

InCELL Hunter™ eXpress HEK 293 BRDT(1,2) Bromodomain Assay

Catalog Number: 96-0046E1

Lot Number: See Vial

Contents: 1.2 x 10⁶ cells per vial in 0.1 mL

Background

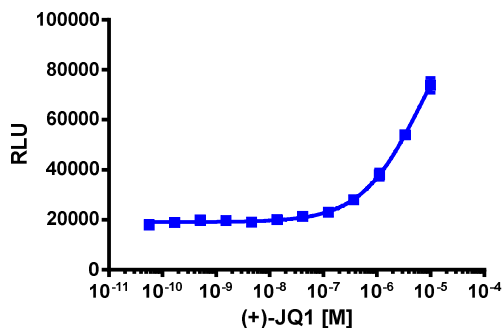
The InCELL Hunter technology measures compound binding to an intracellular target by detecting changes in protein stability. In this assay, the target is fused to the enhanced ProLabel®(ePL) enzyme fragment and expressed in the selected cell background. The amount of fusion protein is quantified through 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:	BRDT(1,2)	Amino Acid Range: (aa 28 - 376)
Description:	Bromodomain containing testis-specific, 1st & 2nd bromodomains	
Accession Number:	NM_001726	
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% CO₂. 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 7 Reagent
Cell Incubation Time (Hours):	48
Compound Incubation Time (minutes):	360
Compound Incubation Temperature (°C):	37
EC₅₀ for Compound (nM):	6811
Signal:Background at E_{max}:	4.1
1:5 EA Dilution Required?	No

Additional Ligand Information

Control Compound: (+)-JQ1

Vendor: Eurofins DiscoverX® (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 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
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com