

PathHunter[®] CHO-K1 NUC-EA Parental Cell Line Starter Pack

Product Code: 493-0004

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 |
|---------|--|
| 93-0004 | PathHunter [®] CHO-K1 NUC-EA Parental Cell Line |

Starter Pack Contents

| Product Use | Description | DiscoverX [®] Part # | Configuration |
|-----------------|--|-------------------------------|---------------|
| Cell Culture | AssayComplete [™] Cell Culture Kit-107 | 92-3107G | Kit |
| Cell Thawing | AssayComplete [™] Thawing Reagent T2 | 92-4102TR | 100 mL |
| Cell Freezing | AssayComplete [™] Freezing Reagent F2 | 92-5102FR | 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. For additional information, please visit www.discoverx.com.