

TECHNICAL MANUAL

Lumit[®] IL-4 (Human) Immunoassay

Instructions for Use of Products
W6060, W6061 and W6062

Lumit[®] IL-4 (Human) Immunoassay

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 | 3 |
| 3. Before You Begin | 4 |
| 4. Direct (No-Transfer) Protocol for Cultured Cells | 6 |
| 4.A. Cell Plating and Treatment..... | 6 |
| 4.B. Preparing Human IL-4 Standard Dilution | 7 |
| 4.C. Adding 5X Anti-hIL-4 Antibody Mixture to Assay Wells..... | 8 |
| 4.D. Adding Lumit [®] Detection Reagent B to Assay Wells | 9 |
| 5. Optional Sample Transfer Protocol | 10 |
| 5.A. Cell Plating and Treatment..... | 10 |
| 5.B. Preparing Human IL-4 Standard Dilutions..... | 11 |
| 5.C. Adding 2X Anti-hIL-4 Antibody Mixture to Sample Wells..... | 12 |
| 5.D. Adding Lumit [®] Detection Reagent B to Sample Wells..... | 13 |
| 6. Calculating Results | 14 |
| 7. Representative Data | 14 |
| 8. Troubleshooting | 16 |
| 9. Appendix..... | 17 |
| 9.A. Assessing Cross-Reactivity of Lumit [®] Cytokine Immunoassays | 17 |
| 9.B. Multiplexing Assays with Lumit [®] IL-4 (Human) Immunoassay..... | 18 |
| 9.C. References | 19 |
| 9.D. Related Products..... | 19 |
| 10. Summary of Changes | 21 |

1. Description

The Lumit[®] IL-4 (Human) Immunoassay^(a,b) is a homogeneous, bioluminescent assay for detecting interleukin-4 (IL-4) released from cells without the need for sample transfers or wash steps. Produced by a subset of immune cells, including T helper type 2 (Th2) cells, basophils, eosinophils and mast cells, IL-4 is an important immunoregulatory cytokine involved in antibody production, hematopoiesis, inflammation and effector T cell responses (1). IL-4 has a major role in asthma and allergic inflammation (2) and a more complex involvement in cancer biology (3,4), with potential relevance to disease treatment. It has also been reported to function in immunity and higher functions of the brain, including memory and learning (5). Hence, measuring IL-4 production in cell culture models is an essential activity in multiple areas of basic research and drug discovery.

The Lumit[®] IL-4 (Human) Immunoassay has been developed for use with cell culture samples. Lumit[®] reagents can be dispensed directly into microplate wells containing cells and culture medium. Alternatively, medium from cell wells can be transferred to a separate plate for analysis. Assay performance with additional sample types must be determined by the user.

The Lumit[®] IL-4 (Human) Immunoassay is based on NanoLuc[®] Binary Technology (NanoBiT[®]). NanoBiT[®] is a luminescent structural complementation system designed for biomolecular interaction studies (6,7). The system is composed of two subunits, Large BiT (LgBiT; 18kDa) and Small BiT (SmBiT; 11 amino acid peptide), that have been optimized for stability and minimal spontaneous association. In this assay, a sample is incubated with a pair of anti-human IL-4 monoclonal antibodies covalently labeled with SmBiT or LgBiT. When the labeled antibodies recognize and bind to released IL-4, the complementary LgBiTs and SmBiTs are brought into proximity, thereby reconstituting the NanoBiT[®] enzyme and generating luminescence in the presence of the Lumit[®] substrate. Luminescence generated is directly proportional to the amount of analyte present in the sample.

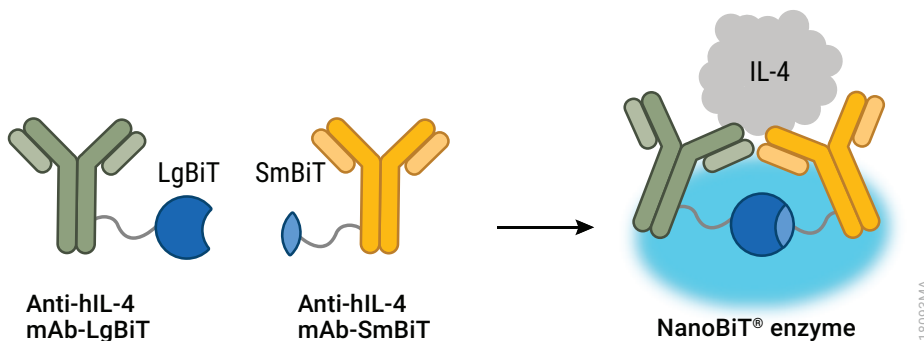


Figure 1. Assay principle. Primary monoclonal antibodies to human IL-4 are labeled with SmBiT and LgBiT. In the presence of IL-4, SmBiT and LgBiT are brought into close proximity, forming the NanoBiT[®] enzyme. When Lumit[®] Detection Reagent B is added, a bright luminescent signal is generated.

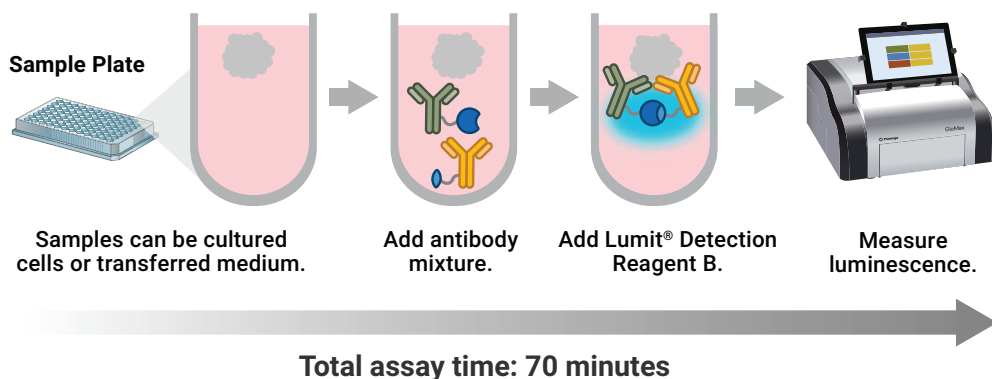


Figure 2. Assay protocol. The Lumit® IL-4 (Human) Immunoassay is performed directly on cells in culture or on medium transferred from the cell culture plate to a new assay plate. The Lumit® Immunoassay protocol does not require wash steps and is complete in 70 minutes.

2. Product Components and Storage Conditions

| PRODUCT | SIZE | CAT. # |
|--|-------------------|--------------|
| Lumit® IL-4 (Human) Immunoassay | 100 assays | W6060 |

Sufficient for 100 assays in 96-well plates; volumes can be adjusted for alternate plate sizes. Includes:

- 30µl Anti-hIL-4 mAb-SmBiT, 500X
- 30µl Anti-hIL-4 mAb-LgBiT, 500X
- 25µl Human IL-4 Standard
- 160µl Lumit® Detection Substrate B
- 3.2ml Lumit® Detection Buffer B

| PRODUCT | SIZE | CAT. # |
|--|---------------------|--------------|
| Lumit® IL-4 (Human) Immunoassay | 1,000 assays | W6061 |

Sufficient for 1,000 assays in 96-well plates; volumes can be adjusted for alternate plate sizes. Includes:

- 300µl Anti-hIL-4 mAb-SmBiT, 500X
- 300µl Anti-hIL-4 mAb-LgBiT, 500X
- 25µl Human IL-4 Standard
- 1.25ml Lumit® Detection Substrate B
- 25ml Lumit® Detection Buffer B

2. Product Components and Storage Conditions (continued)

| PRODUCT | SIZE | CAT. # |
|--|-----------------------|--------------|
| Lumit® IL-4 (Human) Immunoassay | 5 × 100 assays | W6062 |

Sufficient for 500 assays in 96-well plates; volumes can be adjusted for alternate plate sizes. Includes:

- 5 × 30µl Anti-hIL-4 mAb-SmBiT, 500X
- 5 × 30µl Anti-hIL-4 mAb-LgBiT, 500X
- 25µl Human IL-4 Standard
- 5 × 160µl Lumit® Detection Substrate B
- 5 × 3.2ml Lumit® Detection Buffer B

Storage Conditions: Store complete kit at less than –65°C upon receipt. Alternatively, store the Human IL-4 Standard at less than –65°C and all other components at –30°C to –10°C. After thawing, store Human IL-4 Standard at +2°C to +10°C for up to 1 month. If storing the Human IL-4 Standard for more than 1 month after thawing, dispense into aliquots and store at less than –65°C. After thawing, store the Anti-hIL-4 mAb-SmBiT and Anti-hIL-4 mAb-LgBiT at –30°C to –10°C. After thawing, store Lumit® Detection Buffer B at room temperature. Store Lumit® Detection Substrate B protected from light.

3. Before You Begin

There are two protocols for measuring human IL-4. The direct protocol typically achieves higher sensitivity than the transfer protocol and requires less sample manipulation. The optional sample transfer protocol provides flexibility for same well sampling during treatment exposure time courses and split-sample analysis for assessment of multiple cytokine levels from the same sample.

Direct (No-Transfer) Protocol for Cultured Cells (Section 4): Measure human IL-4 directly in cell culture wells. Add 20µl of a 5X antibody mixture to 80µl of cells or IL-4 standard dilutions in culture medium and incubate for 60 minutes. Following incubation, add 25µl of Lumit® Detection Reagent B and record luminescence.

Optional Sample Transfer Protocol (Section 5): Measure human IL-4 in medium samples transferred from treated cell wells. Transfer 50µl of culture medium from cell wells to a separate assay plate. Add 50µl of a 2X antibody mixture to 50µl of transferred sample or standard dilutions and incubate for 60 minutes. Following incubation, add 25µl of Lumit® Detection Reagent B and record luminescence.

Note: Assay volumes are scalable and can be adjusted based on sample sizes. The protocols in Sections 4 and 5 list common volumes for 96- and 384-well plates. Other volumes may be used, maintaining the recommended antibody and detection reagent final concentrations. We recommend using standard tissue culture medium supplemented with 5–10% fetal bovine serum (FBS). Lesser concentrations of FBS may produce higher background (nonanalyte-mediated antibody pairing) and higher replicate variability. The use of medium without phenol red may increase assay sensitivity and reduce inner-filter effects from luminescence quenching.

Reagent Preparation and Storage

Prepare the Human IL-4 Standard dilution series (Section 4.B or 5.B), the Anti-hIL-4 Antibody Mixture (Section 4.C or 5.C) and Lumit[®] Detection Reagent B (Section 4.C or 5.C) on the day of use. Do **not** reuse Human IL-4 Standard dilutions, the Anti-hIL-4 Antibody Mixture or the Lumit[®] Detection Reagent B.



Use personal protective equipment and follow your institution's safety guidelines and disposal requirements when working with biohazardous materials such as cells and cell culture reagents.

Plate Map

| | IL-4 Standard Curve (pg/ml)* | | Test Samples | | | | | | | | | |
|----------|---------------------------------|--------|--------------|---|---|---|---|---|---|----|----|----|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | 25,000 | 25,000 | | | | | | | | | | |
| B | 7,500 | 7,500 | | | | | | | | | | |
| C | 2,250 | 2,250 | | | | | | | | | | |
| D | 675 | 675 | | | | | | | | | | |
| E | 203 | 203 | | | | | | | | | | |
| F | 60.8 | 60.8 | | | | | | | | | | |
| G | 18.2 | 18.2 | | | | | | | | | | |
| H | 0 | 0 | | | | | | | | | | |

***Note:** The indicated Human IL-4 Standard dilution series listed above and described in the subsequent protocol are only a recommended dilution series.

Materials to Be Supplied by the User

- cells (human-derived)
- culture medium (e.g., RPMI 1640, GIBCO[®] Cat.# 22400-089 + 10% heat-inactivated fetal bovine serum, GIBCO[®] Cat.# A3840001)
- white, multiwell tissue culture plates (solid white or white with clear bottom) compatible with a luminometer (e.g., 96-well Corning[®] Cat.# 3917)
- multichannel pipette or automated pipetting station
- dilution tubes or multichamber, dilution reservoir (e.g., Dilux[®] D-1002)
- reagent reservoir trays (e.g., Midwest Scientific Cat.# RR25)
- plate shaker for mixing multiwell plates
- luminometer capable of reading multiwell plates (e.g., GloMax[®] Discover System, Cat.# GM3000)

4. Direct (No-Transfer) Protocol for Cultured Cells

This protocol describes how to detect IL-4 released directly in assay wells containing cells and culture medium. For quantitation purposes, a standard curve is generated using an IL-4 standard diluted in culture medium.

4.A. Cell Plating and Treatment

1. Plate cells into a 96- or 384-well, white (or white with clear bottom) tissue culture plate. Leave columns 1 and 2 empty. Allow cells to attach if using adherent cells.

Note: While the broad dynamic range of the assay offers considerable flexibility for cell number, the optimal number of cells dispensed per well for a specific cell model should be empirically determined. Ensure the maximum level of cytokine released does not exceed the linear range of the detection chemistry. Within that constraint, cell number can be increased to meet the detection requirements of low-level cytokine production. See example cell numbers used for human Th2 and PBMC in Figures 5 and 6.

2. Treat cells by adding a volume of test agent to each well such that the total volume is as follows:

96-well plate: 80µl per well.

384-well plate: 20µl per well.

For example, if 60µl of cells are plated per well in a 96-well plate, add 20µl of 4X treatment agent in culture medium. Cells are typically treated for 6–24 hours or longer, depending on stimuli, to observe release of significant levels of IL-4.

Optional: If manually dispensing into the 384-well assay format, after various additions, briefly centrifuge the microplate (700–900rpm for approximately 10 seconds) followed by a brief mix with a plate shaker to ensure proper mixing.

4.B. Preparing Human IL-4 Standard Dilution

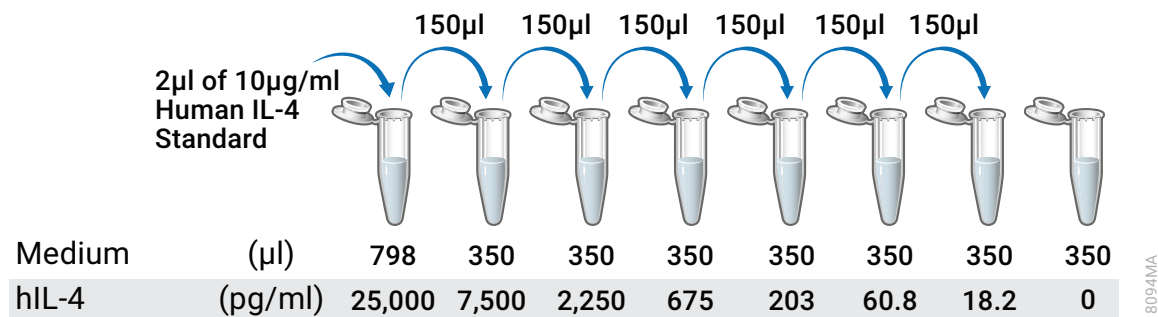


Figure 3. Human IL-4 dilution series.

Shortly before completing cell treatments, prepare IL-4 dilutions in the identical culture medium used for cell samples.

Note: If using multiple cell models each requiring different culture medium, separate standard dilution series must be generated in each medium used in the study.

1. Thaw the Human IL-4 Standard (approximately 15 minutes at room temperature) immediately before use.
2. Briefly centrifuge the tube before opening, then mix by pipetting.
3. Prepare an initial concentration of 25,000pg/ml human IL-4 by diluting Human IL-4 Standard (10µg/ml) 1:400 in pre-warmed cell culture medium. For example, prepare 800µl of 25,000pg/ml human IL-4 by adding 2µl of the Human IL-4 Standard stock to 798µl of culture medium (see Figure 3).
4. Set up seven tubes (or seven chambers in a dilution reservoir) with 350µl of culture medium in each.
5. Prepare 3.33-fold serial dilutions of standard. Transfer 150µl from the 25,000pg/ml initial human IL-4 dilution (Step 3) to 350µl of culture medium for the second dilution. Mix by pipetting and repeat five times to generate seven standard dilutions with a range of 25,000pg/ml to 18.2pg/ml. The last well or chamber should contain only culture medium as the background control.

Note: Change pipette tips between each dilution step and use aerosol filter tips to avoid analyte carryover.

6. After the cell treatment is complete, add the standard dilutions and background control in duplicate to two columns in the assay plate (see the plate map in Section 3).

96-well plate: Dispense 80µl per well.

384-well plate: Dispense 20µl per well.

Notes:

- a. Unused Human IL-4 Standard (10µg/ml) can be stored at +2°C to +10°C for 1 month. If storing the standard for more than 1 month, dispense into aliquots and store at less than -65°C. Avoid multiple freeze-thaw cycles.
- b. We recommend incorporating Human IL-4 Standard controls on each assay plate for normalization.
- c. Additional Lumit® Human IL-4 Standard (10µg/ml) is available for standalone purchase (Cat. # W125A-C).

4.C. Adding 5X Anti-hIL-4 Antibody Mixture to Assay Wells

Note: If using multiple cell models each requiring different culture medium, separate 5X Anti-hIL-4 antibody mixtures must be generated in each medium used in the study.

1. Remove the Anti-hIL-4 antibodies from storage immediately before use. Thaw if necessary.
Note: Remove Lumit[®] Detection Buffer B from storage at the same time and equilibrate to room temperature if not already thawed. Use a water bath to accelerate buffer warming as necessary. If using the 25ml buffer in Cat.# W6061, you may need to initiate buffer warming further in advance of its use in Section 4.D.
2. Briefly centrifuge the Anti-hIL-4 antibody tubes before opening, then mix by pipetting.
3. Immediately prior to use, prepare a 5X antibody mixture by diluting both antibodies 1:100 into a single volume of prewarmed culture medium. Pipet to mix the antibody solution. To assay a complete 96- or 384-well plate, including some reagent volume for pipetting loss, prepare the 5X antibody mixture as follows:

| Reagent | Volume |
|----------------------|---------|
| culture medium | 2,352μl |
| Anti-hIL-4 mAb-SmBiT | 24μl |
| Anti-hIL-4 mAb-LgBiT | 24μl |

4. Add the 5X Anti-hIL-4 antibody mixture to wells containing cultured cells or standard dilutions, carefully avoiding cross contamination between wells by changing pipette tips if moving from high to low analyte levels.
96-well plate: Dispense 20μl/well of 5X Anti-hIL-4 antibody mixture to 80μl/well of cells or IL-4 standard dilutions.
384-well plate: Dispense 5μl/well of 5X Anti-hIL-4 antibody mixture to 20μl/well of cells or IL-4 standard dilutions.
5. Briefly mix with a plate shaker (e.g., 10 seconds at 250–350rpm).
6. Incubate for 45 minutes at 37°C in a humidified 5% CO₂ incubator.

4.D. Adding Lumit[®] Detection Reagent B to Assay Wells

While cells are incubating with the Anti-hIL-4 antibody mixture (Section 4.C), prepare the Lumit[®] Detection Reagent B.

1. Equilibrate the required volume of Lumit[®] Detection Buffer B to room temperature.
2. Remove the Lumit[®] Detection Substrate B from storage and mix. If Lumit[®] Detection Substrate B has collected in the cap or on the sides of the tube, briefly centrifuge.
3. Prepare a 1:20 dilution of Lumit[®] Detection Substrate B in room temperature Lumit[®] Detection Buffer B to create enough volume of Lumit[®] Detection Reagent B for the number of wells to be assayed. To assay a 96- or 384-well assay plate, including some reagent volume for pipetting loss, prepare 5X Lumit[®] Detection Reagent B as follows:

| Reagent | Volume |
|--|---------|
| Lumit [®] Detection Buffer B | 3,040µl |
| Lumit [®] Detection Substrate B | 160µl |

Notes:

- a. The 1,000 assay size Lumit[®] IL-4 (Human) Immunoassay (Cat.# W6061) contains 25ml of Lumit[®] Detection Buffer B and 1.25ml of Lumit[®] Detection Substrate B. There is sufficient overfill to prepare Lumit[®] Detection Reagent B for analyzing 5 or 10 plates at one time. If Cat.# W6061 is used for assaying 10 plates individually, mix 2,375µl of Lumit[®] Detection Buffer B + 125µl of Lumit[®] Detection Substrate B for each plate.
 - b. Once reconstituted, the Lumit[®] Detection Reagent B will lose 10% activity in approximately 3 hours at 20°C. At 4°C, the reconstituted reagent will lose 10% activity in approximately 7 hours.
4. After the incubation in Section 4.C, Step 6 is complete, equilibrate the assay plate with cells to room temperature for 15 minutes.
 5. Add room temperature 5X Lumit[®] Detection Reagent B to each assay well of the plate.
96-well plate: Dispense 25µl per well.
384-well plate: Dispense 6.25µl per well.
 6. Briefly mix with a plate shaker (e.g., 10 seconds at 300–500rpm).
 7. Incubate at room temperature for 3–5 minutes.
 8. Read luminescence.

Note: Assay signal is stable with a half-life of approximately 2 hours, compatible with batch processing of multiple assay plates. We recommend incorporating Human IL-4 Standard controls on each assay plate for normalization.

5. Optional Sample Transfer Protocol

This protocol describes transfer of sample medium from treated cell wells into a separate assay plate for subsequent cytokine detection, leaving the cells and remaining medium for additional uses. For quantitation purposes, a dilution series of Human IL-4 Standard prepared in culture medium is used to generate a standard curve.

5.A. Cell Plating and Treatment

1. Plate cells into a 96- or 384-well tissue culture plate. Leave columns 1 and 2 empty. Allow cells to attach if using adherent cells.

Note: While the broad dynamic range of the assay offers considerable flexibility for cell number, the optimal number of cells dispensed per well for a specific cell model should be empirically determined. Ensure that the maximum level of cytokine released does not exceed the linear range of the detection chemistry. Within that constraint, cell number can be increased to meet the detection requirements of low-level cytokine production. See example cell numbers used for human Th2 and PBMC in Figures 5 and 6.

2. Treat cells by adding a volume of test agent to each well. The final treatment volume is flexible. Typical volumes are:

96-well plate: 100–200µl per well

384-well plate: 25–50µl per well

Optional: If manually dispensing into the 384-well assay format, after various additions, briefly centrifuge the microplate (700–900rpm for approximately 10 seconds) then briefly mix with a plate shaker to ensure proper mixing.

3. After cell treatment is complete, transfer cell medium from each well to the corresponding wells of a separate white assay plate:

96-well plate: Transfer 50µl per well.

384-well plate: Transfer 12.5µl per well.

Notes:

- a. If lower sample volumes are transferred, dilute with culture medium to 50µl for 96-well format and 12.5µl for 384-well format, subsequently accounting for sample dilution when calculating actual cytokine concentration released in cell wells.
- b. For enhanced 384-well assay sensitivity, if needed, you can transfer 20µl of sample (and standards) to assay wells, then add 5µl/well of 5X Anti-hIL-4 antibody mixture (prepared in Section 4.C).

5.B. Preparing Human IL-4 Standard Dilutions

Shortly before completing cell treatments, prepare IL-4 dilutions in the identical culture medium used for cell samples.

Note: If using multiple cell models, each requiring different culture medium, separate standard dilution series must be generated in each medium used.

1. Thaw the Human IL-4 Standard (approximately 15 minutes) immediately before use.
2. Briefly centrifuge the tube before opening, then mix by pipetting.
3. Prepare an initial concentration of 25,000pg/ml human IL-4 by diluting Human IL-4 Standard (10µg/ml) 1:400 in prewarmed cell culture medium. For example, prepare 800µl of 25,000pg/ml human IL-4 by adding 2µl of Human IL-4 Standard stock to 798µl of culture medium (Figure 3).
4. Set up seven tubes (or seven chambers in a dilution reservoir) with 350µl of culture medium in each.
5. Proceed with 3.33-fold serial dilutions of standard. Transfer 150µl from the 25,000pg/ml stock to 350µl of culture medium for the second dilution. Mix and repeat five times to generate seven standard dilutions with a range of 25,000pg/ml to 18.2pg/ml. The last well or chamber should contain only culture medium as the background control.

Note: Change pipette tips between each dilution step and use aerosol filter tips to avoid analyte carryover.

6. After transferring the culture medium from the treated cell wells to a separate assay plate, add the standard dilutions and background control in duplicate to two columns in the transfer plate (see the plate map in Section 3).

96-well plate: Dispense 50µl per well.

384-well plate: Dispense 12.5µl per well.

Notes:

- a. Unused Human IL-4 Standard (10µg/ml) can be stored at +2°C to +10°C for 1 month. If storing the standard for more than 1 month, dispense into aliquots and store at less than -65°C. Avoid multiple freeze-thaw cycles.
- b. We recommend incorporating Human IL-4 Standard controls on each assay plate for normalization.
- c. Additional Lumit® Human IL-4 Standard (10µg/ml) is available for standalone purchase (Cat. # W125A-C).

5.C. Adding 2X Anti-hIL-4 Antibody Mixture to Sample Wells

Note: If using multiple cell models each requiring different culture medium, separate 2X Anti-hIL-4 antibody mixtures must be generated in each medium used in the study.

1. Remove the Anti-hIL-4 antibodies from storage immediately before use. Thaw if necessary.
Note: Remove Lumit[®] Detection Buffer B from storage at the same time and equilibrate to room temperature if not already thawed. Use a water bath to accelerate buffer warming as necessary. If using the 25ml buffer in Cat.# W6061, you may need to initiate buffer warming further in advance of its use in Section 5.D.
2. Briefly centrifuge the Anti-hIL-4 antibody tubes before opening, then mix by pipetting.
3. Immediately prior to use, prepare a 2X antibody mixture by diluting both antibodies 1:250 into a single volume of prewarmed cell culture medium. Pipet to mix the antibody solution. To assay a complete 96- or 384-well plate, including some reagent volume for pipetting loss, prepare the 2X antibody mixture as follows:

| Reagent | Volume |
|----------------------|---------|
| culture medium | 5,952µl |
| Anti-hIL-4 mAb-SmBiT | 24µl |
| Anti-hIL-4 mAb-LgBiT | 24µl |

4. Add the 2X Anti-hIL-4 antibody mixture to transferred culture medium (samples) or IL-4 standard dilutions, carefully avoiding cross contamination between wells by changing pipette tips if moving from high to low analyte levels.

96-well plate: Dispense 50µl/well of 2X Anti-hIL-4 antibody mixture to 50µl/well of medium or standard dilutions.

384-well plate: Dispense 12.5µl/well of 2X Anti-hIL-4 antibody mixture to 12.5µl/well of medium or standard dilutions.

5. Briefly mix with a plate shaker (e.g., 10 seconds at 250–350rpm).
6. Incubate for 60 minutes at room temperature.

Note: To incubate at room temperature, a HEPES-containing medium will provide best results. Without HEPES, the buffering capacity outside of a CO₂ incubator is limited. The plates can also be incubated at 37°C in a CO₂ incubator if subsequently equilibrated to room temperature prior to addition of detection reagent.

5.D. Adding Lumit[®] Detection Reagent B to Sample Wells

While the samples and standard dilutions are incubating with the Anti-hIL-4 antibody mixture (Section 5.C), prepare the Lumit[®] Detection Reagent B.

1. Equilibrate the required volume of Lumit[®] Detection Buffer B to room temperature.
2. Remove the Lumit[®] Detection Substrate B from storage and mix. If the Lumit[®] Detection Substrate B has collected in the cap or on the sides of the tube, briefly centrifuge.
3. Prepare a 1:20 dilution of Lumit[®] Detection Substrate B into room temperature Lumit[®] Detection Buffer B to create enough volume of Lumit[®] Detection Reagent B for the number of wells to be assayed. For a 96- or 384-well assay plate, including some excess reagent volume, prepare 5X Lumit[®] Detection Reagent B as follows:

| Reagent | Volume |
|--|---------|
| Lumit [®] Detection Buffer B | 3,040µl |
| Lumit [®] Detection Substrate B | 160µl |

Notes:

- a. The 1,000 assay size of Lumit[®] IL-4 (Human) Immunoassay (Cat.# W6061) contains 25ml of Lumit[®] Detection Buffer B and 1.25ml of Lumit[®] Detection Substrate B. There is sufficient overfill to prepare Lumit[®] Detection Reagent B for analyzing 5 or 10 plates at once. If Cat.# W6061 is used for assaying 10 plates individually, mix 2,375µl of Lumit[®] Detection Buffer B + 125µl of Lumit[®] Detection Substrate B for each plate.
 - b. Once reconstituted, the Lumit[®] Detection Reagent B loses 10% activity in approximately 3 hours at 20°C. At 4°C the reconstituted reagent loses 10% activity in approximately 7 hours.
4. After the incubation in Section 5.C, Step 6, is complete, add room temperature Lumit[®] Detection Reagent B to each assay well of the plate:
96-well plate: Dispense 25µl per well.
384-well plate: Dispense 6.25µl per well.
 5. Briefly mix with a plate shaker (e.g., 10 seconds at 300–500rpm).
 6. Incubate at room temperature for 3–5 minutes.
 7. Read luminescence.

Note: Assay signal is stable with a half-life of approximately 2 hours, compatible with batch processing of multiple assay plates.

6. Calculating Results

Create a standard curve for the known cytokine concentrations using software (e.g., GraphPad® Prism) capable of nonlinear regression analysis or cubic spline curve fitting.

Subsequently, interpolate the concentration of cytokine in various cell samples. The broad dynamic range of the Lumit® standard curve closely approaches linearity and is well-suited for second- or third-order polynomial regression curve fitting, as well as cubic spline curve fitting. Four-parameter logistic (4PL) curve fitting is also commonly used, but may not be ideal since the broad, linear dynamic range for the Lumit® standard curve is not well-suited for sigmoidal curve fitting (8).

Alternatively, while somewhat less accurate, a Log-Log plot of average RLU (background-subtracted) vs. cytokine standard concentrations can be fit with the Power trendline in Microsoft Excel® (see Section 7) and subsequently used for interpolation of the concentration of cytokine release in various cell samples.

7. Representative Data

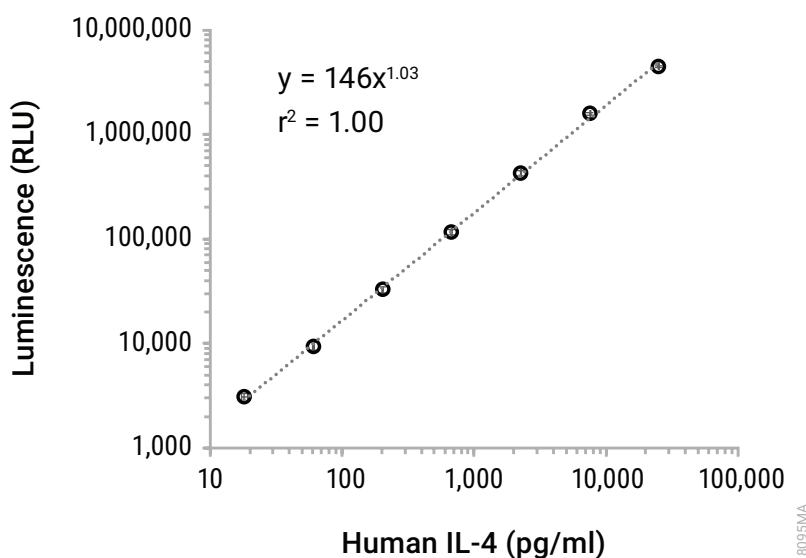


Figure 4. Standard curve for the Lumit® IL-4 (Human) Immunoassay. This is a representative standard curve and should not be used for calculation of unknowns. Generate a standard curve on each assay plate to interpolate the concentration of cytokine in experimental samples. **Note:** The plotted luminescence values were determined by subtracting background RLU.

Table 1. Assay Precision. Three samples of known concentrations of human IL-4 were tested 20 times on two plates to assess assay precision. A standard curve was used on each assay plate to interpolate the IL-4 quantities in each well by using cubic spline fitting with GraphPad® Prism software.

| Assay Precision | | | |
|-----------------------------|--------|--------|--------|
| Number of Replicates | 40 | 40 | 40 |
| Expected (pg/ml) | 20,000 | 2,000 | 200 |
| Mean (pg/ml) | 18,632 | 1,865 | 167 |
| Standard Deviation | 620 | 101 | 13 |
| Percent CV | 3.3 | 5.4 | 7.6 |
| Average Percent of Expected | 93.2 | 93.2 | 83.7 |
| Percent Range | 84–105 | 83–112 | 73–108 |

Note: The minimal detectable dose (MDD), determined at two standard deviations above background, was determined to be 4.5 ± 0.3 pg/ml of hIL-4.

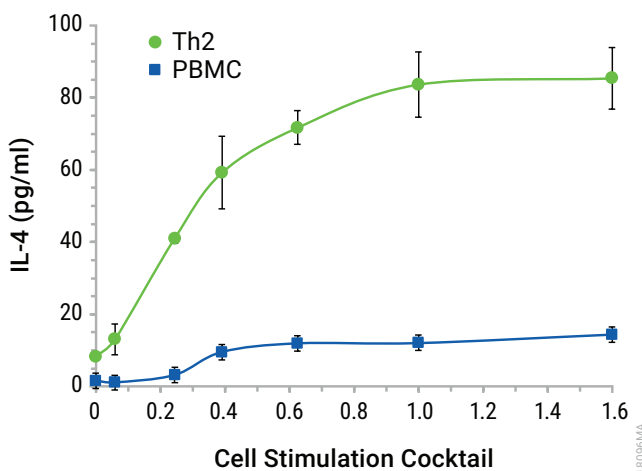


Figure 5. Lumit® detection of IL-4 released from human Th2 and PBMC. Human T helper type 2 cells (Th2; differentiated from CD4+ T cells) and peripheral blood mononuclear cells (PBMC; BioIVT preparation pooled from four donors) were plated in 96-well plate format at 30,000 cells/well in RPMI 1640 (for Th2 cells, supplemented with 0.5mM β -ME and 1mM pyruvate) + 10% heat-inactivated FBS. The next day, cells were treated in triplicate for 24 hours with increasing concentrations of Cell Stimulation Cocktail (CSC; Invitrogen Cat. # 00-4970-93). 1X CSC is a mixture of 81nM phorbol 12-myristate 13-acetate (PMA) and 1.34 μ M ionomycin. Following treatment, the 5X Anti-hIL-4 antibody mixture was dispensed to the cell wells and incubated for 1 hour before Lumit® Detection Reagent B was added. Luminescence readings were analyzed using cubic spline interpolation (GraphPad® Prism 8) against a standard curve to determine levels of IL-4 released.

8. 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 |
|--|---|
| No signal from treated cells | Make sure that human cells are used with the Lumit [®] IL-4 (Human) Immunoassay. The Lumit [®] IL-4 (Human) Immunoassay will not effectively detect IL-4 from other species. |
| Human IL-4 standard curve is not linear | The 3.33-fold dilutions should be carefully created without carryover from a higher concentration. We recommend changing aerosol filter pipette tips after each dilution step to prevent carryover. The sensitivity of the assay and the broad linear range (>3 logs) means that any carryover will disrupt the linear range. Also, make sure that no IL-4 contaminates the background control. |
| The relative light units (RLU) for the standard curve are low and/or variable. | Warm culture medium, Human IL-4 Standard, and Lumit [®] Detection Buffer B to room temperature before use. There may be some variation in RLU due to culture conditions, temperature, etc., but as long as the standard curve is run on the same plate as the test samples under the same conditions, released IL-4 can be accurately quantitated. |
| Assay background is significantly higher than previous experience | On rare occasion, assay background can be influenced by the particular lot of FBS used in the culture medium. Test assay performance using an alternative lot of FBS. In addition, ensure that the antibody mixture used in the assay is not prepared longer before use than necessary. |

9. Appendix

9.A. Assessing Cross-Reactivity of Lumit® Cytokine Immunoassays

Table 2. Testing Immunoassay Cross-Reactivity. The specificity of Lumit® cytokine immunoassays were assessed against a panel of their respective recombinant standards. Each Lumit® cytokine immunoassay was applied to 10ng/ml of the indicated cytokine standards in triplicate assay wells, in addition to wells containing only medium for background determination. Background luminescence was subtracted prior to normalization of average assay signals with each cytokine to that of the intended target of the Lumit® cytokine immunoassay. For the Lumit® IL-4 (Human) Immunoassay, signals obtained from the tested cytokines were less than or equal to 0.06% of the signal from human IL-4.

| Standard (10ng/ml) | Lumit® Cytokine Immunoassay (Percent Assay Control) | | | | | | |
|-----------------------|---|-------|-------|-------|--------|--------|--------|
| | hIL-1β | hIL-2 | hIL-4 | hIL-6 | hIL-10 | hIFN-γ | hTNF-α |
| hIL-1β | 100 | 0.01 | 0.02 | 0 | 0 | 0.02 | 0 |
| hIL-2 | 0 | 100 | 0.02 | 0 | 0 | 0.01 | 0 |
| hIL-4 | 0.06 | 0.01 | 100 | 0 | 0.02 | 0.01 | 0 |
| hIL-6 | 0.07 | 0.02 | 0.06 | 100 | 0.02 | 0.03 | 0 |
| hIL-10 | 0.18 | 0.03 | 0.03 | 0.03 | 100 | 0.04 | 0 |
| hIFN-γ | 0.10 | 0.03 | 0.03 | 0.03 | 0.02 | 100 | 0.03 |
| hTNF-α | 0.01 | 0.01 | 0.02 | 0.02 | 0.01 | 0.03 | 100 |

9.B. Multiplexing Assays with Lumit[®] IL-4 (Human) Immunoassay

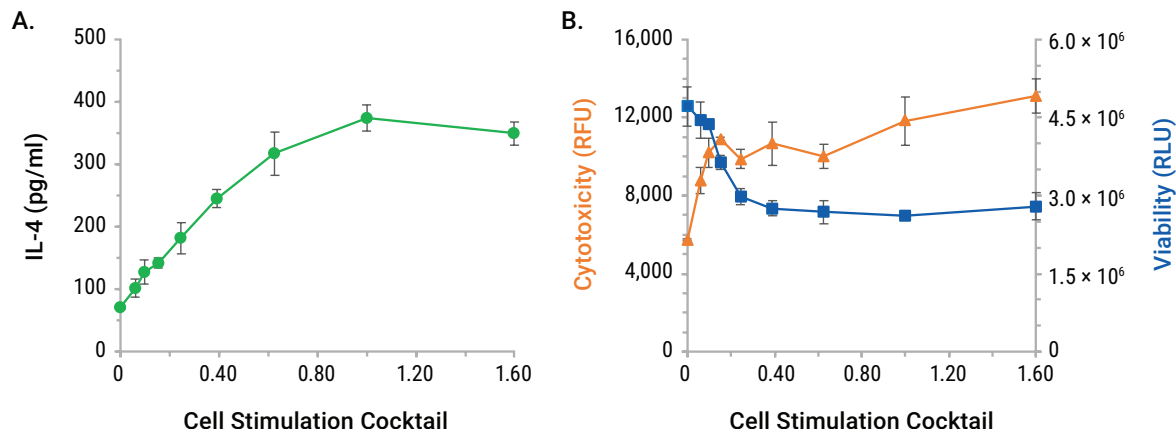


Figure 6. Same-well determination of IL-4 release, cytotoxicity and viability for treated human Th2 cells. Human T helper type 2 (Th2) cells, prepared by differentiated CD4⁺ T cells, were plated in RPMI 1640 (supplemented with 0.5mM β-ME and 1mM pyruvate) + 10% heat-inactivated FBS at 30,000 cells/well in an all-white 96-well plate, then primed overnight with 5ng/ml recombinant human IL-2. Subsequently, cells were treated in a total volume of 70μl with increasing concentrations of Cell Stimulation Cocktail (CSC) for 24 hours. Following treatment, 10μl of 10X CellTox[™] Green Reagent (prepared in Assay Buffer as described in *CellTox[™] Green Cytotoxicity Assay Technical Manual #TM375*) and 20μl of 5X Anti-hIL-4 antibody mixture (Section 4) were added to each well. The assay plate was incubated for 45 minutes at 37°C in a humidified 5% CO₂ incubator prior to measuring fluorescence (RFU) indicating cytotoxicity. The assay plate was allowed to equilibrate to room temperature for 15 minutes before adding 25μl of Lumit[®] Detection Reagent B to assess human IL-4 release. Lastly, 100μl of the lytic CellTiter-Glo[®] Reagent was added to each well (as described in the *CellTiter-Glo[®] Luminescent Cell Viability Assay Technical Bulletin #TB288*) and the plate was shaken for 2 minutes at 350rpm, then incubated for 10 minutes before luminescence (RLU) signals were measured. Luminescence reflects cell viability.

Panel A. Dose-dependent release of human IL-4 was observed in response to CSC treatment. **Panel B.** Some degree of increased cytotoxicity and decreased cell viability was observed in response to 24-hour treatment with increasing CSC.

Note: Maximal cell death provoked by adding a reference cytotoxic agent (50μg/ml digitonin) produced fluorescence readings (using the CellTox[™] Green Cytotoxicity Assay, Cat.# G8741) of more than 20,000 RFU, while viability signals (shown as luminescence using the CellTiter-Glo[®] Viability Assay, Cat.# G7570) were reduced by more than 98% relative to untreated cells (data not shown).

9.C. References

1. Brown, M.A. and Hural, J. (1997) Functions of IL-4 and control of its expression. *Crit. Rev. Immunol.* **17(1)**, 1–32.
2. Steinke, J.W. and Borish, L. (2001) Th2 cytokines and asthma – Interleukin-4: its role in the pathogenesis of asthma, and targeting it for asthma treatment with interleukin-4 receptor antagonists. *Respir. Res.* **2(2)**, 66–70.
3. Li, Z., Chen, L. and Qin, Z. (2009) Paradoxical roles of IL-4 in tumor immunity. *Cell. Mol. Immunol.* **6(6)**, 415–22.
4. Suzuki, A. *et al.* (2015) Targeting of IL-4 and IL-13 receptors for cancer therapy. *Cytokine* **75(1)**, 79–88.
5. Gadani, S.P. *et al.* (2012) IL-4 in the brain: A cytokine to remember. *J. Immunol.* **189(9)**, 4213–9.
6. Dixon, A.S. *et al.* (2016) NanoLuc complementation reporter optimized for accurate measurement of protein interactions in cells. *ACS Chem. Biol.* **11**, 400–8.
7. Hwang, B.B. *et al.* (2020) A homogeneous bioluminescent immunoassay to probe cellular signaling pathway regulation. *Commun. Biol.* **3**, 8.
8. Herman, R.A., Scherer, P.N. and Shan, G. (2008) Evaluation of logistic and polynomial models for fitting sandwich-ELISA calibration curves. *J. Immunol. Methods* **339**, 245–58.

9.D. Related Products

Lumit® Immunoassays

| Product | Size | Cat.# |
|--|------------|---------|
| Lumit® Human IL-4 Standard | 25µl | W125A-C |
| Lumit® FcRn Binding Immunoassay | 100 assays | W1151 |
| Lumit® HMGB1 (Human/Mouse) Immunoassay | 100 assays | W6110 |
| Lumit® Human IL-1β Immunoassay | 100 assays | W6010 |
| Lumit® IFN-γ (Human) Immunoassay | 100 assays | W6040 |
| Lumit® IL-2 (Human) Immunoassay | 100 assays | W6020 |
| Lumit® IL-6 (Human) Immunoassay | 100 assays | W6030 |
| Lumit® IL-10 (Human) Immunoassay | 100 assays | W6070 |
| Lumit® Mouse IL-1β Immunoassay | 100 assays | W7010 |
| Lumit® TNF-α (Human) Immunoassay | 100 assays | W6050 |

Additional sizes available.

9.D. Related Products (continued)

Lumit® Immunoassay Reagents

| Product | Size | Cat. # |
|---|------------|--------|
| Lumit® Detection Reagent B* | 100 assays | VB4050 |
| Lumit® Immunoassay Cellular Systems–Starter Kit | 200 assays | W1220 |
| Lumit® Immunoassay Labeling Kit* | 1 each | VB2500 |

*Additional sizes available.

Inflammation Assays

| Product | Size | Cat. # |
|---------------------------------------|------------|--------|
| Caspase-Glo® 1 Inflammasome Assay | 10ml | G9951 |
| RealTime-Glo™ Extracellular ATP Assay | 200 assays | GA5010 |

Additional sizes available.

Cell Viability Assays

| Product | Size | Cat. # |
|---|---------------|--------|
| CellTiter-Glo® 2.0 Cell Viability Assay | 10ml | G9241 |
| CellTiter-Glo® Cell Viability Assay | 10ml | G7570 |
| RealTime-Glo™ MT Cell Viability Assay | 100 reactions | G9711 |
| CellTiter-Fluor™ Cell Viability Assay | 10ml | G6080 |

Additional sizes available.

Cytotoxicity Assays

| Product | Size | Cat. # |
|-----------------------------------|------|--------|
| LDH-Glo™ Cytotoxicity Assay | 10ml | J2380 |
| CytoTox-Glo™ Cytotoxicity Assay | 10ml | G9290 |
| CellTox™ Green Cytotoxicity Assay | 10ml | G8741 |

Additional sizes available.

Energy Metabolism and Oxidative Stress Assays

| Product | Size | Cat. # |
|--|------|--------|
| NAD/NADH-Glo™ Assay | 10ml | G9071 |
| NADP/NADPH-Glo™ Assay | 10ml | G9081 |
| ROS-Glo™ H ₂ O ₂ Assay | 10ml | G8820 |
| GSH/GSSG-Glo™ Assay | 10ml | V6611 |
| Lactate-Glo™ Assay | 5ml | J5021 |
| Glucose-Glo™ Assay | 5ml | J6021 |
| Glutamine/Glutamate-Glo™ Assay | 5ml | J8021 |
| Glycerol-Glo™ Assay | 5ml | J3150 |
| Triglyceride-Glo™ Assay | 5ml | J3160 |
| Cholesterol/Cholesterol Ester-Glo™ Assay | 5ml | J3190 |

Additional sizes available.

Apoptosis Assays

| Product | Size | Cat. # |
|--|------------|--------|
| Caspase-Glo® 3/7 Assay System | 2.5ml | G8090 |
| Caspase-Glo® 8 Assay System | 2.5ml | G8200 |
| Caspase-Glo® 9 Assay System | 2.5ml | G8210 |
| RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay | 100 assays | JA1011 |

Additional sizes available.

10. Summary of Changes

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

1. Added Human IL-4 Standard to Sections 4.B, 5.B and 9.D.
2. Updated a third party trademark attribution.
3. Updated Lumit to a registered trademark.
4. Updated the fonts.



^(a)U.S. Pat. Nos. 9,797,889; 9,797,890; 10,107,800; 10,648,971; and other patents and patents pending.

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

© 2022–2025 Promega Corporation. All Rights Reserved.

Caspase-Glo, CellTiter-Glo, Lumit, NanoBiT and NanoLuc are registered trademarks of Promega Corporation. CellTiter-Fluor, CellTox, CytoTox-Glo, Ester-Glo, Glucose-Glo, Glutamine/Glutamate-Glo, Glycerol-Glo, GSH/GSSG-Glo, Lactate-Glo, LDH-Glo, NAD/NADH-Glo, NADP/NADPH-Glo, RealTime-Glo, ROS-Glo and Triglyceride-Glo are trademarks of Promega Corporation.

Corning is a registered trademark of Corning, Inc. Dilux is a registered trademark of Dilux, L.L.C. Excel and Microsoft are registered trademarks of Microsoft Corporation. GIBCO is a registered trademark of Life Technologies Corporation. GraphPad Prism is a registered trademark of GraphPad Software, 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.