

PathHunter® eXpress β -Catenin Nuclear Translocation Assay

Catalog Number: 93-0743E3

Lot Number: See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

PathHunter® Nuclear Translocation cells are engineered to co-express two proteins: a) Enzyme Donor (ED) tagged target protein; b) an Enzyme Acceptor (EA), which is localized 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: β -Catenin

Accession #: NM_001904

Description: Catenin (cadherin-associated protein), beta 1

ED Tag: PL

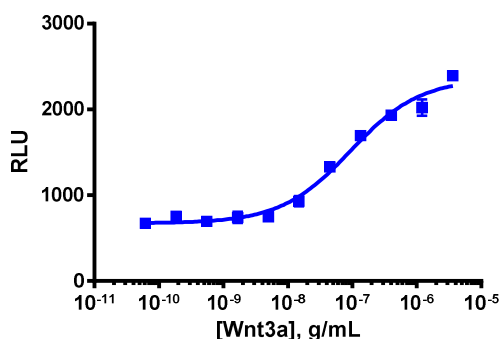
Protein Species: Human

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% CO₂. 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 22 Reagent

Control Compound: Wnt3a protein

Cell Incubation Time (Hours): 24

Compound Incubation Time (minutes): 360

Compound Incubation Temperature (°C): 37

EC₅₀ for Compound Stimulation: 93.6

Signal:Background at Compound E_{max}: 3.6

Additional Ligand Information

Control Compound: Wnt3a protein

Vendor: Eurofins DiscoverX® (Catalog No. 92-1100)

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 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

Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com

General product information: www.discoverx.com