

## PathHunter® U2OS JAK3 Activity Assay

**Catalog Number:** 93-0796C3 **Lot Number:** See Vial  
**Contents:** 2 vials, 4 x 10<sup>6</sup> cells per vial in 1 mL

### Background

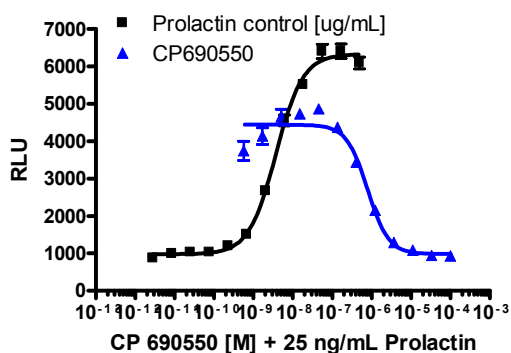
PathHunter Cytosolic Tyrosine Kinase (CTK) Activity cell lines are engineered to express the catalytic domain from CTK of interest fused to the cytokine receptor binding domain (CRBD) of a JAK kinase. The cells co-express an Enzyme Acceptor (EA) tagged SH2 domain and a ProLink™ (PK) tagged Cytokine receptor. The CRBD is designed to bind the CTK catalytic domain to the receptor-PK. Activation of the cytokine receptor-PK induces receptor phosphorylation by the CTK catalytic domain. This leads to SH2-EA recruitment, 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

**CTK Kinase Domain:** JAK3 [Accession #: NM\_000215] Amino Acid Range: (aa 800 - 1124)  
**Description:** Janus kinase 3  
**CRBD:** JAK2 [Accession #: BC039695] Amino Acid Range: (aa 1 - 255)  
**CTK Receptor** PRLR [Accession #: NM\_000949]  
**Species:** Human  
**SH2 Domain:** PLCG1  
**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 384-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 pre-incubated with inhibitor, followed by stimulation with the 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.



**Cell Number/Well:** 5000  
**Control Inhibitor:** CP-690550  
**Inhibitor Pre-incubation Time (minutes):** 60  
**Inhibitor Pre-incubation Temperature (°C):** RT  
**IC<sub>50</sub> for the Inhibitor (nM):** 792.6  
**Signal:Background of the Inhibitor:** 5.2  
**Control Agonist:** Prolactin  
**Agonist Incubation Time (minutes):** 150  
**Agonist Incubation Temperature (°C):** RT

### Passage Stability

This cell line has been confirmed to be stable through 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® Detection Kit	93-0001
Cell Culture	AssayComplete™ Cell Culture Kit-103	92-3103G
Cell Plating	AssayComplete™ Cell Plating 16 Reagent	93-0563R16A
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	250	92-0029
AssayComplete™ G418	500	92-0030

### Additional Ligand Information

**Control Agonist:** Prolactin

**Vendor:** DiscoverX® (Catalog No. 92-1144)

**Control Inhibitor** CP-690550

**Vendor:** Selleck (Catalog No. S5001)

For order placement or technical support, please call 1.866.448.4864 (North America) or +44.121.260.6142 (Europe) or e-mail [info@discoverx.com](mailto:info@discoverx.com). For additional information, please visit [discoverx.com](http://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](https://discoverx.com/license) for a list of products that are governed by limited use label license terms and relevant patent and trademark information.