

PathHunter® CHO-K1 NPR1 Functional Assay

Catalog Number: 93-0804C2 **Lot Number:** See Vial
Contents: 2 vials, 1 x 10⁶ cells per vial in 1 mL

Background

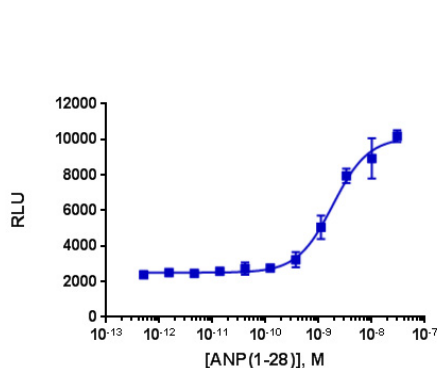
PathHunter NPR cell line express wild-type NPR1 receptors and are designed to detect increases in cGMP levels in response to agonist stimulation of the receptor. This cell line is designed to be used in conjunction with the HitHunter® cGMP Detection Kit.

Product Information

Target Protein: NPR1
Accession Number: NM_000906.3
Description: Natriuretic peptide receptor A
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/5% CO₂ to allow the cells to attach and grow. Cells were then stimulated with a control agonist, using the assay conditions described below. Following stimulation, signal was detected using the HitHunter cGMP 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 Agonist: ANP(1-28)

Ligand Incubation Time (minutes): 30

Ligand Incubation Temperature (°C): 37

EC₅₀ for compound stimulation (nM): 1.9

Signal:Background at agonist E_{max}: 4.3

Cells were plated in a 384-well plate and assayed in suspension mode. Cells were stimulated with a control agonist using the assay conditions described below. Following stimulation, signal was detected using the HitHunter cGMP Detection Kit according to the recommended protocol. Please refer to page 2 for the recommended assay reagents, detection reagents and control compounds.

Passage Stability

This cell line has been confirmed to be stable through 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	HitHunter [®] cGMP Assay	90-0039
Cell Culture	AssayComplete™ Cell Culture Kit-107	92-3107G
Cell Plating	AssayComplete™ Cell Plating 43 Reagent	93-0563R43A
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	Not Applicable	Not Applicable
AssayComplete™ G418	800	92-0030

Additional Ligand Information

Control Agonist: ANP(1-28)

Vendor: DiscoverX[®] (Catalog No. 92-1337)

For order placement or technical support, please call 1.866.448.4864 (North America) or +44.121.260.6142 (Europe) or e-mail info@discoverx.com. For additional information, please visit 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.