

## KILR<sup>®</sup> 4T1 Cell Pool Starter Pack

**Product Code:** 497-1028P041

### 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
97-1028P041	KILR <sup>®</sup> 4T1 Cell Pool

### Starter Pack Contents

Product Use	Description	DiscoverX <sup>®</sup> Part #	Configuration
Cell Culture	AssayComplete <sup>™</sup> Cell Culture Kit-101	92-3101G	Kit
Cell Plating	AssayComplete <sup>™</sup> Cell Plating 39 Reagent	93-0563R39A	100 mL
Cell Thawing	AssayComplete <sup>™</sup> Thawing Reagent T6	92-4106TR	100 mL
Cell Freezing	AssayComplete <sup>™</sup> Freezing Reagent F5	92-5105FR	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).