

TECHNICAL MANUAL

ICOS Bioassay, Core Kit

Instructions for Use of Products
JA6801 and JA6805

ICOS Bioassay, Core Kit

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	9
3. Before You Begin	10
3.A. Materials to Be Supplied by the User	11
4. Assay Protocol for FcγRIIb-Dependent or FcγRIIb-Independent Antibodies	11
4.A. Preparing Cell Plating Medium, Assay Buffer, Bio-Glo-NL™ Reagent and Antibody Samples	13
4.B. Plate Layout Design	14
4.C. Plating FcγRIIb aAPC/CHO-K1 Cells	15
4.D. Preparing Antibody Serial Dilutions	16
4.E. Preparing ICOS Effector Cells	17
4.F. Adding ICOS Effector Cells and Antibody to Assay Plates	18
4.G. Preparing and Adding Bio-Glo-NL™ Reagent	18
4.H. Data Analysis	19
5. Troubleshooting	20
6. References	21
7. Appendix	22
7.A. Representative Assay Results	22
7.B. Related Products	22
8. Summary of Changes	25

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. T cells play a central role in cell-mediated immunity against pathogens; however, T cells also contribute to the pathogenesis and exacerbation of autoimmune disorders. Costimulatory T cell receptors have been identified, which upon activation, can enhance T cell-mediated immune responses. Activating these receptors with agonist monoclonal antibodies has emerged as a strategy to enhance T cell responses in cancer or during persistent infections.

Optimal activation of T cells is initiated by engagement of the T cell antigen receptor (TCR)/CD3 complex and activation of a costimulatory receptor such as Inducible T Cell Co-Stimulator (ICOS, CD278; 1–3). ICOS binds to its ligand ICOSL (B7RP-1, CD275), which is constitutively expressed on B cells, monocytes, and dendritic cells, and can be induced on endothelial and epithelial cells during inflammation (4–7). ICOS costimulation induces the production of effector T cell cytokines such as interferon (IFN)- γ , interleukin-4 (IL-4) and IL-10 (1–3).

Quantitative functional bioassays are needed in the development of biologic drugs designed to activate ICOS. Current methods rely on primary human T cells and antigen-presenting cells (APCs), and measurement of functional endpoints such as cell proliferation, cell surface marker expression and cytokine production. These assays are laborious and highly variable due to their reliance on donor cells, complex assay protocols and unqualified assay reagents. Current methods are, as a result, difficult to establish in a quality-controlled setting.

The ICOS Bioassay, Core Kit^(a-e) (Cat.# JA6801, JA6805), is a bioluminescent cell-based assay that overcomes the limitations of existing assays. It can be used to measure the potency and stability of antibodies and other biologics that activate ICOS (8). The bioassay consists of the following genetically engineered cell lines:

- **ICOS Effector Cells:** Jurkat T cells expressing ICOS and endogenous TCR/CD3, and a NanoLuc[®] (NL) luciferase reporter driven by ICOS and TCR/CD3 pathway-dependent response elements.

The ICOS Bioassay should be performed with one of the following genetically engineered cell lines:

- **Fc γ RIIb aAPC/CHO-K1 Cells (Cat.# JA9331, JA9335):** CHO-K1 cells engineered to express a proprietary TCR-activating protein as well as human Fc γ RIIb. Fc γ RIIb serves as a crosslinking receptor that binds to antibody Fc domains and clusters agonist antibodies for maximal agonist activity.
- **aAPC/CHO-K1 Cells (Cat.# JA9441, JA9445):** CHO-K1 cell engineered to express an engineered cell-surface protein designed to activate the TCR complex in an antigen-independent manner. Can be substituted for use with Fc γ RIIb-independent antibodies or other biologics that do not require crosslinking for optimal activity.

The ICOS Effector Cells, Fc γ RIIb aAPC/CHO-K1 Cells and aAPC/CHO-K1 Cells are provided in thaw-and-use format, eliminating the need for cell culture and propagation.

In addition to the ICOS Bioassay Core Kit, Fc γ RIIb aAPC/CHO-K1 and aAPC/CHO-K1 Cells, we offer Control Ab, Anti-ICOS (Cat.# K1241), for use as a positive control for assay optimization and routine quality control. When screening for ICOS agonist antibodies, we recommend performing the assay with Fc γ RIIb aAPC/CHO-K1 Cells in parallel with aAPC/CHO-K1 Cells in order to determine the role of Fc γ RIIb-mediated crosslinking in enhancing the effect of agonist antibodies.

Inducing the ICOS Effector Cells with an ICOS agonist antibody results in increased promoter-mediated luminescence (Figure 1). The bioluminescent signal is quantified using the Bio-Glo-NL™ Luciferase Assay System^(a,c,d) (Cat.# J3081) and a standard luminometer such as the GloMax® Discover System (see Related Products, Section 7.B).

The ICOS Bioassay reflects the mechanism of action (MOA) of biologics designed to activate ICOS. Specifically, ICOS-mediated luminescence activation is increased following the addition of an ICOS activating antibody but not following addition of anti-4-1BB or anti-OX40 agonist antibodies (Figure 2). The bioassay is prequalified following International Council for Harmonisation of Technical Requirements for Pharmaceuticals for Human Use (ICH) guidelines and shows the precision, accuracy and linearity required for routine use in potency and stability studies (Table 1 and Figure 3). The bioassay is performed in a two-day timeframe, and the 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 antibody samples; Figure 5), indicating potential for further development into a neutralizing antibody bioassay.

It is increasingly common during drug development to analyze potential therapeutic antibodies for Antibody-Dependent Cell-Mediated Cytotoxicity (ADCC) activity. Another application of the ICOS Bioassay is the ability to measure ADCC activity of anti-ICOS antibodies by combining ADCC Reporter Bioassay Effector cells, available separately (Cat. #G7010), with ICOS Effector Cells used as target cells in this application (Figure 6).

Blockade of ICOS is a separate immunotherapy strategy to reduce autoimmunity and alloimmunity. The ICOS Bioassay is not designed to detect blocking activity of ICOS antibodies. We offer separately an ICOS Blockade Bioassay (Cat.# JA6001, JA6005), which can be used for screening and potency testing of ICOS antagonists.

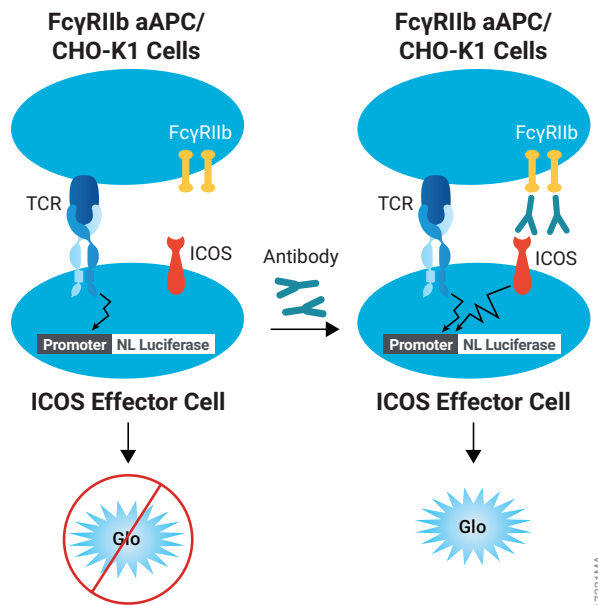


Figure 1. Representation of the ICOS Bioassay. The ICOS Bioassay for FcγRIIb-dependent agonist antibodies consists of two cell lines, ICOS Effector Cells and FcγRIIb aAPC/CHO-K1 Cells. In the absence of agonist antibody, ICOS is not activated and the luminescence signal is low. In the presence of FcγRIIb aAPC/CHO-K1 Cells, anti-ICOS antibody can be crosslinked, thereby inducing ICOS pathway-activating luminescence in a dose-dependent manner.

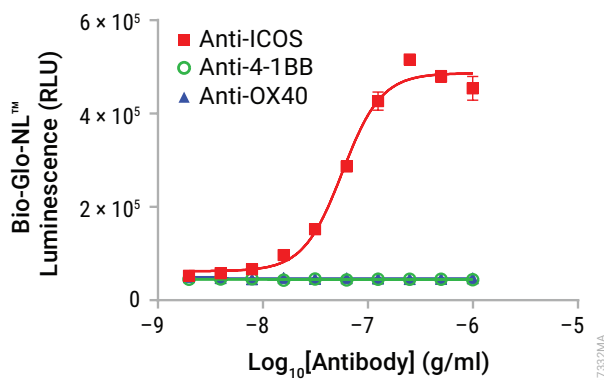


Figure 2. The ICOS Bioassay reflects the mechanism of action (MOA) and shows specificity for antibodies designed to activate ICOS. ICOS Effector Cells were incubated with FcγRIIb aAPC/CHO-K1 Cells in the presence of serial titrations of agonist antibodies as indicated. After a 6-hour induction, Bio-Glo-NL™ 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.

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

Parameter	Results	
Accuracy	% Expected Relative Potency	% Recovery
	50	46.6
	70	69.0
	100	102.8
	140	143.1
	200	211.7
Repeatability (% CV)	100% (Reference)	2.7
Intermediate Precision (% CV)		5.8
Linearity (r^2)		0.9996
Linearity ($y = mx + b$)		$y = 1.094x - 7.835$
<p>A 50–200% theoretical potency series of Control Ab, Anti-ICOS, was analyzed in triplicate in three independent experiments performed on three days by two analysts using the ICOS Bioassay for FcγRIIb-dependent antibodies. Bio-Glo-NL™ 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.</p>		

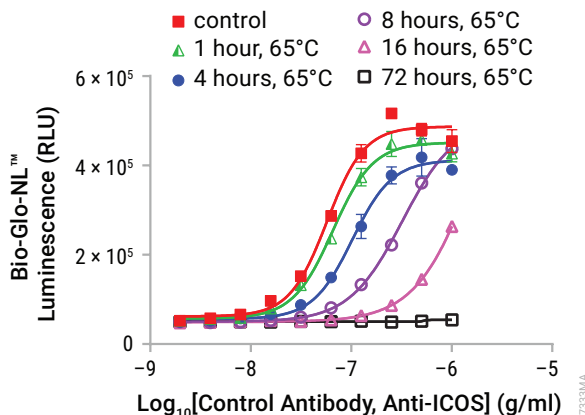


Figure 3. The ICOS Bioassay is stability-indicating. Samples of Control Ab, Anti-ICOS, were maintained at 4°C (control) or heat-treated at the indicated times and temperatures, then analyzed using the ICOS Bioassay for FcγRIIb-dependent antibodies. Bio-Glo-NL™ Reagent was added and luminescence quantified 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.

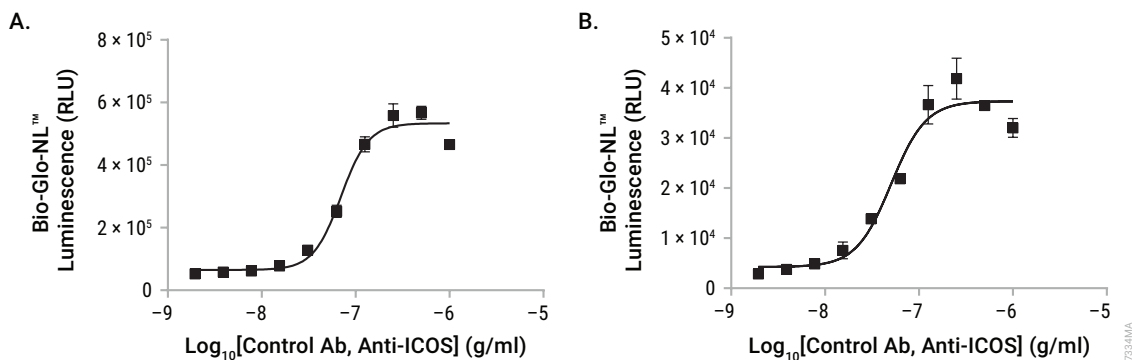


Figure 4. The ICOS Bioassay is amenable to 384-well plate format. Panel A. The ICOS Bioassay for FcγRIIb-dependent antibodies was performed in 96-well plates as described in this technical manual with a titration of Control Ab, Anti-ICOS. **Panel B.** The ICOS Bioassay for FcγRIIb-dependent antibodies was performed in 384-well format as briefly described here. FcγRIIb aAPC/CHO-K1 cells were harvested and 8×10^3 cells/15μl/well plated 16–24 hours prior to assay, in a 384-well white assay plate (e.g., Corning® Cat.# 3570). On the day of the assay, 5μl of 5X serially diluted Control Ab, Anti-ICOS, was added, followed by the addition of 2×10^4 /5μl/well of ICOS Effector Cells. After a 6-hour assay incubation at 37°C, 5% CO₂, 25μl of Bio-Glo-NL™ Reagent was added per well and luminescence was quantified using the GloMax® Discover System. Data were fitted to four-parameter logistic curves using GraphPad Prism® software. The EC₅₀ values were 7.0 and 5.1ng/ml for the 96-well and 384-well format, respectively, and the fold induction was 10.2 and 13.1 for 96-well and 384-well format, respectively. Data were generated using thaw-and-use cells.

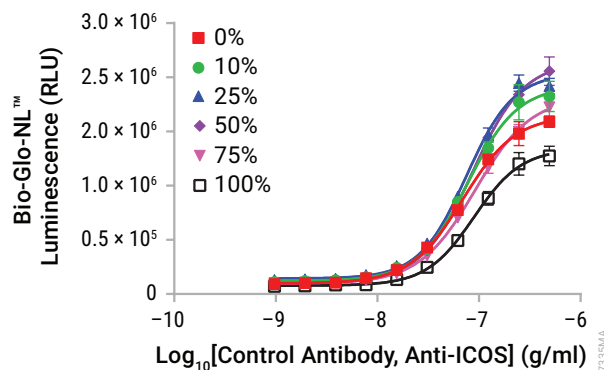


Figure 5. The ICOS Bioassay is tolerant to human serum. Following the ICOS Bioassay protocol for FcγRIIb- dependent antibodies, Control Ab, Anti-ICOS (Cat.# K1241), was analyzed in the absence or presence of pooled normal human serum (0–100% in the antibody sample). After the 6-hour assay induction at 37°C, Bio-Glo-NL™ 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. The ICOS Bioassay is tolerant to human serum with this serum pool. A different human serum pool showed similar effects on the assay (data not shown).

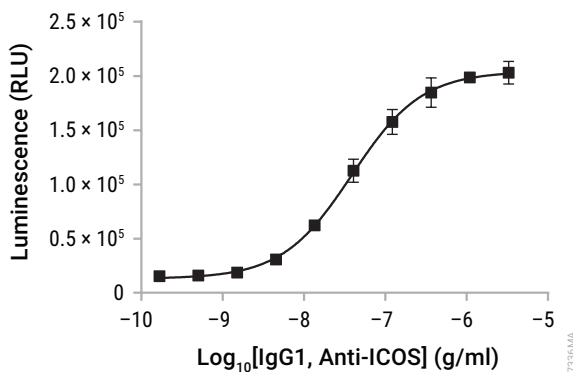


Figure 6. The ICOS Bioassay can be used to measure ADCC activity. A 2:1 ratio of ICOS Effector Cells (used as target cells in this application) and ADCC Bioassay Effector Cells (Cat.# G7010), were incubated for 6 hours in the presence of a research-grade anti-ICOS antibody (human IgG1 isotype). For reference, we use 3.3×10^{-6} g/ml as a starting concentration (1X) and a threefold serial dilution when testing hulgG1 anti-ICOS. 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.

2. Product Components and Storage Conditions

PRODUCT	SIZE	CAT.#
ICOS Bioassay, Core Kit	1 each	JA6801

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 ICOS Effector Cells
- 36ml RPMI 1640 Medium
- 4ml Fetal Bovine Serum
- 1 vial Bio-Glo-NL™ Luciferase Assay Substrate
- 10ml Bio-Glo-NL™ Luciferase Assay Buffer

PRODUCT	SIZE	CAT.#
ICOS Bioassay, Core Kit 5X	1 each	JA6805

Not for Medical Diagnostic Use. Each kit contains sufficient reagents for 600 assays using the inner 60 wells of ten 96-well plates. Includes:

- 5 vials ICOS Effector Cells
- 5 × 36ml RPMI 1640 Medium
- 5 × 4ml Fetal Bovine Serum
- 5 vials Bio-Glo-NL™ Luciferase Assay Substrate
- 5 × 10ml Bio-Glo-NL™ Luciferase Assay Buffer

Note: The ICOS Bioassay components and ICOS Effector Cells are shipped separately because of different temperature requirements. ICOS Effector Cells are shipped on dry ice. Bio-Glo-NL™ Luciferase Assay Substrate, Bio-Glo-NL™ Luciferase Assay Buffer and Fetal Bovine Serum are shipped on dry ice, separately from the cells. 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) for long-term storage. Do not store cell vials submerged in liquid nitrogen. **Do not** store cell vials at –80°C because this will negatively impact cell viability and cell performance.
- Store Bio-Glo-NL™ Luciferase Assay Substrate, Bio-Glo-NL™ Luciferase Assay Buffer and Fetal Bovine Serum at –30°C to –10°C. The Bio-Glo-NL™ Luciferase Assay Substrate remains liquid and does not freeze. Avoid multiple freeze-thaw cycles of the serum.
- Store RPMI 1640 at 4°C, protected from fluorescent light.

Available Separately

PRODUCT	SIZE	CAT.#
FcγRIIb aAPC/CHO-K1 Cells	1 each	J9331
FcγRIIb aAPC/CHO-K1 Cells 5X	1 each	J9335
aAPC/CHO-K1 Cells	1 each	J9441
aAPC/CHO-K1 Cells 5X	1 each	J9445

Not for Medical Diagnostic Use.

3. Before You Begin

Please read through 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.



Note: The ICOS Bioassay, uses the Bio-Glo-NL™ Luciferase Assay System (Cat.# J3081, J3082, J3083), for detection. **Do not** use the Bio-Glo™ Luciferase Assay System (Cat.# G7940, G7941).

The ICOS Bioassay is intended to be used with user-provided antibodies or other biologics designed to activate ICOS. Control Ab, Anti-ICOS (Cat.# K1241), FcγRIIb aAPC/CHO-K1 Cells (Cat.# JA9331, JA9335), and aAPC/CHO-K1 Cells (Cat.# JA9441, JA9445), are available separately for use in assay optimization and routine quality control. We strongly recommend using FcγRIIb aAPC/CHO-K1 Cells and Control Ab, Anti-ICOS, as a positive control in the first few assays to gain familiarity with the assay. Data generated using FcγRIIb aAPC/CHO-K1 Cells with Control Ab, Anti-ICOS, are shown in Figures 2–5 and Section 7.A, Representative Assay Results.

The ICOS 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. Do not overmix or overwarm the cell reagents.

The ICOS Bioassay produces a bioluminescent signal and requires a sensitive luminescence plate reader. Bioassay development and performance data included in this Technical Manual were generated using the GloMax® Discover System (see Section 7.B, Related Products). An integration time of 0.5 second/well was used for all readings. The bioassay is compatible with most other plate-reading luminometers, though relative luminescence unit (RLU) readings will vary with the sensitivity and settings of each instrument. If using a reader with adjustable gain, we recommend a high-gain setting. The use of different instruments and gain adjustment will affect the magnitude of the raw data but should not affect the measured relative potency of test samples.

3.A. Materials to Be Supplied by the User

- user-defined anti-ICOS antibodies or other biologics samples
- solid-white, flat-bottom 96-well assay plates (e.g., Corning® Cat.# 3917) or 384-well assay plates (e.g., Corning® Cat.# 3570) for plating and reading luminescence
- sterile clear V-bottom 96-well plate with lid (e.g., Costar® Cat.# 3896) 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®/Costar® Cat.# 4870)
- 37°C, 5% CO₂ incubator
- 37°C water bath
- sensitive plate reader with glow luminescence measuring capability or luminometer (e.g., GloMax® Discover System or equivalent system)
- **optional:** Control Ab, Anti-ICOS (Cat.# K1241; This antibody is FcγRIIb-dependent and requires the use of FcγRIIb aAPC/CHO-K1 Cells for optimal signal.)
- **optional:** FcγRIIb aAPC/CHO-K1 Cells (Cat.# JA9331, JA9335; for FcγRIIb-dependent antibodies)
- **optional:** aAPC/CHO-K1 Cells (Cat.# JA9441, JA9445; for FcγRIIb-independent antibodies)

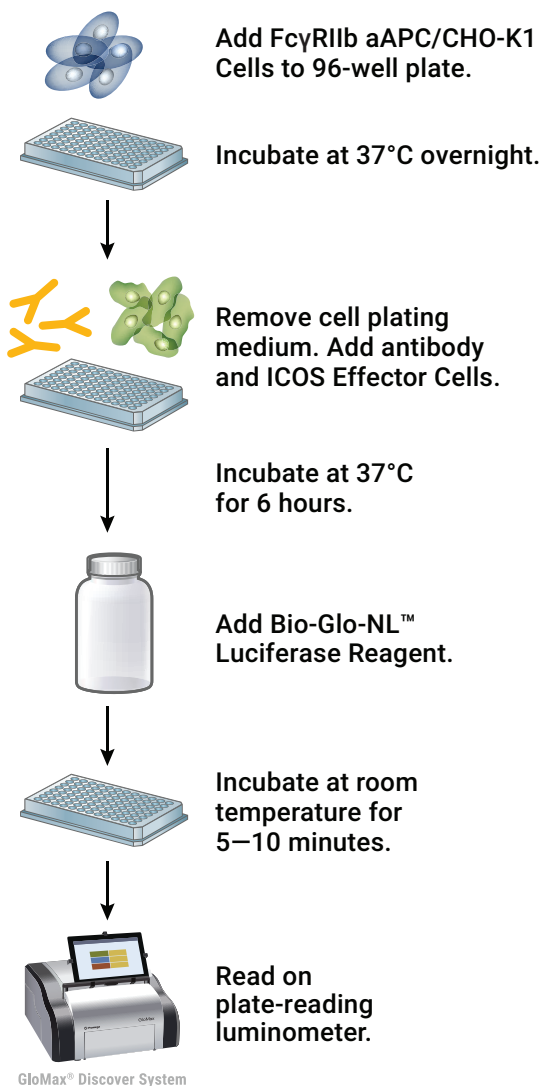
4. Assay Protocol for FcγRIIb-Dependent or FcγRIIb-Independent Antibodies

This assay protocol requires two engineered cell lines: ICOS Effector Cells, included with this assay, and either FcγRIIb aAPC/CHO-K1 Cells (Cat.# JA9331), for FcγRIIb-dependent antibodies, or aAPC/CHO-K1 Cells (Cat.# JA9441), for FcγRIIb-independent antibodies. These two cell lines are available separately as listed in Section 3.A. Section 4.A describes preparation of cell plating medium containing Ham's F12 Medium, which is provided with FcγRIIb aAPC/CHO-K1 Cells and aAPC/CHO-K1 Cells.

The procedure below illustrates the use of the ICOS Bioassay to test two FcγRIIb-dependent antibody samples against a reference sample in a single assay (Figure 7). 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.

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 1 µg/ml as a starting concentration (1X) and twofold serial dilution when testing Control Ab, Anti-ICOS, using the FcγRIIb-dependent method.

For testing antibodies not dependent on FcγRIIb-mediated crosslinking, use aAPC/CHO-K1 Cells instead of FcγRIIb aAPC/CHO-K1 Cells in this protocol. The aAPC/CHO-K1 Cells and FcγRIIb aAPC/CHO-K1 Cells are thawed and plated using the same protocol. When using the aAPC/CHO-K1 Cells, we use the same 1 µg/ml as a starting concentration (1X) of Control Ab, Anti-ICOS, and a twofold serial dilution series.



17337/NA

Figure 7. Schematic protocol for the ICOS Bioassay.

4.A. Preparing Cell Plating Medium, Assay Buffer, Bio-Glo-NL™ Reagent and Antibody Samples

1. **FcγRIIb aAPC/CHO-K1 Cells Plating Medium:** On the day before the assay, prepare an appropriate amount of cell plating medium (90% Ham's F-12/10% FBS). Thaw the FBS overnight at 4°C or in a 37°C water bath on the day of use. Add 1.5ml of FBS to 13ml of Ham's F-12 medium. Mix well and warm to 37°C before use. For reference, 14.5ml of cell plating medium is sufficient to thaw and plate 1 vial of FcγRIIb aAPC/CHO-K1 Cells. If multiple vials will be thawed, adjust the amount of cell plating medium accordingly. Warm the remaining Ham's F12 Medium to 37°C. Store the remaining FBS at 4°C for use in preparing the assay buffer on the day of the assay.
2. **Assay Buffer:** On the day of the assay, prepare 20ml of assay buffer (99% RPMI 1640/1% FBS). Add 0.2ml of FBS to 19.8ml RPMI 1640 Medium. Mix well and warm to 37°C before use. Warm the remaining RPMI 1640 Medium to 37°C.

Note: The recommended assay buffer contains 1% FBS. This concentration of FBS works well for the Control Ab, Anti-ICOS that we tested. If you experience assay performance issues when using this assay buffer, we recommend testing different serum concentrations in the range of 0.5–10%.

3. **Bio-Glo-NL™ Reagent:** For reference, 10ml of Bio-Glo-NL™ Reagent is sufficient to assay 120 wells in a 96-well assay format. The Bio-Glo-NL™ Luciferase Assay Substrate should always be stored at –20°C. Thaw the Bio-Glo-NL™ Luciferase Assay Buffer at room temperature (do not exceed 25°C) during the 6-hour induction period of the assay. We recommend preparing the reconstituted Bio-Glo-NL™ Luciferase Reagent immediately before use. For instructions on use of the Bio-Glo-NL™ Luciferase Assay System, please refer to the *Bio-Glo-NL™ Luciferase Assay System Quick Protocol #FB227*.



Note: The ICOS Bioassay is compatible only with Bio-Glo-NL™ Luciferase Assay System (Cat.# J3081, J3082, J3083). **Do not** use the Bio-Glo™ Luciferase Assay System (Cat.# G7940, G7941) with the ICOS Bioassay.

4. **Test and Reference Samples:** Using assay buffer as the diluent, prepare stock starting dilutions (dilu1, 2X final concentration) of two test antibodies (300µl each) and one reference antibody (600µl) 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-ICOS (Cat.# K1241), as a reference antibody in your assay, prepare a 50µg/ml working stock of Control Ab, Anti-ICOS, by adding 2µl of Control Ab, Anti-ICOS stock (1.0mg/ml) to 38µl of assay buffer. Prepare a 600µl starting dilution of 2µg/ml of Control Ab, Anti-ICOS (dilu1, 2X final concentration) by adding 24µl of Control Ab, Anti-ICOS working stock (50µg/ml) to 576µl of assay buffer.
- b. The highest concentration of Control Ab, Anti-ICOS, 2µg/ml, is appropriate for both FcγRIIb-dependent and FcγRIIb-independent antibodies.

4.B. Plate Layout Design

For the protocol described here, use the plate layout illustrated in Figure 8 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 8. Example plate layout showing nonclustered sample locations of test antibody and reference antibody dilution series, and wells containing assay buffer (denoted by “B”) alone.

4.C. Plating FcγRIIb aAPC/CHO-K1 Cells

Thaw-and-use FcγRIIb aAPC/CHO-K1 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 manipulation is required or recommended. We recommend that you thaw and dilute a maximum of two vials of thaw-and-use cells at any one time.

Follow the instructions below when using aAPC/CHO-K1 Cells for FcγRIIb-independent antibodies.

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

Perform the following steps in a sterile cell culture hood.

1. On the day before performing the assay, prepare 14.5ml of cell plating medium (90% Ham's F12/10% FBS) as described in Section 4.A.
2. Remove one vial of FcγRIIb aAPC/CHO-K1 Cells from storage at -140°C and transfer to the bench on dry ice. Thaw the cells in a 37°C water bath until just thawed (about 2 minutes). While thawing, gently agitate and visually inspect the vial. Do not invert.
3. Gently mix the cell suspension by pipetting, then transfer 0.5ml cells to the tube containing 14.5ml of cell plating medium. Mix well by gently inverting the tube 1–2 times.
4. Transfer the suspension to a sterile reagent reservoir. Using a multichannel pipette, immediately dispense 100μl of the cell suspension to each of the inner 60 wells of two 96-well white flat-bottom assay plates.
5. Add 100μl of prewarmed (37°C) Ham's F12 Medium to each of the outside wells of the assay plates.
6. Place lids on the assay plates and incubate in a 37°C , 5% CO_2 incubator overnight (18–22 hours).

4.D. Preparing Antibody Serial Dilutions

The instructions described here are for preparation of a single stock of twofold serial dilutions of a single antibody for analysis in triplicate (150µl of each dilution provides a sufficient volume for analysis in triplicate). Alternatively, you can prepare three independent stocks of serial dilutions to generate triplicate samples. To prepare twofold serial dilutions, you will need 600µl of reference antibody at 2X the highest antibody concentration in your dose-response curve. You will need 300µl of each test antibody at 2X 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 Ab, Anti-ICOS (Cat.# K1241), as a control in the assay, follow the instructions below to prepare twofold serial dilutions. A twofold serial dilution for test antibodies is listed as an example below as well.

1. On the day of the assay, prepare an appropriate amount of assay buffer as described in Section 4.A.
2. To a sterile clear V-bottom 96-well plate, add 300µl of reference antibody starting dilution (dilu1, 2X final concentration) to wells A11 and B11 (see Figure 9).
3. Add 300µl of test antibodies 1 and 2 starting dilution (dilu1, 2X final concentration) to wells E11 and G11, respectively (see Figure 9).
4. Add 150µl of assay buffer to other wells in these four rows, from column 10 to column 2.
5. Transfer 150µl of the antibody starting dilutions from column 11 into column 10. Mix well by pipetting. Avoid creating bubbles.
6. Repeat equivalent twofold serial dilutions across the columns from right to left through column 3. Do not dilute into column 2.

Note: Wells A2, B2, E2 and G2 contain 150µl of assay buffer without antibody as a negative control.

7. Cover the antibody dilution plate with a lid and keep at ambient temperature (22–25°C) while preparing the ICOS 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 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 9. Example plate layout showing antibody serial dilutions.

4.E. Preparing ICOS Effector Cells

The thaw-and-use ICOS 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 manipulation is required or recommended. We recommend that you thaw and dilute a maximum of two vials of thaw-and-use cells at any one time.

1. Label a sterile 15ml conical tube "Effector Cells". Add 6ml of prewarmed (37°C) assay buffer to the 15ml conical tube.
2. Remove one vial of ICOS Effector Cells from storage at -140°C and transfer to the bench on dry ice. Thaw the cells in a 37°C water bath until just thawed (about 2 minutes). While thawing, gently agitate and visually inspect the vial.
3. Gently mix the cell suspension by pipetting, then transfer 0.5ml cells to the 15ml conical tube containing 6ml of assay buffer. Mix well by gently inverting the tube.

4.F. Adding ICOS Effector Cells and Antibody to Assay Plates

1. Take the 96-well assay plates containing FcγRIIb aAPC/CHO-K1 Cells out of the incubator. 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. Alternatively, remove 95µl of medium from each of the wells using a manual multichannel pipette.
2. Using a multichannel pipette, add 40µl of the appropriate antibody dilution (Figure 9) to the assay plates according to the plate layout in Figure 8.
3. Mix the ICOS Effector Cells by inverting the tube and transfer the suspension to a sterile reagent reservoir. Using a multichannel pipette, immediately dispense 40µl of the cell suspension to each of the inner 60 wells of the assay plates. Gently swirl the assay plates to ensure mixing of the ICOS Effector Cells and antibody.
4. Add 80µl of assay buffer to each of the outside wells of the assay plates.
5. Cover the assay plates with lids and incubate in a 37°C, 5% CO₂ incubator for 6 hours.

Note: The 6-hour assay time was optimized using the Control Ab, Anti-ICOS. We recommend optimizing assay time (5–24 hours) with your own antibody or other biologic samples.

4.G. Preparing and Adding Bio-Glo-NL™ Reagent

We recommend preparing the Bio-Glo-NL™ Reagent immediately before use. Ensure that the Bio-Glo-NL™ Luciferase Assay Buffer is equilibrated to room temperature (do not exceed 25°C) before reconstituting the reagent. **Do not** store the reconstituted reagent. Once reconstituted, the reagent will lose 10% activity in approximately 8 hours at room temperature.



Note: The ICOS Bioassay is compatible only with Bio-Glo-NL™ Luciferase Assay System (Cat.# J3081, J3082, J3083).

Do not use Bio-Glo™ Luciferase Assay System (Cat.# G7940, G7941) with the ICOS Bioassay.

1. Remove the Bio-Glo-NL™ Luciferase Assay Substrate from –20°C storage and mix by pipetting. Briefly centrifuge the tubes if the substrate has collected in the cap or on the sides of the tubes.
2. Prepare the desired amount of reconstituted Bio-Glo-NL™ Reagent by combining one volume of substrate with 50 volumes of buffer. For example, if the experiment requires 10ml of reagent, add 200µl of substrate to 10ml of buffer. Ten milliliters (10ml) of Bio-Glo-NL™ Reagent is sufficient for 120 wells (two assay plates, using the inner 60 wells of each plate).
3. Remove assay plates from the incubator after the incubation period and equilibrate to room temperature for 10–15 minutes.
4. Using a manual multichannel pipette, add 80µl of Bio-Glo-NL™ Reagent to the inner 60 wells of the assay plates, taking care not to create bubbles.
5. Add 80µl of Bio-Glo-NL™ Reagent to wells B1, D1 and F1 of each assay plate to measure background signal.
6. Wait 5–10 minutes, then measure the luminescence in a GloMax® Discover System or a plate reader with glow-type luminescence reading capabilities. The luminescence intensity will decay gradually, with a signal half-life of approximately 120 minutes at room temperature.

Note: Varying the Bio-Glo-NL™ Reagent incubation time will affect the raw relative light unit (RLU) values but should not significantly change the EC₅₀ value and maximum fold induction.

4.H. Data Analysis

1. Determine the plate background by calculating the average RLU from wells B1, D1 and F1.
2. Calculate fold induction =
$$\frac{\text{RLU (induced-background)}}{\text{RLU (no antibody control-background)}}$$
3. Graph data as RLU versus Log_{10} [antibody] and fold induction versus Log_{10} [antibody]. Fit curves and determine the EC_{50} value of antibody response using appropriate curve fitting software (such as GraphPad Prism® software).

5. 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>Ensure that you are using Bio-Glo-NL™ Reagent, which is designed for NanoLuc® luciferase reporter bioassays. The ICOS Bioassay is not compatible with Bio-Glo™ Reagent, which is designed for firefly luciferase reporter bioassays.</p> <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>Some models of luminometers with low sensitivity should be avoided. If using a reader with an adjustable gain, we recommend a high-gain setting.</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>Low activity of Bio-Glo-NL™ Reagent leads to low RLU. Store and handle the Bio-Glo-NL™ Reagent according to the instructions. For best results, prepare immediately before use.</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 with the ICOS Bioassay may vary from the EC₅₀ value obtained using other methods such as primary T cell-based assays.</p> <p>The assay is sensitive to the concentration of FBS in assay buffer. Optimize the FBS concentration from 0.5%-10% in assay buffer if assay performance is not ideal.</p> <p>Optimize the assay incubation time within a range of 5–24 hours.</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>

6. References

1. Hutloff, A. *et al.* (1999) ICOS is an inducible T-cell co-stimulator structurally and functionally related to CD28. *Nature* **397**, 263–6.
2. Yoshinaga, S. *et al.* (1999) T-cell co-stimulation through B7RP-1 and ICOS. *Nature* **402**, 827–32.
3. Coyle, A. *et al.* (2000) The CD28-related molecule ICOS is required for effective T cell-dependent immune responses. *Immunity* **13**, 95–105.
4. Wang, S. *et al.* (2000) Costimulation of T cells by B7-H2, a B7-like molecule that binds ICOS. *Blood* **96**, 2808–13.
5. Aicher, A. *et al.* (2000) Characterization of human inducible costimulatory ligand expression and function. *J. Immunol.* **164**, 4689–96.
6. Khayyamian, S. *et al.* (2002) ICOSLigand expressed on human endothelial cells, costimulates Th1 and Th2 cytokine secretion by memory CD4+ T cells. *Proc. Natl. Acad. Sci. USA* **99**, 6198–203.
7. Wahl, P. *et al.* (2002) Renal tubular epithelial expression of the costimulatory molecule B7RP-1 (inducible costimulatory ligand). *J. Am. Soc. Nephrol.* **13**, 1517–26.
8. Grailer, J. *et al.* (2019) Quantitative cell-based bioassays to advance immunotherapy programs targeting immune checkpoint receptors. *Cancer Res.* **79**, 4113.

7. Appendix

7.A. Representative Assay Results

The following data were generated using the ICOS Bioassay, and Control Ab, Anti-ICOS (Figure 10).

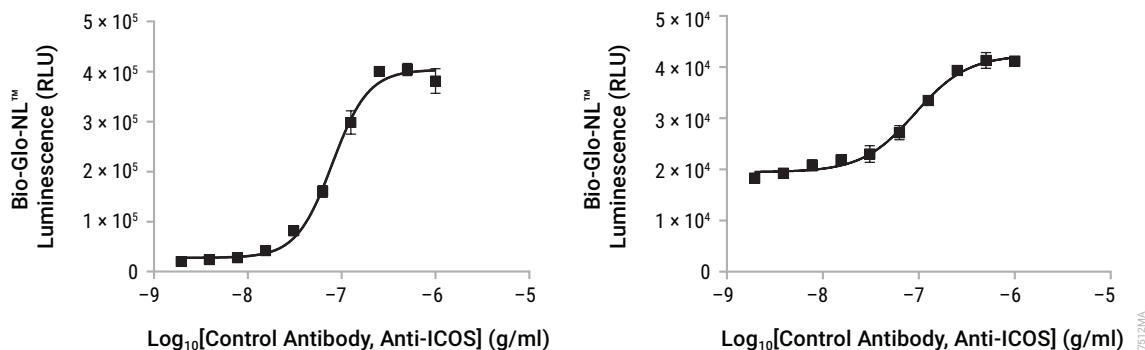


Figure 10. The ICOS Bioassay measures the activity of Control Ab, Anti-ICOS. FcγRIIb aAPC/CHO-K1 Cells (**Panel A**) or aAPC/CHO-K1 Cells (**Panel B**) were added to a 96-well assay plate 20–24 hours prior to the assay. On the day of assay, ICOS Effector Cells and a titration of Control Ab, Anti-ICOS (Cat.# K1241), were added. After a 6-hour induction at 37°C, Bio-Glo-NL™ Reagent was added and luminescence measured using the GloMax® Discover System. Data were fitted to a four-parameter logistic curve using GraphPad Prism® software. The EC₅₀ was 79.6ng/ml (**Panel A**) and 91.9ng/ml (**Panel B**), and the fold induction was 20 and 2.3 when using FcγRIIb aAPC/CHO-K1 Cells and aAPC/CHO-K1 Cells, respectively.

7.B. 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
FcγRIIa-H ADCP Reporter Bioassay, Complete Kit**	1 each	G9901
FcγRIIa-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

*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
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 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

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

7.B. Related Products (continued)

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.

Control Antibodies and Proteins

Product	Size	Cat.#
Control Ab, Anti-4-1BB	50µg	K1161
Control Ab, Anti-CD-20	5 µg	GA1130
Control Ab, Anti-OX40	50µg	K1191
Control Ab, Anti-CD40	50µg	K1181
Control Ab, Anti-CTLA-4	100µg	JA1020
Control Ab, Anti-LAG-3	100µg	K1150
Control Ab, Anti-PD-1	100µg	J1201
Control Ab, Anti-TIGIT	100µg	J2051
Control Ab, Anti-TIM-3	100µg	K1210
Recombinant VEGF ligand	10µg	J2371

Detection Reagent

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

Not for Medical Diagnostic Use.

Luminometers

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

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

Note: Additional Fc Effector, Immune Checkpoint, T Cell Activation, Cytokine, Macrophage, Primary Cell and Target Cell Killing Bioassays are available. To view and order Promega Bioassay products visit:

www.promega.com/products/reporter-bioassays/ 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/

8. Summary of Changes

The following changes were made to the 5/25 revision of this document:

1. In Section 3, text about the label was revised.
2. A patent statement was updated.



^(e)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-NL™ 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. Nos. 8,557,970, 8,669,103, 9,777,311, 9,840,730, 9,951,373, 10,233,485, 10,633,690, 10,774,364, 10,844,422, 11,365,436, 11,661,623, 11,667,950; European Pat. Nos. 2456864, 2635595, 2990478, 3181687, 3409764; Japanese Pat. Nos. 6038649, 6155424, 6227615, 6374420, 6539689; and other patents and patents pending.

^(c)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.

^(d)U.S. Pat. No. 8,809,529, European Pat. No. 2635582, Japanese Pat. No. 5889910 and other patents and patents pending.

^(e)Patent Pending.

© 2021, 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.

Corning and Costar are registered trademarks of Corning, Inc. GraphPad Prism is a registered trademark of GraphPad Software, Inc. JMP is a registered trademark of SAS Institute, Inc.

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.