

## Recombinant Human Neuregulin-1 (Heregulin-β2)

Neuregulin-1 is a glycoprotein that acts on EGFR family receptors, particularly ERBB receptor activation, increasing their tyrosine phosphorylation. Plays a critical role in growth and development, heart structure and function, and cardiomyocyte differentiation.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1091</b>
<b>Quantity Supplied:</b>	1 vial(s), 50 µg per vial

### 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>	7055 g/mol
<b>Source:</b>	Recombinant
<b>Purity:</b>	≥ 96% by RP-HPLC
<b>Endotoxin Level:</b>	< 0.1 ng/µg (<1EU/µg).
<b>Formulation:</b>	N/A
<b>Storage:</b>	≤ -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>	100 µg/mL
<b>Reconstitution Volume:</b>	0.5 mL
<b>Reconstitution Solvent:</b>	H <sub>2</sub> O

***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)