product release assays for biologic drugs and vaccines. No other commercial use is allowed. “Commercial use” means any and all uses of this product by recipient for monetary or other consideration, including providing a service, information or data to unaffiliated third parties, and resale of this product for any use. Recipient has no right to propagate, modify, derivatize, genetically engineer or otherwise create variations of the cells or genes stably transfected within the cells. In addition, recipient must use Bio-Glo™ Luciferase Assay System purchased from Promega Corporation for all luminescence assays using this product or contact Promega to obtain a license for use of this product with reagents other than Promega’s. PROMEGA MAKES NO REPRESENTATIONS OR WARRANTIES OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING AS TO MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE WITH REGARDS TO THIS PRODUCT. The terms of this agreement shall be governed under the laws of the State of Wisconsin, USA.

<sup>(b)</sup>U.S. Pat. No. 8,008,006.

<sup>(c)</sup>Product cannot be used for proficiency testing.

<sup>(d)</sup>Licensed from Lonza Cologne GmbH under U.S. Pat. Nos. 7,700,357 and 8,003,389, European Pat. Nos. 1607484 and 1741778 and other pending and issued patents.

<sup>(e)</sup>Patents Pending.

© 2019–2024 Promega Corporation. All Rights Reserved.

GloMax is a registered trademark of Promega Corporation. Bio-Glo is a trademark of Promega Corporation.

Corning is a registered trademark of Corning, Inc. Echo is a registered trademark of Labcyte, Inc. GraphPad Prism is a registered trademark of GraphPad Software, Inc. JMP is a registered trademark of SAS Institute, Inc. Mantis is a registered trademark of Formulatrix, Inc.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.