

## InCELL Hunter™ eXpress HEK 293 BRD2(1) Bromodomain Assay

**Catalog Number:** 96-0004E1

**Lot Number:** See Vial

**Contents:** 1.2 x 10<sup>6</sup> cells per vial in 0.1 mL

### Background

The 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 protein is quantified through the addition of cell lysis buffer and a complementary enzyme acceptor fragment (EA). The two enzyme fragments complement to form a functional enzyme that is capable of hydrolyzing substrate and generating 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:** BRD2(1) Amino Acid Range: (aa 74 - 188)

**Description:** Bromodomain containing 2, 1st bromodomain

**Accession Number:** NM\_005104.3

**Target Species:** Human

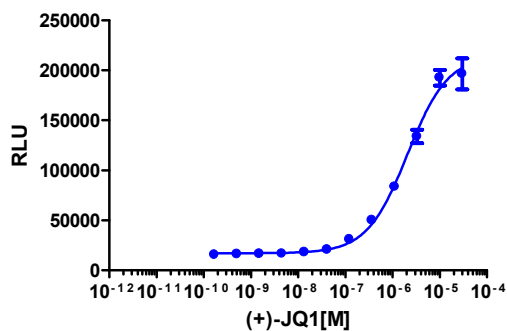
**ProLink™ Tag:** ePL **ProLink™ Tag located at (terminus):** C-Terminus

**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:** (+)-JQ1

**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> for Compound (nM):** 2071

**Signal:Background at E<sub>max</sub>:** 12.4

**1:5 EA Dilution Required?** No

### Additional Ligand Information

**Control Compound:** (+)-JQ1

**Vendor:** Eurofins DiscoverX<sup>®</sup> (Catalog No. 92-1149)

### 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)