

## PathHunter® U2OS BCMA Internalization Cell Line

**Catalog Number:** 93-1145C3

**Lot Number:** See Vial

**Contents:** 2 x 10<sup>6</sup> cells per vial in 1 mL

### Background

PathHunter Receptor Internalization cell lines are engineered to co-express the ProLink™ (PK) tagged receptor, and an Enzyme Acceptor (EA) tag localized to the endosomes. Receptor antibody binding to the receptor-PK fusion protein results in receptor internalization and colocalization with EA-tagged endosomes, forcing complementation of the two β-galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

### Product Information

**Target Receptor:** BCMA

**Description:** Tumor necrosis factor receptor superfamily member 17

**Accession Number:** NM\_001192.2

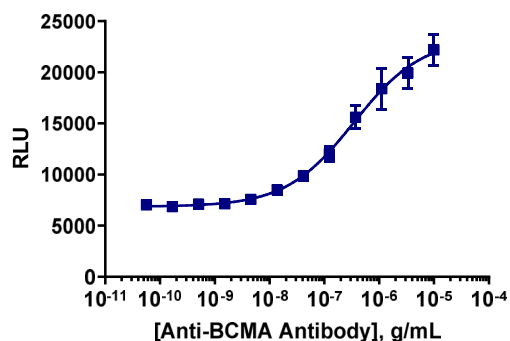
**Target Species:** Human

**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

Cells were plated in a 96-well plate and incubated overnight at 37°C and 5% CO<sub>2</sub> to allow the cells to attach and grow. Cells were then stimulated with a control agonist, 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>	4
<b>Control Antibody:</b>	Anti-BCMA Antibody
<b>Assay Incubation Time (minutes):</b>	960
<b>Assay Incubation Temperature (°C):</b>	37
<b>EC<sub>50</sub> for Agonist Stimulation (ng/mL):</b>	346
<b>Signal:Background at Agonist E<sub>max</sub>:</b>	3.2
<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® Flash Detection Kit	93-0247
Cell Culture	AssayComplete™ Cell Culture Kit-103	92-3103G
Cell Plating	AssayComplete™ Cell Plating 5 Reagent	93-0563R5A
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	Not Applicable	Not Applicable
AssayComplete™ Hygromycin B	250	92-0029
AssayComplete™ G418	500	92-0030

### Additional Ligand Information

**Control Agonist:** Anti-BCMA Antibody

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1356)

### 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.