

PathHunter® U2OS PD-L1/TCR Activator Ligand Cell Line

Catalog Number: 93-1154C3 **Lot Number:** See Vial
Contents: 2 vials, 1.2 x 10⁶ cells per vial in 1 mL

Background

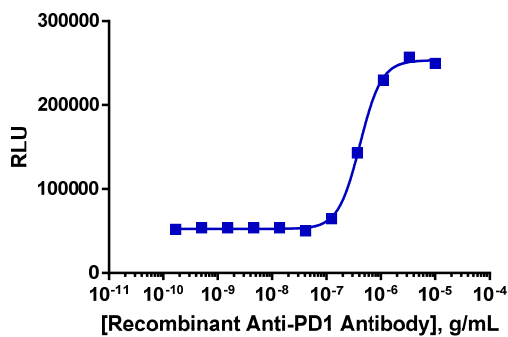
Ligand / TCR Activator Cell Lines are companion products to the Pathway Checkpoint Reporter Assays, which have been engineered to exogenously co-express a cell surface TCR activator, to stimulate TCR signaling, and a Checkpoint Receptor ligand, to counteract TCR activation. When co-cultured with its cognate pathway reporter assay cells, the ligand(s) expressed on the Ligand / TCR Activator Cell Line will attenuate TCR activation in the pathway reporter cell line. Addition of checkpoint receptor antagonists to block TCR inhibition results in increased TCR activation and increased ED-tagged reporter protein expression that is detected by the addition of exogenous Enzyme Acceptor (EA), and buffer to force complementation of the ED and EA enzyme fragments, and results in functional enzyme that hydrolyzes substrate to generate a chemiluminescent signal

Product Information

Target Protein : PD-L1
Accession #: NM_014143.3
Description: Programmed cell death 1 ligand 1
Species: Human
Cell Type: U2OS
Storage: Short term (<24 h): Store at -80°C; Long term (>24 h): Store in vapor phase of liquid nitrogen.

Functional Performance

Pathway Reporter assay cells were seeded in a 96-well plate and then pre-incubated with a control antibody, followed by co-incubation with the ligand cell line 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.



Pathway Reporter Assay: PathHunter Jurkat PD-1 Pathway Reporter Cell Line
Signaling Assay Cell Number/Well: 20000
Control Antibody: Recombinant Anti-PD1 Antibody [NAT105]
Antibody Pre-Incubation Time (minutes): 60
Antibody Pre-Incubation Temperature (°C): 37
IC₅₀ for Antibody (ng/mL): 411.6
Signal:Background at agonist E_{max}: 4.8
Ligand Cell Number/Well: 10000
Receptor:Ligand Cell Ratio: 2:1
Ligand Incubation Time (minutes): 300
Ligand Incubation Temperature (°C): 37

Passage Stability

This cell line has been confirmed to be stable through a minimum of 10 passages with no significant drop in expression levels.

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-103	92-3103G
Cell Plating	AssayComplete™ Cell Plating 0 Reagent	93-0563R0A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T3	92-4103TR
Cell Freezing	AssayComplete™ Freezing Reagent F3	92-5103FR

*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	Not Applicable	Not Applicable
AssayComplete™ G418	500	92-0030

To be used with: PathHunter Jurkat PD-1 Pathway Reporter Cell Line (93-1153C19)

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: 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 for a list of products that are governed by limited use label license terms and relevant patent and trademark information.