

## InCELL Hunter™ eXpress BCL2L1-BAX Protein Binding Assay

**Catalog Number:** 96-0009E3

**Lot Number:**

See Vial

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

### Background

InCELL Hunter assay detects degradation of a partner protein that is ordinarily stabilized by its natural interaction with the target protein. The cell lines are engineered to co-express an untagged target protein, and an Enzyme Donor (ED) tagged partner protein. Either protein may be full length or truncated. Addition of inhibitors disrupting the interaction of the two proteins 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:** BCL2L1 [Accession #: NM\_138578.1] Amino Acid Range: (aa 1 - 233)

**Target Description:** BCL2-like 1

**Partner Protein:** BAX [Accession #: NM\_138761] Amino Acid Range: (aa 52 - 72)

**Partner Description:** BCL2-associated X protein

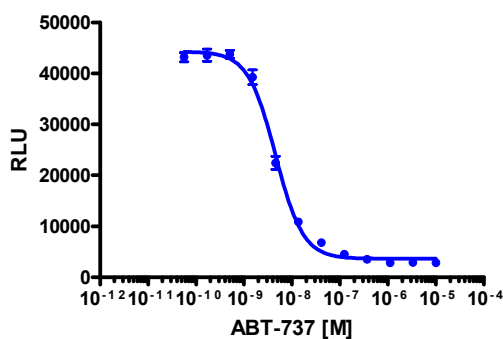
**Target Species:** Human

**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% 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:** ABT-737

**Cell Plating Reagent:** AssayComplete™ Cell Plating 16 Reagent

**Cell Incubation Time (Hours):** 24

**Compound Incubation Time (minutes):** 180

**Compound Incubation Temperature (°C):** 37

**EC<sub>50</sub> of Compound (nM):** 4.6

**Signal:Background at E<sub>max</sub>:** 15.2

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

### Additional Ligand Information

**Control Compound:** ABT-737

**Vendor:** Eurofins DiscoverX® (Catalog No. 92-1151)

### 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)