

PathHunter® HVEM Ligand Cell Line (U2OS) Starter Pack

Catalog Number: 493-1185C3

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 |
|-----------|--|
| 93-1185C3 | PathHunter® HVEM Ligand Cell Line (U2OS) |

Starter Pack Contents

| Product Use | Description | DiscoverX® Part # | Configuration |
|-----------------|--|-------------------|---------------|
| Cell Culture | AssayComplete™ Cell Culture Kit-103 | 92-3103G | Kit |
| Cell Plating | Not included with Ligand cell lines | N/A | 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)

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com