

# PathHunter® GLP1 (7-37) Bioassay Kit

For chemiluminescent detection of GLP1 (7-37) activity

***Simple Solutions for Complex Biology***



# Table of Contents

Overview .....	1
Technology Principle: PathHunter® GLP1 (7-37) Bioassay .....	1
Materials Provided.....	2
Storage Conditions.....	2
Additional Materials Required .....	3
Protocol Schematic .....	4
Quickstart Guide .....	4
<b>Detailed Protocols</b>	
Cell Thawing Method .....	5
Assay Protocol .....	6
<b>Supplemental Information .....</b>	<b>9-10</b>
Representative Plate Maps for Agonist/Inhibitor Curves .....	9
Instrument Compatibility .....	10
<b>Troubleshooting Guide.....</b>	<b>11</b>
<b>Limited Use License Agreement.....</b>	<b>12</b>

**Read the entire product insert before beginning the assay**

**For additional information or Technical Support, contact [support@discoverx.com](mailto:support@discoverx.com) or visit [www.discoverx.com](http://www.discoverx.com).**

## Overview

### Intended Use

PathHunter® GLP1 (7-37) Bioassay kits provide a robust, highly sensitive and easy-to-use, cell-based functional assay to study GLP1 (7-37) potency and neutralizing antibodies. The Bioassay kits contain all the reagents needed for a complete assay including cells, detection reagents, cell plating reagent, positive control agonist, and assay plates. The pre-validated, frozen cells have been manufactured for single use and are provided in a ready-to-assay format that saves time and adds convenience.

### Technology Principle: PathHunter® GLP1 (7-37) Bioassay

The PathHunter GLP1 (7-37) Bioassay monitors GPCR activity by detecting the interaction of  $\beta$ -Arrestin with the activated GPCR using  $\beta$ -galactosidase ( $\beta$ -gal) enzyme fragment complementation. In this system, the GPCR of interest is fused in frame with the small, 42 amino acid fragment of  $\beta$ -gal called ProLink™ and co-expressed in cells stably expressing a fusion protein of  $\beta$ -Arrestin and the larger, N-terminal deletion mutant of  $\beta$ -gal (called enzyme acceptor or EA). Activation of the GPCR stimulates binding of  $\beta$ -Arrestin to the ProLink-tagged GPCR and forces complementation of the two enzyme fragments of  $\beta$ -galactosidase, resulting in the formation of an active  $\beta$ -gal enzyme. This action leads to an increase in enzyme activity that can be measured using chemiluminescent PathHunter Bioassay Detection Reagents. Because arrestin recruitment occurs independent of G-protein coupling, these assays provide a direct, universal platform for measuring receptor activation, thus providing an ideal tool to characterize functional biologics.

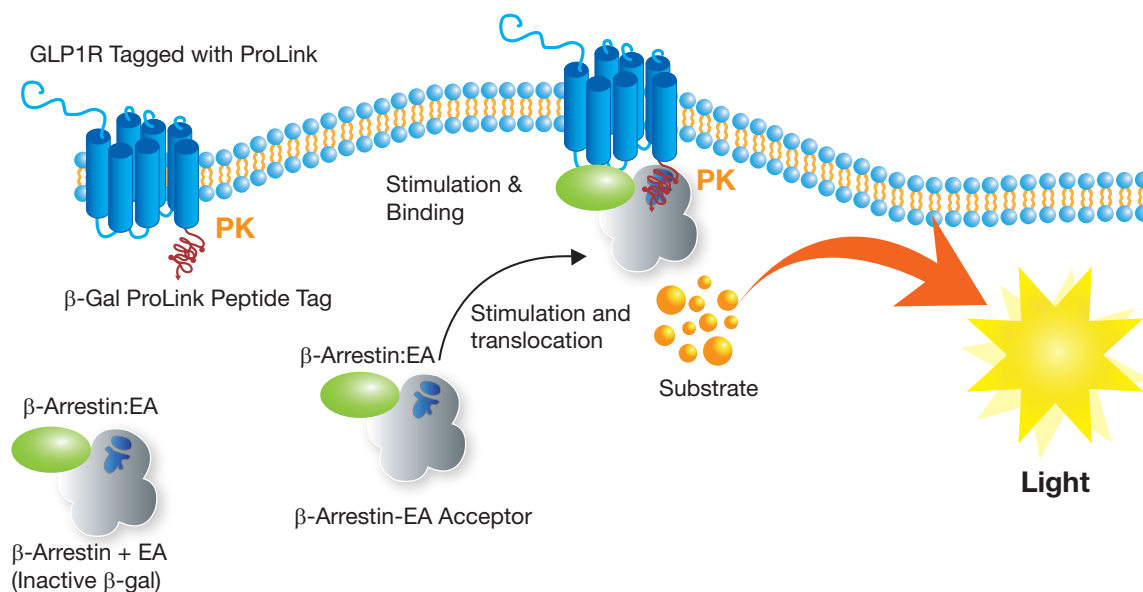


Figure 1. PathHunter GLP1 (7-37) Bioassay Principle

## Materials Provided

List of Components	(93-0300Y2-00027)	(93-0300Y2-00028)
Description Kit Size	Contents	Contents
PathHunter CHO-K1 GLP1R Bioassay Cells	2 vials	10 vials
PathHunter Bioassay Detection Kit	200 dp	1,000 dp
• Detection Reagent 1	2 mL	10 mL
• Detection Reagent 2	8 mL	40 mL
AssayComplete Cell Plating Reagent 0	1 x 100 mL	3 x 100 mL
Protein Dilution Buffer	1 x 50 mL	2 x 50 mL
Control Agonist (Exendin-4)	1 vial	1 vial
96-well Tissue Culture Treated Plates	2 plates	10 plates

## Storage Conditions

### PathHunter CHO-K1 GLP1R Bioassay Cells

Cells are shipped on dry ice and should arrive in a frozen state. To ensure maximum cell viability, thaw the vials as soon as possible upon receipt.

If continued storage of the frozen vials is necessary, store as follows:

- Short term (2 weeks or less): Store vials at  $-80^{\circ}\text{C}$  immediately upon arrival. (DO NOT store at  $-80^{\circ}\text{C}$  for more than 2 weeks).
- Long term (greater than 2 weeks): Vials should ONLY be stored in the vapor phase of liquid nitrogen ( $\text{LN}_2$ ).

**Safety Warning:** A face shield, gloves and lab coat should be worn at all times when handling frozen vials. The manufacturer of the cryovial recommends storing the vials in the vapor phase above the liquid  $\text{N}_2$ . Upon thawing, if  $\text{LN}_2$  is present in the cryovial, it rapidly converts back to its gas phase which can result in the explosion of the vial upon its removal from the liquid nitrogen tank.

### PathHunter Bioassay Detection Kit

Store at  $-20^{\circ}\text{C}$ . Once thawed, the detection reagents can be kept at  $4^{\circ}\text{C}$  for up to 4 days. For longer storage (up to the expiration date listed in the kit certificate of analysis), the reagent should be aliquoted and stored at  $-20^{\circ}\text{C}$  until needed. Avoid multiple freeze-thaw cycles.

To make aliquots suitable for testing one assay plate each, 1mL of Detection Reagent 1 per aliquot should be dispensed and frozen down. 4mL of Detection Reagent 2 per aliquot should be dispensed and frozen down separately. Do not mix the two reagents during aliquoting.

### AssayComplete Cell Plating Reagent 0

Once thawed, the Cell Plating Reagent can be stored at  $4^{\circ}\text{C}$  for up to 4 weeks. For longer storage (up to the expiration date listed in the kit certificate of analysis), the reagent should be aliquoted and stored at  $-20^{\circ}\text{C}$  until needed. Avoid multiple freeze-thaw cycles. To make aliquots suitable for testing one assay plate each, 30mL of reagent per aliquot should be dispensed and frozen down.

### 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 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, 10mL of reagent per aliquot should be dispensed and frozen down. This amount may vary depending on your stock sample concentrations and should be adjusted accordingly.

### Recombinant Human Exendin-4 Control Agonist

Store at -20°C until ready to use (up to the expiration date listed in the kit certificate of analysis). Centrifuge the vial prior to opening to maximize recovery.

### 96-well Tissue Culture Treated Plates

Store at room temperature.

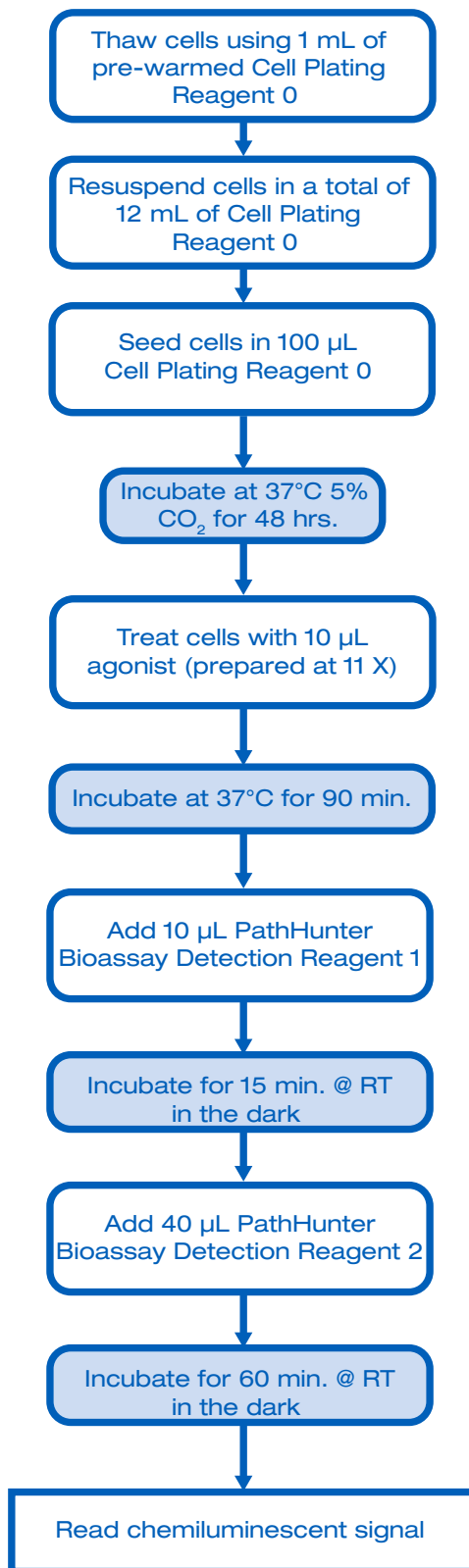
## Additional Materials Required

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

Equipment
Single and multichannel micro-pipettors and pipette tips
Multimode or luminescence plate reader
V-Bottom 96-well compound dilution plates (DiscoverX 92-0011 or similar)
Disposable Reagent Reservoir (Thermo Scientific, Cat. #8094 or similar)

## Protocol Schematic

**Quick-Start Procedure:** In a 96-well tissue culture treated plate perform the following:



## Detailed Protocols

### Day 1: PathHunter CHO-K1 GLP1R Bioassay cell suspension preparation: \_\_\_\_\_

The following protocol is for thawing and plating frozen PathHunter CHO-K1 GLP1R Bioassay cells from cryovials.

1. Before thawing the cells, ensure that all the necessary materials required are set up in the tissue culture hood. This includes:
  - a. One 100 mL reagent reservoir
  - b. One 15 mL conical tube
  - c. A micropipettor (P1000) set to dispense 1 mL.
  - d. A multichannel pipette and tips set to dispense 100  $\mu$ L.
  - e. A bottle of Cell Plating Reagent 0 (CP0, pre-warmed in a 37°C water bath for 15min.)
  - f. A white-walled, clear-bottom 96-well assay plate.
2. Dispense 12 mL of cell plating reagent into the 15 mL conical tube
3. Remove the cryovial from liquid nitrogen and immediately place in dry ice.

**DO NOT** use heated water bath to thaw the vial. Wipe down the cryovial quickly with 70% EtOH, and bring it into the tissue culture hood right away. **DO NOT** touch the sides or bottom of the vial to avoid thawing of the cell pellet through body heat.

4. Immediately add 1 mL (using P1000) of pre-warmed Cell Plating Reagent from the 15 mL conical tube to the cryovial, thawing the cell pellet. The medium should be added slowly along the side of the wall of cryovial tube. Mix the cells gently by pipetting up and down several times to break up any clumps. Transfer the cell suspension to the conical tube containing the remaining 11 mL of Cell plating reagent. Remove any medium/suspension left in the tube to ensure complete recovery of all the cells from the vial.
5. Mix the tube by inversion to ensure the cells are properly mixed in the medium without creating any froth in the suspension and immediately pour the suspension into the 100 mL reservoir.
6. Add 100  $\mu$ L of cells (10,000 cells/well) to each well of the 96 well assay plate using the multichannel pipette.
7. Let the assay plate sit for 15 min. at room temperature, then gently place the assay plate in a tissue culture incubator set to 37°C, 5% CO<sub>2</sub>. Allow the cells to recover for 48 hours before continuing.

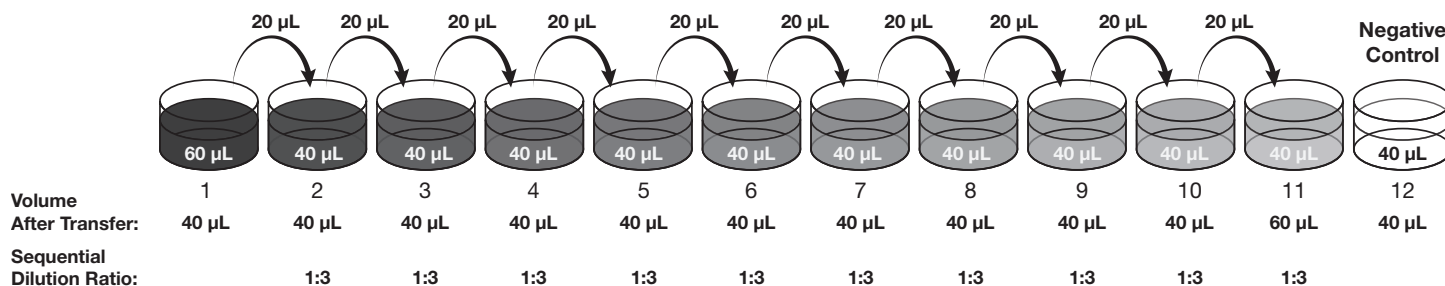


### Day 3: Ligand Preparation

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. Therefore, standard curves of control can typically be prepared in neat serum or plasma and added directly to cells without further dilution. For the best results, the optimized minimum required dilution (MRD) of crude samples should be empirically determined.

#### Method 1

A 1:3 serial dilution. This method is simple to run and covers a broad concentration range, making it ideal for initial assay optimization. The volumes listed below are designed for running samples from one dose response curve in duplicate on the assay plate (as pictured in the assay plate map on page 11).



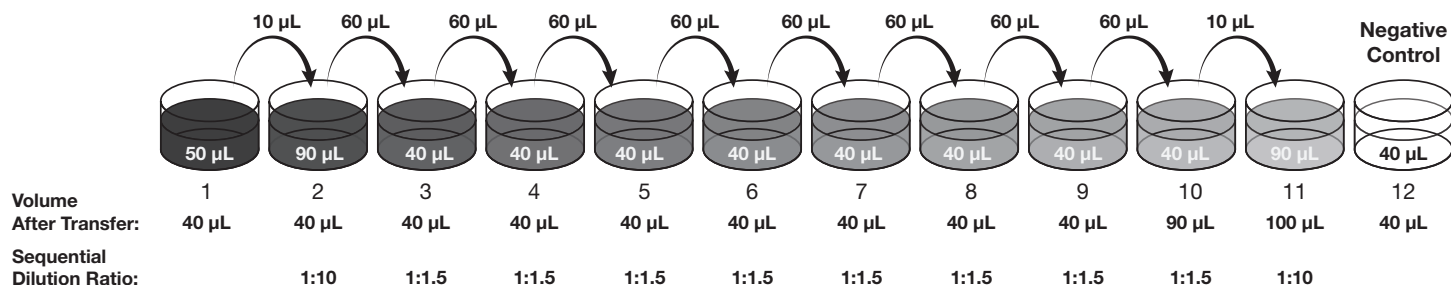
- Prepare the reference agonist (DiscoverX Exendin-4) dose response curve, which will serve as a positive control in this assay. Agonist is prepared at 11x the desired final concentration as it will be diluted by adding to the 100 µL of media present in the assay plate.
  - Add 40 µL of Protein Dilution Buffer (PDB) to columns 2-12, row A of a master dilution plate (e.g. a V-bottom polypropylene 96-well dilution plate, DiscoverX 92-0011 or similar).
  - Add 500 µL of PDB to the exendin-4 vial containing 0.5 mg of lyophilized powder to make a 1 mg/mL stock solution.
  - Add 57.4 µL of PDB to column one, row A of the master dilution plate. Add 2.6 µL of the 1 mg/mL exendin-4 stock to this well. Mix thoroughly by pipetting up and down several times. This results in an 44 µg/mL solution (11x the final 4 µg/mL curve top).
  - Using a clean tip, transfer 20 µL from column 1 into column 2 and mix by pipetting up and down several times. Replace the pipette tip, and transfer 20 µL from column 2 into column 3. Mix by pipetting up and down several times. Repeat this process until column 11 is reached, resulting in an eleven point, 1:3 dilution series. No ligand is transferred to column 12 as this will serve as a negative control.
- Prepare GLP1 (7-37) reference curve.
  - Add 40 µL of Protein Dilution Buffer (PDB) to columns 2-12 in a new row of the master dilution plate.
  - Add 60 µL of GLP1 (7-37) prepared at 11x the desired final concentration to column 1 of this row on the master dilution plate.
  - Using a clean tip, transfer 20 µL from column 1 into column 2 and mix by pipetting up and down several times. Replace the pipette tip, and transfer 20 µL from column 2 into column 3. Mix thoroughly by pipetting up and down several times. Repeat this process until column 11 is reached, resulting in an eleven point, 1:3 dilution series. Do not transfer ligand into column 12 as this will serve as a negative control.

3. Prepare biosimilar curves. Biosimilars are prepared at 11x the desired final concentration.
  - a. Add 40  $\mu\text{L}$  of Protein Dilution Buffer (PDB) to columns 2-12 in a new row of the master dilution plate.
  - b. Add 60  $\mu\text{L}$  of the highest concentration of biosimilar to column 1 of this row on the master dilution plate.
  - c. Using a clean tip, transfer 20  $\mu\text{L}$  from column 1 into column 2 and mix by pipetting up and down several times. Replace the pipette tip, and transfer 20  $\mu\text{L}$  from column 2 into column 3. Mix by pipetting up and down several times. Repeat this process until column 11 is reached, resulting in an eleven point, 1:3 dilution series. Do not transfer ligand into column 12 as this will serve as a negative control.
4. Remove the assay plate from the 37°C, 5% CO<sub>2</sub> incubator and bring into the tissue culture hood.
5. Add 10  $\mu\text{L}$  from each well of the master dilution plate to the appropriate wells of the assay plate. Refer to the example plate map on page 11.
6. Return the assay plate to the 37°C, 5% CO<sub>2</sub> incubator and incubate for 90 minutes before proceeding.

## Method 2

This method covers a smaller concentration range in order to generate more data points within the linear range of the dose response curve, and is suitable for testing ligands with known pharmacology in this assay platform. The volumes listed below are designed for running samples from one dose response curve in duplicate on the assay plate (as pictured in the assay plate map on page 11).

1. Prepare the reference agonist (DiscoverRx Exendin-4) dose response curve, which will serve as a positive control in this assay. Agonist is prepared at 11x the desired final concentration as it will be diluted by adding to the 100  $\mu\text{L}$  of media present in the assay plate.
  - a. Add 90  $\mu\text{L}$  of Protein Dilution Buffer (PDB) to columns 2 and 11, row A of a master dilution plate (e.g. a V-bottom 96-well dilution plate, DiscoverRx 92-0011 or similar). Then add 40  $\mu\text{L}$  of Protein Dilution Buffer (PDB) to the remaining wells in the same row.
  - b. Add 500  $\mu\text{L}$  of PDB to the exendin-4 vial containing 0.5 mg of lyophilized powder to make a 1 mg/mL stock solution.
  - c. Add 47.8  $\mu\text{L}$  of PDB to column one, row A of the master dilution plate. Add 2.2  $\mu\text{L}$  of the 1 mg/mL stock to this well. Mix thoroughly by pipetting up and down several times. This results in an 44  $\mu\text{g}/\text{mL}$  solution (11x the final 4  $\mu\text{g}/\text{mL}$  curve top).
  - d. Using a clean tip, transfer 10  $\mu\text{L}$  from column 1 into column 2 and mix by pipetting up and down several times.
  - e. Replace the pipette tip, and transfer 60  $\mu\text{L}$  from column 2 into column 3. Mix by pipetting up and down several times. Repeat this process until column 10 is reached.



- f. Using a clean tip, transfer 10 µL from column 10 into column 11 and mix by pipetting up and down several times.
2. Prepare GLP1 (7-37) reference curve.
    - a. Add 90 µL of Protein Dilution Buffer (PDB) to columns 2 and 11 in a new row on the master dilution plate. Then add 40 µL of Protein Dilution Buffer (PDB) to the remaining wells in the same row.
    - b. Prepare 50 µL of GLP1 (7-37) at 11x the desired final concentration and add to column 1 of this row on the master dilution plate.
    - c. Using a clean tip, transfer 10 µL from column 1 into column 2 and mix by pipetting up and down several times.
    - d. Replace the pipette tip, and transfer 60 µL from column 2 into column 3. Mix by pipetting up and down several times. Repeat this process until column 10 is reached.
    - e. Using a clean tip, transfer 10 µL from column 10 into column 11 and mix by pipetting up and down several times.
  3. Prepare biosimilar curves. Biosimilars are prepared at 11x the desired final concentration.
    - a. Add 90 µL of Protein Dilution Buffer (PDB) to columns 2 and 11 in a new row on the master dilution plate. Then add 40 µL of Protein Dilution Buffer (PDB) to the remaining wells in the same row.
    - b. Add 45 µL of the highest concentration of biosimilar to column 1 of a new row on the master dilution plate.
    - c. Using a clean tip, transfer 10 µL from column 1 into column 2 and mix by pipetting up and down several times.
    - d. Replace the pipette tip, and transfer 60 µL from column 2 into column 3. Mix by pipetting up and down several times. Repeat this process until column 10 is reached.
    - e. Using a clean tip, transfer 10 µL from column 10 into column 11. Mix by pipetting up and down several times.
  4. Remove the assay plate from the 37°C, 5% CO<sub>2</sub> incubator and bring it into the tissue culture hood.
  5. Add 10 µL from each well of the master dilution plate to the appropriate wells of the assay plate. Refer to the example plate map on page 11.
  6. Return the assay plate to the 37°C, 5% CO<sub>2</sub> incubator and incubate for 90 minutes before proceeding.

### Day 3: Detection

---

1. Using a multichannel pipette, add 10 µL of Detection Reagent 1 to each well of the assay plate.
2. Incubate the plate at room temperature for 15 minutes in the dark.
3. Using a multichannel pipette add 40 µL of 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 sample on a standard luminescence plate reader at 0.1 to 1 second/well for PMT readers or 5-10 seconds for imager.

**Note:** For crude biologic samples, gently removing the liquid from all wells and replacing with 100 µL of Cell Plating Reagent before the addition of the detection reagents can result in higher signal. Additional Cell Plating Reagent is necessary for this method.

Representative Plate Maps for Agonist/Inhibitor Curve

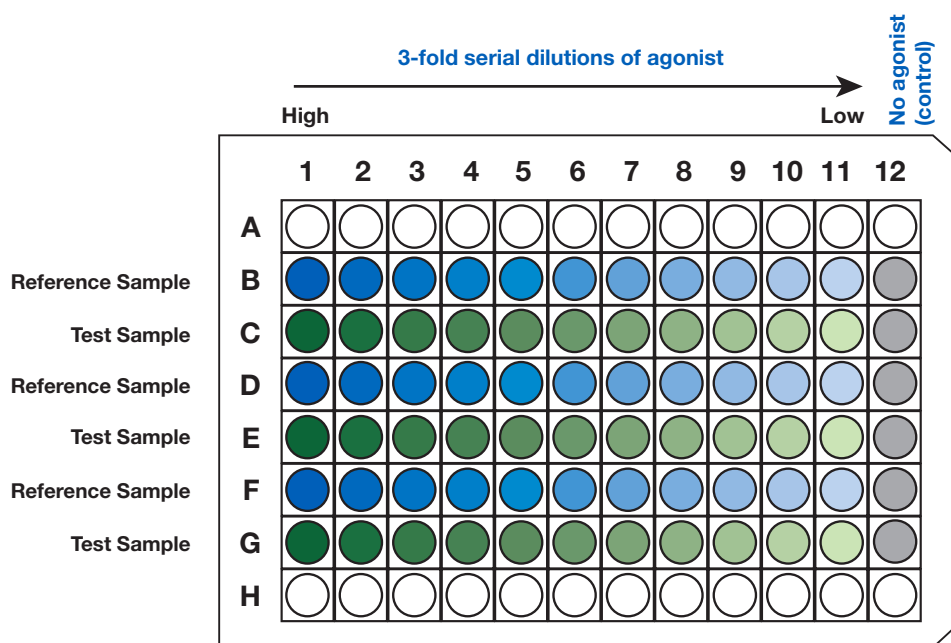


Figure 1. Serial dilution method 1. This plate map shows 11-point dose curves with three data points at each concentration for one reference and one test sample per plate.

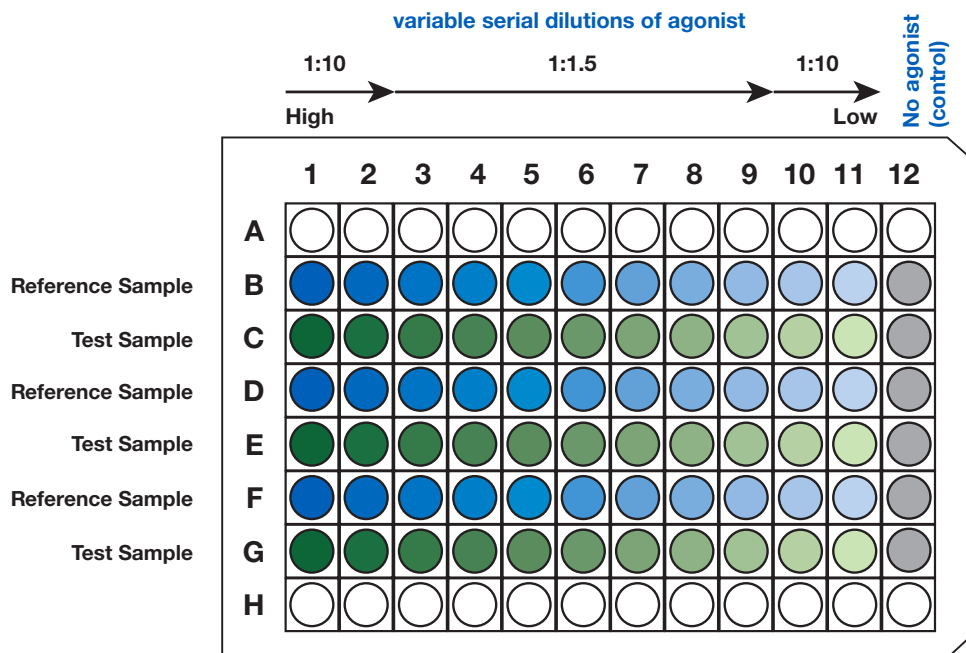


Figure 2. Serial dilution method 2. This plate map shows a 11 point dose curve with 3 data points at each concentration for one reference and one test sample per plate, with a variable serial dilution scheme.

## Related Products

Description	Ordering Information
PathHunter Bioassay Detection Reagents	<a href="http://www.discoverx.com/detectionreagents">www.discoverx.com/detectionreagents</a>
Cell Culture Kits, Reagents & Consumables	<a href="http://www.discoverx.com/cell-culture-kits-reagents-consumables">www.discoverx.com/cell-culture-kits-reagents-consumables</a>
AssayComplete Cell Plating Reagents	<a href="http://www.discoverx.com/cellplatingreagents">www.discoverx.com/cellplatingreagents</a>
Control Ligands	<a href="http://www.discoverx.com/controlligands">www.discoverx.com/controlligands</a>

## Instrument Compatibility Chart

Assay	Instrument	Read-Out
All PathHunter® assays	<b>COMPATIBLE WITH ANY LUMINOMETER</b> <b>BMG:</b> PheraStar, Cytostar, LumiStar	Luminescence
HitHunter® cAMP HitHunter® cGMP	<b>Perkin Elmer:</b> TopCount, Victor 2 or V, Fusion, LumiCount, Envision, Micro-beta (Trilux), Viewlux, Northstar  <b>GE:</b> LEAD seeker, Farcyte  <b>Molecular Devices:</b> FLIPR, SpectraMax M3/M4/M5/M5e, FlexStation 3, SpectraMax L  <b>Tecan:</b> Ultra, Evolution  <b>Turner BioSystems:</b> Modulus Microplate  <b>Caliper</b> LabChip 3000 & EZ Reader  <b>Berthold Technologies:</b> Mithras LB940, CentroLIApc  <b>Hamamatsu</b> FDS6000, FDSS/RayCatcher  <b>Thermo Scientific:</b> Luminoskan Ascent  <b>Biotek:</b> Synergy 2	

\*For other instruments not listed here, please contact technical support at [support@discoverx.com](mailto:support@discoverx.com) to ensure compatibility.

## Troubleshooting Guide

Problem	Cause	Solution
<b>No Response</b>	Improper thawing procedure	Refer to thawing instructions on page 4 of this user manual.
	Improper ligand used or improper ligand incubation time	See datasheet for recommended ligand and assay conditions
	Improper preparation of ligand (agonist or antagonist)	Refer to vendor specific datasheet to ensure proper handling, dilution and storage of ligand
	Improper time course for induction	Optimize time course of induction with agonist and antagonist
<b>Decreased Response</b>	Cells are not adherent and exhibit incorrect morphology	Confirm adherence of cells using microscope
<b>Low or No Signal</b>	Improper preparation of detection reagents	Detection reagents should ideally be prepared just prior to use and are sensitive to light
	Problem with microplate reader	Microplate reader should be in luminescence mode. Read at 0.1-1 sec/well.
<b>Experimental S:B does not match datasheet value</b>	Incorrect incubation temperature	Confirm assay conditions Check and repeat assay at correct incubation temperature as indicated on the assay datasheet
	Improper preparation of ligand (agonist or antagonist)	Some ligands are difficult to handle. Confirm the final concentration of ligands
<b>EC<sub>50</sub> is right-shifted</b>	Improper ligand handling or storage	Check ligand handling requirements
	Difference in agonist binding affinity	Confirm that the ligand used is comparable to the ligand in the Product Insert
	Problems with plate type and compound stability	Hydrophobic compounds should be tested for solubility and may be diluted in Protein Dilution Buffer
		Non-binding surface plates may be necessary for hydrophobic compounds

For additional information or technical support, please contact technical support at [support@discoverx.com](mailto:support@discoverx.com).

## Limited Use License Agreement

**A.** This product and/or its use is covered by U.S. patents #6,342,345 B1, #7,135,325, B2, #8,101,373 B2 and/or related foreign patents and pending applications and trade secrets that are either owned by or licensed to DiscoverX corporation. The cells and detection reagents (collectively "Materials") purchased from DiscoverX are expressly restricted in their use. DiscoverX has developed a Cell-Based assay ("Assay") that employs genetically modified cells and vectors (collectively, the "Cells"), and related detection reagents (the "Reagents") (collectively referred to as "Materials"). By purchasing and using the Materials, the Purchaser agrees to comply with the following terms and conditions of this label license and recognizes and agrees to such restrictions:

1. Purchaser will not further propagate the Cells included in the Assay kit.
2. Purchaser is permitted to use the Cells only for use in the Assay and in connection with Reagents purchased from DiscoverX Corporation or its authorized distributor.
3. The Materials are not transferable and will be used only at the site for which they were purchased. Transfer to another site owned by Purchaser will be permitted only upon written request by Purchaser followed by subsequent written approval by DiscoverX.
4. The Reagents contain or are based upon the proprietary and valuable know-how developed by DiscoverX, and the Reagents have been optimized by DiscoverX to function more effectively with the Cells in performing the Assay. Purchaser will not analyze or reverse engineer the Materials nor have them analyzed on Purchaser's behalf.
5. In performing the Assay, Purchaser will use only Reagents supplied by DiscoverX or an authorized DiscoverX distributor for the Materials.
6. Purchaser will not use the Cells with any other reagents or substrates, other than the Reagents that are provided by or purchased from DiscoverX or an authorized DiscoverX distributor, in connection with the Materials.
7. The number of Assays performed will not exceed the authorized number for which Materials were purchased.

**B.** The use of this cell line may require third party licenses. Purchaser shall have the sole responsibility for obtaining any such licenses prior to use of the cell lines.

If the purchaser has any further questions regarding the rights conferred with purchase of the Materials, please contact:

Licensing Department  
DiscoverX Corporation  
42501 Albrae Street  
Fremont, CA 94538 USA  
t | 1.510.979.1415 x 104  
e | info@discoverx.com