



TECHNICAL MANUAL

RANKL Bioassay, Propagation Model

Instructions for Use of Product
J3102

RANKL Bioassay, Propagation Model

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

1. Description	2
2. Product Components and Storage Conditions	8
3. Before You Begin	8
4. Preparing RANKL Bioassay Cells	10
4.A. Cell Thawing and Initial Cell Culture	10
4.B. Cell Maintenance and Propagation	10
4.C. Cell Freezing and Banking	11
5. Stimulation Protocol	12
5.A. Plate Layout Design	13
5.B. Day One: Preparing and Plating RANKL Bioassay Cells	14
5.C. Day Two: Assay Day with Addition of Test and Reference Samples	14
5.D. Preparing Serial Dilutions	15
5.E. RANKL Stimulation Assay	16
5.F. Adding Bio-Glo™ Reagent	16
5.G. Data Analysis	16
6. Inhibition Protocol	17
6.A. RANKL Inhibition Assay	17
6.B. Adding Bio-Glo™ Reagent	18
6.C. Data Analysis	18
7. Troubleshooting	19
8. References	20
9. Appendix	20
9.A. Composition of Buffers and Solutions	20
9.B. Related Products	21
10. Summary of Changes	24

1. Description

Receptor activator of nuclear factor- κ B (NF- κ B; RANK/TRANCE receptor/TNFRSF11A) is a member of the tumor necrosis factor receptor (TNFR) family. Binding of its ligand (RANKL) to the receptor regulates osteoclast formation, activation and survival in bone modeling and remodeling, and several other pathologic conditions characterized by increased bone turnover (1).

The osteoclastogenesis signaling pathway is activated by osteoblasts producing RANKL, which binds to and activates the RANK receptor on osteoclast precursors. The adapter protein TRAF6 is recruited to the RANK receptor and activates NF- κ B, which leads to its translocation to the nucleus. This increases the expression of c-FOS which, together with NFATc1, increases the transcription of osteoclastogenic genes (2). Osteoprotegerin (OPG) binds to and inhibits RANKL. In cells with excess RANKL or insufficient OPG, upregulated RANKL/RANK signaling leads to superfluous osteoclast formation and bone resorption, causing pathologic bone loss and destruction (3).

In osteosarcoma, in addition to cancer-induced bone destruction, RANKL is also involved in tumorigenesis and metastasis. RANKL inhibition significantly delays mammary tumor formation in carcinogen and hormone-induced breast cancer mouse models (4, 5). Denosumab is a human IgG2 monoclonal antibody that targets and inhibits RANKL. It is used to treat osteoporosis and hypercalcemia, as well as bone cancer and other cancer-associated bone disease. Denosumab blocks the RANK-RANKL interaction, inhibiting osteoclast formation, function and survival; thus, it decreases bone resorption and interrupts cancer-induced bone destruction (6).

The RANKL Bioassay, Propagation Model^(a,b) (Cat.# J3102) is a bioluminescent cell-based assay designed to measure RANK/RANKL pathway stimulation or inhibition. The RANKL Bioassay Cells are provided in a Cell Propagation Model (CPM) format, which includes cryopreserved cells that can be thawed, propagated and banked for long-term use (also offered in a thaw-and-use format; Cat.# JA2701, JA2705).

The RANKL Bioassay is comprised of a human cell line engineered to express the RANK receptor and a luciferase reporter driven by a response element (RE). When RANKL binds the RANK receptor, transduced intracellular signals activate the RE, resulting in luminescence (Figure 1). The bioluminescent signal is detected and quantified using the Bio-Glo™ Luciferase Assay System (Cat.# G7940, G7941) and a standard luminometer, such as the GloMax® Discover System (see Related Products, Section 9.B).

1. Description (continued)

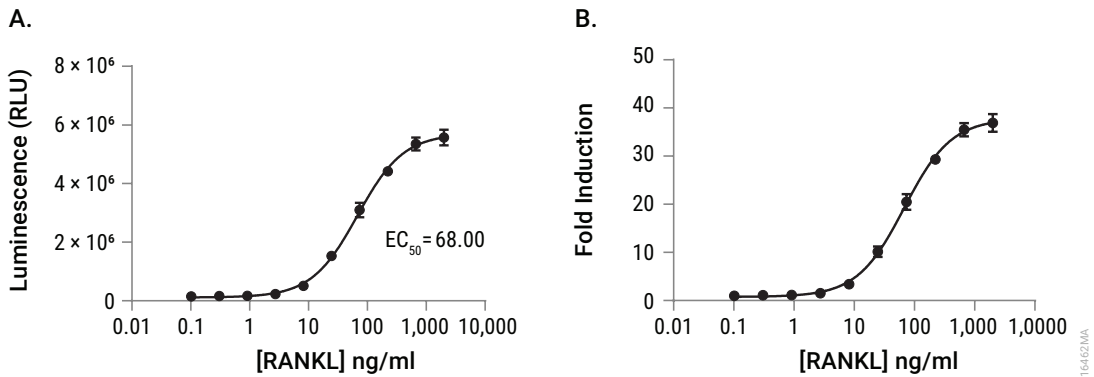


Figure 2. The RANKL Bioassay responds to recombinant RANKL. RANKL Bioassay Cells were grown and prepared as described in this protocol, and incubated with serial dilutions of recombinant RANKL. After a 6-hour incubation, Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were fitted to a four-parameter logistic curve using GraphPad Prism® software. **Panel A** shows raw luminescence measurements. **Panel B** displays the calculated fold induction. Data were generated using CPM cells.

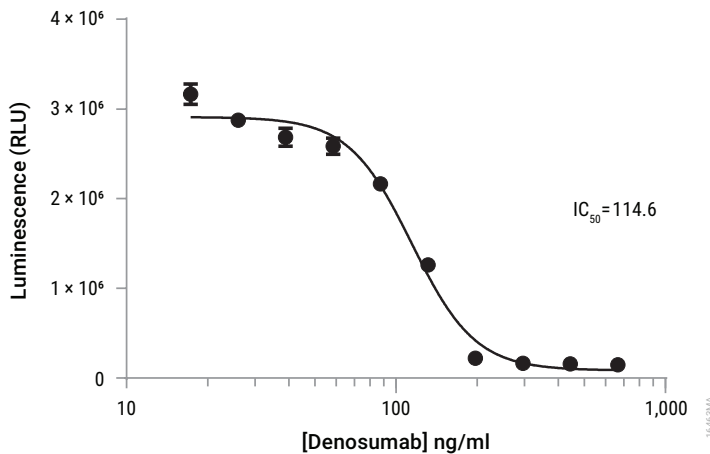


Figure 3. The RANKL Bioassay responds to denosumab. RANKL Bioassay Cells were grown and prepared as described in this protocol, and incubated with serial dilutions of denosumab (anti-RANKL) antibody and 100 ng/ml RANKL. After a 6-hour incubation, Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were fitted to a four-parameter logistic curve using GraphPad Prism® software. Data were generated using CPM cells.

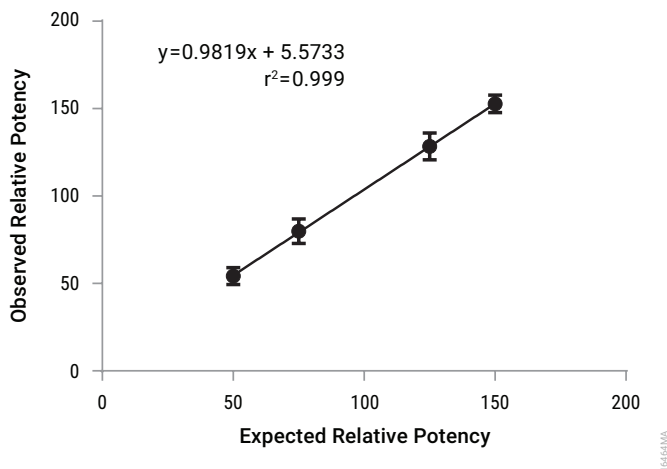


Figure 4. The RANKL Bioassay shows precision, accuracy, and linearity. A 50–150% theoretical potency series of RANKL was analyzed in triplicate in three independent experiments performed on three days by two analysts using the RANKL Bioassay (for a total of six independent experiments). Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Linearity and r^2 values were determined using GraphPad Prism® software. Data were generated using thaw-and-use cells.

1. Description (continued)

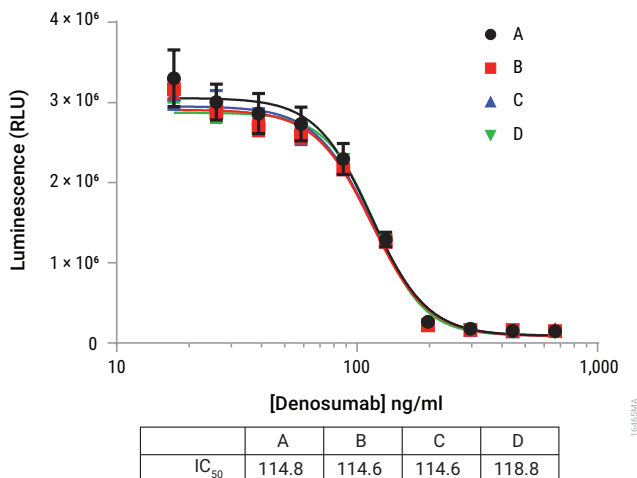


Figure 5. The RANKL Bioassay demonstrates repeatability. Four separate serial dilution series of denosumab were analyzed on four individual assay plates using the RANKL Bioassay. Bio-Glo™ Reagent was added, and luminescence was quantified using the GloMax® Discover System. Data were fitted to a four-parameter logistic curve using GraphPad Prism® software. Data were generated using thaw-and-use cells.

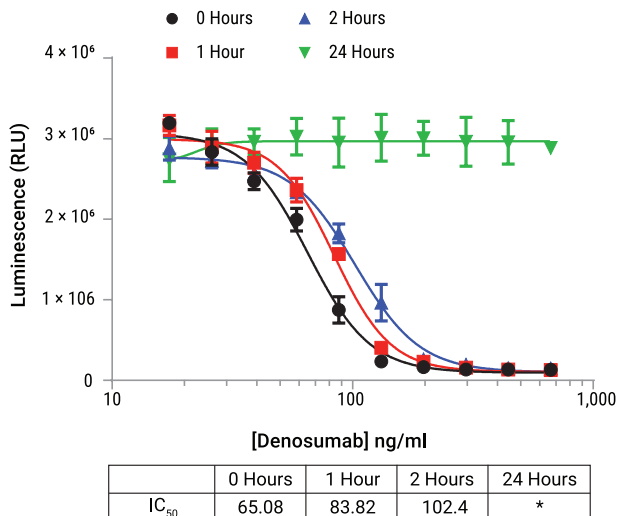


Figure 6. The RANKL Bioassay indicates stability. Denosumab was heated at 65°C for four different periods of time and used in the inhibition assay. Prolonged heating induced inactivated denosumab based on time heated, which is reflected by the RANKL assay. Data were generated using thaw-and-use cells. *Data for 24-hour treatment were not calculable.

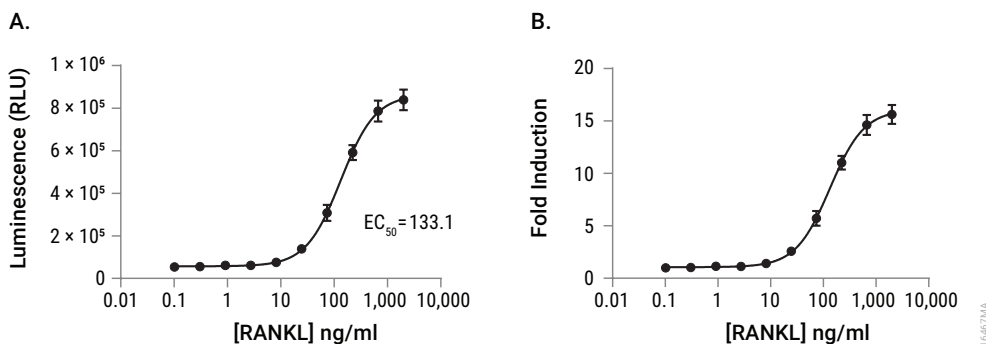


Figure 7. The RANKL Bioassay is amenable to 384-well plate format. The RANKL Bioassay was tested in 384-well format. RANKL Bioassay Cells were prepared and dispensed as 30 μ l/well. Serial threefold dilutions of recombinant human RANKL were prepared and added to cells after seeding overnight and removing of the seeding media. After 6 hours of stimulation, 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. **Panel A** shows raw luminescence measurements. **Panel B** displays the calculated fold induction. Data were generated using thaw-and-use cells.



2. Product Components and Storage Conditions

PRODUCT	SIZE	CAT. #
RANKL Bioassay, Propagation Model	1 each	J3102

Not for Medical Diagnostic Use. Includes:

- 2 vials RANKL Bioassay Cells, 1.0×10^7 cells/ml (1ml per vial)

Note: Thaw and propagate one vial to create frozen cell banks before use in an assay. Reserve the second vial for future use.

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

3. Before You Begin

Please read 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 the Certificate of Analysis.

Cell thawing, propagation and banking should be performed exactly as described in Section 4. Cell seeding and propagation densities have been optimized to ensure stable cell growth, which is reflected in a steady cell doubling rate, to achieve optimal, consistent performance.

The RANKL Bioassay is intended to be used with user-provided biologics designed to activate or inhibit the RANK/RANKL signaling pathway. The recommended cell plating density, induction time and assay buffer components described in Section 5 were established using research-grade recombinant human RANKL. You may need to adjust the parameters provided here and optimize assay conditions for other biologic samples. Data generated using these reagents is shown above in Figure 2.

The RANKL Bioassay produces a bioluminescent signal and requires a sensitive luminometer or luminescence plate reader for the detection of luminescence. Bioassay development and performance data included in this Technical Manual were generated using the GloMax[®] Discover System. An integration time of 0.5 second/well was used for all readings. The bioassay is compatible with most other plate-reading luminometers; however, relative luminescence unit readings may vary due to the sensitivity and settings of each instrument. The use of different instruments should not affect the measured relative potency of test samples.

Materials to Be Supplied by the User

Composition of Buffers and Solutions is provided in Section 9.A.

Reagents

- user-defined biologics samples
- RPMI 1640 (e.g., Sigma Cat.# R7388-1L)
- McCoy's 5A (e.g., GIBCO® Cat.# 16600-082)
- fetal bovine serum (e.g., HyClone Cat.# SH30070)
- hygromycin B (e.g., Invitrogen Cat.# 10687-010)
- Antibiotic G-418 Sulfate (e.g., Cat.# V8091)
- D-PBS (e.g., Invitrogen Cat.# 14190144)
- Accutase® or equivalent (e.g., Innovative Cell Technologies Cat.# AT104)
- trypan blue solution (e.g., Sigma Cat.# T8154)
- Bio-Glo™ Luciferase Assay System (Cat.# G7940 or G7941)
- DMSO (e.g., Sigma Cat.# D2650-5X5ML)
- **optional:** recombinant human RANKL (Biolegend Cat.# 591106)

Supplies and Equipment

- white, flat-bottom 96-well assay plates (e.g., Corning® Cat.# 3917)
- sterile clear 96-well plate with lid (e.g., Corning® Cat.# 3896 or Falcon Cat.# 353077) for preparing sample dilutions
- sterile 12-well reagent reservoir (e.g., Axygen Cat.# RES-MW12-LP-SI) for preparing sample dilutions
- pipettes (single-channel and 12-channel)
- sterile 15ml and 50ml conical tubes
- T75 tissue culture flask
- sterile reagent reservoirs
- humidified 37°C, 5% CO₂ incubator
- 37°C water bath
- plate reader that measures glow luminescence or luminometer (e.g., GloMax® Discover System)

4. Preparing RANKL Bioassay Cells

4.A. Cell Thawing and Initial Cell Culture

Note: RANKL Bioassay Cells are grown as **adherent** cultures.



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

1. Prepare 60ml of thaw medium (see Section 9.A) and prewarm to 37°C. This medium will be used for culturing the cells immediately after thawing.
2. Transfer 8ml of thaw medium into a 15ml conical tube.
3. Remove one vial of RANKL Bioassay Cells from storage at -140°C and thaw in a 37°C water bath with gentle agitation (no inversion) until just thawed (typically 2–3 minutes).
4. Spray the vial with 70% ethanol and transfer to a cell culture hood.
5. Transfer all of the cells to the 15ml conical tube containing 8ml of prewarmed thaw medium.
6. Centrifuge at 150 × *g* for 5 minutes.
7. Carefully aspirate the medium and resuspend the cell pellet in 40ml of prewarmed thaw medium in a 50ml conical tube.
8. Count cells with trypan blue and determine cell number and viability.
9. Transfer the cell suspension evenly into two T75 flasks. Place the flasks horizontally in a humidified 37°C, 5% CO₂ incubator and incubate for 2 days.

4.B. Cell Maintenance and Propagation

Note: For cell maintenance and propagation starting from the second cell passage, use growth medium containing selection antibiotic (see Section 9.A), and monitor cell viability and doubling rate during propagation. Cell growth rate will stabilize by approximately 7–10 days after thawing. At this time, cell viability is typically >95% and the average cell doubling rate is approximately 21 hours. Passage number should be recorded for each passage. Cells are expected to retain their functionality for up to 35 passages.

1. On the day of cell passage, visualize the cells under a microscope and estimate confluency.
2. Aspirate the growth medium and wash the cells with D-PBS.
3. Add 2ml of Accutase® to each T75 flask and rock the flask several times to mix and coat the cell surface.
4. Incubate at 37°C until cells begin to detach (approximately 3–5 minutes).

5. Add 8ml of prewarmed growth medium (see Section 9.A) and resuspend the cells.
6. Sample and count by Trypan blue exclusion.
7. Add fresh growth medium and transfer the cells to a new flask. Mix gently.
8. Recommended density for passaging cells is as follows:
 - a. For 2-day culture: 2×10^4 cells/cm²
 - b. For 3-day culture: 1×10^4 cells/cm²

Note: We recommend using the following media volumes for routine cell propagation: 14ml for a T75 flask, 28ml for a T150 flask and 42ml for a T225 flask. Scale according to the surface area of the flask.

9. Place the flasks horizontally in a humidified 37°C, 5% CO₂ incubator.

4.C. Cell Freezing and Banking

Note: We recommend making master and working cell banks at the earliest possible passage.

1. On the day of cell freezing, prepare new cell freeze medium (see Section 9.A) and keep on ice.
2. Harvest the cells as described in Section 4.B.
3. Remove a sample for cell counting by trypan blue staining. Calculate the volume of freeze medium needed based on desired cell freezing density. We recommend a freezing density range of 2×10^6 – 2×10^7 cells/ml.
4. Transfer the cells to 50ml sterile conical tubes or larger sized centrifuge tubes, and centrifuge at $150 \times g$ for 10–15 minutes.
5. Gently aspirate the supernatant, being careful not to disturb the cell pellet.
6. Carefully resuspend the cell pellet in ice-cold freeze medium to the desired final cell density. Combine the cell suspensions into a single tube and dispense 1ml into cryovials.
7. Freeze using a controlled-rate freezer (or use an insulated Mr. Frosty® or a Styrofoam® type of cell freezing container at –80°C overnight).
8. Transfer to –140°C or below for long-term storage.

5. Stimulation Protocol

The RANKL Bioassay can be used to test RANKL and RANKL-blocking antibodies. This stimulation protocol illustrates the use of the RANKL Bioassay to examine two test samples against a reference sample in a single assay run. Each test and reference sample is run in triplicate, in a ten-point dilution series, in a single 96-well assay plate using the inner 60 wells (see Figure 9). Other experimental and plate layouts are possible but may require further optimization. The inhibition protocol (Section 6) illustrates the use of the RANKL Bioassay to block RANKL activity.

Notes:

- When preparing test and reference samples, 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 0–2 μ g/ml final concentration of recombinant human RANKL (Biolegend Cat.# 591106) as a sample range, with serial threefold dilutions to achieve full dose curves as a ten-point series. Concentration ranges and dilution schemes may need to be optimized for your samples.
- While maintaining the RANKL Bioassay Cells in culture, follow the recommended cell seeding density during routine propagation. Changes in cell culture volume or seeding density could affect subsequent assay performance. Only use cells in this assay after the doubling rate has stabilized during propagation. Use actively growing, healthy cells harvested as part of a routine 2- or 3-day passage; culture viability should be >95% prior to use in the RANKL Bioassay.

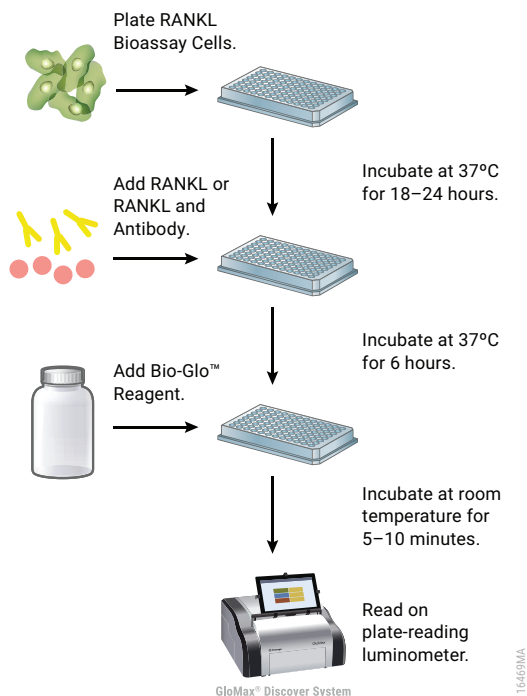


Figure 8. RANKL Bioassay schematic protocol.

5.A. Plate Layout Design

For the protocol described here, use the plate layout illustrated in Figure 9 as a guide. The protocol describes serial replicate dilutions (n = 3) of test and reference samples 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 drug	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Replicate 1
C	B	no drug	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Replicate 1
D	B	no drug	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Replicate 2
E	B	no drug	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Replicate 2
F	B	no drug	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Reference Replicate 3
G	B	no drug	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1	B	Test Replicate 3
H	B	B	B	B	B	B	B	B	B	B	B	B	Assay Buffer (B)

Figure 9. Example plate layout. This suggested layout shows nonclustered locations for three replicates of each test and reference sample dilution series (dilu1–dilu9) and wells containing assay buffer (denoted by “B”) alone.

5.B. Day One: Preparing and Plating RANKL Bioassay Cells

1. Prepare 50ml of seeding media as described in Section 9.A and warm to 37°C before use.
2. Aspirate the growth medium and wash the cells with D-PBS.
3. Add 2ml of Accutase® to a T75 flask and rock the flask several times to mix and coat the cell surface.
4. Incubate at 37°C until cells begin to lift off (approximately 3–5 minutes).
5. Add 8ml of seeding media (see Section 9.A) and resuspend the cells.
6. Sample and count by trypan blue exclusion.
7. Based on the number of samples and plates, estimate the number of cells required and include 50–100% extra to account for loss during centrifugations. For each assay plate, a minimum of 1.5×10^6 cells are required (2.5×10^4 cells/well \times 60 wells).
8. Place the cells into 50ml centrifuge tubes and centrifuge at $150 \times g$ for 5–10 minutes.
9. Remove the supernatant. Resuspend the cells in seeding media to an estimated 2×10^6 cells/ml and count again by Trypan blue exclusion.
10. Adjust to 2.5×10^5 cells/ml using additional seeding media.
11. Dispense 100 μ l/well (2.5×10^4 cells/well) using a multichannel pipette into the inner 60 wells of two solid white 96-well plates. Add 100 μ l/well of assay buffer to the outer 36 wells.
12. Incubate overnight (18–24 hours) at 37°C, 5% CO₂.

5.C. Day Two: Assay Day with Addition of Test and Reference Samples

Preparing Reagents for the Assay Day

1. **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 in a refrigerator 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 the Bio-Glo™ Luciferase Assay Buffer into the amber bottle containing the Bio-Glo™ Luciferase Assay Substrate and mix by inversion until the Substrate is thoroughly dissolved. Equilibrate and store the reconstituted Bio-Glo™ Reagent at ambient temperature (22–25°C) protected from light before adding to assay plates.

If you are using a large (100ml) size of Bio-Glo™ Luciferase Assay System, dispense the reconstituted Bio-Glo™ Reagent into 10ml aliquots and store at –20°C for up to 6 weeks. Avoid repeated freeze-thaw cycles. On the day of the assay, thaw the appropriate amount of reconstituted Bio-Glo™ Reagent in a room temperature water bath for at least 1–2 hours before use. Approximate stability of Bio-Glo™ Reagent after reconstitution is 18% loss of luminescence after 24 hours at ambient temperature and 12% loss of luminescence after 5 days at 4°C.
2. **Assay Buffer:** Use RPMI 1640 without serum as assay buffer. Warm to 37°C prior to use. For reference, 30ml of assay buffer is typically sufficient for 120 wells in a 96-well assay format using the inner 60 wells.

- Test and Reference Samples:** Prepare starting dilutions (denoted as dilu1) of test and reference samples (see Figure 10). Using assay buffer as the diluent, prepare a minimum of 800µl of reference sample starting dilution and a minimum of 400µl of each test sample starting dilution in 1.5ml tubes. Store the tubes containing starting dilutions appropriately before making serial dilutions.

5.D. Preparing Serial Dilutions

The instructions described here are for preparation of a single stock of threefold serial dilutions of a single sample for analysis in triplicate (250µl of each dilution provides a sufficient volume). Alternatively, you can prepare three independent stocks of serial dilutions to generate triplicate samples.

Note: For RANKL stimulation using recombinant human RANKL as your reference sample (Biolegend Cat.# 591106), we recommend starting with a 1X concentration of 2µg/ml and performing serial threefold dilutions. When using other reference sources of RANKL, the starting concentration may need to be adjusted. To perform the dilutions as described here, you will need a total of four dilution reservoirs, two for the test samples and two for the reference samples.

- Add 375µl of each sample starting dilution (dilu1) to column 11 of a sterile 12-well dilution reservoir (see Figure 10).
- Add 250µl of assay buffer to the other wells, from column 10 to column 2.
- Transfer 125µl of the sample starting dilution from column 11 into column 10. Mix well by pipetting. Avoid creating bubbles.
- Repeat equivalent threefold serial dilutions across the columns from right to left until you reach column 3. Remove 125µl from column 3 so all wells have 250µl volume. Do not dilute into column 2.
- Create the same dilution series for the remaining samples.

Recommended Reservoir Layout for Sample Dilutions													
	1	2	3	4	5	6	7	8	9	10	11	12	
A		no drug	dilu9	dilu8	dilu7	dilu6	dilu5	dilu4	dilu3	dilu2	dilu1		Sample

Figure 10. Example reservoir layout showing sample serial dilutions. Well 2 contains 250µl of assay buffer without sample.

5.E. RANKL Stimulation Assay

1. 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, use a multichannel pipette or aspirator to remove seeding media (~100µl) from each well.
2. Dispense 75µl of each sample to the preplated cells according to the plate layout in Figure 9.
3. Cover each assay plate with a lid and incubate in a humidified 37°C, 5% CO₂ incubator for 6 hours.
4. After the 6-hour incubation is over, proceed to Section 5.F.

5.F. Adding Bio-Glo™ Reagent

Note: Prepare Bio-Glo™ Reagent as described in Section 5.C. Bio-Glo™ Reagent should be at ambient temperature (22–25°C) when added to assay plates.

1. Remove the assay plates from the incubator, remove the plate lid and equilibrate to ambient temperature for 10–15 minutes.
2. Using a multichannel pipette, add 75µl of Bio-Glo™ Reagent to the inner 60 wells of the assay plates, taking care to not create bubbles.
3. Add 75µl of Bio-Glo™ Reagent to wells B1, C1 and D1 of each assay plate to measure the background signal.
4. Incubate at ambient temperature for 5–10 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.G. Data Analysis

1. Measure plate background by calculating the average RLU from wells B1, C1 and D1.

2. Calculate fold induction =
$$\frac{\text{RLU (sample-background)}}{\text{RLU (no drug control-background)}}$$

Note: When calculating fold induction, if the no-drug control sample RLU is at least 100X the plate background RLU, there is no need to subtract plate background from sample RLU.

3. Graph data as RLU versus log₁₀[sample] and fold induction versus log₁₀[sample]. Fit curves and determine the EC₅₀ value of RANKL response using appropriate curve fitting software (such as GraphPad Prism®).

6. Inhibition Protocol

The RANKL Bioassay Cells can be used to measure inhibition of RANK/RANKL signaling using a blocking antibody such as denosumab, which targets RANKL. A preliminary stimulation experiment with RANKL is necessary to determine the EC_{80} concentration, which is used during an inhibition assay. The protocol may be modified for other blocking antibodies, depending on their mechanism of action.

6.A. RANKL Inhibition Assay

1. Prepare and plate RANKL Bioassay Cells as described in Sections 5.A and 5.B (see Figure 9). Incubate overnight in a 37°C, 5% CO₂ humidified incubator.
2. In a separate clear 96-well plate, prepare serial dilutions of denosumab in prewarmed assay buffer to a 2X final concentration. We recommend a final concentration of denosumab of 0–660ng/ml as serial 10-point 1.5-fold dilutions. Prepare at least 125µl of denosumab solution for each sample.
3. Prepare a RANKL solution at 2X the concentration of the predetermined EC_{80} response concentration in prewarmed assay buffer. Prepare at least 125µl of RANKL solution for each sample.
4. Combine an equal volume of RANKL solution with each antibody dilution. Mix by pipetting.
Note: Each component is now at 1X its final concentration.
5. Remove the 96-well assay plates containing the overnight preplated RANKL Bioassay Cells from the incubator.
6. 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, use a multichannel pipette or aspirator to remove all of the seeding media (~100µl) from each well of the assay plate.
7. Add 75µl of RANKL and denosumab dilutions from the dilution plate in Step 4 into each replicate well of the assay plate.
8. Cover each assay plate with a lid and incubate in a 37°C, 5% CO₂ humidified incubator for 6 hours.
9. After the 6-hour incubation is completed, proceed to Section 6.B.

6.B. Adding Bio-Glo™ Reagent

Note: Prepare Bio-Glo™ Reagent as described in Section 5.C. Bio-Glo™ Reagent should be at ambient temperature (22–25°C) when added to assay plates.

1. Remove the assay plates from the incubator, remove the plate lid and equilibrate to ambient temperature for 10–15 minutes.
2. Using a multichannel pipette, add 75µl of Bio-Glo™ Reagent to the inner 60 wells of the assay plates, taking care to not create bubbles.
3. Add 75µl of Bio-Glo™ Reagent to wells B1, C1 and D1 of each assay plate to measure the background signal.
4. Incubate at ambient temperature for 5–10 minutes.
Note: Varying the incubation time will affect the raw RLU values but should not significantly change the IC₅₀ value and fold induction.
5. Measure luminescence using a luminometer or luminescence plate reader.

6.C. Data Analysis

1. Measure plate background by calculating the average RLU from wells B1, C1 and D1.

2. Calculate fold induction =
$$\frac{\text{RLU (sample-background)}}{\text{RLU (no drug control-background)}}$$

Note: When calculating fold induction, if the no-drug control sample RLU is at least 100X the plate background RLU, there is no need to subtract plate background from sample RLU.

3. Graph data as RLU versus log₁₀[sample] and fold induction versus log₁₀[sample]. Fit curves and determine the IC₅₀ value of RANKL inhibition response using appropriate curve fitting software (such as GraphPad Prism® software).

7. Troubleshooting

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

Symptoms	Causes and Comments
Low luminescence measurements (RLU readout)	<p>Choose an instrument designed for plate-reading luminescence detection. Instruments designed primarily for fluorescence detection are not recommended. Luminometers measure and report luminescence as relative values, and actual RLU numbers will vary between instruments.</p> <p>Insufficient cells per well can lead to low RLU. Handle and plate the cells according to the instructions to ensure a sufficient number of viable cells per well.</p> <p>Low cell viability can lead to low luminescence readout and variability in assay performance.</p> <p>Low activity of Bio-Glo™ Reagent leads to low RLU. Store and handle the Bio-Glo™ Reagent according to the instructions.</p>
Assay performance is variable	<p>Ensure that incubation times are consistent between assays.</p> <p>Ensure the Preparing and Plating protocol (see Section 5.B) is strictly followed.</p> <p>Cells must be treated the same way prior to assay for each assay. Variability in cell growth rates and preculture plating densities will result in variable assay results.</p> <p>Ensure RANKL is prepared and stored properly with carrier protein. Follow manufacturer's protocol for initial rehydration of cytokine. Single use frozen aliquots are recommended for each assay.</p> <p>RANKL lot-to-lot activity differences may be observed. Consult cytokine provider for details.</p>
Weak assay response (low fold induction)	<p>RANKL frozen single-use aliquot has lost biological activity. Follow manufacturer's recommendation for storage and stability.</p> <p>If untreated control RLU is less than 100X above plate reader background RLU, subtract plate reader background RLU from all samples prior to calculating fold induction.</p>



8. References

1. Boyce, BF. *et al.* (2007) Biology of RANK, RANKL, and osteoprotegerin. *Arthritis Res. Ther.* **9**, S1.
2. Walsh, MC. *et al.* (2014) Biology of the RANKL–RANK–OPG system in immunity, bone, and beyond. *Fron. Immunol.* **5(511)**, 1–11.
3. Sigl, V. *et al.* (2016) RANKL/RANK: from bone loss to the prevention of breast cancer. *Open Biol.* **6(11)**, 160230.
4. Renema, N. *et al.* (2016) RANK–RANKL signalling in cancer. *Biosci Rep.* **36(4)**, e00366.
5. Infante, M. *et al.* (2019) RANKL/RANK/OPG system beyond bone remodeling: involvement in breast cancer and clinical perspectives. *J Exp Clin Cancer Res.* **38(12)**.
6. Dubois, EA. (2011) Denosumab. *Br J Clin Pharmacol.* **71(6)**, 804–6.

9. Appendix

9.A. Composition of Buffers and Solutions

thaw medium

90% McCoy's 5A

10% fetal bovine serum

Prepare and use within 2 weeks, stored at 4°C.

freeze medium

85% McCoy's 5A

10% fetal bovine serum

5% DMSO

Prepare immediately before use. Maintain at 4°C during use.

growth medium

90% McCoy's 5A

10% fetal bovine serum

600µg/ml Antibiotic G-418 Sulfate

200µg/ml hygromycin B

Prepare and use within 2 weeks, stored at 4°C.

seeding medium

99% RPMI 1640

1% fetal bovine serum

assay buffer

100% RPMI 1640

9.B. Related Products

Cytokine and Growth Factor Bioassays

Product	Size	Cat.#
RANKL Bioassay	1 each	JA2701
RANKL Bioassay 5X	1 each	JA2705
VEGF Bioassay	1 each	GA2001
VEGF Bioassay 5X	1 each	GA2005
VEGF Bioassay, Cell Propagation Model	1 each	GA1082
Recombinant VEGF	10µg	J2371
IL-2 Bioassay	1 each	JA2201
IL-2 Bioassay 5X	1 each	JA2205
IL-2 Bioassay, Propagation Model	1 each	J2952
IL-6 Bioassay	1 each	JA2501
IL-6 Bioassay 5X	1 each	JA2505
IL-6 Bioassay, Propagation Model	1 each	J2992
IL-15 Bioassay	1 each	JA2011
IL-15 Bioassay 5X	1 each	JA2015
IL-15 Bioassay, Propagation Model	1 each	J2962

Not for Medical Diagnostic Use.

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 (IL-2)	1 each	J1651
T Cell Activation Bioassay (IL-2) 5X	1 each	J1655

Not for Medical Diagnostic Use.

9.B. Related Products (continued)

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
Fc γ R11a-H ADCP Reporter Bioassay, Complete Kit**	1 each	G9901
Fc γ R11a-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
Control Ab, Anti-CD40	50 μ g	K1181
CTLA-4 Blockade Bioassay	1 each	JA3001
CTLA-4 Blockade Bioassay 5X	1 each	JA3005
CTLA-4 Blockade Bioassay, Propagation Model	1 each	JA1400
Control Antibody, Anti-CTLA-4	100 μ g	JA1020
LAG-3/MCHII Blockade Bioassay	1 each	JA1111
LAG-3/MHCII Blockade Bioassay 5X	1 each	JA1115
LAG-3/MHCII Blockade Bioassay, Propagation Model	1 each	JA1112
TCR Activating Antigen Stock Solution	500 μ l	K1201
Control Ab, Anti-LAG-3	100 μ g	K1150
PD-1/PD-L1 Blockade Bioassay	1 each	J1250
PD-1/PD-L1 Blockade Bioassay 5X	1 each	J1255
PD-1 Blockade Bioassay, Propagation Model	1 each	J1252
PD-L1 Negative Cells	1 each	J1191
Control Ab, Anti-PD-1	100 μ g	J1201

Immune Checkpoint Bioassays (continued)

Product	Size	Cat.#
OX40 Bioassay	1 each	JA2191
OX40 Bioassay 5X	1 each	JA2195
OX40 Bioassay, Propagation Model	1 each	J2172
Control Antibody, Anti-OX40	50µg	K1191
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-TIGIT	100µg	J2051
4-1BB Bioassay	1 each	JA2351
4-1BB Bioassay 5X	1 each	JA2355
4-1BB Bioassay, Propagation Model	1 each	J2332
Control Ab, Anti-4-1BB	50µg	K1161

Detection Reagents

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

Not for Medical Diagnostic Use.

Detection Instrument

Product	Size	Cat.#
GloMax® Discover System	1 each	GM3000

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/



10. Summary of Changes

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

1. Updated the fonts and cover image.
2. Revised text about the label in Section 3.
3. Removed an expired patent statement.
4. Updated a third party trademark.
5. Made miscellaneous text edits.

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

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

^(b) Patents Pending.

© 2020–2025 Promega Corporation. All Rights Reserved.

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

Accutase is a registered trademark of Innovative Cell Technologies, Inc. Corning is a registered trademark of Corning, Inc. GIBCO is a registered trademark of Life Technologies Corporation. GraphPad Prism is a registered trademark of GraphPad Software, Inc. JMP is a registered trademark of SAS Institute, Inc. Mr. Frosty is a registered trademark of Nalge Nunc International Corporation. Styrofoam is a registered trademark of Dow Chemical Company.

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.