

PathHunter® Jurkat CD47-Presenting Cell Line Starter Pack

Catalog Number: 493-1149C19

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
93-1149C19	PathHunter® Jurkat CD47-Presenting Cell Line

Starter Pack Components

Product Use	Description	DiscoverX® Part #
Cell Culture	AssayComplete™ Cell Culture Kit-101	92-3101G
Cell Plating	Not included with Ligand cell lines	N/A
Cell Thawing	AssayComplete™ Thawing Reagent T6	92-4106TR
Cell Freezing	AssayComplete™ Freezing Reagent F5	92-5105FR
Cell Detachment	AssayComplete™ 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. For additional information, please visit discoverx.com.