

KILR[®] HCT-116 Cell Pool Starter Pack

Catalog Number: 497-1040P047

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
97-1040P047	KILR [®] HCT-116 Cell Pool

Starter Pack Components

Product Use	Description	DiscoverX [®] Part #
Cell Culture	AssayComplete [™] Cell Culture Kit-102	92-3102G
Cell Plating	AssayComplete [™] Cell Plating 0 Reagent	93-0563R0A
Cell Thawing	AssayComplete [™] Thawing Reagent T8	92-4108TR
Cell Freezing	AssayComplete [™] Freezing Reagent F6	92-5106FR
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.