

## PathHunter® LPS Signaling Reporter Cell Line (THP-1)

**Catalog Number:** 93-1187C14

**Lot Number:** See Vial

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

### Background

PathHunter Pathway Reporter cell lines are engineered to express an Enzyme Donor (ED) tagged reporter protein controlled by a pathway-inducible transcriptional response element. Pathway activation via endogenous or exogenous target results in induced expressions 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

**Response Element** NF-Kb

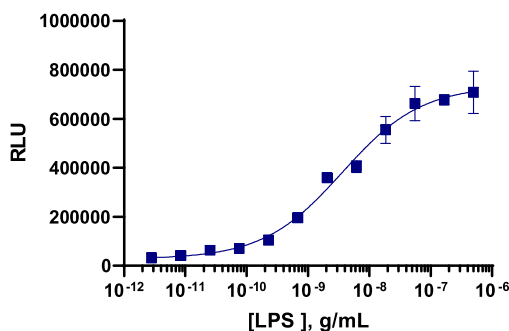
**Reporter ED Tag:** ePL

**Cell Type:** THP-1

**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 96-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>	24
<b>Control Agonist:</b>	LPS
<b>Ligand Incubation Time (minutes):</b>	300
<b>Ligand Incubation Temperature (°C):</b>	37
<b>EC<sub>50</sub> for compound stimulation ng/mL</b>	3.64
<b>Signal:Background at agonist E<sub>max</sub>:</b>	22.3
<b>Detection Incubation Time (hrs.)</b>	1

Cell line growth media should include 55 µM β-mercaptoethanol. ATCC recommends culture seeding at 0.2- 0.4M cells/mL for optimal growth. Cell line may require a few passages after thaw to achieve best results.

### 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 <sup>®</sup> ProLabel <sup>®</sup> /ProLink <sup>™</sup> Detection Kit	93-0812
Cell Culture	AssayComplete <sup>™</sup> Cell Culture Kit-101	92-3101G
Cell Plating	AssayComplete <sup>™</sup> Cell Plating 0 Reagent	93-0563R0A
Cell Detachment	Not Applicable	Not Applicable
Cell Thawing	AssayComplete <sup>™</sup> Thawing Reagent T6	92-4106TR
Cell Freezing	AssayComplete <sup>™</sup> Freezing Reagent F5	92-5105FR

\*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 <sup>™</sup> Puromycin	Not Applicable	Not Applicable
AssayComplete <sup>™</sup> Hygromycin B	150	92-0029
AssayComplete <sup>™</sup> G418	Not Applicable	Not Applicable

### Additional Ligand Information

**Control Agonist:** LPS

**Vendor:** Sigma (Catalog No. L2630) or equivalent

### 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 <https://www.discoverx.com/legal-policies/> for limited-use label license agreement and trademark information for this product.