

PathHunter® IL-2 Signaling Reporter Cell Line (REH)

Catalog Number: 93-1180C055

Lot Number: See Vial

Contents: 3 x 10⁶ 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.

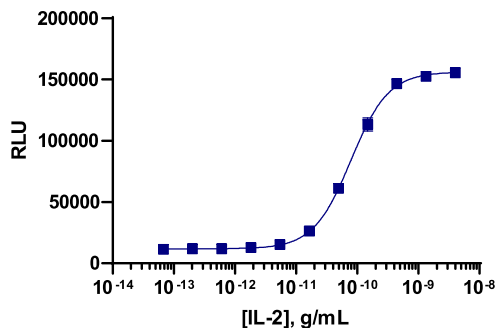
Product Information

Target Receptor	IL2RA	Co-Receptor	IL2RB
Target Accession #	NM_000417.2	Accession #	NM_000878.5
Description:	interleukin 2 receptor subunit alpha	Description:	interleukin 2 receptor subunit beta
Target Species:	Human	Target Protein 3:	IL2RG
Response Element	STAT5	Accession #:	NM_000206.2
Reporter ED Tag:	ePL	Description:	Interleukin 2 receptor, gamma
Cell Type:	REH		

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% CO₂ 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.



Cell Number/Well:	10000
Cell Seeding Time (hours):	0
Control Agonist:	IL-2
Ligand Incubation Time (minutes):	300
Ligand Incubation Temperature (°C):	37
EC₅₀ for compound stimulation ng/mL	0.079
Signal:Background at agonist E_{max}:	14.4
Detection Incubation Time (hrs.)	1

To maintain healthy and responsive REH cells during culture, always pellet the cells, resuspend cells in fresh growth media and use the cell suspension to plate in fresh media. In addition, cultures should be maintained at a density of <1.5 million cells/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 EC₅₀.

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	150	92-0029
AssayComplete [™] G418	500	92-0030
AssayComplete [™] Blastcidin	3	R21001 (Invitrogen)

Additional Ligand Information

Control Agonist: IL-2

Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1253)

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.