

PathHunter® eXpress CXCR7 Activated GPCR Internalization Assay

Catalog Number: 93-0738E3

Lot Number: See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

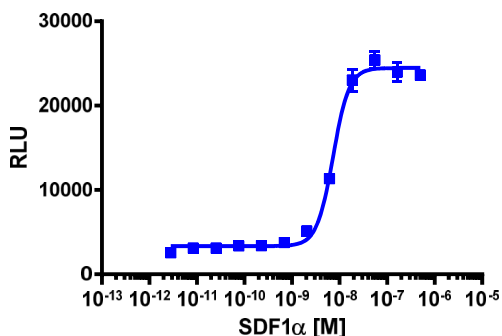
PathHunter eXpress β -Arrestin GPCR cells are engineered to co-express the ProLink™ (PK) tagged GPCR and the Enzyme Acceptor (EA) tagged β -Arrestin. Activation of the GPCR-PK induces β -Arrestin-EA recruitment, forcing complementation of the two β -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal. 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:	CXCR7
Description:	Chemokine (C-X-C motif) receptor 7
Receptor Family:	Chemokine
Coupling:	None
Accession Number:	NM_020311.2
GPCR Species:	Human
β-Arrestin Isoform:	β -Arrestin-2
ProLink™ Tag:	None
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:	Recombinant Human SDF-1 α (CXCL12)
Cell Plating Reagent:	AssayComplete™ Cell Plating 27 Reagent
Cell Incubation Time (Hours):	24
Agonist Incubation Time (Minutes):	240
Agonist Incubation Temperature (°C):	37
EC₅₀ for Agonist Stimulation (nM):	7.3
Signal:Background at Agonist E_{max}:	9.5

Additional Ligand Information

Control Agonist: Recombinant Human SDF-1 α (CXCL12)

Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1011)

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