

## InCELL Hunter™ eXpress CLK1 Kinase Binding Assay

**Catalog Number:** 96-0059E1

**Lot Number:** See Vial

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

### Background

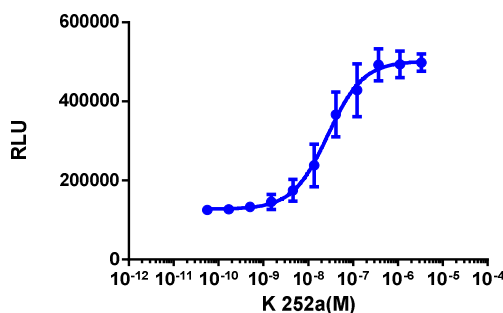
InCELL Hunter technology measures compound binding to an intracellular target by detecting changes in protein stability. In this assay, the target is fused to the enhanced ProLabel®(ePL) enzyme fragment and expressed in the selected cell background. The amount of fusion is quantified through the addition of substrate and complementary enzyme acceptor fragment (EA). The small size, but robust activity of the ePL tag makes it an ideal fusion partner. The small disordered peptide has minimal effects on target half-life, and generates high levels of enzyme activity, resulting in a sensitive measurement of protein levels. 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:** CLK1 [Accession #: NM\_004071.3]  
**Description:** CDC-Like Kinase 1  
**Accession Number:** NM\_004071.3  
**Target Species:** Human  
**ProLink™ Tag:** ePL **ProLink™ Tag located at (terminus):** C-Terminus  
**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 InCELL Hunter 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  
**Control Compound:** K 252a  
**Cell Plating Reagent:** AssayComplete™ Cell Plating 0 Reagent  
**Cell Incubation Time (Hours):** 48  
**Compound Incubation Time (minutes):** 360  
**Compound Incubation Temperature (°C):** 37  
**EC<sub>50</sub> of Compound (nM):** 27.5  
**Signal:Background at E<sub>max</sub>:** 3.9  
**1:5 EA Dilution Required?** Yes

### Additional Ligand Information

**Control Compound:** K 252a

**Vendor:** DiscoverX® (Catalog No. 92-1245)

### 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)