

InCELL Hunter™ Jurkat RIPK3 Kinase Cell Line Starter Pack

Catalog Number: 496-0077C19

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
96-0077C19	InCELL Hunter™ Jurkat RIPK3 Kinase Cell Line

Starter Pack Contents

Product Use	Description	DiscoverX® Part #	Configuration
Cell Culture	AssayComplete™ Cell Culture Kit-101	92-3101G	Kit
Cell Plating	AssayComplete™ Cell Plating 25 Reagent	93-0563R25A	100 mL
Cell Thawing	AssayComplete™ Thawing Reagent T6	92-4106TR	100 mL
Cell Freezing	AssayComplete™ Freezing Reagent F5	92-5105FR	2 x 50 mL

Storage And Shipping Information

Storage: Store at -20°C

Shipping: Frozen on Dry Ice (-70°C)

Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit <https://www.discoverx.com/legal-policies/> for limited-use label license agreement and trademark information for this product.

© 2024 Eurofins DiscoverX, a Eurofins Discovery Company. All Rights Reserved.