



TECHNICAL MANUAL

CD40 Bioassay, Propagation Model

Instructions for use of Product
J2132

CD40 Bioassay, Propagation Model

All technical literature is available at: www.promega.com/protocols/
 Visit the website to verify that you are using the most current version of this Technical Manual.
 Email Promega Technical Services if you have questions on use of this system: techserv@promega.com

1.	Description	2
2.	Product Components and Storage Conditions	8
3.	Before You Begin	9
4.	Preparing CD40 Effector Cells	10
4.A.	Cell Thawing and Initial Cell Culture	10
4.B.	Cell Maintenance and Propagation	11
4.C.	Cell Freezing and Banking	11
5.	Assay Protocol for FcγRIIb-Dependent Antibodies	12
5.A.	Preparing Assay Buffer, Bio-Glo™ Reagent and Antibody Samples	12
5.B.	Plate Layout Design	15
5.C.	Preparing and Plating CD40 Effector Cells	16
5.D.	Preparing Antibody Serial Dilutions	16
5.E.	Preparing FcγRIIb CHO-K1 Cells	18
5.F.	Adding Antibody Samples and FcγRIIb CHO-K1 Cells to Assay Plates	18
5.G.	Adding Bio-Glo™ Reagent	19
5.H.	Data Analysis	19
6.	Assay Protocol for Ligand or FcγRIIb-Independent Antibodies	20
6.A.	Preparing Assay Buffer, Bio-Glo™ Reagent and Antibody Samples	20
6.B.	Plate Layout Design	22
6.C.	Preparing and Plating CD40 Effector Cells (the day before assay)	22
6.D.	Preparing Ligand or Antibody Serial Dilutions (the day of assay)	23
6.E.	Adding Ligand or Antibody Samples to Assay Plates	24
6.F.	Adding Bio-Glo™ Reagent	25
6.G.	Data Analysis	25
7.	Troubleshooting	26
8.	References	27
9.	Appendix	28
9.A.	Representative Assay Results with FcγRIIb-Dependent Antibody	28
9.B.	Representative Assay Results with CD40 Ligand	29
9.C.	Composition of Buffers and Solutions	29
9.D.	Related Products	30
10.	Summary of Changes	33

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 costimulatory 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 agonist monoclonal antibodies has emerged as a novel strategy to enhance anti-tumor immune responses and promote immune-mediated tumor rejection (1,2).

The cell surface molecule CD40 expressed by B cells, dendritic cells and monocytes is a member of the tumor necrosis factor receptor superfamily. CD40 ligand (CD154) is the primary ligand for CD40 and is expressed by activated T cells, which are critical regulators of cellular and humoral immunity. Signaling via CD40 triggers activation of antigen-presenting cells (APC). Agonist CD40 antibodies were found to mimic the signal of CD40 ligand and were capable of substituting for the function of CD4+ helper T cells in murine models of T-cell-mediated immunity. Therefore, agonist CD40 antibodies can rescue the function of APC in tumor-bearing hosts and restore effective immune responses against tumor antigens. Subsequent data from multiple preclinical models has demonstrated synergistic enhancement from combining CD40 agonists with cytotoxics, especially chemotherapy (2,3).

Current methods used to measure the activity of biologic drugs targeting CD40 rely on primary human T cells and measurement of functional endpoints such as cell proliferation, cell surface marker expression, and interferon gamma (IFN γ) 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 unqualified assay reagents. As a result, these assays are difficult to establish in a drug-development setting.

The CD40 Bioassay, Propagation Model^(a-e) (Cat.# J2132), 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 and other biologics that can bind and activate CD40 (4,5). The assay consists of CD40 Effector cells that express human CD40. CD40 Effector Cells are also genetically engineered with a luciferase reporter driven by a response element that can respond to CD40 ligand/agonist antibody stimulation. The CD40 Effector Cells are provided in Cell Propagation Model (CPM) format, which includes cryopreserved cells that can be thawed, propagated and banked for long-term use.

Based on the antibody properties to be tested, the CD40 Bioassay can be conducted in either single-cell systems or with Fc γ RIIb CHO-K1 Cells (Cat.# J2232) to determine the agonist antibodies that can activate CD40 in a Fc γ R-dependent manner. Induction of the CD40 Effector Cells with CD40 ligand or an agonist CD40 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 9.D, Related Products).

The CD40 Bioassay, Propagation Model, reflects the mechanism of action (MOA) of biologics designed to activate the CD40 receptor. Specifically, CD40-mediated luminescence is detected following the addition of CD40 agonist antibodies or CD40 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 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 50% human serum (in antibody samples) with some loss in fold induction (Figure 5), indicating potential for further development into a neutralizing antibody bioassay.

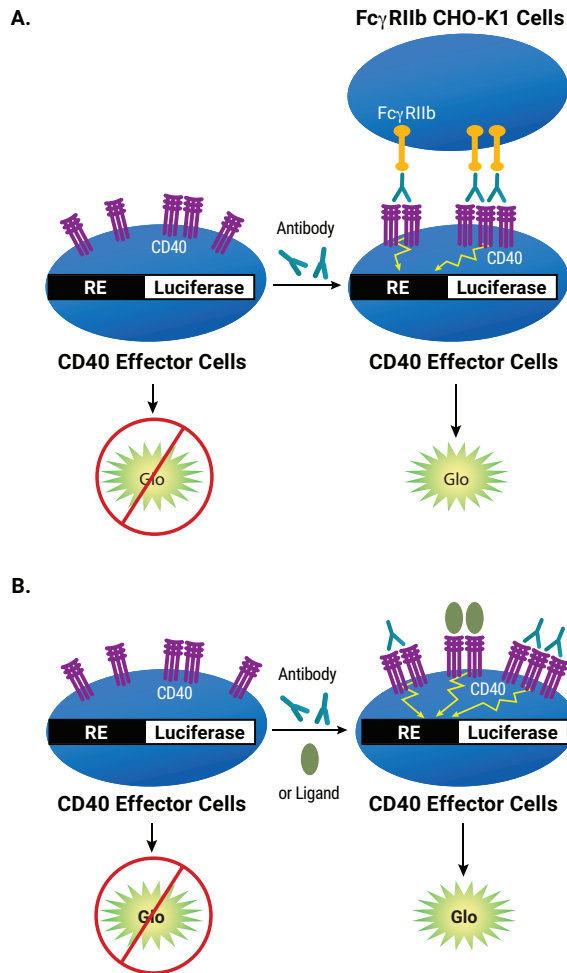


Figure 1. Representation of the CD40 Bioassay. Panel A. Assay with Fc γ RIIb-dependent antibody. The bioassay consists of two engineered cell lines, CD40 Effector Cells and Fc γ RIIb CHO-K1 Cells. In the presence of Fc γ RIIb CHO-K1 Cells, the anti-CD40 antibody is cross-linked, thereby inducing CD40 pathway-activated luminescence, which can be detected in a dose-dependent manner by addition of Bio-Glo™ Reagent and quantitation with a luminometer. **Panel B.** Assay with Fc γ RIIb-independent antibody or ligand. The bioassay consists of one engineered cell line, CD40 Effector Cells. In the absence of agonist antibody or CD40 ligand, the CD40 receptor is not activated and luminescence signal is low. The addition of CD40 ligand or antibody induces the CD40 pathway-activated luminescence, which can be detected in a dose-dependent manner by addition of Bio-Glo™ Reagent and quantified with a luminometer.

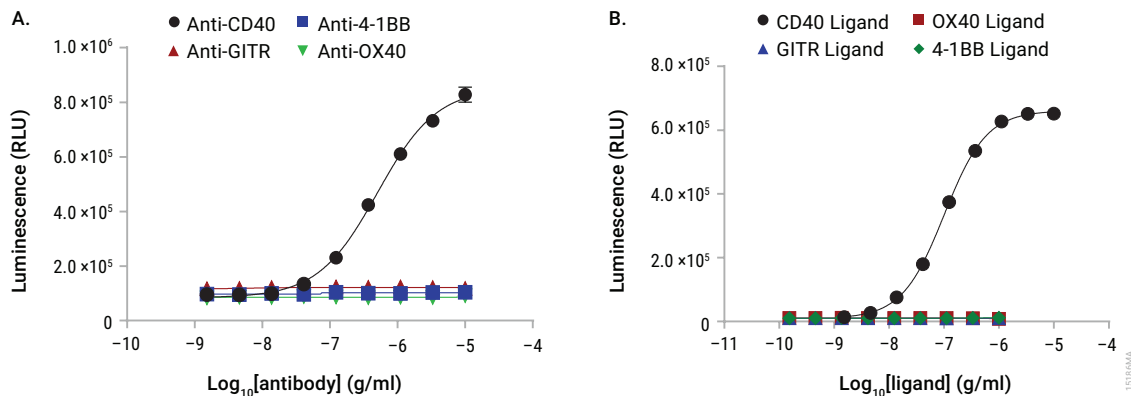


Figure 2. The CD40 Bioassay reflects the mechanism of action (MOA) and specificity of biologics designed to activate CD40 receptor. Panel A. CD40 Effector Cells were induced respectively with a serial titration of anti-CD40 antibody, anti-4-1BB antibody, anti-OX40 antibody or anti-GITR antibody, as indicated, in the presence of FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255). Panel B. CD40 Effector Cells were induced with a serial titration of CD40 Ligand, 4-1BB Ligand, OX40 Ligand or GITR Ligand, as indicated. For both panels, Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were fitted to a four-parameter logistic curve using GraphPad Prism® software. Data were generated using thaw-and-use cells.

Table 1. The CD40 Bioassay Shows Precision, Accuracy and Linearity.

Parameter	Results	
	% Expected Relative Potency	% Recovery
Accuracy	50	98.7
	70	99.8
	140	104.0
	200	100.8
Repeatability (% CV)	100% (Reference)	5.1
Intermediate Precision (% CV)		7.0
Linearity (r ²)		0.997
Linearity (y = mx + b)		y = 1.028x - 0.0284

A 50–200% theoretical potency series of Control Ab, Anti-CD40 (Cat.# K1181), was analyzed in triplicate in three independent experiments performed on three days by two analysts. 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.

1. Description (continued)

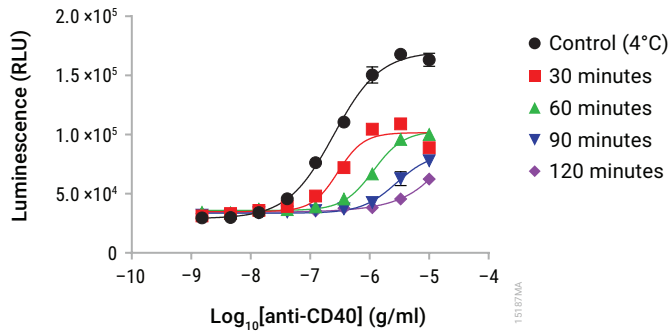


Figure 3. The CD40 Bioassay is stability-indicating. Samples of Control Ab, Anti-CD40 (Cat.# K1181), were maintained at 4°C (control) or heat-treated at 65°C for the indicated times, and then analyzed using the CD40 Bioassay with FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255). Bio-Glo™ Reagent was added, and luminescence quantified using the GloMax® Discover System. Data were fitted to a four-parameter logistic curve using GraphPad Prism® software. Data were generated using thaw-and-use cells.

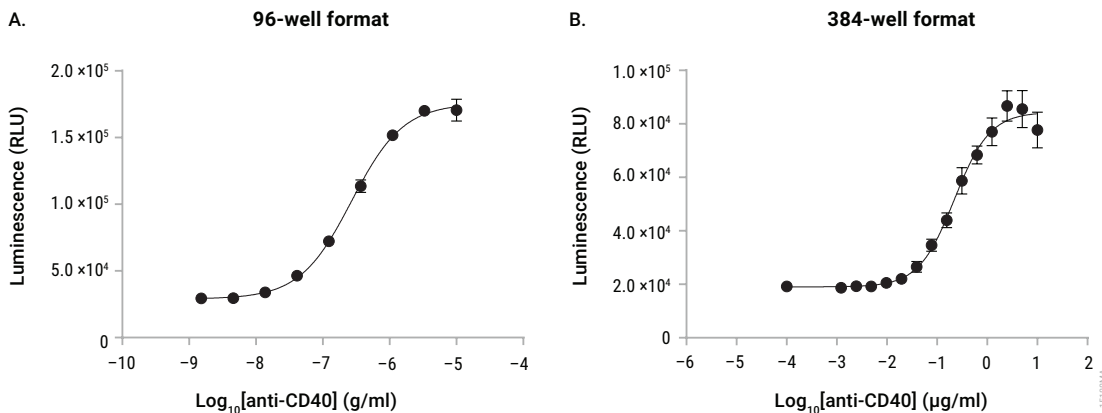


Figure 4. The CD40 Bioassay is amenable to 384-well plate format and compatible with laboratory automation.

Panel A. The CD40 Bioassay was performed in 96-well plates as described in this technical manual using Control Ab, Anti-CD40 (Cat.# K1181). **Panel B.** The CD40 Bioassay was performed in 384-well format using a ThermoFisher Multidrop™ Combi nL Reagent Dispenser to dispense the cells and assay buffer and Echo® Acoustic liquid handler for antibody handling. On the day before assay, CD40 Effector Cells were plated at 8×10^3 cells/10µl/well. The next day, assay buffer was added to the plate at 5µl/well. Control Ab, Anti-CD40, was dispensed in submicroliter volumes to the assay plate and serially titrated by direct dilution. Finally, FcγRIIb CHO-K1 cells were added at 10×10^3 cells/5µl/well. After a 6-hour incubation, 20µl of Bio-Glo™ Reagent was added and luminescence was quantified using the GloMax® Discover System. Data were fitted to a four-parameter logistic curve using GraphPad Prism® software. The EC₅₀ values were 0.27µg/ml and 0.23µg/ml, and the fold inductions were 6.3 and 4.7 for 96-well and 384-well format, respectively. Data were generated using thaw-and-use cells.

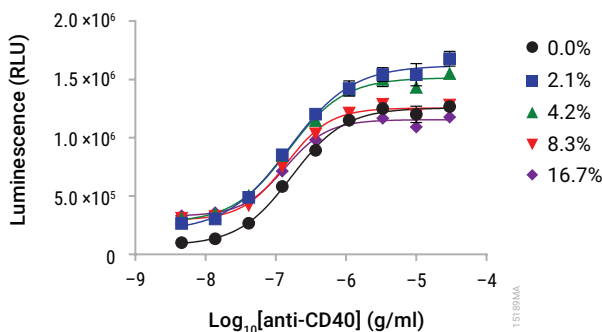


Figure 5. The CD40 Bioassay is tolerant to human serum. Control Ab, Anti-CD40 (Cat.# K1181) was analyzed in the presence of FcγRIIb CHO-K1 Cells and increasing concentrations of pooled normal human serum (0–50% in the antibody sample), resulting in final assay concentration of human serum (0–16.7%). Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were fitted to a four-parameter logistic curve using GraphPad Prism® software. Data were generated using thaw-and-use cells. The CD40 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.#
CD40 Bioassay, Propagation Model	1 each	J2132

Not for Medical Diagnostic Use. Includes:

- 2 vials CD40 Effector Cells (CPM), 7.5×10^6 cells/ml (1.0ml per vial)

Note: Thaw and propagate one vial to create frozen cell banks before use in an assay. The second vial should be reserved for future use.

Storage Conditions: Upon arrival, immediately transfer the cell vials to below -140°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°C because this will decrease cell viability and cell performance.

3. Before You Begin

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

Note the catalog number and lot number from the cell vial box label. This information can be used to download documents for the specified product from the website such as Certificate of Analysis.

The CD40 Bioassay, Propagation Model, is intended to be used with user-provided ligands, antibodies or other biologics designed to activate or block CD40 receptor signal. Control Ab, Anti-CD40 (Cat.# K1181) and FcγRIIb CHO Cells (Cat.# J2232) are available separately for use in assay optimization and routine quality control. We strongly recommend including Control Ab, Anti-CD40, as a positive control in the first few assays to gain familiarity with the assay. Data generated using these reagents are shown in Figures 3–5 and Sections 9.A (antibodies) and 9.B (ligands), Representative Assay Results.

Cell thawing, propagation and banking should be performed exactly as described in Section 4. Cell seeding and propagation densities have been optimized to ensure stable cell growth, which is reflected in a steady cell doubling rate, to achieve optimal, consistent performance. An accurate, reliable and reproducible cell counting method is required for routine cell culturing and optimal bioassay performance. The recommended cell plating densities, induction time and assay buffer components described in Sections 5 and 6 were established using CD40 Ligand and Control Ab, Anti-CD40, respectively. You may need to adjust the parameters provided here and optimize assay conditions for your own antibody or biologic.

The CD40 Bioassay produces a bioluminescent signal and is expected to be compatible with all standard luminometers or luminescence plate readers. Bioassay development and performance data included in this Technical Manual were generated using the GloMax[®] Discover System (see Section 9.D, Related Products). An integration time of 0.5 seconds/well was used for all readings.

Materials to Be Supplied by the User

(Composition of buffers and solutions is provided in Section 9.C.)

Reagents


- Control Ab, Anti-CD40 (Cat.# K1181) available as a standalone reagent or user-defined anti-CD40 antibodies or other biologics
- CD40 Ligand (R&D Systems, Cat.# 6420-CL/CF), optional as control
- RPMI 1640 Medium with L-glutamine and HEPES (e.g., Corning[®] Cat.# 10-041-CV or GIBCO[®] Cat.# 22400)
- McCoy's 5A Medium with L-glutamine (e.g., GIBCO[®] Cat.# 16600)
- fetal bovine serum (e.g., HyClone Cat.# SH30070.03 or GIBCO[®] Cat.# 16000044)
- hygromycin B (e.g., GIBCO[®] Cat.# 10687010)
- DMSO (e.g., Sigma Cat.# D2650)
- DPBS (e.g., GIBCO[®] Cat.# 14190)
- Accutase[®] solution (e.g., Sigma Cat.# A6964)
- Trypan blue solution (e.g., Sigma Cat.# T8154)
- Bio-Glo[™] Luciferase Assay System (Cat.# G7940, G7941)

3. Before You Begin (continued)

Supplies and Equipment

- white, flat-bottom 96-well assay plates (e.g., Corning® Cat.# 3917)
- T75 tissue culture flask (e.g., Corning® Cat.# CLS430641)
- sterile, clear V-bottom 96-well plate with lid (e.g., Costar® 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® Cat.# 4870)
- 37°C, 5% CO₂ incubator
- 37°C water bath
- plate reader with glow luminescence measuring capability or luminometer (e.g., GloMax® Discover System or equivalent)

4. Preparing CD40 Effector Cells

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

4.A. Cell Thawing and Initial Cell Culture

1. Prepare 50ml of initial cell culture medium by adding 5ml of FBS to 45ml of McCoy's 5A medium prewarmed to 37°C. This initial cell culture medium will be used for culturing the cells immediately after thawing.
2. Transfer 9ml of prewarmed initial cell culture medium to a 15ml conical tube.
3. Remove one vial of CD40 Effector Cells from storage at -140°C and thaw in a 37°C water bath with gentle agitation (no inversion) until just thawed (typically 2–3 minutes).
4. Transfer all of the cells (approximately 1ml) to the 15ml conical tube containing 9ml of prewarmed initial cell culture medium.
5. Centrifuge at 180 × *g* for 5 minutes.
6. Carefully aspirate the medium, and resuspend the cell pellet in 30ml of prewarmed initial cell culture medium.
7. Transfer the cell suspension to two T75 tissue culture flasks (15ml per flask), and place the flasks horizontally in a 37°C, 5% CO₂ incubator.
8. Incubate for approximately 48 hours before passaging the cells.

4.B Cell Maintenance and Propagation

Note: For cell maintenance and propagation starting from the second cell passage, use the cell growth medium containing antibiotics (see Section 9.C) and monitor cell viability and doubling rate during propagation. The cell growth rate will stabilize by 7–10 days post-thaw, at which time cell viability is typically >95%, and the average cell doubling rate is ~24 hours. Passage number should be recorded for each passage. In our experience, cells maintain their functionality for up to 25 passages if passaging is performed on a Monday-Wednesday-Friday schedule.

1. On the day of cell passage, aspirate the cell culture medium and wash the cells with DPBS.
2. Add 3ml of Accutase® solution to each T75 flask and place in a 37°C, 5% CO₂ incubator for 5–7 minutes or until the cells round up and detach from the bottom of the flask.
3. Add 7ml of cell culture medium to each T75 flask. Transfer the cell suspension to a sterile 15ml or 50ml conical tube.
4. Count the cells by Trypan blue staining. We suggest seeding the cells at a density of 4×10^4 cells/cm² if passaging every two days (e.g., Monday–Wednesday or Wednesday–Friday) or 1.8×10^4 cells/cm² if passaging every three days (e.g., Friday–Monday).
5. Add an appropriate amount of cell growth media (see Section 9.C) to a new flask.
6. Transfer the appropriate volume of cell suspension to achieve the desired cell seeding density per area.
7. Place the flasks horizontally in a 37°C, 5% CO₂ incubator.

4.C. Cell Freezing and Banking

1. On the day of cell freezing, make fresh cell freezing medium (Section 9.C) and keep on ice.
2. Aspirate the cell culture medium and wash the cells with DPBS.
3. Add 3ml of Accutase® solution to each T75 flask and place in a 37°C, 5% CO₂ incubator for 5–7 minutes or until the cells round up and detach from the bottom of the flask.
4. Add 7ml of cell culture medium to each T75 flask. Transfer the cell suspension to a sterile 15ml or 50ml conical tube.
5. Remove a sample for cell counting by Trypan blue staining. Calculate the volume of cell freezing medium needed based on desired cell freezing densities of 3×10^6 – 1×10^7 cells/ml.
6. Transfer the cell suspension to 50ml sterile conical tubes or larger sized centrifuge tubes, and centrifuge at $180 \times g$, 4°C, for 5–10 minutes.
7. Carefully aspirate the supernatant and avoid disturbing the cell pellet.
8. Gently resuspend the cell pellet in ice-cold cell freezing medium to a final cell density of 3×10^6 – 1×10^7 cells/ml. Combine the cell suspensions into a single tube and dispense into cryovials.
9. Freeze the cells using a controlled-rate freezer (preferred), or a Mr. Frosty® or a Styrofoam® rack in a –80°C freezer overnight. Transfer the vials to at or below –140°C for long-term storage.

5. Assay Protocol for FcγRIIb-Dependent Antibodies

This assay protocol requires two engineered cell lines: CD40 Effector Cells and FcγRIIb CHO-K1 Cells. We provide FcγRIIb CHO-K1 Cells in thaw and use (Cat.# JA2151, JA2155) and cell propagation model format (Cat.# JA2132). See *FcγRIIb CHO-K1 Cells, Propagation Model Technical Manual*, #TM569, for cell growth, thawing and banking. Either cell format may be used in this assay.

The procedure below illustrates the use of the CD40 Bioassay to test two FcγRIIb-dependent antibody samples against a reference sample in a single assay run. Each test and reference antibody is run in triplicate, in a ten-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 full 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-CD40.

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

1. **Cell Recovery Medium:** On the day before the assay, prepare 30ml of cell recovery medium (90% RPMI 1640/10% 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 3ml of FBS to 27ml 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-CD40, 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.

If you are using a large (100ml) size of Bio-Glo™ Luciferase Assay System, dispense the reconstituted Bio-Glo™ Reagent into 10ml aliquots and store at –20°C for up to six weeks. Avoid repeated freeze-thaw cycles. On the day of the assay, thaw an appropriate amount of reconstituted Bio-Glo™ Reagent in a room temperature water bath for at least 1–2 hours before use. Bio-Glo™ Reagent can be stored at room temperature after reconstitution with ~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 (180 μ l each) and one reference antibody (360 μ l) in 1.5ml tubes. Store the tubes containing antibody starting dilutions appropriately before making antibody serial dilutions.

Notes:

- a. If you are using Control Ab, Anti-CD40 (Cat.# K1181), as a reference antibody in your assay, prepare 360 μ l of starting dilution with 30 μ g/ml anti-CD40 antibody (dilu1, 3X final concentration) by adding 10.8 μ l of anti-CD40 stock (1.0mg/ml) to 349.2 μ l of assay buffer. Store the antibody starting dilution on ice until ready to use in the assay.
- b. To streamline assay setup, prepare antibody serial dilutions prior to harvesting and plating cells.

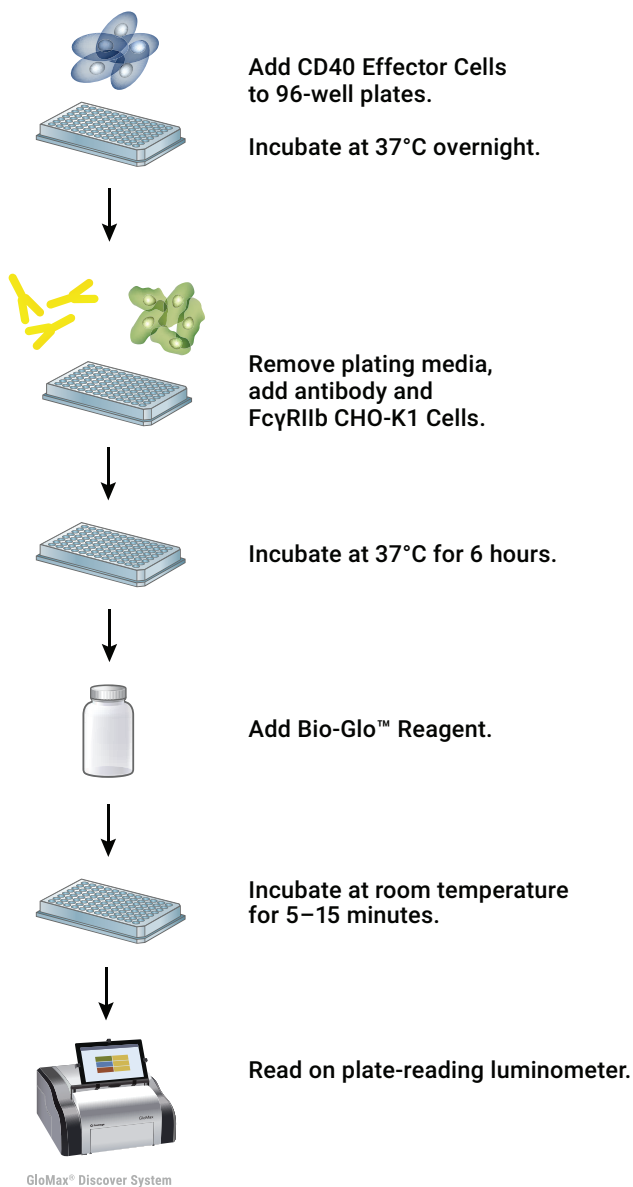


Figure 6. Schematic protocol for FcγRIIb-dependent CD40 Bioassay.

5.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 ten-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 sample locations of test and reference antibody dilution series and wells containing assay buffer (denoted by "B") alone.

5.C. Preparing and Plating CD40 Effector Cells

While maintaining the CD40 Effector Cells, follow the recommended cell seeding density. Changes in cell culture volume or seeding density will affect cell growth rate and assay performance. Do not allow the cells to reach 100% confluence. Only use the cells in the assay after the cell doubling rate has stabilized during propagation.

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

1. We recommend passaging the cells two days or three days before performing the assay as described in Section 4 to ensure optimal and consistent assay performance.
2. On the day before performing the assay, prepare fresh CD40 Effector cell recovery medium (RPMI 1640/10% FBS).
3. Aspirate the cell culture medium from the CD40 Effector Cells and wash with DPBS.
4. Add 3ml of Accutase® solution to each T75 flask and place in a 37°C, 5% CO₂ incubator for 5–7 minutes or until the cells round up and detach from the bottom of the flask.
5. Add 7ml of initial cell culture medium (McCoy's 5A Medium + 10% FBS) to each T75 flask. Transfer the cell suspension to a sterile 15ml or 50ml conical tube.
6. Gently mix and count the CD40 Effector Cells by Trypan blue staining.
7. Pellet the cells by centrifugation at 180 × *g* for 5–10 minutes.
8. Gently resuspend the cell pellet in cell recovery medium to achieve a concentration of 1.5 × 10⁵ cells/ml.
9. Transfer the CD40 Effector Cells to a sterile reagent reservoir. Using a multichannel pipette, dispense 100µl (1.5 × 10⁴ cells) of CD40 Effector Cells into the inner 60 wells of the assay plates.
10. Add 75µl of cell recovery medium to the outside wells of the 96-well assay plates.
11. Place lids on the assay plates and incubate in a 37°C, 5% CO₂ incubator overnight (18–22 hours).

5.D. Preparing Antibody Serial Dilutions

The instructions described here are for preparation of 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 360µl 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-CD40 (Cat.# K1181), 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 5.A.
2. 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 11).
3. Add 180µl of test antibodies 1 and 2 starting dilution (dilu1, 3X final concentration) to wells E11 and G11, respectively (see Figure 11).

4. Add 120µl of assay buffer to other wells in these four rows, from column 10 to column 2.
5. Transfer 60µl of the antibody starting dilutions from column 11 into column 10. Mix well by pipetting. Avoid creating bubbles.
6. Repeat equivalent threefold serial dilutions across the columns from right to left through column 3. Do not dilute into column 2.
7. Cover the assay plate with a lid and keep at ambient temperature (22–25°C) while preparing the CD40 Effector Cells.

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 Ligand	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		Reference Ligand
B		no Ligand	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 8. Example plate layout showing antibody serial dilutions. Wells A2, B2, E2 and G2 contain 120µl of assay buffer without antibody as a negative control.

5.E. Preparing FcγRIIb CHO-K1 Cells

While maintaining the FcγRIIb CHO-K1 Cells, it is important to follow the recommended cell seeding density. Changes in cell culture volume or seeding density will affect cell growth rate and assay performance. Only use the cells in the assay after the cell doubling rate has stabilized during propagation.

1. Passage the cells two days or three days before performing the assay. To ensure optimal and consistent assay performance, maintain the cell density, upon harvest, in the range of $0.5\text{--}1.0 \times 10^5$ cells/ml and cell viability at greater than 90%.
2. Count the FcγRIIb CHO-K1 Cells by Trypan blue staining, and calculate the cell density and viability.
3. Transfer an appropriate amount of FcγRIIb CHO-K1 Cells from the culture vessel to a 50ml conical tube or larger sized centrifuge tube.
4. Centrifuge the cells at $180 \times g$ for 5–10 minutes at ambient temperature and resuspend the pellet in assay buffer to generate a targeted cell density of 8.0×10^5 cells/ml.
5. You will need at least 8ml of FcγRIIb CHO-K1 Cells to fill 120 assay wells, or the inner 60 wells of two assay plates.

5.F. Adding Antibody Samples and FcγRIIb CHO-K1 Cells to Assay Plates

1. Take the 96-well assay plates containing CD40 Effector Cells out of the incubator. Using a manual multichannel pipette, carefully remove all medium from the inner 60 wells of each plate.
2. Using an electronic multichannel pipette, add 25μl of the appropriate antibody titration to the assay plates according to the plate layout in Figure 8.
3. Transfer the FcγRIIb CHO-K1 Cells prepared in Section 5.E to a sterile reagent reservoir. Using a multichannel pipette, dispense 50μl (4.0×10^4 cells) of FcγRIIb CHO-K1 Cells into the wells containing antibody or ligand.
4. Cover the assay plates with lids and incubate in a 37°C, 5% CO₂ incubator for 6 hours.

5.G. Adding Bio-Glo™ Reagent

Note: Bio-Glo™ Reagent should be at ambient temperature (22–25°C) when added to assay plates.

1. Take the assay plates out from the incubator and equilibrate to ambient temperature for 10–15 minutes.
2. Using a manual multichannel pipette, add 75µl of Bio-Glo™ Reagent to the inner 60 wells of the assay plates, taking care not to create bubbles.
3. Add 75µl 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 EC₅₀ value and fold induction.

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

5.H. 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 Log₁₀ [antibody] and fold induction versus Log₁₀ [antibody]. Fit curves and determine the EC₅₀ value of antibody response using appropriate curve fitting software (such as GraphPad Prism® software).

6. Assay Protocol for Ligand or FcγRIIb-Independent Antibodies

This assay protocol illustrates the use of the CD40 Bioassay to test two ligand or antibody samples against a reference sample in a single assay run. Each test and reference antibody is run in triplicate, in a ten-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 and reference 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 10µg/ml as a starting concentration (1X) and threefold dilution when testing a CD40 ligand.

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

1. **Cell Recovery Medium:** On the day before the assay, prepare 30ml of cell recovery medium (90% RPMI 1640/10% 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 3ml of FBS to 27ml 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 recommended CD40 Ligand (CD40L).

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.

If you are using a large (100ml) size of Bio-Glo™ Luciferase Assay System, dispense the reconstituted Bio-Glo™ Reagent into 10ml aliquots and store at –20°C for up to six weeks. Avoid repeated freeze-thaw cycles. On the day of the assay, thaw an appropriate amount of reconstituted Bio-Glo™ Reagent in a room temperature water bath for at least 1–2 hours before use. 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 or ligands (180µl each) and one reference ligand or reference antibody (360µl) in 1.5ml tubes. Store the tubes containing ligand or antibody starting dilutions appropriately before making antibody serial dilutions.

Note: If you are using a CD40 Ligand (CD40L) as a reference ligand in your assay, prepare 360µl starting dilution with 30µg/ml CD40L (dilu1, 3X final concentration). Store the antibody starting dilution on ice until ready to use in the assay.

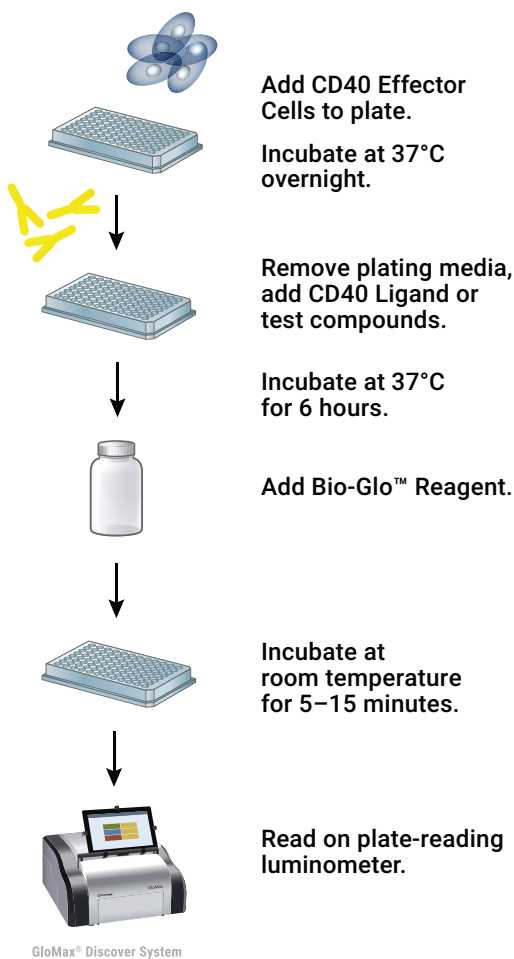


Figure 9. Schematic protocol for CD40 Ligand or FcγRIIb-independent CD40 Bioassay.

6.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 antibodies to generate two ten-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 10. Example plate layout showing non-clustered sample locations of test and reference antibody dilution series and wells containing assay buffer (denoted by "B") alone.

6.C. Preparing and Plating CD40 Effector Cells (the day before assay)

While maintaining the CD40 Effector Cells, follow the recommended cell seeding density. Changes in cell culture volume or seeding density will affect cell growth rate and assay performance. Do not allow the cells to reach 100% confluence. Only use the cells in the assay after the cell doubling rate has stabilized during propagation.

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

1. We recommend passaging the cells two days or three days before performing the assay as described in Section 4 to ensure optimal and consistent assay performance.
2. On the day before performing the assay, prepare fresh CD40 Effector Cell recovery medium (RPMI 1640/10% FBS).

3. Aspirate the cell culture medium from the CD40 Effector Cells and wash with DPBS.
4. Add 3ml of Accutase solution to each T75 flask and place in a 37°C, 5% CO₂ incubator for 5–7 minutes or until the cells round up and detach from the bottom of the flask.
5. Add 7ml of initial cell culture medium (i.e., 90% McCoys medium + 10% FBS) to each T75 flask. Transfer the cell suspension to a sterile 15ml or 50ml conical tube.
6. Gently mix and count the CD40 Effector Cells by Trypan blue staining.
7. Pellet the cells by centrifugation at 180 × *g* for 5–10 minutes.
8. Gently resuspend the cell pellet in cell recovery medium to achieve a concentration of 1.5 × 10⁵ cells/ml.
9. Transfer the CD40 Effector Cells to a sterile reagent reservoir. Using a multichannel pipette, dispense 100µl (1.5 × 10⁴ cells) of CD40 Effector Cells into the inner 60 wells of the assay plates.
10. Add 75µl of cell recovery medium to the outside wells of the 96-well assay plates.
11. Place lids on the assay plates and incubate in a 37°C, 5% CO₂ incubator overnight (18–22 hours).

6.D. Preparing Ligand or Antibody Serial Dilutions (the day of assay)

The instructions described here are for preparation of a single stock of threefold serial dilutions of a ligand for analysis in triplicate (120µl of each 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 threefold serial dilutions for CD40 ligand, you will need 360µl of reference ligand at 3X the highest concentration in your dose-response curve. To prepare threefold serial dilutions for test antibodies, you will need 180µl 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 Control Ligand, CD40L, as a control in the assay, follow the instructions below to prepare threefold serial dilutions.

1. On the day of assay, prepare an appropriate amount of assay buffer as described in Section 5.A.
2. To a sterile, clear V-bottom 96-well plate, add 180µl of reference ligand starting dilution (dilu1, 3X final concentration) to wells A11 and B11 (see Figure 8).
3. Add 180µl of test antibodies 1 and 2 starting dilution (dilu1, 3X final concentration) to wells E11 and G11, respectively (see Figure 8).
4. Add 120µl of assay buffer to other wells in these four rows, from column 10 to column 2.
5. Transfer 60µl of the ligand or antibody starting dilutions from column 11 into column 10. Mix well by pipetting. Avoid creating bubbles.
6. Repeat equivalent threefold serial dilutions across the columns from right to left through column 3. Do not dilute into column 2.
7. Cover the assay plate with a lid and keep at ambient temperature (22–25°C) while preparing the CD40 Effector Cell plates.

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 Ligand	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		Reference Ligand
B		no Ligand	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. Wells A2, B2, E2 and G2 contain 120µl of assay buffer without antibody as a negative control.

6.E. Adding Ligand or Antibody Samples to Assay Plates

1. Remove assay plates from incubator and carefully remove all cell recovery medium from the inner 60 wells of each assay plate.
2. Using an electronic multichannel pipette, add 25µl of the appropriate antibody or ligand titration to the assay plates according to the plate layout in Figure 10.
3. Add additional 50µl of assay buffer to all assay wells for a final volume of 75µl per well.
4. Cover the assay plates with lids and incubate in a 37°C, 5% CO₂ incubator for 6 hours.

6.F. Adding Bio-Glo™ Reagent

Note: Bio-Glo™ Reagent should be at ambient temperature (22–25°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 75µl of Bio-Glo™ Reagent to the inner 60 wells of the assay plates, taking care not to create bubbles.
3. Add 75µl 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 EC₅₀ value and fold induction.

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

6.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 Log₁₀ [antibody] and fold induction versus Log₁₀ [antibody]. Fit curves and determine the EC₅₀ value of antibody response using appropriate curve fitting software (such as GraphPad Prism® software).

7. Troubleshooting

For questions not addressed here, please contact your local Promega Branch Office or Distributor. Contact information available at: www.promega.com. Email: techserv@promega.com

Symptoms	Causes and Comments
Low luminescence measurements (RLU readout)	<p>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.</p> <p>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.</p> <p>If performing the assay for the first time, we recommend that you use the FcγRIIb cells since your Ab of your interest may be dependent on crosslinking by FcγRIIb. In the case of ligands, crosslinking by an antibody may be necessary.</p>
Variability in assay performance	<p>Low activity of Bio-Glo™ Reagent leads to low RLU. Store and handle the Bio-Glo™ Reagent according to the instructions. Assay performance can be affected by variations in cell growth conditions including plating and harvest density, centrifuge times and speeds, and freezing and/or DMSO exposure times during cell banking.</p> <p>Poor cell viability and variations in doubling time may affect assay performance. Ensure consistent cell growth by handling the cells exactly according to the instructions. Avoid one-day cell passages whenever possible, especially with the CD40 Effector Cells. Ensure you are using high-quality cell culture reagents (especially serum) and plasticware for maintaining cells in culture. Ensure accurate and consistent cell counting methods.</p>
Weak assay response (low fold induction)	<p>Optimize the concentration range of your test sample(s) to achieve a full dose response with complete upper and lower asymptotes. The EC₅₀ value obtained in the CD40 Bioassay may vary from the EC₅₀ value obtained using other methods.</p> <p>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.</p>

Symptoms

Weak assay response (low fold induction)

Causes and Comments

If performing the assay for the first time, we recommend trying the Fc γ RIIb cells since your Ab of interest may be dependent on crosslinking by Fc γ RIIb. (See Figure 1 for details on the use of Fc γ RIIb.) If you are using ligands, crosslinking by an antibody may be necessary.

8. References

1. Mahoney, K.M. *et al.* (2015) Combination cancer immunotherapy and new immunomodulatory targets. *Nat. Rev. Drug Disc.* **14**, 561–84.
2. Vonderheide, R.H. *et al.* (2013) CD40 immunotherapy for pancreatic cancer. *Cancer Immunol. Immunother.* **62(5)** 949–54.
3. Melero, I. *et al.* (2015) Evolving synergistic combinations of target immunotherapies to combat cancer. *Nat. Rev. Cancer* **15**, 457–72.
4. Chester, C. *et al.* (2016) 4-1BB agonism: adding the accelerator to cancer immunotherapy. *Cancer Immunol Immunother.* **65(10)**, 1243–8.
5. Garvin, D. *et al.* (2018) Measurement of Fc10/19-mediated ADCC and CDC activity of anti-TNF α and anti-VEGF therapeutic antibodies using reporter-based bioassays and engineered TNF α + and VEGF+ target cells. *American Association for Cancer Research (AACR) Annual Meeting*, Poster abstract# 5628.

9. Appendix

9.A. Representative Assay Results with FcγRIIb-Dependent Antibody

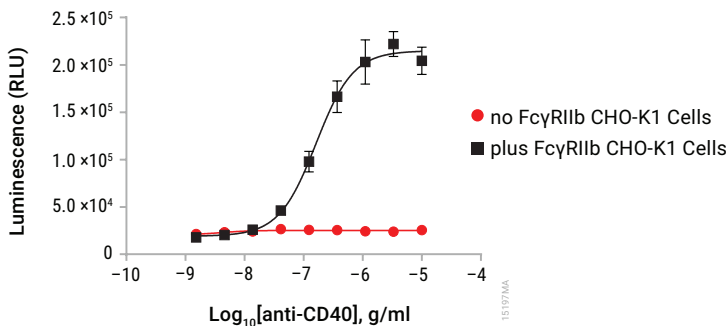


Figure 12. The CD40 Bioassay measures the activity of Control Ab, Anti-CD40. The day before assay, CD40 Effector Cells were plated and allowed to attach overnight. On the assay day, a titration of Control Ab, Anti-CD40, was added followed by the addition of FcγRIIb CHO-K1 Cells. After a 4-hour induction 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 EC₅₀ response determined was 0.163µg/ml, and the fold induction was 11.7.

9.B. Representative Assay Results with CD40 Ligand

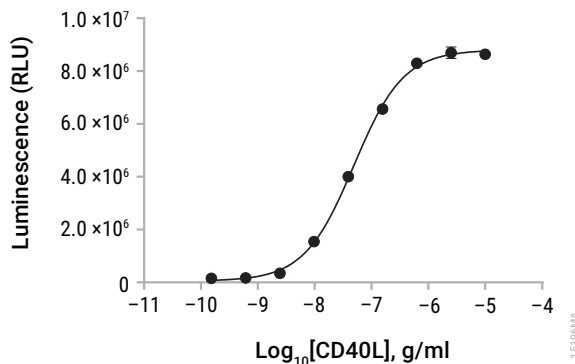


Figure 13. The CD40 Bioassay measures the activity of CD40 Ligand. The day before assay, CD40 Effector Cells were plated in a 96-well plate at 25,000 cells/well. On the day of assay, a serial dilution of CD40L was added to assay plates. After a 4-hour induction at 37°C, Bio-Glo™ Luciferase Assay Reagent was added and luminescence was determined using a GloMax® Discover System. Four-parameter logistic curve analysis was performed with GraphPad Prism® software. The EC₅₀ response determined was 49ng/ml, and the fold induction was 63-fold.

9.C. Composition of Buffers and Solutions

initial cell culture medium for CD40

Effector Cells

- 90% McCoy's 5A with L-glutamine
- 10% FBS

cell growth medium for CD40 Effector Cells

- 90% McCoy's 5A with L-glutamine
- 10% FBS
- 200µg/ml hygromycin B

cell freezing medium for CD40 Effector Cells

- 85% McCoy's 5A with L-glutamine
- 10% FBS
- 5% DMSO

cell recovery medium for CD40 Effector Cells

- 90% RPMI 1640 with L-glutamine
- 10% FBS assay buffer

assay buffer

- 99% RPMI 1640 with L-glutamine
- 1% FBS

9.D. 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 and sizes are available.

Immune Checkpoint Bioassays

Product	Size	Cat.#
4-1BB Bioassay	1 each	JA2351
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

Product	Size	Cat.#
PD-1/PD-L1 Blockade Bioassay	1 each	J1250
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 and sizes are available.

Macrophage-Directed Bioassays

Product	Size	Cat.#
SIRP α /CD47 Blockade Bioassay	1 each	JA6011
SIRP α /CD47 Blockade Bioassay, Fc-Dependent	1 each	JA4801
TLR Bioassay	1 each	JA9011
ADCP Reporter Bioassay (THP-1)	1 each	JA9411

Not for Medical Diagnostic Use. Additional kit formats and sizes 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 (TCR $\alpha\beta$ -KO, CD4+)	1 each	GA1172
T Cell Activation Bioassay (TCR $\alpha\beta$ -KO, CD8+)	1 each	GA1162
T Cell Activation Bioassay (TCR $\alpha\beta$ -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.

9.D. 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.

Detection Instruments

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 and Cytokine, Macrophage and Target Cell Killing Bioassays are available. To view and order Promega Bioassay products visit:

www.promega.com/products/reporter-bioassays/ or email: EarlyAccess@promega.com. For information on custom bioassay development and services visit the Promega Tailored R&D Solutions website:

www.promega.com/custom-solutions/tailored-solutions/

10. Summary of Changes

The following changes were made to the 6/25 revision of this document:

1. Removed an expired patent statement and updated another patent statement.
2. Updated text about the label in Section 3.
3. Made miscellaneous text edits.

©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 modify, derivatize, genetically engineer or otherwise create variations of the cells or genes stably transfected within the cells except that recipient may propagate and store the cells for long-term use. 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.

©U.S. Pat. No. 8,008,006.

©Licensed from Lonza Cologne GmbH under U.S. Pat. Nos. 7,700,357, 8,192,990 and 8,003,389, European Pat. Nos. 1297119, 1522587, 1607484 and 1741778 and other pending and issued patents.

©Patent Pending.

© 2019–2025 Promega Corporation. All Rights Reserved.

GloMax and Lumit are registered trademarks of Promega Corporation. Bio-Glo and Bio-Glo-NL are trademarks of Promega Corporation.

Accutase is a registered trademark of Innovative Cell Technologies, Inc. Corning and Costar are registered trademarks of Corning, Inc. Echo is a registered trademark of Labcyte, Inc. GIBCO is a registered trademark of Life Technologies Corporation. Mr. Frosty is a registered trademark of Nalge Nunc International Corporation. Styrofoam is a registered trademark of Dow Chemical Company. GraphPad Prism is a registered trademark of GraphPad Software, Inc. JMP is a registered trademark of SAS Institute, Inc. Multidrop is a trademark of Thermo Fisher Scientific, Ltd.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our website 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.