

## PathHunter® U2OS ErbB2/ErbB4 Dimerization Cell Line

**Catalog Number:** 93-0960C3

**Lot Number:** See Vial

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

### Product Information

**Target Protein 1:** ErbB2

**Target Protein 2:** ErbB4

**Amino Acid Range:** 1 - 686

**Amino Acid Range:** 1 - 686

**Accession #:** NM\_004448.2

**Accession #:** NM\_005235.2

**Description:** v-erb-b2 erythroblastic leukemia viral oncogene homolog 2

**Description:** v-erb-b2 erythroblastic leukemia viral oncogene homolog 4

**Target Tag 1:** PK1

**Target Tag 2:** EA

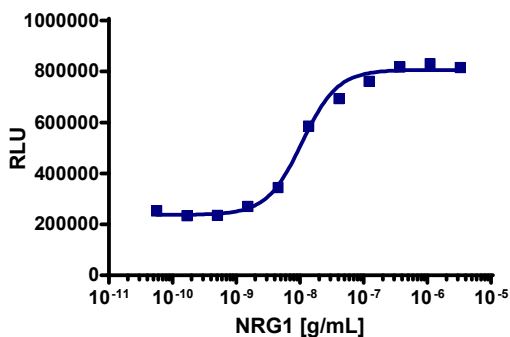
**Target Species:** Human

**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 seeded in a 96-well plate, incubated at 37°C / 5% CO<sub>2</sub> followed by stimulation with a control ligand, as defined in the assay conditions below. After stimulation, assay 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

**Cell Seeding Time (Hours):** 0

**Control Compound:** Heregulin-β1

**Compound Incubation Time (minutes):** Overnight

**Compound Incubation Temperature (°C):** 37

**EC<sub>50</sub> for Compound Stimulation (ng/mL):** 10.9

**Signal:Background at Compound E<sub>max</sub>:** 3.5

Optimal ligand incubation time is 16hrs.

### 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 <sup>®</sup> 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      |
| Ligand Dilution | AssayComplete™ Protein Dilution Buffer      | 92-0023M       |

\*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 Compound:** Heregulin-β1

**Vendor:** Peprotech (Catalog No. 100-03)

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.