

TECHNICAL MANUAL

OX40 Bioassay

Instructions for use of Products
JA2191 and JA2195

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

OX40 Bioassay

All technical literature is available at: 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

1. Description.....	2
2. Product Components and Storage Conditions	8
3. Before You Begin.....	10
4. Assay Protocol for FcγRIIb-Dependent Antibodies.....	11
4.A. Preparing Assay Buffer, Bio-Glo™ Reagent and Antibody Samples.....	12
4.B. Plate Layout Design	14
4.C. Preparing and Plating FcγRIIb CHO-K1 Cells	15
4.D. Preparing and Plating OX40 Effector Cells.....	16
4.E. Preparing Antibody Serial Dilutions and Setting Up Assay	17
4.F. Adding Bio-Glo™ Reagent	19
4.G. Data Analysis	19
5. Assay Protocol for Ligand or FcγRIIb-Independent Antibodies	20
5.A. Preparing Assay Buffer, Bio-Glo™ Reagent and Antibody Samples.....	20
5.B. Plate Layout Design	22
5.C. Preparing and Plating OX40 Effector Cells.....	23
5.D. Preparing Ligand or Antibody Serial Dilutions and Setting Up Assay	24
5.E. Adding Bio-Glo™ Reagent	26
5.F. Data Analysis	26
6. Troubleshooting.....	27
7. References.....	28
8. Appendix.....	29
8.A. Representative Assay Results with FcγRIIb-Dependent Antibody.....	29
8.B. Representative Assay Results with OX40 Ligand	30
8.C. Related Products	30
9. 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 tumor necrosis factor receptor family-related protein (GITR), OX40, 4-1BB, 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).

OX40 (CD134/TNFRSF4), a member of the tumor necrosis factor (TNF) receptor superfamily, is a costimulatory receptor expressed primarily on activated T cells, and on neutrophils and natural killer (NK) cells to a lesser extent (4). When present on the cell surface, OX40 interacts with OX40 ligand (OX40L), and induces subsequent cell proliferation, survival and production of cytokines, particularly in T cells (5,6).

Current methods used to measure the activity of biologic drugs targeting OX40 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 quality controlled drug-development setting.

The OX40 Bioassay^(a-f) (Cat.# JA2191, JA2195), is a bioluminescent cell-based assay that overcomes the limitations of existing assays. It can be used to measure the potency and stability of ligands or agonist antibodies that can bind and activate OX40 (5,7). The assay consists of a genetically engineered Jurkat T cell line that expresses human OX40 and a luciferase reporter driven by a response element that can respond to OX40 ligand/agonist antibody stimulation. The OX40 Effector Cells 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 OX40 Bioassay should be conducted with Fc γ RIIb CHO-K1 Cells (Cat.# JA2251, JA2255) to test whether agonist antibodies activate OX40 in an Fc γ RIIb-dependent manner. Fc γ RIIb 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 costimulatory immune checkpoints, you perform the assay both with and without Fc γ RIIb CHO-K1 cells to ascertain the need for these cells in enhancing the effect of the agonist antibodies raised against the costimulatory immune checkpoint targets.

Induction of the OX40 Effector Cells with an OX40 ligand or agonist antibody results in response element-mediated luminescence (Figure 1). The bioluminescent signal is quantified using the Bio-Glo™ Luciferase Assay System^(f) and a standard luminometer such as the GloMax® Discover System (see Section 8.C, Related Products).

The OX40 Bioassay reflects the mechanism of action (MOA) of biologics designed to activate OX40. Specifically, OX40-mediated luminescence is detected following the addition of OX40 ligand and OX40 agonist antibodies (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, 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 100% human serum (in ligand or antibody samples; Figure 5), indicating potential for further development into a neutralizing antibody bioassay.

1. Description (continued)

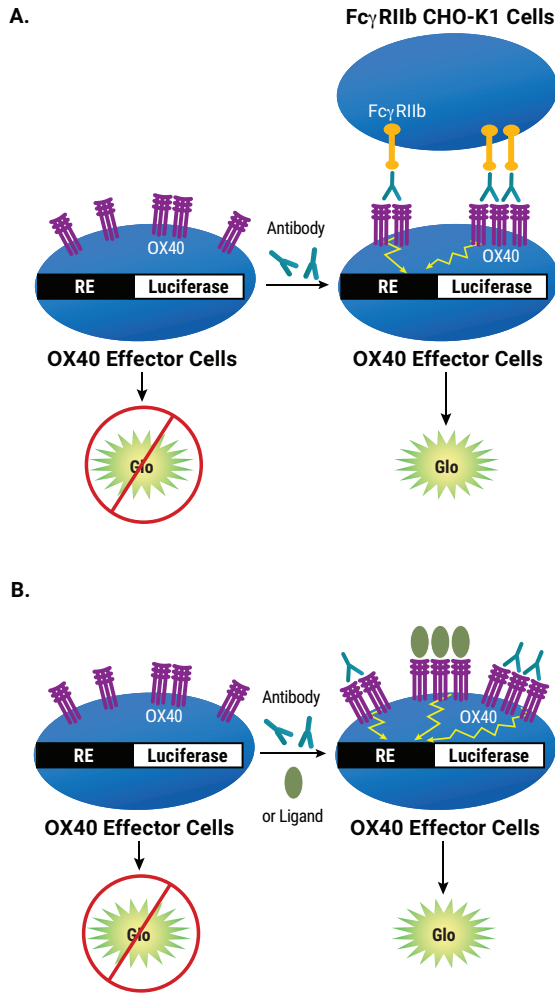


Figure 1. Representation of the OX40 Bioassay. Panel A. Assay with Fc γ RIIb-dependent agonist antibody. The bioassay consists of two engineered cell lines, OX40 Effector Cells and Fc γ RIIb CHO-K1 Cells. In the presence of Fc γ RIIb CHO-K1 Cells, the anti-OX40 antibody can be crosslinked, thereby inducing OX40 pathway-activated luminescence. **Panel B.** Assay with Fc γ RIIb-independent agonist antibody or ligand. The bioassay consists of one engineered cell line, OX40 Effector Cells. In the absence of agonist antibody or OX40 ligand, the OX40 receptor is not activated and the luminescence signal is low. The addition of agonist antibody or OX40 ligand induces the OX40 pathway-activated luminescence, which can be detected in a dose-dependent manner.

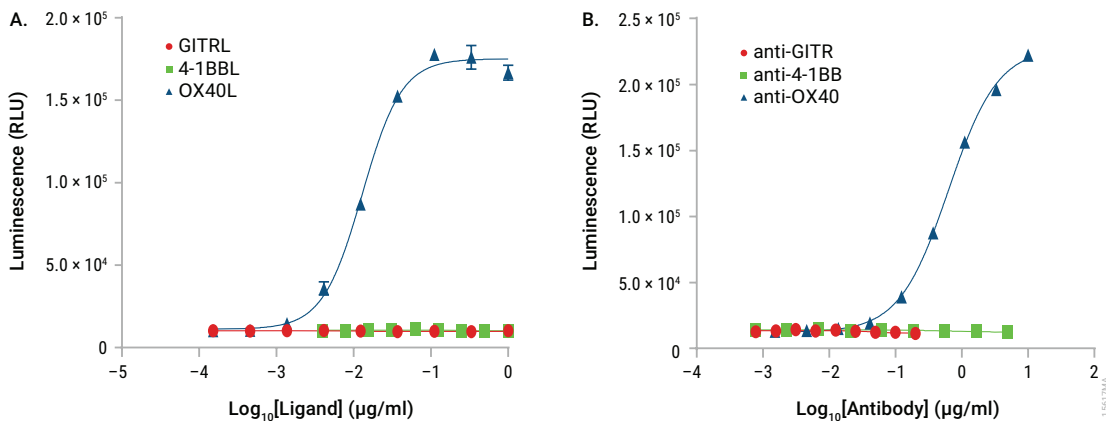


Figure 2. The OX40 Bioassay reflects the mechanism of action (MOA) and shows specificity of biologics designed to activate OX40. Panel A. OX40 Effector Cells were induced with a serial titration of ligands: GITRL crosslinked by anti-HA Ab; 4-1BBL crosslinked by anti-His Ab; or OX40L, as indicated. **Panel B.** OX40 Effector Cells were induced with a serial titration of anti-GITR antibody, anti-4-1BB antibody or anti-OX40 antibody, as indicated, in the presence of FcγRIIb CHO-K1 Cells (Cat. # JA2251, JA2255). After a 5-hour induction, Bio-Glo™ Reagent^(f) was added, and luminescence was quantified using the GloMax® Discover System. Data were generated using thaw-and-use cells and fitted to a four-parameter logistic curve using GraphPad Prism® software.

Table 1. The OX40 Bioassay Shows Precision, Accuracy and Linearity.

Parameter	Results	
	% Expected Relative Potency	% Recovery
Accuracy	50	99.4
	70	100.8
	140	101.2
	200	109.9
Repeatability (% CV)	100% (Reference)	5.4
Intermediate Precision (% CV)		9.0
Linearity (r ²)		0.996
Linearity (y = mx + b)		y = 1.123x – 8.895
A 50–200% theoretical potency series of Control Ab, Anti-OX40, was analyzed in triplicate in three independent experiments performed on three days by two analysts. Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were analyzed and relative potencies were calculated after parallelism determination using JMP® software. Data were generated using thaw-and-use cells.		

1. Description (continued)

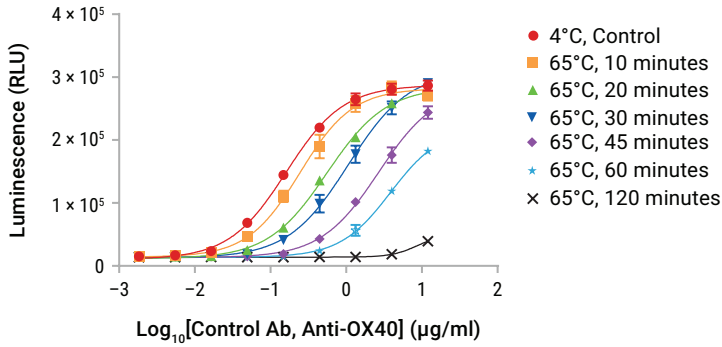


Figure 3. The OX40 Bioassay is stability-indicating. Samples of Control Ab, Anti-OX40 (Cat.# K1191), were maintained at 4°C (control) or heat-treated at 65°C for the indicated times, then analyzed using the OX40 Bioassay with FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255). After a 5-hour induction, Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were generated using thaw-and-use cells and fitted to a four-parameter logistic curve using GraphPad Prism® software.

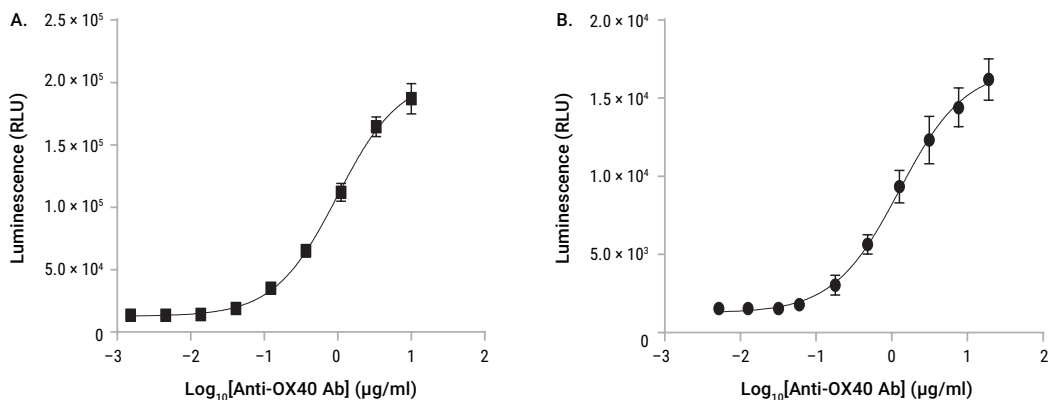


Figure 4. The OX40 Bioassay is amenable to 384-well plate format and compatible with laboratory automation. Panel A. The OX40 Bioassay was performed in 96-well plates as described in this technical manual using anti-OX40 antibody. **Panel B.** The OX40 Bioassay was performed in 384-well format using a Mantis[®] liquid handler to dispense the cells and Echo[®] Acoustic liquid handler for antibody handling. On the day before the assay, FcγRIIb CHO-K1 Cells (Cat. # JA2251, JA2255) were plated at 8×10^3 cells/10µl/well and incubated at 37°C for 5 hours. OX40 Effector Cells were then added to the plate at 1×10^4 cells/10µl/well to co-incubate with FcγRIIb CHO-K1 Cells at 37°C overnight. On the day of the assay, anti-OX40 antibody was serially diluted and added to the plate at 0.2µl/well. After a 5-hour induction, 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 0.99µg/ml and 1.26µg/ml, and the fold inductions were 15 and 11 for 96-well and 384-well format, respectively. Data were generated using thaw-and-use cells.

1. Description (continued)

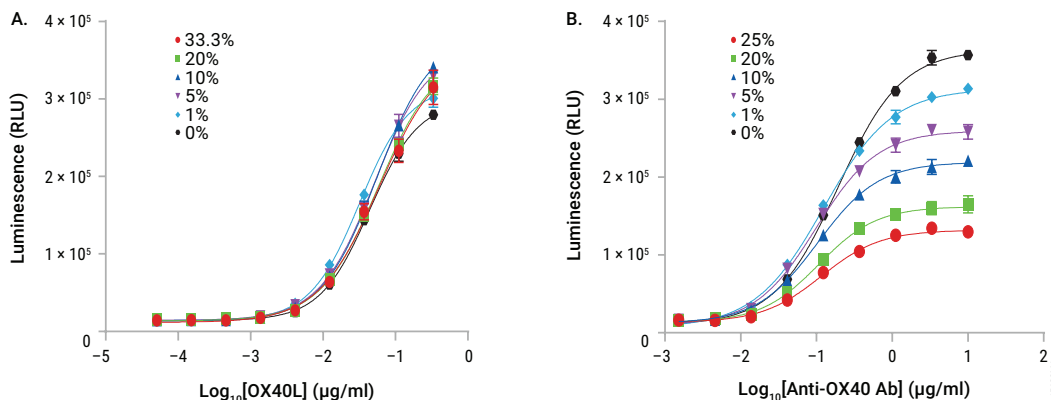


Figure 5. The OX40 Bioassay is tolerant to human serum. **Panel A.** OX40 ligand (OX40L, Biologend Cat.# 555704) was analyzed in the presence of increasing concentrations of pooled normal human serum (0–100% in the ligand sample), resulting in final assay concentration of human serum (0–33.3%). **Panel B.** Anti-OX40 antibody was analyzed in the presence of FcγRIIb 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 (0–25%). After a 5-hour induction at 37°C, 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 OX40 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.#
OX40 Bioassay	1 each	JA2191

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 OX40 Effector Cells (0.5ml)
- 36ml RPMI 1640 Medium
- 4ml Fetal Bovine Serum
- 1 vial Bio-Glo™ Luciferase Assay Substrate (lyophilized)
- 10ml Bio-Glo™ Luciferase Assay Buffer

PRODUCT	SIZE	CAT.#
OX40 Bioassay 5X	1 each	JA2195

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 OX40 Effector Cells (0.5ml)
- 5 × 36ml RPMI 1640 Medium
- 5 × 4ml Fetal Bovine Serum
- 5 vials Bio-Glo™ Luciferase Assay Substrate (lyophilized)
- 5 × 10ml Bio-Glo™ Luciferase Assay Buffer

Note: OX40 Bioassay components are shipped separately because of different temperature requirements. OX40 Effector Cells are shipped on dry ice. 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°C (freezer or liquid nitrogen vapor phase). 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.
- Store Bio-Glo™ Luciferase Assay Substrate, Bio-Glo™ Luciferase Assay Buffer and Fetal Bovine Serum at –30°C to –10°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 –20°C for up to 6 weeks.
- Store RPMI 1640 Medium at 4°C protected from fluorescent light.

Items Available Separately

PRODUCT	SIZE	CAT.#
FcγRIIb CHO-K1 Cells	1 each	JA2251

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 FcγRIIb Cells (0.5ml)
- 36ml RPMI 1640 Medium

PRODUCT	SIZE	CAT.#
FcγRIIb CHO-K1 Cells 5X	1 each	JA2255

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 FcγRIIb Cells (0.5ml per vial)
- 5 × 36ml RPMI 1640 Medium



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 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 OX40 Bioassay is intended to be used with user-provided antibodies or ligands designed to activate OX40. Control Ab, Anti-OX40 (Cat.# K1191) and FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255) are available separately for use in assay optimization and routine quality control. We strongly recommend including OX40 Ligand or Control Ab, Anti-OX40, and 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 OX40 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 appropriate concentration for the assay. The cells are sensitive, and care should be taken to follow cell thawing and plating procedures as described.

The OX40 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[®] Discover System (see Section 8.C, Related Products). An integration time of 0.5 seconds/well was used for all readings.

Materials to Be Supplied by the User

- user-defined anti-OX40 antibodies or other biologics samples
- solid-white, flat-bottom 96-well assay plates (e.g., Corning® Cat.# 3917) for plating and reading luminescence
- 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 125ml media bottles (e.g., Fisher Scientific Cat.# 03-313-901)
- sterile reagent reservoirs (e.g., Costar®/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 system)
- **optional:** solid-white, flat-bottom 384-well assay plates (e.g., Corning® Cat.# 3570)
- **optional:** FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255; We recommend this cell line when using this assay for the first time and/or testing an Ab that may be dependent on crosslinking FcγRIIb.)
- **optional:** Control Ab, Anti-OX40 (Cat.# K1191)
- **optional:** OX40 ligand (OX40L, Biologend Cat.# 555704)

4. Assay Protocol for FcγRIIb-Dependent Antibodies

This assay protocol requires two engineered cell lines: OX40 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.# J2232; see *FcγRIIb CHO-K1 Propagation Model Technical Manual*, #TM569, for details). Either cell format can be used in this assay.

The procedure below illustrates the use of the OX40 Bioassay to test two FcγRIIb-dependent antibody samples against a reference sample in a single assay run using the FcγRIIb CHO-K1 Cells in 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 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-OX40, which is FcγRIIb-dependent and requires the use of FcγRIIb CHO-K1 Cells.

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

1. **Assay Buffer:** On the day before assay, prepare 70ml of assay buffer (95% RPMI 1640/5%FBS) in a 125ml media bottle. Thaw the FBS overnight at 4°C or in a 37°C water bath on the day of use. Add 3.5ml of FBS to 66.5ml of RPMI 1640 medium. Mix well and warm to 37°C before use. For reference, 70ml 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 5% FBS. This concentration of FBS works well for the Control Ab, Anti-OX40, that we tested. If assay performance is not ideal using 5% FBS assay buffer when testing your antibody of interest, try optimizing the FBS concentration from 0.5–10% in assay buffer.

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. When stored at ambient temperature, the reconstituted Bio-Glo™ Reagent will lose ~18% signal after 24 hours.

3. **Test and Reference Samples:** Using assay buffer as the diluent, prepare starting dilutions (dilu1, 4X 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-OX40 (Cat.# K1191), as a reference antibody in your assay, prepare a 400µl starting dilution with 40µg/ml of anti-OX40 antibody (dilu1, 4X final concentration) by adding 16µl of anti-OX40 stock (1,000µg/ml) to 384µl of assay buffer. Store the antibody starting dilution on ice until ready to use in the assay.

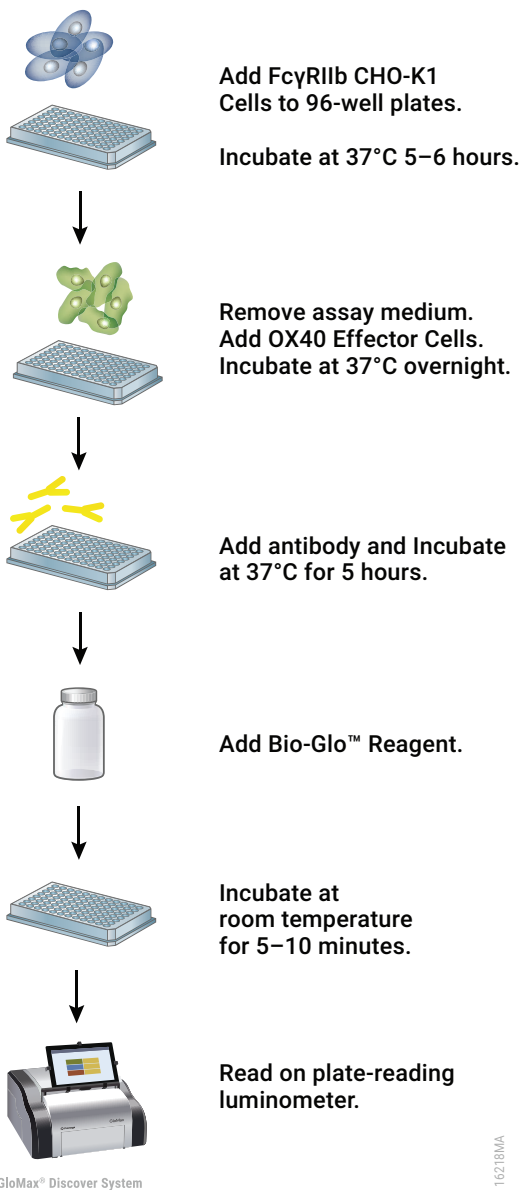


Figure 6. Schematic protocol for OX40 Bioassay with Fc γ RIIb-dependent antibody.

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 alone (denoted by “B”).

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

The thaw-and-use FcγRIIb CHO-K1 Cells (Cat. # JA2251, JA2255) are available separately. These cells 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.

Note: Perform the following steps using aseptic technique in a sterile cell culture hood early in the morning on the day before the assay.

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

1. On the day before the assay, prepare 70ml of fresh assay buffer (RPMI/5% FBS) as described in Section 4.A. Divide the assay buffer into three aliquots, one for FcγRIIb CHO-K1 preparation (30ml), one for OX40 Effector Cells preparation (30ml) and one for antibody serial dilution preparation (10ml). Store two aliquots (30ml and 10ml) at 4°C for later use.
2. Warm the assay buffer (30ml aliquot) in a 37°C water bath for 15 minutes then transfer 14.5ml of assay buffer 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 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 assay buffer. 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 the cell suspension to each of the inner 60-wells of two 96-well, white, flat-bottom assay plates.
6. Add 100µl of assay buffer per well to the outside wells of the assay plates.
7. Cover the assay plates with a lid and incubate the cells in a 37°C, 5% CO₂ incubator for 5–6 hours.

4.D. Preparing and Plating OX40 Effector Cells

Note: The thaw-and-use OX40 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. Warm the assay buffer (30ml aliquot) in a 37°C water bath for 15 minutes and transfer 11.5ml of assay buffer to a 15ml conical tube.
2. Remove one vial of thaw-and-use OX40 Effector Cells from –140°C storage and transfer to the 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 the vial.
3. Gently mix the cell suspension in the vial by pipetting, then transfer 0.5ml of cells to the 15ml conical tube containing 11.5ml of assay buffer. Mix well by gently inverting 1–2 times.
4. Remove the 96-well assay plates containing FcγRIIb CHO-K1 Cells from the incubator. Using a manual multichannel pipette, remove 95μl of medium from each of the wells. Alternatively, invert the assay plate above a sink to remove the medium. Then place the inverted plate on paper towels for 5–10 seconds to drain any remaining medium.
5. Transfer the OX40 Effector Cells prepared in Step 3 to a sterile reagent reservoir. Immediately, using a multichannel pipette, dispense 60μl of cell suspension to each of the inner 60-wells of the plates containing preplated FcγRIIb CHO-K1 Cells.
6. Add 80μl of assay buffer to the outside wells of the 96-well assay plates.
7. Cover the assay plates with lids and incubate in a 37°C, 5% CO₂ incubator overnight.

4.E. Preparing Antibody Serial Dilutions and Setting Up Assay

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 400µl of reference antibody at 4X the highest antibody concentration in your dose-response curve. You will need 180µl of each test antibody at 4X 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-OX40, as a control in the assay, follow the instructions below to prepare threefold serial dilutions. A threefold serial dilution for test antibodies was listed as an example below as well.

1. On the day of assay, warm the assay buffer (10ml aliquot, described in Section 4.C, Step 1) in a 37°C water bath for 15 minutes.
2. To a sterile, clear V-bottom 96-well plate, add 180µl of reference antibody starting dilution (dilu1, 4X final concentration) to wells A11 and B11 (see Figure 8).
3. Add 180µl of test antibodies 1 and 2 starting dilution (dilu1, 4X 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 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. Remove the preplated FcγRIIb CHO-K1 and OX40 Effector Cells from the incubator. Using an electronic multichannel pipette, add 20µl of the appropriate antibody titration to the assay plates according to the plate layout in Figure 7.
8. Cover the plate with a lid and incubate in a 37°C, 5% CO₂ incubator for 5 hours.

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

4.E. Preparing Antibody Serial Dilutions and Setting Up Assay (continued)

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.

4.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 80µl of Bio-Glo™ Reagent to the inner 60 wells of the assay plates, taking care not to create bubbles.
3. Add 80µl 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.

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 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).

5. Assay Protocol for Ligand or FcγRIIb-Independent Antibodies

This assay protocol illustrates the use of the OX40 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 threefold dilution when testing an OX40 ligand.

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

1. **Assay Buffer:** On the day before the assay, prepare 36ml of assay buffer (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 1.8ml of FBS to 34.2ml of RPMI 1640 medium. Mix well and warm to 37°C before use. For reference, 36ml 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 5% FBS. This concentration of FBS works well with the OX40 ligand that we tested.

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. When stored at ambient temperature, the reconstituted Bio-Glo™ Reagent will lose ~18% signal after 24 hours.

3. **Test and Reference Samples:** Using assay buffer as the diluent, prepare starting dilutions (dilu1, 4X final concentration) of two test antibodies (180 μl each) and one reference ligand (400 μl) in 1.5ml tubes. Store the tubes containing ligand or antibody starting dilutions appropriately before making antibody serial dilutions.

Note: If you are using OX40 ligand (OX40L, Biolegend Cat.# 555704) as a reference ligand, prepare a 400 μl starting dilution with 4 μg/ml of OX40L (dilu1, 4X final concentration) by adding 8 μl of OX40L stock (200 μg/ml) to 392 μl of assay buffer. The final (1X) starting concentration is 1 μg/ml of OX40L. Store the ligand starting dilution on ice until ready to use in the assay.

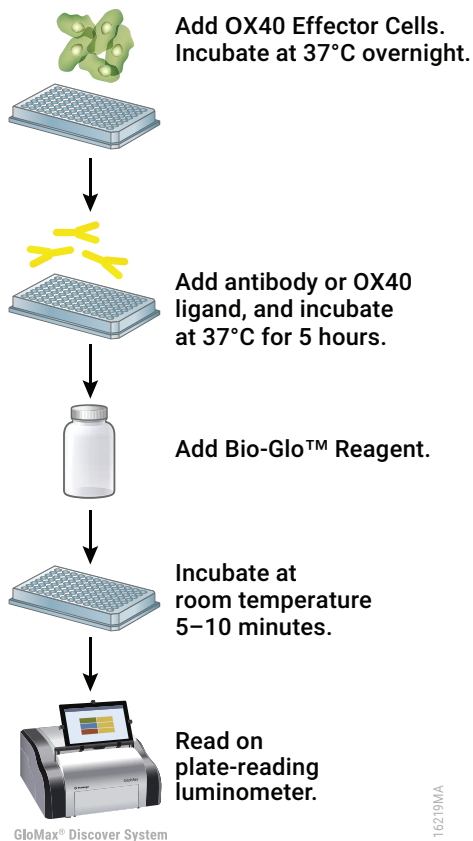


Figure 9. Schematic protocol for OX40 Bioassay with OX40 ligand or FcγRIIb-independent antibody.

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 alone (denoted by “B”).

5.C. Preparing and Plating OX40 Effector Cells

Note: The thaw-and-use OX40 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. On the day before the assay, prepare 36ml of assay buffer (95% RPMI 1640/5% FBS) as described in Section 5.A. Divide the assay buffer into two aliquots, one for OX40 Effector Cells preparation (26ml) and the other for antibody serial dilution preparation (10ml). Store the 10ml aliquot at 4°C for later use.
2. Warm the assay buffer (26ml aliquot) in a 37°C water bath for 15 minutes and transfer 11.5ml to a 15ml conical tube.
3. Remove one vial of thaw-and-use OX40 Effector Cells from –140°C storage and transfer to the 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, then transfer 0.5ml of cells to the 15ml conical tube containing 11.5ml of assay buffer. Mix well by gently inverting 1–2 times.
5. Transfer the OX40 Effector Cell suspension to a sterile reagent reservoir. Immediately, using a multichannel pipette, dispense 60µl of cell suspension to each well of the inner 60 wells of two 96-well, white, flat-bottom assay plates. .
6. Add 80µl of assay buffer to the outside wells of the 96-well assay plates.
7. Cover the assay plates with lids and incubate in a 37°C, 5% CO₂ incubator overnight.

5.D. Preparing Ligand or Antibody Serial Dilutions and Setting Up 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 ligand, you will need 400µl of reference ligand at 4X 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 4X 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 OX40L (see Section 3) as a control in the assay, use the following instructions to prepare threefold serial dilutions. A threefold serial dilution for test antibodies is shown in Figure 11.

1. On the day of the assay, warm the 10ml assay buffer aliquot prepared in Section 5.C., Step 1.
2. To a sterile, clear V-bottom 96-well plate, add 180µl of reference ligand starting dilution (dilu1, 4X final concentration) to wells A11 and B11 (see Figure 11).
3. Add 180µl of test antibodies 1 and 2 starting dilution (dilu1, 4X final concentration) to wells E11 and G11, respectively (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 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. Remove the preplated OX40 Effector Cells from the incubator. Using an electronic multichannel pipette, add 20µl of the OX40L or appropriate antibody titration to the assay plates according to the plate layout in Figure 10.
8. Cover the plate with a lid and incubate in a 37°C, 5% CO₂ incubator for 5 hours.

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 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 reference ligand and antibody serial dilutions.

5.E. 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 80µl of Bio-Glo™ Reagent to the inner 60 wells of the assay plates, taking care not to create bubbles.
3. Add 80µl 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.F. Data Analysis

1. Determine the plate background by calculating the average RLU from wells B1, C1 and D1.
2. Calculate fold induction =
$$\frac{\text{RLU (sample-background)}}{\text{RLU (no drug control-background)}}$$
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. 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> <hr/> <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> <hr/> <p>If performing the assay for the first time, we recommend that you use FcγRIIb CHO-K1 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>Inappropriate cell handling during cell thawing, including long water bath incubation times, can cause low assay performance and high assay variation.</p>
Weak assay response (low fold induction)	<p>Low activity of Bio-Glo™ Reagent leads to low RLU. Store and handle the Bio-Glo™ Reagent according to the instructions.</p> <hr/> <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 OX40 Bioassay may vary from the EC₅₀ value obtained using other methods such as primary T cell-based assays.</p> <hr/> <p>Determine if the antibody used is dependent on crosslinking by testing in the presence of FcγRIIb Cells.</p> <hr/> <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>



7. References

1. Mahoney, K.M. *et al.* (2015) Combination cancer immunotherapy and new immunomodulatory targets. *Nat. Rev. Drug Disc.* **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. Buchan, S.L. *et al.* (2018) The immunobiology of CD27 and OX40 and their potential as targets for cancer immunotherapy. *Blood* **131(1)**, 39–48.
4. Willoughby, J. *et al.* (2017) OX40: Structure and function - What questions remain? *Mol. Immunol.* **83**, 13–22.
5. Croft, M. *et al.* (2009) The significance of OX40 and OX40L to T-cell biology and immune disease. *Immunol. Rev.* **229 (1)**, 173–91.
6. Ishii, N. *et al.* (2010) OX40-OX40 ligand interaction in T-cell-mediated immunity and immunopathology. *Adv Immunol.* **105**, 63–98.
7. 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

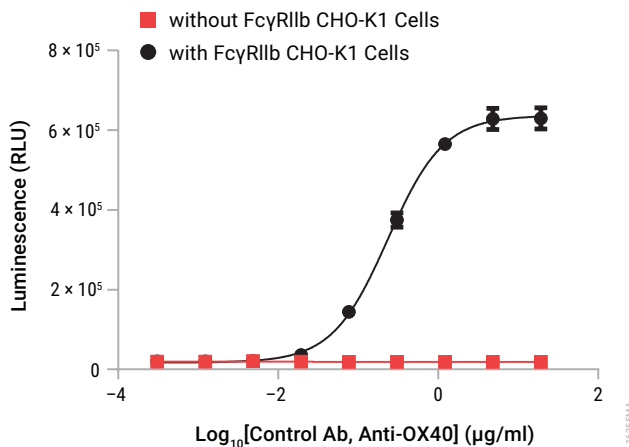


Figure 12. The OX40 Bioassay measures the activity of Control Antibody, Anti-OX40. Thaw-and-use FcγRIIb CHO-K1 Cells (Cat.# JA2251, JA2255) and OX40 Effector Cells (Cat.# JA2191, JA2195) were thawed and plated as instructed, and incubated overnight. The following day, a titration of Control Ab, Anti-OX40, (Cat.# K1191; Section 4.C) was added. After 5 hours of induction at 37°C, Bio-Glo™ 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 ~0.24µg/ml, and the fold induction was ~30.

8.B. Representative Assay Results with OX40 Ligand

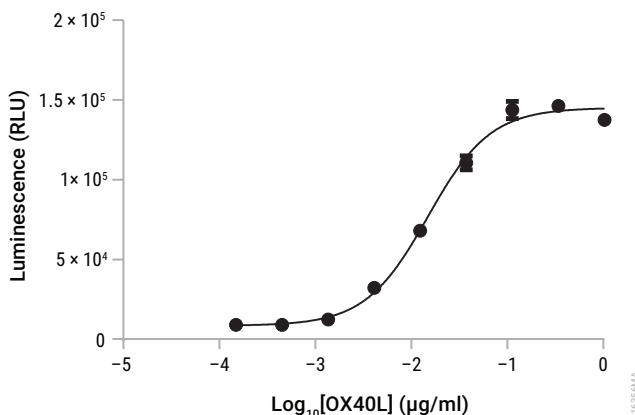


Figure 13. The OX40 Bioassay measures the activity of OX40 Ligand. Thaw-and-use OX40 Effector Cells were plated overnight. The following day, cells were incubated with various concentrations of OX40 ligand (Section 5.C). After 5 hours of induction at 37°C, Bio-Glo™ Reagent was added, and luminescence was determined using a GloMax® Discover System. Four-parameter logistic curve analysis was performed with GraphPad Prism® software. The EC₅₀ response determined was ~0.015µg/ml, and the fold induction was ~15.2.

8.C. Related Products

Fc Effector Bioassays

Product	Size	Cat.#
ADCC Reporter Bioassay, Complete Kit (Raji)*	1 each	G7015
ADCC Reporter Bioassay, Target Kit (Raji)*	1 each	G7016
ADCC Reporter Bioassay, Core Kit*	1 each	G7010
ADCC Reporter Bioassay, F Variant, Core Kit**	1 each	G9790
FcγRIIa-H ADCP Reporter Bioassay, Complete Kit**	1 each	G9901
FcγRIIa-H ADCP Reporter Bioassay, Core Kit**	1 each	G9991

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

**Not for Medical Diagnostic Use.

Additional kit formats are available.

Immune Checkpoint Bioassays

Product	Size	Cat.#
CD40 Bioassay	1 each	JA2151
CD40 Bioassay 5X	1 each	JA2155
CD40 Bioassay, Propagation Model	1 each	J2132
LAG-3/MHCII Blockade Bioassay	1 each	JA1111
LAG-3/MHCII Blockade Bioassay 5X	1 each	JA1115
LAG-3/MHCII Blockade Bioassay, Propagation Model	1 each	JA1112
OX40 Bioassay, Propagation Model	1 each	J2172
4-1BB Bioassay	1 each	JA2351
4-1BB Bioassay 5X	1 each	JA2355
4-1BB Bioassay, Propagation Model	1 each	J2332
PD-1/PD-L1 Blockade Bioassay	1 each	J1250
PD-1/PD-L1 Blockade Bioassay 5X	1 each	J1255
PD-L1 Negative Cells	1 each	J1191
CTLA-4 Blockade Bioassay	1 each	JA3001
CTLA-4 Blockade Bioassay 5X	1 each	JA3005
TIGIT Negative Cells	1 each	J1921
PD-1+TIGIT Combination Bioassay	1 each	J2211
PD-1+TIGIT Combination Bioassay 5X	1 each	J2215
Control Ab, Anti-CTLA-4	100µg	JA1020
Control Ab, Anti-PD-1	100µg	J1201
Control Ab, Anti-TIGIT	100µg	J2051
Control Ab, Anti-4-1BB	50µg	K1161
Control Ab, Anti-OX40	50µg	K1191
Control Ab, Anti-GITR	50µg	K1171
Control Ab, Anti-CD40	50µg	K1181

Not for Medical Diagnostic Use.

Additional kit formats are available.



8.C. Related Products (continued)

T Cell Activation Bioassays

Product	Size	Cat.#
T Cell Activation Bioassay (NFAT)	1 each	J1621
T Cell Activation Bioassay (NFAT) 5X	1 each	J1625
T Cell Activation Bioassay (NFAT), Propagation Model	1 each	J1601
T Cell Activation Bioassay (IL-2)	1 each	J1651
T Cell Activation Bioassay (IL-2) 5X	1 each	J1655
T Cell Activation Bioassay (IL-2), Propagation Model	1 each	J1631

Not for Medical Diagnostic Use.

Cytokine and Growth Factor Bioassays

Product	Size	Cat.#
IL-2 Bioassay	1 each	JA2201
IL-2 Bioassay 5X	1 each	JA2205
IL-2 Bioassay, Propagation Model	1 each	J2952
IL-15 Bioassay	1 each	JA2011
IL-15 Bioassay 5X	1 each	JA2015
IL-15 Bioassay, Propagation Model	1 each	J2962
VEGF Bioassay	1 each	GA2001
VEGF Bioassay 5X	1 each	GA2005
VEGF Bioassay, Cell Propagation Model	1 each	GA1082
Recombinant VEGF	10µg	J2371

Not for Medical Diagnostic Use.

Detection Reagent

Product	Size	Cat.#
Bio-Glo™ Luciferase Assay System	10ml	G7941
	100ml	G7940

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 and Cytokine Bioassays are available from Promega Custom Assay Services. See the Early Access listings at:

www.promega.com/applications/biologics-drug-discovery/functional-bioassays/target-pathway-assays/
 or email: **CAS@promega.com**

9. Summary of Changes

The following changes were made to the 4/22 revision of this document:

1. In Section 2, “long-term storage” was removed from the storage conditions for the cells and storage temperature ranges were updated to match those on the product labels.
2. The cover image was updated.



^(b)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.

^(b)U.S. Pat. No. 8,008,006 and European Pat. No. 1341808.

^(c)Product cannot be used for proficiency testing.

^(d)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.

^(e)Patents Pending.

^(f)U.S. Pat. No. 10,077,244 and other patents.

© 2019, 2022 Promega Corporation. All Rights Reserved.

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

Corning and Costar are registered trademarks 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. 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 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.