

User Manual

PathHunter[®] U2OS IL1RL1/IL1RAP Dimerization Bioassay Kit

For Detection of IL-33-Mediated IL-33 Receptor Heterodimerization

For Bioassay Kits:

93-1067Y3-00079: 2-Plate Kit

93-1067Y3-00080: 10-Plate Kit

This page is intentionally left blank.

Table of Contents

Important: Please read this entire user manual before proceeding with the assay.

Overview.....	1
Assay Principle.....	1
Materials Provided.....	2
Storage Conditions.....	2
Additional Materials Required.....	3
Protocol Schematic.....	4
Detailed Protocol.....	5
Day 1: Bioassay Cell Preparation.....	5
Day 1: Sample Preparation.....	6
Day 1: Ligand-Neutralizing Sample Preparation.....	7
Day 1: Detection.....	8
Typical Results.....	9
Troubleshooting Guide.....	10
Document Revision History.....	11
Limited Use License Agreement.....	12
Contact Information.....	12

Table of Figures

Figure 1. Assay Principle.....	1
Figure 2. Agonist Serial Dilutions.....	6
Figure 3. Representative Assay Plate Map.....	8
Figure 4. Typical Results.....	9

For additional information or Technical Support, see contact information at the bottom of this page.

Overview

The PathHunter U2OS IL1RL1/IL1RAP Dimerization Bioassay Kit provides a robust, highly sensitive, and easy-to-use functional cell-based assay to determine drug potency and detect neutralizing antibodies. The bioassay kit contains all the materials needed for a complete assay, including cryopreserved cells, detection reagents, cell plating reagent, positive control agonist, and assay plates. The ready-to-use cryopreserved cells have been manufactured to ensure assay reproducibility and faster implementation from characterization to lot release. This bioassay has been optimized for a 96-well plate format. The protocol can be optimized further for running the assay in a high-throughput 384-well plate format.

Assay Principle

The PathHunter U2OS IL1RL1/IL1RAP Dimerization Bioassay Kit utilizes Enzyme Fragment Complementation (EFC) technology, where the β -galactosidase (β -gal) enzyme is split into two fragments, a small 42 amino acid fragment called ProLink™ (PK), and the larger, N-terminal deletion mutant of β -gal called Enzyme Acceptor (EA). Independently, these fragments have no β -gal activity; however, when forced to complement through protein-protein interactions, they form an active β -gal enzyme.

This bioassay is an application of Eurofins DiscoverX's dimerization platform. The assay is designed to detect IL-33-induced functional dimerization of interleukin 1 receptor-like 1 (IL1RL1) and interleukin-1 receptor accessory protein (IL1RAP). The cells have been engineered to co-express one receptor subunit fused to EA, and a second dimer partner fused to PK. Binding of an agonist, such as IL-33, to one receptor subunit induces it to interact with its dimer partner, forcing complementation of the two enzyme fragments. This results in the formation of a functional enzyme that hydrolyzes a substrate to generate a chemiluminescent signal.

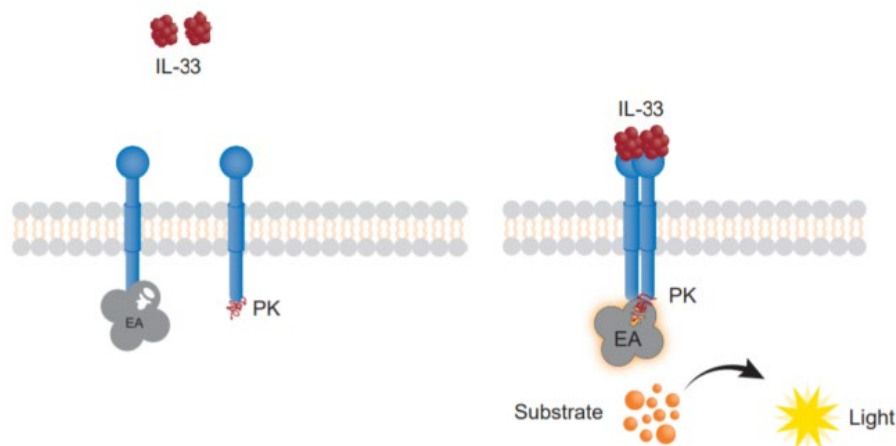


Figure 1. Assay Principle: IL-33-mediated stimulation of one of the tagged receptor units facilitates dimerization with its dimer partner. Dimerization results in EA-PK complementation and forms an active β -gal enzyme that hydrolyzes a substrate to produce a chemiluminescent signal.

Materials Provided

List of Components	93-1067Y3-00079	93-1067Y3-00080
PathHunter U2OS IL1RL1/IL1RAP Dimerization Bioassay Cells (1.2 x 10 ⁶ cells in 1 mL per vial)	2	10
AssayComplete™ Cell Plating 0 Reagent (100 mL per bottle)	1	3
Control Agonist: Recombinant Human IL-33 (10 µg per vial)	1	2
AssayComplete Protein Dilution Buffer (50 mL per bottle)	1	2
PathHunter Bioassay Detection Kit		
Detection Reagent 1 (Bottle)	1 x 3 mL	1 x 15 mL
Detection Reagent 2 (Bottle)	1 x 12 mL	1 x 60 mL
96-Well White, Clear Flat-Bottom, TC-Treated, Sterile Plates with Lid	2	10

Storage Conditions

PathHunter U2OS IL1RL1/IL1RAP Dimerization Bioassay Cells

Cells are shipped on dry ice and should arrive in a frozen state. To ensure maximum cell viability, store the vials of bioassay cells in the vapor phase of liquid nitrogen as soon as possible upon receipt. Please contact technical support immediately, if the cells received were already thawed.

- Short-term (24 hours or less): Store vials at -80°C immediately upon arrival. (DO NOT store at -80°C for longer than 24 hours).
- Long-term (greater than 24 hours): Vials should ONLY be stored in the vapor phase of liquid nitrogen.



Safety Warning: A face shield, gloves and lab coat should be worn at all times when handling frozen vials. Use tongs to remove cryovials from liquid nitrogen storage, and place the vials immediately on dry ice in a covered container. Wait at least 1 minute for any liquid nitrogen that may be present inside the vial to evaporate. Do not touch the bottom of the vials at any time to avoid inadvertent thawing of the cells.

AssayComplete™ Cell Plating 0 Reagent (CP0)

Upon receipt, store at -20°C. Once thawed, the Cell Plating Reagent can be stored at 4°C for up to 4 weeks. For longer storage (up to the expiration date listed on the kit's Certificate of Analysis), the reagent should be aliquoted and stored at -20°C until needed. Avoid multiple freeze-thaw cycles.

To make aliquots suitable for testing one assay plate each, 25 mL of reagent per aliquot can be dispensed and frozen down.

PathHunter® U2OS IL1RL1/IL1RAP Dimerization Bioassay Kit User Manual

Recombinant Human IL-33 (Control Agonist)

Upon receipt, store at -20°C until ready to use (up to the date listed in the kit's Certificate of Analysis). Centrifuge the vial before opening to maximize recovery. Once prepared, the excess stock solution can be stored as suitable aliquots (e.g. 30 µL) at -20°C until needed. Do not freeze/thaw more than twice.

AssayComplete™ Protein Dilution Buffer

Once thawed, the Protein Dilution Buffer can be stored at 4°C for up to 4 weeks. For longer storage (up to the expiration date listed in the kit's Certificate of Analysis), the reagent should be aliquoted and stored at -20°C until needed. Avoid multiple freeze-thaw cycles.

PathHunter Bioassay Detection Kit

Upon receipt, store at -20°C. Once thawed, the detection reagents can be kept at 4°C for up to 4 days. For long-term storage (up to the expiration date listed on the kit's Certificate of Analysis), the reagent should be aliquoted and stored at -20°C until needed. Avoid multiple freeze-thaw cycles.

For the ten-plate kit: If not using all the plates at the same time, we recommend making five aliquots for each of the two detection reagents. Each aliquot will be adequate for two assay plates. Make five aliquots of 3 mL each for Detection Reagent 1, and five aliquots of 12 mL each for Detection Reagent 2. Sufficient volumes for each reagent component have been provided in the kit to make these aliquots.

96-Well Tissue Culture-Treated Plates

Upon receipt, store at room temperature.

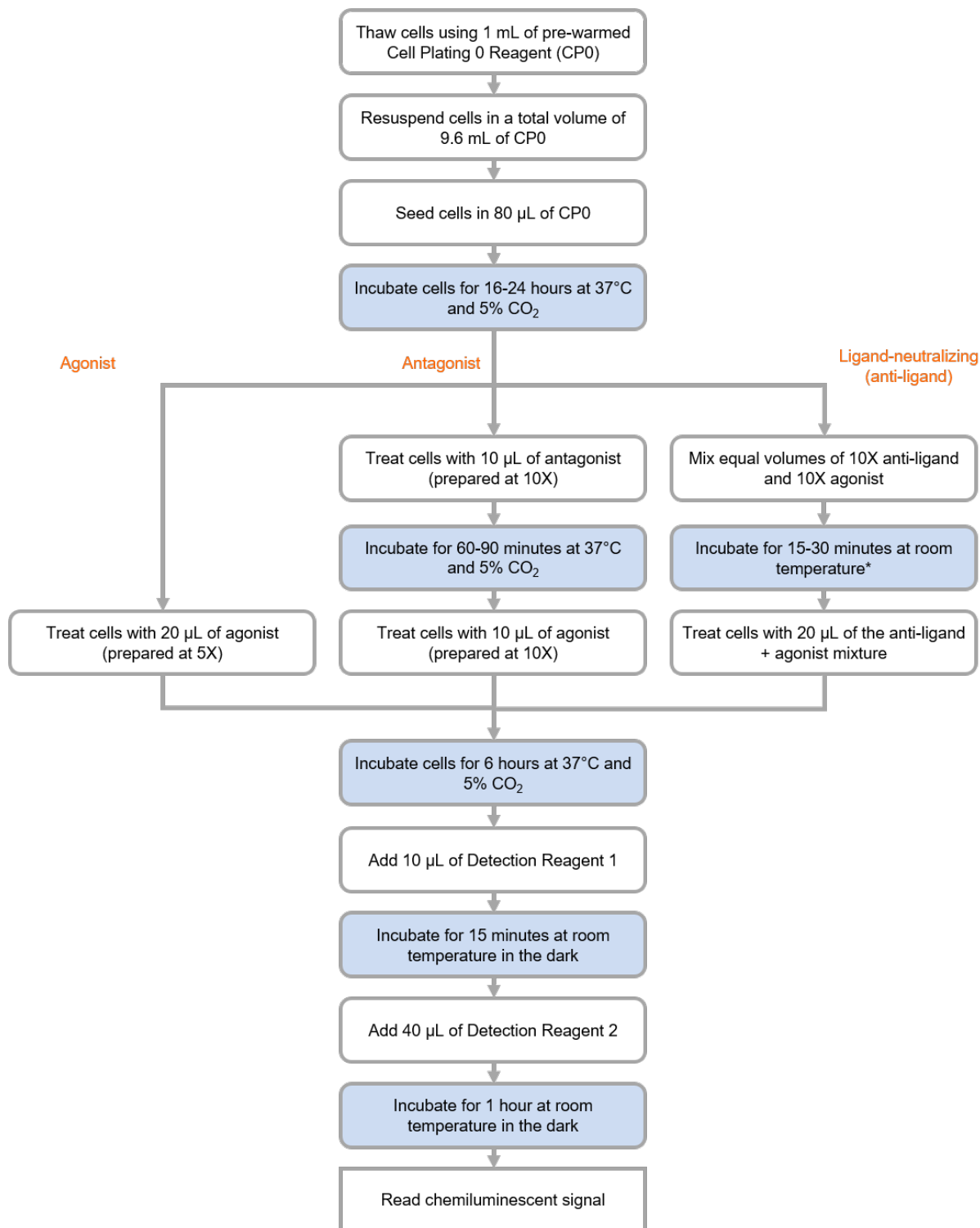
Additional Materials Required

The following equipment and additional materials are required to perform these assays:

Material	Ordering Information
96-Well Green, V-Bottom, Untreated, Non-Sterile Dilution Plates	92-0011
Multimode or luminescence plate reader	Refer to the Instrument Compatibility Chart at discoverx.com/instrument-compatibility
Sterile disposable reagent reservoir	Thermo Fisher Scientific, Cat. No. 8094 or similar
Single and multichannel micropipettes and pipette tips (10 µL-1000 µL)	
Polypropylene tubes (50 mL and 15 mL)	
Microcentrifuge tubes (1.5 mL)	
Tissue culture disposable pipettes (1 mL-25 mL) and tissue culture flasks (T25 and T75 flasks, etc.)	

Protocol Schematic

Quick-start Procedure: In a 96-well plate, perform the following steps:



*Room temperature refers to a range of 23-25°C

Detailed Protocol

This user manual provides a protocol for determining potency in a 96-well format. To run this bioassay in a 384-well format, or to optimize the assay for the detection of neutralizing antibodies, contact Technical Support at DRX_SupportUS@eurofinsUS.com.

Day 1: Bioassay Cell Preparation

The following protocol is for thawing and plating cryopreserved U2OS IL1RL1/IL1RAP bioassay cells from cryovials.

1. Before thawing the cells, ensure that all the materials required are set up in the tissue culture hood. These include:
 - a. One sterile 25 mL reagent reservoir
 - b. One sterile 15 mL conical tube
 - c. A micropipette (P1000) set to dispense 1 mL
 - d. A multichannel pipette and tips set to dispense 80 µL
 - e. A bottle of AssayComplete™ Cell Plating 0 Reagent (CP0), pre-warmed in a 37°C water bath for 15 minutes
 - f. A 96-Well White, Clear Flat-Bottom, Tissue Culture-Treated Sterile Assay Plate (provided with the kit)
 - g. A 96-Well Green, V-Bottom, Untreated, Non-Sterile Dilution Plate, to be used as the master dilution plate
2. Dispense 9.6 mL of CP0 into the 15 mL conical tube.
3. Remove the cryovial from the liquid nitrogen vapor and immediately place it in dry ice.



DO NOT use a heated water bath to thaw the vial. Hold the cryovials by the cap; **DO NOT** touch the sides or bottom of the vial to avoid thawing of the cell pellet. Wipe down the outside surface of the cryovial quickly with 70% ethanol, and immediately bring it into the tissue culture hood.

4. Add 1 mL of pre-warmed CP0 from the 15 mL conical tube into the cryovial, to thaw the cell pellet. The reagent should be added slowly along the wall of the cryovial tube. Slowly pipet the cell suspension up and down, several times, to uniformly resuspend the cells.
5. Transfer the cell suspension to the conical tube containing the remaining 8.6 mL of CP0. Remove any remaining suspension from the cryovial to ensure maximum recovery of all the cells.
6. Replace the cap on the conical tube and gently invert it several times to ensure that the cells are uniformly resuspended in CP0, without creating any froth in the suspension. Immediately pour the suspension into the sterile 25 mL reagent reservoir.
7. Transfer 80 µL of the cell suspension to each well of the 96-well assay plate using a multichannel pipette. Replace the lid and leave the plate at room temperature for 15 minutes, to allow the cells to settle uniformly in the well and minimize potential edge effects.
8. Incubate the assay plate at 37°C and 5% CO₂ for 16-24 hours before proceeding with the assay.

Day 1: Sample Preparation

The following protocol is designed for testing purified biologics. The PathHunter assays can also be run in the presence of high levels of serum or plasma without significantly impacting assay performance. In some assays, standard curves of the control can be prepared in neat serum or plasma, and added directly to cells without further dilution. For best results, the optimized minimum required dilution of crude samples should be empirically determined.

A 1:3 serial dilution for the control agonist, IL-33, has been used in this protocol, as shown in **Figure 2. Agonist Serial Dilutions**. The volumes listed below are designed for running samples from one dose-response curve in triplicate on the assay plate (Refer to **Figure 3: Representative Assay Plate Map**).

1. Add 80 µL of the AssayComplete™ Protein Dilution Buffer (PDB) to Wells A2 to A12 of the master dilution plate.
2. Prepare the control agonist IL-33 dose-response curve:
IL-33 will serve as a positive control in this assay. The agonist is prepared at 5X the desired final concentration as it will be diluted by adding 20 µL to 80 µL of the medium present in the assay plate.
 - 2.1. Add 120 µL of the control agonist prepared at 5X to Well A1 of the master dilution plate. Using a clean pipette tip, transfer 40 µL from Well A1 to Well A2, and mix thoroughly by pipetting up and down several times.
 - 2.2. Replace the pipette tip and transfer 40 µL from Well A2 to A3. Repeat this process until Well A11 is reached, resulting in an 11-point, 1:3 dilution series. No sample is transferred to Well A12 as this is the negative control.
3. Prepare test samples in a similar manner in additional rows of the master dilution plate.
4. Remove the assay plate from the incubator and place it in the tissue culture hood.
5. Transfer 20 µL from each well of the agonist dilution series from the master dilution plate to the appropriate wells of the assay plate.
6. Incubate the assay plate in a humidified incubator at 37°C and 5% CO₂ for 6 hours.

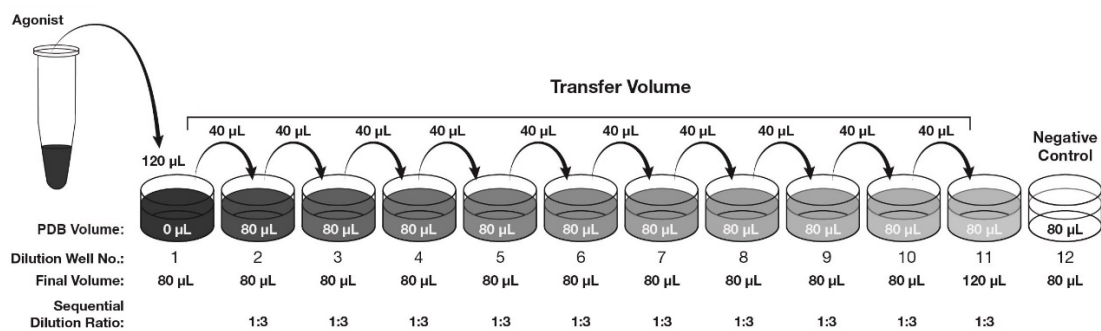


Figure 2. Agonist Serial Dilutions: Make eleven 3-fold serial dilutions of the agonist in a master dilution plate.

7. Prepare the antagonist dilution series using the following protocol. If antagonist samples are not being tested, then skip **Steps 7-12** and proceed to **Day 1: Ligand-Neutralizing Sample Preparation** or **Day 1: Detection**.

- 7.1. The antagonist is prepared at 10X the desired final concentration, in 11-point dilution series, similar to that described for IL-33 in [Steps 1 and 2](#). However, an appropriate top dose and dilution scheme (e.g. a 1:3 dilution scheme) must be determined empirically. No antagonist is added to any wells in Column 12, as these are the negative control wells.
- 7.2. Remove the assay plate from the incubator and place it in the tissue culture hood.
- 7.3. Transfer 10 µL from each well of the antagonist dilution series from the master dilution plate to the appropriate wells of the assay plate using a multichannel pipette.
- 7.4. Incubate the assay plate at 37°C and 5% CO₂ for 60-90 minutes in a humidified incubator. The optimal incubation time for your antagonist must be determined empirically.
8. Prepare the agonist challenge for the antagonist dilution series at 10X the desired final concentration of the agonist (i.e. at 10X EC₈₀ of the agonist).
9. Remove the assay plate from the incubator and place it in the tissue culture hood.
10. Transfer the 10X agonist challenge into a 25 mL reagent reservoir.
11. Transfer 10 µL of the 10X agonist challenge prepared in [Step 8](#), to the appropriate wells (i.e. wells with antagonist dilution series) of the assay plate using a multichannel pipette.
12. Incubate the assay plate at 37°C and 5% CO₂ in a humidified incubator for 6 hours.

Day 1: Ligand-Neutralizing Sample Preparation

The following protocol is used for quantitation of anti-ligand samples (e.g. antibodies or Fc-fusion proteins, etc.). The ligand is typically pre-incubated with the neutralizing sample for 15-30 minutes. However, the optimal incubation time should be determined empirically before addition to the assay.

1. Prepare the anti-ligand antibody dilution series at 10X the desired final concentration in a new row of the master dilution plate, using a similar protocol as described in the [Day 1: Sample Preparation section](#).
2. Prepare the agonist challenge at 10X the desired final concentration, similar to that described in [Step 8 in Sample Preparation](#).
3. For each dose-response curve, add 40 µL of the agonist challenge to Wells 1-11 of the appropriate row in a pre-mixing plate (e.g. a V-bottom polypropylene 96-well plate). Transfer 40 µL of the 10X anti-ligand antibody dilution series to the appropriate row of the pre-mixing plate already containing 40 µL of the agonist challenge. Mix by gently pipetting up and down several times.
4. Add 80 µL of PDB to Column 12 (negative control wells) of the pre-mixing plate.
5. Incubate the plate at room temperature for 15-30 minutes. The optimal pre-incubation time for each antibody should be determined empirically.
6. Transfer 20 µL of the 10X anti-ligand antibody and 10X agonist mixture to the appropriate wells of the assay plate.
7. Incubate the assay plate at 37°C and 5% CO₂ in a humidified incubator for 6 hours.

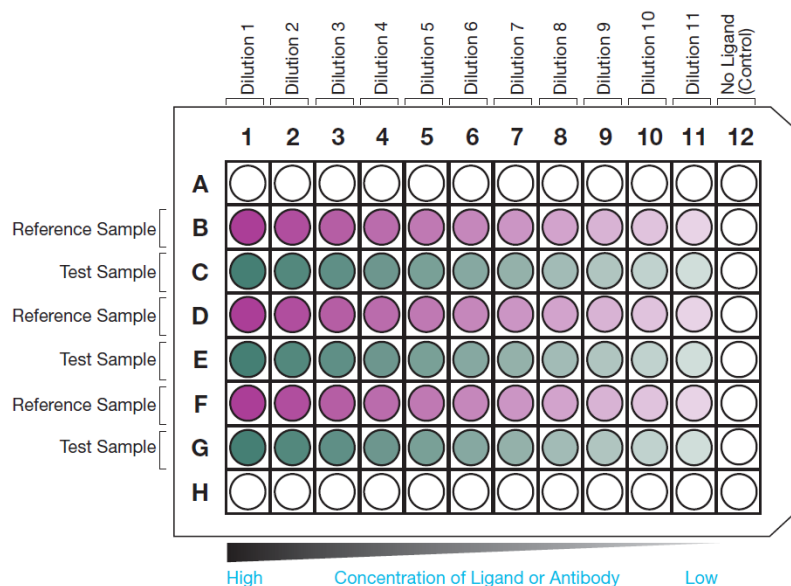


Figure 3. Representative Assay Plate Map: This plate map shows two interdigitated 11-point dose curves, with three replicates per dose point, for a test and reference sample tested using the same dilution scheme.

Day 1: Detection

The following steps for adding detection reagents are applicable when testing purified biologics. For crude biologic samples, gently removing the liquid from all wells of the assay plate and replacing it with 100 μ L of the Cell Plating Reagent before the addition of detection reagents can result in a higher signal. Additional Cell Plating Reagent will be required for this method.

1. Add 10 μ L of the Detection Reagent 1 to each well of the assay plate.
2. Incubate the plate at room temperature (23-25°C) for 15 minutes in the dark.



Room temperature refers to a range of 23-25°C



Detection Reagents are light sensitive, hence incubation in the dark is necessary.

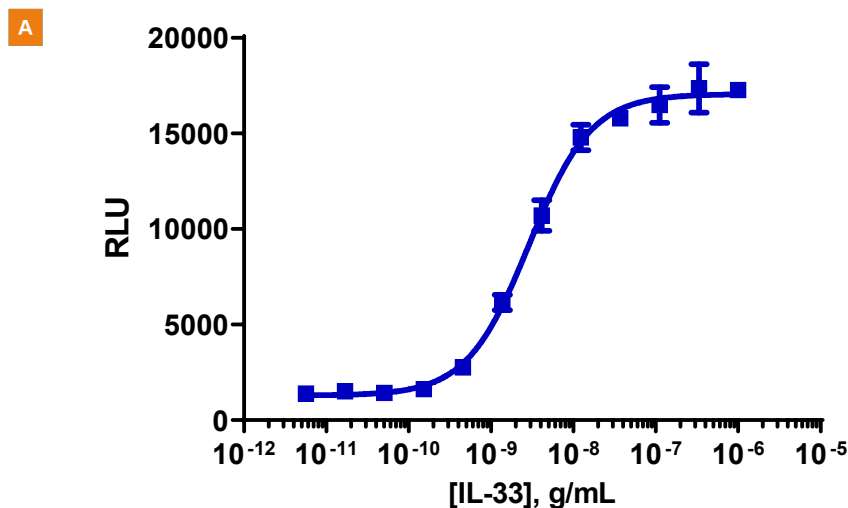
3. Add 40 μ L of the Detection Reagent 2 to each well of the assay plate.
4. Incubate the plate at room temperature for 1 hour in the dark.
5. Read the sample on a standard luminescence plate reader at 0.1 to 1 second/well for photomultiplier tube (PMT) readers or 5-10 seconds for an imager. To determine instrument compatibility, visit discoverx.com/instrument-compatibility.

Data analysis can be performed using any statistical analysis software such as GraphPad Prism, SoftMax Pro, Gen5, or Microsoft Excel.

Typical Results

The following graph is an example of a typical dose-response curve for the IL1RL1/IL1RAP Bioassay generated using the protocol outlined in this manual. The data shows potent, dose-dependent dimerization of the IL1RL1 and IL1RAP receptors when treated with the ligand IL-33.

The plate was read on the EnVision® Multimode Plate Reader and data analysis was conducted using GraphPad Prism.



B

EC ₅₀ (ng/mL)	2.9
S/B	12.4

Figure 4. Typical Results: Representative **A**, dose-response curve and **B**, the EC₅₀ and assay window for IL-33-mediated dimerization of IL1RL1 and IL1RAP receptors, as measured in this bioassay.

Troubleshooting Guide

Problem	Potential Cause	Proposed Solution
No response	Incorrect thawing procedure	Refer to the thawing instructions in the Bioassay Cell Preparation section of this user manual.
	Incorrect ligand used or improper ligand incubation time	Refer to the Certificate of Analysis for recommended ligand and assay conditions.
	Incorrect preparation of the ligand (agonist or antagonist) Sub-optimal time course for induction	Refer to the vendor-specific datasheet to ensure proper handling, dilution, and storage of the ligand. Optimize time course of induction with agonist and antagonist.
Decreased or no response	Cells are not adherent and exhibit incorrect morphology	Confirm adherence of cells using a microscope.
	Problem with microplate reader	The microplate reader should be in luminescence mode. Read at 0.1-1 seconds/well.
Low or no signal	Incorrect preparation of detection reagents	Detection reagents are sensitive to light and should ideally be prepared just prior to use.
	Problem with microplate reader	The microplate reader should be in luminescence mode. Read at 0.1-1 second/well.
Experimental S/B does not match the value noted in the Certificate of Analysis provided	Incorrect incubation temperature	Confirm assay conditions. Check and repeat the assay at the correct incubation temperature, as indicated on the assay datasheet.
	Incorrect preparation of ligand (agonist or antagonist)	Some ligands are difficult to handle. Confirm the final concentration of ligands.
EC ₅₀ is right-shifted	Improper ligand handling or storage	Ensure that the ligands are stored and incubated at the proper temperature.
	Difference in agonist binding affinity	Confirm that the ligand used is comparable to the ligand in the Certificate of Analysis.
	Problems with dynamic range or dilutions	Changing tips during dilution can help in avoiding carryover.
High variability between replicates	Instrument calibration	Ensure that the dispensing equipment is properly calibrated, and proper pipetting technique is used.

For questions on using this product, please contact Technical Support at [1.866.448.4864](tel:1.866.448.4864) or DRX_SupportUS@eurofinsUS.com

Document Revision History

Revision Number	Date Released	Revision Details
0	December 2015	<ul style="list-style-type: none">• New document
1	February 2021	<ul style="list-style-type: none">• Protocol updated to reflect steps for agonist, antagonist, and ligand-neutralization assay formats• Branding updates

Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit discoverx.com/license for limited use label license agreement, and trademark information for this product.

Contact Information

Eurofins DiscoverX

42501 Albrae Street, Fremont, CA 94538

Web: discoverx.com/bioassays

Phone: 1.866.448.4864

Technical Support: DRX_SupportUS@eurofinsUS.com