

## PathHunter<sup>®</sup> eXpress FGFR4- $\beta$ -Klotho Functional Assay

**Catalog Number:** 93-1060E3

**Lot Number:**

See Vial

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

### Background

PathHunter RTK Functional cell lines are engineered to co-express a ProLink™ (PK) tagged RTK, an Enzyme Acceptor (EA) tagged SH2 domain, and, in some cases, an untagged co-receptor. Activation of the RTK-PK induces receptor dimerization, leading to SH2-EA recruitment, and forcing complementation of the two  $\beta$ -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme 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 RTK:** FGFR4

**Co-receptor:**  $\beta$ -Klotho

**RTK Accession #:** NM\_002011

**Accession #:** NM\_175737.3

**Description:** Fibroblast growth factor receptor 4

**Description:** KLB

**Kinase Species:** Human

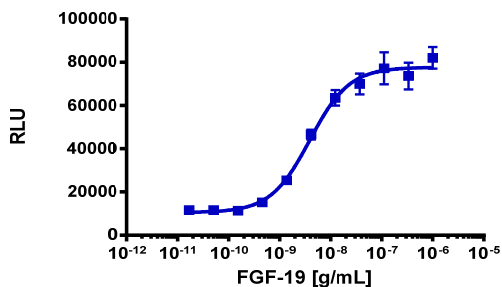
**SH2 Domain:** PLCG1

**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 stimulated with a control agonist, using the assay conditions described below. Following stimulation, signal was detected according to the recommended protocol. Please refer below for information on control compounds.



**Cell Number/Well:**

10000

**Control Agonist:**

FGF-19

**Cell Plating Reagent:**

AssayComplete™ Cell Plating 0 Reagent

**Cell Incubation Time (Hours):**

4

**Agonist Incubation Time (minutes):**

960

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

37

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

3.9

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

7.0

### Additional Information

**Control Agonist:** FGF-19

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1301)

### 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)