

## PathHunter® TREM1-DAP12 NFAT Signaling Reporter Cell Line (Jurkat)

**Catalog Number:** 93-1167C19

**Lot Number:** See Vial

**Contents:** 3 x 10<sup>6</sup> cells per vial in 1 mL

### Background

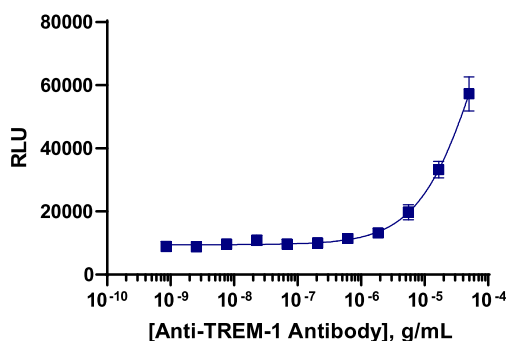
PathHunter Pathway Reporter cell lines are engineered to express an Enzyme Donor (ED) tagged reporter protein controlled by a pathway-inducible transcriptional response element. Pathway activation via endogenous or exogenous target results in induced expressions of the ED-tagged protein. Addition of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

<b>Target Receptor</b>	TREM1	<b>Co-Receptor</b>	DAP12
<b>Target Accession #</b>	NM_018643.3	<b>Accession #</b>	NM_003332.3
<b>Description:</b>	Triggering receptor expressed on myeloid cells 1	<b>Description:</b>	TYRO protein tyrosine kinase-binding protein
<b>Target Species:</b>	Human		
<b>Response Element</b>	NFAT		
<b>Reporter ED Tag:</b>	ePL		
<b>Cell Type:</b>	Jurkat		
<b>Storage:</b>	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>	10000
<b>Cell Seeding Time (hours):</b>	0
<b>Control Agonist:</b>	Anti-TREM1 Antibody
<b>Ligand Incubation Time (minutes):</b>	Overnight
<b>Ligand Incubation Temperature (°C):</b>	37
<b>EC<sub>50</sub> for compound stimulation</b>	NA
<b>Signal:Background at agonist E<sub>max</sub>:</b>	6.7
<b>Detection Incubation Time (hrs.)</b>	1

### 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® ProLabel®/ProLink™ Detection Kit	93-0812
Cell Culture	AssayComplete™ Cell Culture Kit-101	92-3101G
Cell Plating	AssayComplete™ Cell Plating 0 Reagent	93-0563R0A
Cell Detachment	Not Applicable	Not Applicable
Cell Thawing	AssayComplete™ Thawing Reagent T6	92-4106TR
Cell Freezing	AssayComplete™ 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™ Puromycin	0.25	92-0028
AssayComplete™ Hygromycin B	250	92-0029
AssayComplete™ G418	500	92-0030
AssayComplete™ Blasticidin	Not Applicable	R21001 (Invitrogen)

### Additional Ligand Information

**Control Agonist:** Anti-TREM1 Antibody

**Vendor:** R&D Systems (Catalog No. MAB-1278-100) 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.