

## SPRINTer™ BRAF(V600E) Protein Turnover Biosensor Cell Line (A375) Starter

Catalog Number: 491-1008C057

### 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   |
|-------------|---|
| 91-1008C057 | SPRINTer™ BRAF(V600E) Protein Turnover Biosensor Cell Line (A375) |

### Starter Pack Contents

| Product Use     | Description                            | DiscoverX® Part # | Configuration |
|-----------------|--|-------------------|---------------|
| Cell Culture    | AssayComplete™ Cell Culture Kit-105    | 92-3105G          | Kit           |
| Cell Plating    | AssayComplete™ Cell Plating 0 Reagent  | 93-0563R0A        | 100 mL        |
| Cell Thawing    | AssayComplete™ Thawing Reagent T1      | 92-4101TR         | 100 mL        |
| Cell Freezing   | AssayComplete™ Freezing Reagent F1     | 92-5101FR         | 2 x 50 mL     |
| Cell Detachment | AssayComplete™ 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)