

PathHunter[®] EDG5 Activated GPCR Internalization U2OS Cell Line Starter Pack

Product Code: 493-0735C3

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
93-0735C3	PathHunter [®] EDG5 Activated GPCR Internalization U2OS 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 28 Reagent	93-0563R28A	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. For additional information, please visit www.discoverx.com.