

## PathHunter® eXpress HEK 293 IκB Degradation Assay

**Catalog Number:** 93-0538E1

**Lot Number:**

See Vial

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

### Background

PathHunter Degradation cells 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. 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

**Target Protein:** IκB

**Accession #:** NM\_020529

**Description:** nuclear factor of kappa light polypeptide gene enhancer in B-cells inhibitor, alpha

**Target ED Tag:** PL

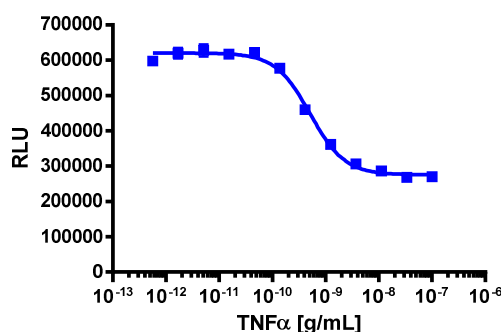
**Target Species:** Human

**Cell Type:** HEK 293

**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 7 Reagent

**Control Compound:**

TNFα (human)

**Cell Incubation Time (Hours):**

24

**Compound Incubation Time (minutes):**

60

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

37

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

0.51

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

2.2

### Additional Ligand Information

**Control Compound:** TNF $\alpha$  (human)

**Vendor:** Eurofins DiscoverX<sup>®</sup> (Catalog No. 92-1097)

### 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)