

PathHunter® CHO-K1 RAR α Protein Interaction Cell Line

Catalog Number: 93-0325C2

Lot Number: See Vial

Contents: 1 x 10⁶ cells per vial in 1 mL

Background

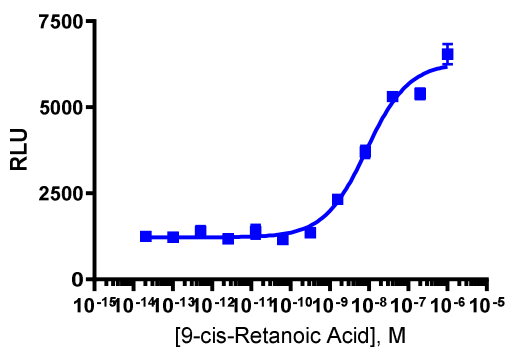
PathHunter NHR Protein Interaction cell lines are engineered to co-express a ProLink™ (PK) tagged NHR and an Enzyme Acceptor (EA) tagged SRC protein. Activation of the receptor results in SRC-EA recruitment, forcing complementation of the two β -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

Product Information

Target NHR:	RAR α
Description:	Retinoic acid receptor, alpha
Accession Number:	NM_000964
SRC Protein:	TRAP220
Target Species:	Human
Cell Type:	CHO-K1
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:	10000
Control Compound:	9- <i>cis</i> -Retinoic acid
Compound Incubation Time (minutes):	360
Compound Incubation Temperature (°C):	37
EC₅₀ for Compound Stimulation (nM):	8.5
Signal:Background at Compound E_{max}:	5.2

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-107	92-3107G
Cell Plating	AssayComplete™ Cell Plating 13 Reagent	93-0563R13A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T2	92-4102TR
Cell Freezing	AssayComplete™ Freezing Reagent F2	92-5102FR

*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	250	92-0029
AssayComplete™ G418	600	92-0030

Additional Ligand Information

Control Agonist: 9-*cis*-Retinoic acid

Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1136)

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.