

## PathHunter® eXpress TORC2 Nuclear Translocation Assay

**Catalog Number:** 93-0813E3

**Lot Number:** See Vial

**Contents:** 1 x 10<sup>6</sup> cells per vial in 0.1 mL

### Background

PathHunter® Nuclear Translocation cells are engineered to co-express two fusion proteins: a) Enzyme Donor (ED) tagged target protein; b) an Enzyme Acceptor (EA) tagged TAZ domain, derived from the CBP/P300 transcription factor, that localizes to the nucleus. Depending on the assay, activation of the signaling pathway can either a) induce translocation of the ED-tagged target protein into the nucleus, which will force complementation of the two enzyme fragments, and result in the formation of a functional enzyme that will hydrolyze substrate and generate a chemiluminescent signal; or b) induce the ED-tagged protein to vacate the nucleus, resulting in a decrease of functional enzyme and a subsequent decrease of chemiluminescent signal. Some nuclear translocation assays will also co-express an untagged secondary protein involved in the pathway of interest. 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

**Translocating Protein:** TORC2

**Accession #:** NM\_181715

**Description:** CREB regulated transcription coactivator 2

**ED Tag:** PK1

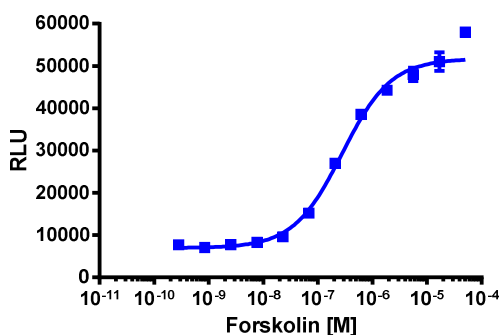
**Protein Species:** Human

**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 at 37°C and 5% CO<sub>2</sub>. 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 19 Reagent

**Control Compound:** Forskolin

**Cell Incubation Time (Hours):** 24

**Compound Incubation Time (minutes):** 180

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

**EC<sub>50</sub> for Compound Stimulation:** 278

**Signal:Background at Compound E<sub>max</sub>:** 7.5

### Additional Ligand Information

**Control Compound:** Forskolin

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-0005)

### 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.

**Ordering:** +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)

**Technical support:** +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)

**General product information:** [www.discoverx.com](https://www.discoverx.com)