

## PathHunter<sup>®</sup> U2OS FcgRIa Clustering Cell Line Starter Pack

Catalog Number: 493-1138C3

### Product Information

The Cell Line Starter Pack contains functionally tested reagents necessary to propagate and maintain the cell line referenced below and run the assay. 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
93-1138C3	PathHunter <sup>®</sup> U2OS FcgRIa Clustering Cell Line

### Starter Pack Components

Product Use	Description	DiscoverX <sup>®</sup> Part #
Cell Culture	AssayComplete <sup>™</sup> Cell Culture Kit-103	92-3103G
Cell Plating	AssayComplete <sup>™</sup> Cell Plating 0 Reagent	93-0563R0A
Cell Thawing	AssayComplete <sup>™</sup> Thawing Reagent T3	92-4103TR
Cell Freezing	AssayComplete <sup>™</sup> Freezing Reagent F3	92-5103FR
Cell Detachment	AssayComplete <sup>™</sup> Cell Detachment Reagent	92-0009

### 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)

