

PathHunter® eXpress NMBR Total GPCR Internalization Assay

Catalog Number: 93-0939E3

Lot Number:

See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

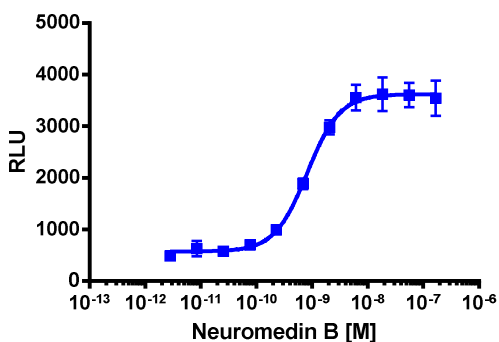
PathHunter eXpress Total GPCR Internalization cells are engineered to co-express the ProLink™ (PK) tagged GPCR, and an Enzyme Acceptor (EA) tag localized to the endosomes. Activation of the GPCR-PK induces internalization of the receptor in EA-tagged endosomes, forcing complementation of the two β-galactosidase enzyme fragments (EA and PK). The resulting functional enzyme 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 GPCR:	NMBR
Description:	Neuromedin B Receptor
Receptor Family:	Bombesin
Coupling:	Gq
Accession Number:	NM_002511.2
GPCR 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 stimulated with a control agonist, using the assay conditions described below. Following stimulation, signal was detected according to the recommended protocol. Please refer below for information on control compounds.



Cell Number/Well:

10000

Control Agonist:

Neuromedin B (porcine)

Cell Plating Reagent:

AssayComplete™ Cell Plating 5 Reagent

Cell Incubation Time (Hours):

24

Agonist Incubation Time (minutes):

180

Agonist Incubation Temperature (°C):

37

EC₅₀ for Agonist Stimulation (nM):

0.8

Signal:Background at Agonist E_{max}:

6.1

Additional Ligand Information

Control Agonist: Neuromedin B (porcine)

Vendor: Eurofins DiscoverX® (Catalog No. 92-1233)

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