

PathHunter® eXpress IL7R/CRLF2 Dimerization Assay

Catalog Number: 93-1019E13

Lot Number:

See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

The PathHunter® Dimerization assay detects ligand induced dimerization of two subunits of a receptor-dimer pair. The cells have been engineered to co-express one receptor subunit fused to Enzyme Donor (ED), and a second dimer partner fused to Enzyme Acceptor (EA). Cytoplasmic tail may have been deleted from one or both receptors. Binding of an agonist to one receptor subunit induces it to interact with its dimer partner, forcing complementation of the two enzyme fragments. This results in the formation of a functional enzyme that hydrolyzes a substrate to generate a chemiluminescent signal. These cells have been modified to prevent long term propagation and expansion using a proprietary compound that has no apparent effect on assay performance.

Product Information

Target Protein 1: IL7R

Target Protein 2: CRLF2

Amino Acid Range: 1 - 270

Amino Acid Range: 1 - 259

Accession #: NM_002185.3

Accession #: NM_022148.3

Description: Interleukin 7 receptor

Description: cytokine receptor-like factor 2

Target Tag 1: PK1

Target Tag 2: EA

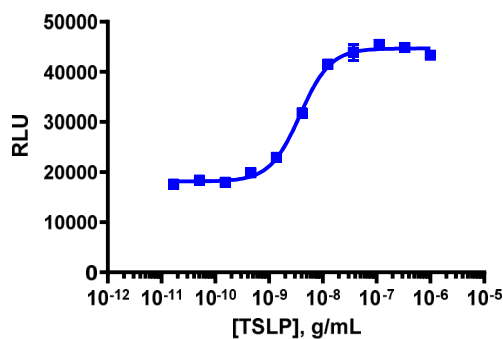
Target Species: Human

Cell Type: DLD1

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/5% CO₂ for the Cell Seeding Time indicated below. Cells were then stimulated with a control compound using the assay conditions described below. Following stimulation, signal was detected using the PathHunter Detection Reagents provided in the kit according to the recommended protocol. For a detailed protocol, please refer to the user manual.



| | |
|---|--|
| Cell Number/Well: | 10000 |
| Cell Plating Reagent: | AssayComplete™ Cell Plating 25 Reagent |
| Cell Seeding Time (Hours): | 4 |
| Control Compound: | TSLP |
| Compound Incubation Time (minutes): | 960 |
| Compound Incubation Temperature (°C): | 37 |
| EC₅₀ for Compound Stimulation ng/mL | 3.8 |
| Signal:Background at Compound E_{max}: | 2.6 |

Additional Ligand Information

Control Compound: TSLP

Vendor: Eurofins DiscoverX® (Catalog No. 92-1276)

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.

