

## PathHunter<sup>®</sup> Jurkat SIRPαSignaling Assay Starter Pack

**Catalog Number:** 493-1135C19

### 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
93-1135C19	PathHunter <sup>®</sup> Jurkat SIRPαSignaling Assay

### Starter Pack Components

Product Use	Description	DiscoverX <sup>®</sup> Part #
Cell Culture	AssayComplete <sup>™</sup> Cell Culture Kit-101	92-3101G
Cell Plating	AssayComplete <sup>™</sup> Cell Plating 0 Reagent	93-0563R0A
Cell Thawing	AssayComplete <sup>™</sup> Thawing Reagent T6	92-4106TR
Cell Freezing	AssayComplete <sup>™</sup> Freezing Reagent F5	92-5105FR
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)

For order placement or technical support, please call 1.866.448.4864 (North America) or +44.121.260.6142 (Europe) or e-mail [info@discoverx.com](mailto:info@discoverx.com). For additional information, please visit [discoverx.com](http://discoverx.com).