

## PathHunter® A549 IκB Degradation Cell Line

**Catalog Number:** 93-0538C15

**Lot Number:** See Vial

**Contents:** 2 x 10<sup>6</sup> cells per vial in 1 mL

### Background

PathHunter Degradation cell lines are engineered to express an Enzyme Donor (ED) tagged target protein. Activation of the pathway results in degradation of the ED-tagged protein. Addition of exogenous Enzyme Acceptor (EA), and buffer, lyses the cell and forces complementation of the ED and EA enzyme fragments. This results in the formation of a functional enzyme that hydrolyzes substrate to generate a chemiluminescent signal.

### Product Information

**Target Protein:** IκB

**Accession #:** NM\_020529

**Description:** nuclear factor of kappa light polypeptide gene enhancer in B-cells inhibitor, alpha

**Target ED Tag:** PL

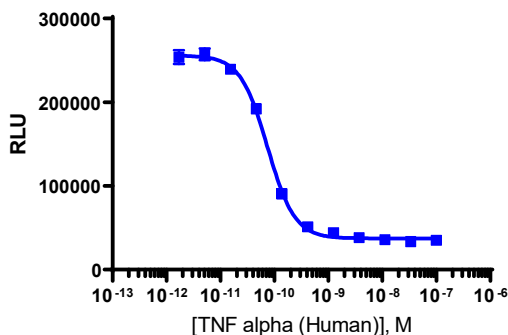
**Target Species:** Human

**Cell Type:** A549

**Storage:** Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

### Functional Performance

Cells were plated in a 384-well plate and incubated at 37°C and 5% CO<sub>2</sub> to allow the cells to attach and grow. Cells were then stimulated with a control compound, using the assay conditions described below. Following stimulation, signal was detected using the PathHunter Detection Kit according to the recommended protocol. Please refer to page 2 for recommended assay reagents, detection reagents, and control compounds.



<b>Cell Number/Well:</b>	5000
<b>Cell Seeding Time (hours):</b>	48
<b>Control Agonist:</b>	TNFα (human)
<b>Ligand Incubation Time (minutes):</b>	120
<b>Ligand Incubation Temperature (°C):</b>	RT
<b>EC<sub>50</sub> for compound stimulation (nM):</b>	0.07
<b>Signal:Background at agonist E<sub>max</sub>:</b>	7.2

To demonstrate functional response of these cells it is recommended to use 11 or more 4-fold serial dilutions of the reference ligand starting at a top concentration of 100 ng/mL.

### Passage Stability

This cell line has been confirmed to be stable through a minimum of 10 passages with no significant drop in assay window or change in EC<sub>50</sub>.

### Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

### Required Materials

The following additional materials are required but not provided:

Product Use*	Product Description	Catalog Number
Detection	PathHunter® ED Bioassay Detection Kit	93-1043
Cell Culture	AssayComplete™ Cell Culture Kit-105	92-3105G
Cell Plating	AssayComplete™ Cell Plating 38 Reagent	93-0563R38A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T1	92-4101TR
Cell Freezing	AssayComplete™ Freezing Reagent F1	92-5101FR

\*Please inquire about our cell line-specific AssayComplete Starter Packs to get you started with your cell culture needs.

### Required Antibiotics

Antibiotic Name	Concentration (µg/mL)	Catalog Number
AssayComplete™ Puromycin	Not Applicable	Not Applicable
AssayComplete™ Hygromycin B	Not Applicable	Not Applicable
AssayComplete™ G418	500	92-0030

### Additional Ligand Information

**Control Agonist:** TNFα (human)

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1097)

**Ordering:** +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)

**Technical support:** +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)

**General product information:** [www.discoverx.com](http://www.discoverx.com)

### 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](https://discoverx.com/license) for a list of products that are governed by limited use label license terms and relevant patent and trademark information.