

## Recombinant Human Heregulin- $\beta$ 1

Heregulin  $\beta$ -1 is a signaling protein for the ErbB2/ErbB4 receptor heterodimer on the cardiac muscle cells, playing an important role in the heart structure and function by inducing the receptor phosphorylation and cardiomyocyte differentiation. Proteolytic processing of the extracellular domain of the transmembrane NRG-1 isoforms release soluble growth factors

### Materials Provided

<b>Catalog Number:</b>	<b>92-1031</b>
<b>Quantity Supplied:</b>	1 vial(s), 10 $\mu$ g per vial
<b>Reconstitution Buffer:</b>	1 mL

### Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

### Product Information

<b>Molecular Weight:</b>	7500 g/mol
<b>Source:</b>	E. coli
<b>Purity:</b>	$\geq$ 98% by SDS-PAGE and HPLC
<b>Endotoxin Level:</b>	< 0.1 ng/ $\mu$ g (<1EU/ $\mu$ g).
<b>Formulation:</b>	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
<b>Storage:</b>	$\leq$ -20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

<b>Stock Concentration:</b>	50 $\mu$ g/mL
<b>Reconstitution Volume:</b>	0.200 mL
<b>Reconstitution Solvent:</b>	Reconstitution Buffer

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)