

## PathHunter<sup>®</sup> U2OS mPD-L1 Ligand Cell Line Starter Pack

**Catalog Number:** 493-1159C3

### Product Information

The Cell Line Starter Pack contains pre-validated, functionally tested reagents necessary to propagate and maintain the cell line referenced below. These reagents are especially developed to provide optimal cell culture environment and the highest possible assay performance.

### Intended Use

The Starter Pack is recommended to be used with the following cell line:

Part #	Description
93-1159C3	PathHunter <sup>®</sup> U2OS mPD-L1 Ligand Cell Line

### Starter Pack Contents

Product Use	Description	DiscoverX <sup>®</sup> Part #	Configuration
Cell Culture	AssayComplete <sup>™</sup> Cell Culture Kit-103	92-3103G	Kit
Cell Plating	Not included with Ligand cell lines	N/A	100 mL
Cell Thawing	AssayComplete <sup>™</sup> Thawing Reagent T3	92-4103TR	100 mL
Cell Freezing	AssayComplete <sup>™</sup> Freezing Reagent F3	92-5103FR	2 x 50 mL
Cell Detachment	AssayComplete <sup>™</sup> Cell Detachment Reagent	92-0009	50 mL

### Storage And Shipping Information

**Storage:** Store at -20°C

**Shipping:** Frozen on Dry Ice ( -70°C)

**Ordering:** +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
**Technical support:** +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
**General product information:** [www.discoverx.com](http://www.discoverx.com)