

PathHunter[®] HEK 293 EDG3 β -Arrestin Cell Line Starter Pack

Product Code: 493-0217C1

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
93-0217C1	PathHunter [®] HEK 293 EDG3 β -Arrestin Cell Line

Starter Pack Contents

Product Use	Description	DiscoverX [®] Part #	Configuration
Cell Culture	AssayComplete [™] Cell Culture Kit-105	92-3105G	Kit
Cell Plating	AssayComplete [™] Cell Plating 7 Reagent	93-0563R7A	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)

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.