

PathHunter® TSLP Signaling Reporter Cell Line (REH) Starter Pack

Catalog Number: 493-1182C055

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 |
|-------------|---|
| 93-1182C055 | PathHunter® TSLP Signaling Reporter Cell Line (REH) |

Starter Pack Contents

| Product Use | Description | DiscoverX® Part # | Configuration |
|---------------|---------------------------------------|-------------------|---------------|
| Cell Culture | AssayComplete™ Cell Culture Kit-101 | 92-3101G | Kit |
| Cell Plating | AssayComplete™ Cell Plating 0 Reagent | 93-0563R0A | 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.