

## Certificate Of Analysis

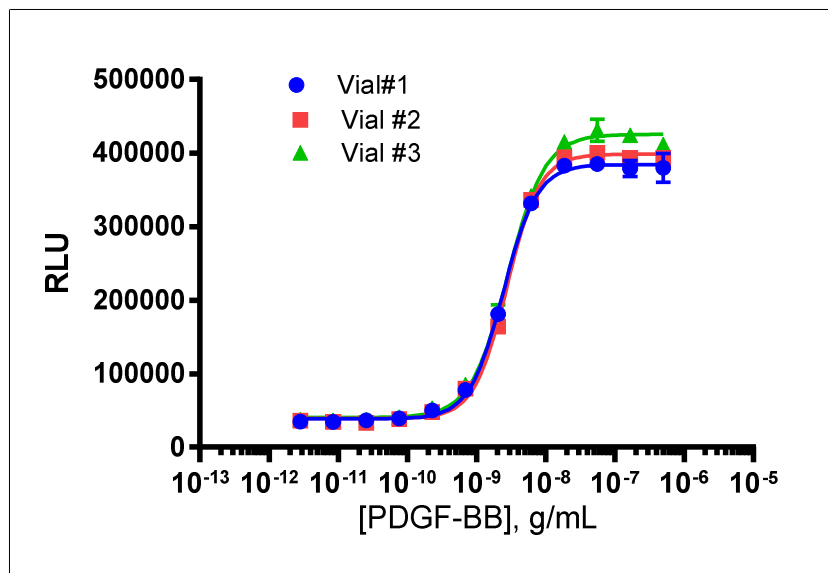
### Background

PathHunter<sup>®</sup> RTK Bioassay Cells are engineered to co-express a ProLink<sup>™</sup> (PK) tagged RTK, an Enzyme Acceptor (EA) tagged SH2 domain, and, in some cases, an untagged co-receptor. Activation of the RTK-PK induces receptor dimerization, leading to receptor phosphorylation, and SH2-EA recruitment, and forcing complementation of the two  $\beta$ -galactosidase enzyme fragments (EA and PK). The resulting functional enzyme hydrolyzes substrate to generate a chemiluminescent signal.

|                             |  |
|-----------------------------|--|
| <b>Product Name</b>         | <b>PathHunter<sup>®</sup> U2OS PDGFRB Bioassay Cells</b> |
| Cryovial Label              | U2OS PDGFRB Bioassay Cells                               |
| Bioassay Catalog #          | 93-0493Y3  |
| Bioassay Manufactured Lot # | 20B1008  |
| Passage # @ Freezing        | 3  |

| <b>Assay Information</b>     |   |
|------------------------------|---|
| Target                       | PDGFRB  |
| Target Accession Number      | NM_002609   |
| Target Description           | Platelet-derived growth factor receptor beta                                |
| Target 2                     | N/A   |
| Target 2 Accession Number    | N/A   |
| Target 2 Description         | N/A   |
| SH2 Domain                   | PLCG1   |
| Species                      | Human   |
| Cell                         | U2OS  |
| CP Reagent                   | AssayComplete <sup>™</sup> Cell Plating 22 Reagent (DiscoverX, 93-0563R22A) |
| Ligand                       | Recombinant Human PDGF-BB (DiscoverX, 92-1110)                              |
| Ligand Diluent               | AssayComplete <sup>™</sup> Protein Dilution Buffer                          |
| Detection Kit                | PathHunter <sup>®</sup> Bioassay Detection Kit (DiscoverX, 93-0933)         |
| Cell Number/Well             | 10,000  |
| Cell Seeding Time (hours)    | 24  |
| Ligand Inc Time (minutes)    | 180   |
| Agonist Inc Temperature (°C) | RT  |

| Cell Density Information                      |                               |
|---|-------------------------------|
| Cell Number (millions)                        | 1.2                           |
| Fill Volume per Vial (mL)                     | 0.1                           |
| Cell Viability                                |                               |
| Viability at Initial Thaw (%)                 | 98                            |
| Recovery After 48 Hours (%)                   | 216                           |
| Mycoplasma and Sterility                      |                               |
| Mycoplasma Test                               | Passed                        |
| Sterility Test                                | Passed                        |
| Functional Performance (3 manufactured vials) |                               |
| S:B Ratio                                     | Vial 1 10.9                   |
|   | Vial 2 10.9                   |
|   | Vial 3 11.0                   |
| EC <sub>50</sub> (g/mL)                       | Vial 1 2.4 x 10 <sup>-9</sup> |
|   | Vial 2 2.8 x 10 <sup>-9</sup> |
|   | Vial 3 2.8 x 10 <sup>-9</sup> |



| <b>Shipping and Storage Information</b> |   |
|---|---|
| Shipping Conditions                     | Dry Ice   |
| Storage Conditions                      | Short term (<24 hours): -80°C; Long term (>24 hours): Vapor phase of liquid nitrogen. |
| Manufacturing Date                      | February 2020   |
| Expiration Date                         | August 2023   |

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: \_\_\_\_\_