



DiscoverX

User Manual

PathHunter[®] Human Growth Hormone
Bioassay Kit

For Chemiluminescent Detection of hGH Activity

Table of Contents

Overview	1
Technology Principle	1
Materials Provided	2
Storage Conditions.....	2
Additional Materials Required	3
Protocol Schematic.....	4
Detailed Protocols.....	5
Day 1: PathHunter Bioassay Cell Preparation	5
Day 1: Sample Preparation	6
Day 2: Detection	6
Representative Plate Map.....	7
Typical Results.....	7
Troubleshooting Guide.....	8
Limited Use License Agreement.....	8



Please read this entire user manual before proceeding with the assay.
For additional information or Technical Support, see contact information below.

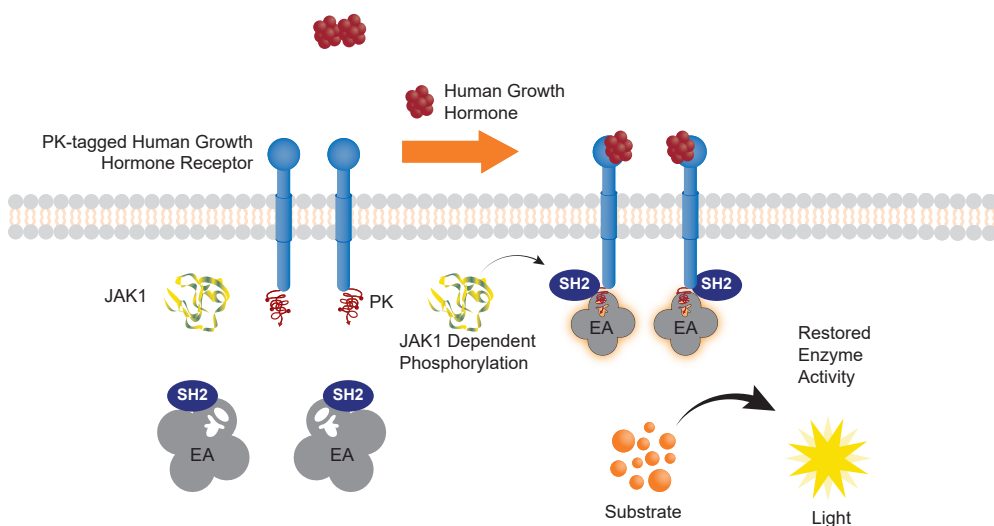
Overview

PathHunter® Human Growth Hormone (hGH) Bioassay kits provide a robust, highly sensitive and easy-to-use, cell-based functional assay to study hGH 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, dilution buffer, and assay plates. The qualified, 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

These assays utilize Enzyme Fragment Complementation (EFC) technology, where the β -galactosidase (β -gal) enzyme is split into two fragments, ProLink™ (PK) and 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.

In the PathHunter assay approach for cytokine receptors, the ProLink tag is fused to the C-terminus of the receptor. The EA is fused to a phosphotyrosine SH2-domain containing protein that is able to bind the activated cytokine. Ligand-induced activation of the receptor results in receptor phosphorylation by a cytosolic tyrosine kinase such as JAK1. The SH2-EA fusion protein binds the phosphorylated receptor, forcing complementation of PK and EA to form an active β -gal enzyme. β -gal enzymatic activity is quantitatively measured using a chemiluminescent substrate in the PathHunter Bioassay Detection Kit (Figure 1).



Materials Provided

List of Components	93-0756Y3-00023	93-0756Y3-00024
PathHunter U2OS GHR-JAK1 Bioassay Cells	2 vials	10 vials
PathHunter Bioassay Detection Kit		
Detection Reagent 1 (mL)	2	10
Detection Reagent 2 (mL)	8	40
AssayComplete™ Cell Plating Reagent 4	1 X 100 mL	3 X 100 mL
Assay Complete Dilution Buffer B5	1 X 100 mL	1 X 100 mL
Control Agonist (hGH)	1 vial	1 vial
96-well Clear-Bottom TC Treated, Sterile Plates w/lid	2 plates	10 plates

*Excess reagent may be left in the bottle, after using the volumes needed to run all the plates

Storage Conditions

PathHunter U2OS GHR-JAK1 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 (24 hours or less): Store vials at -80°C immediately upon arrival. (DO NOT store at -80°C for more 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. The manufacturer of the cryovial recommends storing the vials in the vapor phase above the liquid nitrogen. Upon thawing, if liquid nitrogen 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°C . Once thawed, the detection reagents can be kept at 4°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°C until needed. Avoid multiple freeze-thaws.

For the ten-plate kit, if all the plates will not be used 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 2.3 mL each for Detection Reagent 1, and five aliquots of 9.2 mL each for Detection Reagent 2. Sufficient reagent volumes are provided in the kit to make these aliquots.

If reagents will be used for a single plate, then the remaining Detection Reagents can be frozen. The detection reagents can be thawed and frozen for a total of three times without loss in performance.

AssayComplete Cell Plating Reagent 4 (CP4)

Once thawed, the CP4 can be stored at 4°C for up to 4 weeks. For longer storage (up to the expiration date listed on 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, 30 mL of reagent per aliquot can be dispensed and frozen down.

Assay Complete Dilution Buffer 5 (DB-B5)

Once thawed, the Dilution Buffer can be stored at 4°C for up to 4 weeks. For longer storage (up to the expiration date listed on 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, 10 mL of the Dilution Buffer per aliquot should be dispensed and frozen down. This amount may vary depending on your stock sample concentrations and should be adjusted accordingly.

PathHunter® Human Growth Hormone Bioassay Kit User Manual

Recombinant Human Growth Hormone 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. When ready to use, reconstitute to a concentration of 1 mg/mL by adding 100 µL of sterile deionized water to the vial containing 100 µg of lyophilized powder. Reconstituted ligand is stable for 1 week at 2-8°C. For longer storage (up to the expiration date listed on the kit Certificate of Analysis), it is recommended to store in working aliquots at -20 to -80°C.

96-Well Tissue Culture Treated Plates

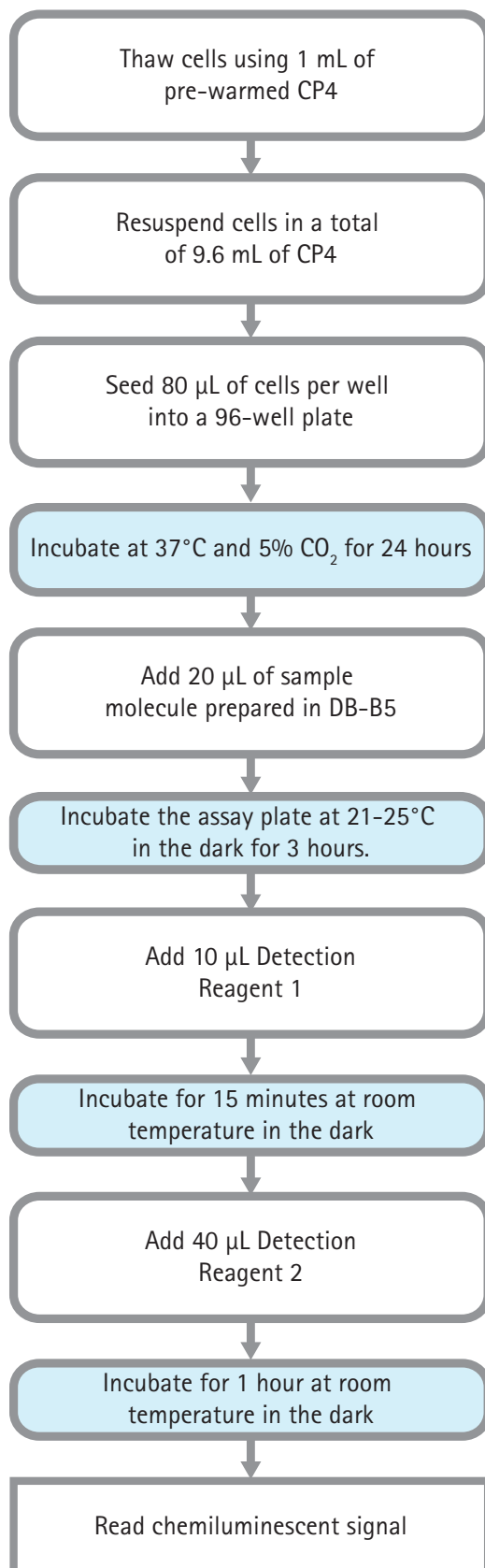
Store at room temperature.

Additional Materials Required

Material	Ordering Information
V-Bottom 96-well ligand dilution plates	92-0011
Multimode or luminescence plate reader	Refer to Instrument Compatibility Chart at discoverx.com/instrument-compatibility
Single and multichannel micro-pipettors and pipette tips	
Disposable reagent reservoir	Thermo Fisher Scientific, Cat. No. 8094 or similar

Protocol Schematic

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



Detailed Protocol

Day 1: PathHunter Bioassay Cell Preparation

The following protocol is for thawing and plating frozen PathHunter U2OS GHR-JAK1 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 25 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 80 μ L.
 - e. A bottle of Cell Plating Reagent 4 (CP4, pre-warmed in a 37°C water bath for 15 minutes).
 - f. A white-walled, Clear-bottom 96-well assay plate (DiscoverX Cat. No. 92-0027 or similar).
2. Dispense 9.6 mL of CP4 into the 15 mL conical tube.
3. Remove the cryovial from liquid nitrogen and immediately place in dry ice.

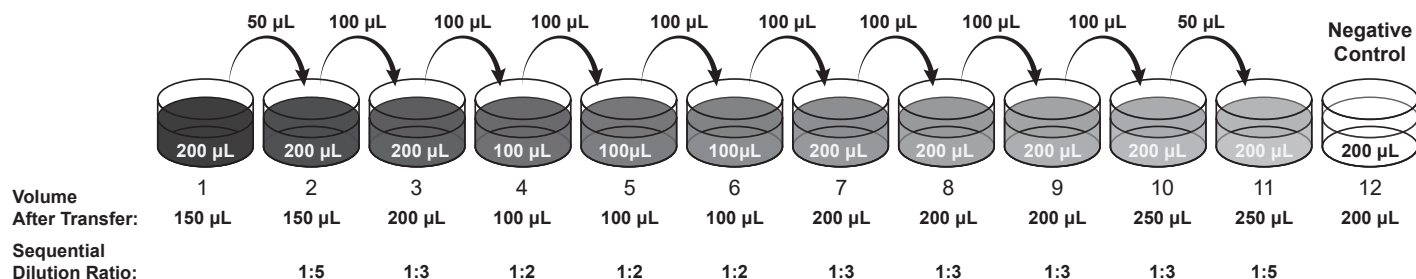


DO NOT use heated water bath to thaw the vial. Wipe down the cryovial quickly with 70% EtOH, and bring it into the tissue culture hood right away. DO NOT touch the sides or bottom of the vial to avoid thawing of the cell pellet through body heat.

4. Thaw the pellet by immediately adding 1 mL (using P1000) of pre-warmed CP4 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.
5. Transfer the cell suspension to the conical tube containing the remaining 8.6 mL of CP4. Remove any medium/ suspension left in the tube to ensure complete recovery of all the cells from the vial.
6. Mix the tube by gentle 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 25 mL reservoir.
7. Add 80 μ L of cells to each well of the 96-well assay plate using the multichannel pipette. Let the plate sit for 15 minutes at room temperature to allow the cells to settle and reduce potential edge effects.
8. Place the plate in a tissue culture incubator (37°C, 5% CO₂) for 24 hours before adding ligand.

Day 1: Sample Preparation

The following protocol is designed for testing purified biologics. 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.

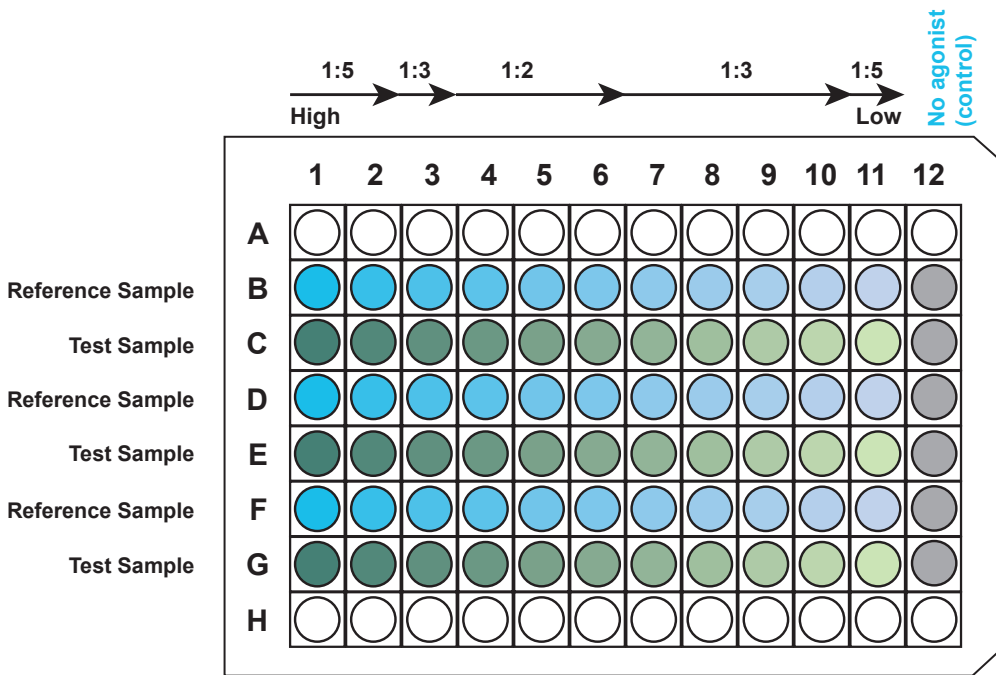


1. Prepare the agonist human growth hormone (hGH) dose response curve. Agonist is prepared at 5X the desired final concentration. Top dose for hGH is 10 µg/mL at final assay concentration.
 - a. Add 200 µL of DB-B5 to wells A2, A3 and A7 to A12 of a master dilution plate. Add 100 µL of DB-B5 to wells A4 to A6.
 - b. Add 100 µL of reconstitution solvent (deionized water) to 100 µg of hGH lyophilized powder to make a stock concentration of 1 mg/mL.
 - c. Add 10 µL of hGH from the stock (1 mg/mL) in 190 µL of DB-B5 in well A1. This creates a working stock of 50 µg/mL of hGH (5X the final concentration).
 - d. Using a clean tip, transfer 50 µL from well A1 into well A2 for a 1:5 dilution, and mix by pipetting up and down several times. Replace the pipette tip, and transfer 100 µL from well A2 into well A3 for a 1:3 dilution. Mix by pipetting up and down several times.
 - e. Transfer 100 µL from A3 to A4 for a 1:2 dilution, and mix by pipetting up and down. Repeat the process till well A6 is reached. Continue transferring 100 µL from A6 to A7 for a 1:3 dilution and repeat process till well A10 is reached. Finally transfer 50 µL from A10 to A11 for a 1:5 dilution. No ligand is transferred to A12 as this will serve as a negative control.
2. Add 20 µL from the hGH curve on the master dilution plate to the appropriate wells of the assay plate.
3. Incubate the assay plate at 21–25°C in the dark for 3 hours

Day 2: Detection

1. Using a multichannel pipette, add 10 µL of Bioassay 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 Bioassay 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 or photomultiplier tube (PMT) readers or 5–10 seconds for imager.

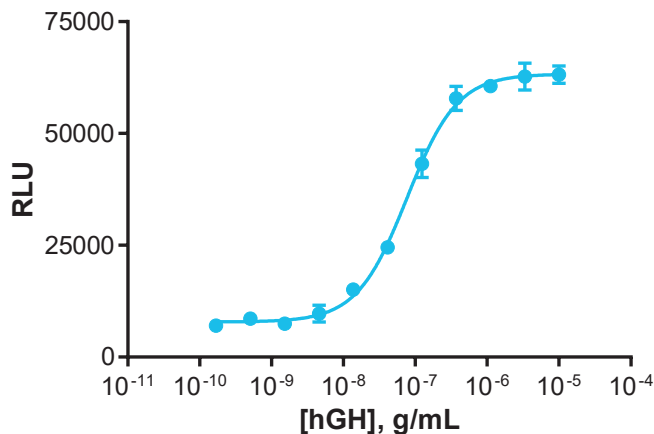
Representative Plate Map



This plate map shows 11-point dose curves with 3 data points at each concentration for one reference and one test sample per plate.

Typical Results

Following graph is a representative dose-response curve with hGH. Using the protocol outlined in this user manual, the data show a potent, dose-dependent phosphorylation of the receptor by the cytosolic tyrosine kinase JAK1 in U2OS cells when treated with a dose-response of hGH.



EC ₅₀ (ng/mL)	76.7
S/B	8.0

Troubleshooting Guide

Problem	Cause	Solution
No response	Improper thawing procedure	Refer to thawing instructions in this user manual. Thawing process can have a significant effect on cell viability.
	Improper ligand used or improper ligand incubation time	See certificate of analysis for recommended ligand and assay conditions.
	Improper preparation of ligand (agonist or antagonist)	Refer to 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.
Decreased response	Cells are not adherent and exhibit incorrect morphology	Confirm adherence of cells using microscope.
Low or no signal	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 seconds/well.
Experimental S:B does not match datasheet value	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.
EC ₅₀ is right-shifted	Improper ligand handling or storage	Make sure 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.
High variability between replicates	Contamination from tips	Changing tips during serial dilutions can help to avoid carryover.
	Instrument calibration	Ensure dispensing equipment is properly calibrated, and proper pipetting technique is used.
		Non-binding surface plates may be necessary for hydrophobic compounds.

For additional information or technical support, please use the contact information below.

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 www.discoverx.com/license for limited use label license and trademark information for this product.