

## SPRINTer<sup>®</sup> HCT-116 c-Myc Protein Turnover Biosensor Cell Line Starter Pack

**Catalog Number:** 491-1003C047

### 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
91-1003C047	SPRINTer <sup>®</sup> HCT-116 c-Myc Protein Turnover Biosensor Cell Line

### 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 0 Reagent	93-0563R0A	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)

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
 Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
 General product information: [www.discoverx.com](http://www.discoverx.com)