

KILR[®] SK-MEL-5 Cell Pool Starter Pack

Product Code: 497-1033P030

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
97-1033P030	KILR [®] SK-MEL-5 Cell Pool

Starter Pack Components

Product Use	Description	DiscoverX [®] Part #
Cell Culture	AssayComplete [™] Cell Culture Kit-103	92-3103G
Cell Plating	AssayComplete [™] Cell Plating 39 Reagent	93-0563R39A
Cell Thawing	AssayComplete [™] Thawing Reagent T3	92-4103TR
Cell Freezing	AssayComplete [™] Freezing Reagent F3	92-5103FR
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 www.discoverx.com.

