

PathHunter® eXpress CCR5 Activated GPCR Internalization Assay

Catalog Number: 93-0653E3 **Lot Number:** See Vial

Contents: 1 x 10⁶ cells per vial in 0.1 mL

Background

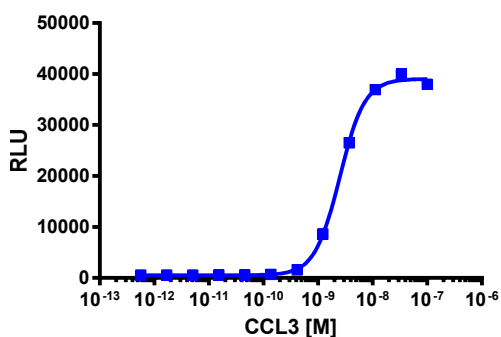
PathHunter eXpress Activated GPCR Internalization cells are engineered to co-express an untagged GPCR, Enzyme Acceptor (EA) tagged β -Arrestin, and a ProLink™ (PK) tag localized to the endosomes. Activation of the GPCR induces β -Arrestin recruitment, followed by internalization of the Receptor/Arrestin-EA complex in PK-tagged endosomes. This forces complementation of the two β -galactosidase enzyme fragments (EA and PK), forming functional enzyme that 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: CCR5
Description: Chemokine (C-C motif) receptor 5
Receptor Family: Chemokine
Coupling: Gi/Go
Accession Number: BC038398
GPCR Species: Human
 β -Arrestin Isoform: β -Arrestin-2
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 MIP-1 α (CCL3)
Cell Plating Reagent: AssayComplete™ Cell Plating 0 Reagent
Cell Incubation Time (Hours): 48
Agonist Incubation Time (minutes): 180
Agonist Incubation Temperature (°C): 37
EC₅₀ for Agonist Stimulation (nM): 2.5
Signal:Background at Agonist E_{max}: 70.7

Additional Ligand Information

Control Agonist: Recombinant Human MIP-1 α (CCL3)

Vendor: Eurofins DiscoverX[®] (Catalog No. 92-1002)

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 <https://www.discoverx.com/legal-policies/> for limited-use label license agreement and trademark information for this product.