

## PathHunter® U2OS CDC25A Degradation Cell Line

**Catalog Number:** 93-0765C3 **Lot Number:** See Vial

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

### Background

PathHunter Degradation cell lines are engineered to express an Enzyme Donor (ED) tagged target protein. Activation of the pathway results in degradation 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 Protein:** CDC25A

**Accession #:** NM\_001789.2

**Description:** Cell division cycle 25 homolog A

**Target ED Tag:** PK2

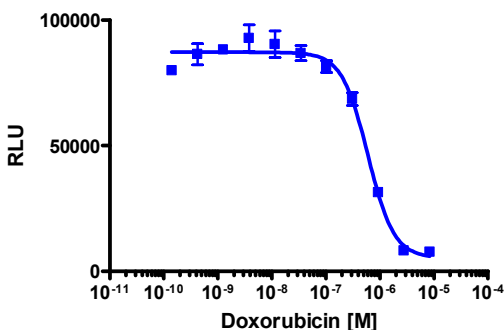
**Target Species:** Human

**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 384-well plate and incubated at 37°C and 5% CO<sub>2</sub> 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:** 2500  
**Cell Seeding Time (hours):** 24  
**Control Agonist:** Doxorubicin HCl  
**Ligand Incubation Time (minutes):** 900  
**Ligand Incubation Temperature (°C):** 37  
**EC<sub>50</sub> for compound stimulation (nM):** 598  
**Signal:Background at agonist E<sub>max</sub>:** 11.2

### 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® ProLabel®/ProLink™ Detection Kit | 93-0812        |
| 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 | Not Applicable        | Not Applicable |
| AssayComplete™ G418         | 500                   | 92-0030        |

### Additional Ligand Information

**Control Agonist:** Doxorubicin HCl

**Vendor:** DiscoverX® (Catalog No. 92-1101)

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.