

PathHunter® HEK 293 AR Nuclear Translocation Cell Line

Catalog Number: 93-0177C1

Lot Number: See Vial

Contents: 1 x 10⁶ cells per vial in 1 mL

Background

PathHunter NHR Nuclear Translocation cell lines are engineered to co-express a ProLabel® (PL) tagged NHR, and Enzyme Acceptor (EA) that is localized in the cell nucleus. Activation of the receptor induces its translocation to the nucleus, forcing complementation of the two β -galactosidase enzyme fragments (EA and PL). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

Product Information

Target NHR AR

Description: Androgen receptor

Accession Number: NM_000044

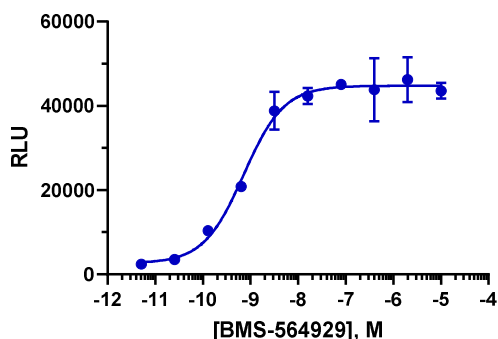
Target Species: Human

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 384-well plate and incubated overnight at 37°C and 5% CO₂ to allow the cells to attach and grow. Cells were then stimulated with a control compound, using the assay conditions described below. Following stimulation, signal was detected using the PathHunter Detection Kit according to the recommended protocol. Please refer to page 2 for recommended assay reagents, detection reagents, and control compounds



Cell Number/Well: 5000

Control Compound: BMS-564929

Compound Incubation Time (minutes): 360

Compound Incubation Temperature (°C): 37

EC₅₀ for Compound Stimulation (nM): 0.72

Signal:Background at Compound E_{max}: 17.2

Passage Stability

This cell line has been confirmed to be stable through a minimum of 10 passages with no significant drop in assay window or change in EC₅₀.

Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

Required Materials

The following additional materials are required but not provided:

Product Use*	Product Description	Catalog Number
Detection	PathHunter® Detection Kit	93-0001
Cell Culture	AssayComplete™ Cell Culture Kit-105	92-3105G
Cell Plating	AssayComplete™ Cell Plating 15 Reagent	93-0563R15A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T1	92-4101TR
Cell Freezing	AssayComplete™ Freezing Reagent F1	92-5101FR

*Please inquire about our cell line-specific AssayComplete Starter Packs to get you started with your cell culture needs.

Required Antibiotics

Antibiotic Name	Concentration (µg/mL)	Catalog Number
AssayComplete™ Puromycin	Not Applicable	Not Applicable
AssayComplete™ Hygromycin B	200	92-0029
AssayComplete™ G418	500	92-0030

Additional Ligand Information

Control Agonist: BMS-564929

Vendor: Eurofins DiscoverX® (Catalog No. 92-1363)

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

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.