

PathHunter® Jurkat mPD-1 Signaling Assay Starter Pack

Catalog Number: 493-1158C19

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
93-1158C19	PathHunter® Jurkat mPD-1 Signaling Assay

Starter Pack Contents

Product Use	Description	DiscoverX® Part #	Configuration
Cell Culture	AssayComplete™ Cell Culture Kit-101	92-3101G	Kit
Cell Plating	AssayComplete™ Cell Plating 0 Reagent	93-0563R0A	100 mL
Cell Thawing	AssayComplete™ Thawing Reagent T6	92-4106TR	100 mL
Cell Freezing	AssayComplete™ Freezing Reagent F5	92-5105FR	2 x 50 mL
Cell Detachment	AssayComplete™ 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
 Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
 General product information: www.discoverx.com