

## PathHunter® U2OS TGFBR1/ACVR2B Dimerization Cell Line Starter Pack

**Product Code:** 493-1046C3

### 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   |
|-----------|---|
| 93-1046C3 | PathHunter® U2OS TGFBR1/ACVR2B Dimerization Cell Line |

### Starter Pack Contents

| Product Use     | Description                            | DiscoverX® Part # | Configuration |
|-----------------|--|-------------------|---------------|
| Cell Culture    | AssayComplete™ Cell Culture Kit-103    | 92-3103G          | Kit           |
| Cell Plating    | AssayComplete™ Cell Plating 5 Reagent  | 93-0563R5A        | 100 mL        |
| Cell Thawing    | AssayComplete™ Thawing Reagent T3      | 92-4103TR         | 100 mL        |
| Cell Freezing   | AssayComplete™ Freezing Reagent F3     | 92-5103FR         | 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)

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).