

**KILR<sup>®</sup> SKBR3 Cell Pool Starter Pack****Product Code:** 497-1002P018**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
97-1002P018	KILR <sup>®</sup> SKBR3 Cell Pool

**Starter Pack Contents**

Product Use	Description	DiscoverX <sup>®</sup> Part #	Configuration
Cell Culture	AssayComplete <sup>™</sup> Cell Culture Kit-102	92-3102G	Kit
Cell Plating	AssayComplete <sup>™</sup> Cell Plating 39 Reagent	93-0563R39A	100 mL
Cell Thawing	AssayComplete <sup>™</sup> Thawing Reagent T8	92-4108TR	100 mL
Cell Freezing	AssayComplete <sup>™</sup> Freezing Reagent F6	92-5106FR	2 x 50 mL
Cell Detachment	AssayComplete <sup>™</sup> Cell Detachment Reagent	92-0009	50 mL

**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 [www.discoverx.com](http://www.discoverx.com).