

Certificate Of Analysis

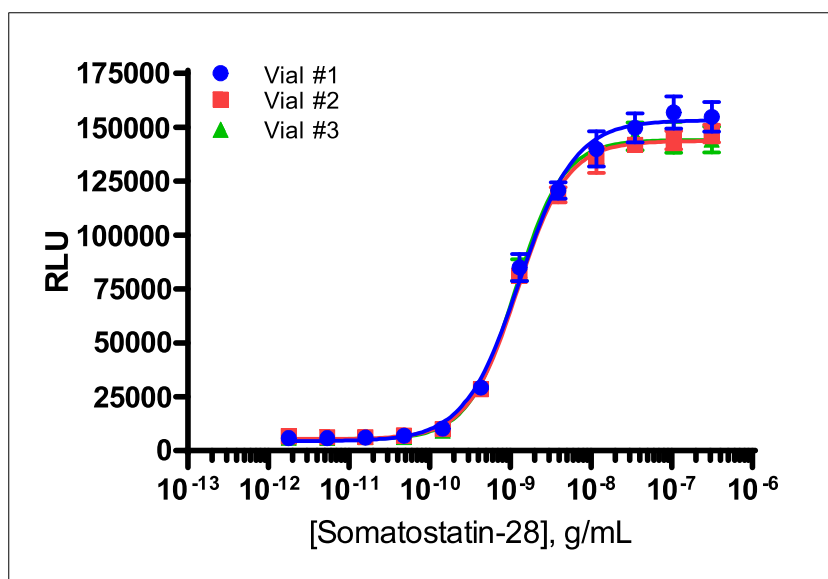
Background

PathHunter[®] β -Arrestin GPCR Bioassay 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.

| | |
|-----------------------------|---|
| Product Name | PathHunter[®] CHO-K1 SSTR2 Bioassay |
| Cryovial Label | CHO-K1 SSTR2 Beta-Arrestin Bioassay Cells |
| Bioassay Catalog # | 93-0181Y2 |
| Bioassay Manufactured Lot # | 21K2812 |
| Passage # @ Freezing | 3 |

| Assay Information | |
|------------------------------|---|
| Target 1 | SSTR2 |
| Target 1 Accession Number | AY236542 |
| Target 1 Description | Somatostatin receptor 2 |
| β -Arrestin Isoform | β -Arrestin-2 |
| Target Species | Human |
| Cell | CHO-K1 |
| CP Reagent | AssayComplete [™] Cell Plating 0 Reagent (DiscoverX, 93-0563R0A) |
| Ligand | Somatostatin 28 (DiscoverX, 92-1068) |
| Ligand Diluent | Protein Dilution Buffer |
| Detection Kit | PathHunter [®] Bioassay Detection Kit (DiscoverX, 93-0933) |
| Cell Number/Well | 10,000 |
| Cell Seeding Time (hours) | 48 |
| Ligand Inc Time (minutes) | 90 |
| Agonist Inc Temperature (°C) | 37 |

| Cell Density Information | |
|---|-------------------------------|
| Cell Number (millions) | 1.2 |
| Fill Volume per Vial (mL) | 0.1 |
| Cell Viability | |
| Viability at Initial Thaw (%) | 95 |
| Recovery After 24 Hours (%) | 219 |
| Mycoplasma and Sterility | |
| Mycoplasma Test | Passed |
| Sterility Test | Passed |
| Functional Performance (3 manufactured vials) | |
| S:B Ratio | Vial 1 27.2 |
| | Vial 2 22.7 |
| | Vial 3 23.6 |
| EC ₅₀ (g/mL) | Vial 1 1.3 x 10 ⁻⁹ |
| | Vial 2 1.2 x 10 ⁻⁹ |
| | Vial 3 1.2 x 10 ⁻⁹ |



| Shipping and Storage Information | |
|---|---|
| Shipping Conditions | Dry Ice |
| Storage Conditions | Short term (<24 hours): -80°C; Long term (>24 hours): Vapor phase of liquid nitrogen. |
| Manufacturing Date | October 2021 |
| Expiration Date | October 2024 |

Shelf life of over 3 years has been established for DiscoverX cell lines and Assay-Ready Cells in general, when stored in the vapor phase of liquid nitrogen.

Documented by / Date: _____

Approved by / Date: _____