

**KILR<sup>®</sup> Hut78 Cell Pool Starter Pack****Product Code:** 497-1024P037**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
97-1024P037	KILR <sup>®</sup> Hut78 Cell Pool

**Starter Pack Contents**

Product Use	Description	DiscoverX <sup>®</sup> Part #	Configuration
Cell Culture	AssayComplete™ Cell Culture Kit-115	92-3115G	Kit
Cell Plating	AssayComplete™ Cell Plating 39 Reagent	93-0563R39A	100 mL
Cell Thawing	AssayComplete™ Thawing Reagent T10	92-4110TR	100 mL
Cell Freezing	AssayComplete™ Freezing Reagent F7	92-5107FR	2 x 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).