

## Recombinant Rat Beta-NGF

Beta-nerve growth factor (NGF) is a prototypical growth factor that functions in sympathetic and sensory neuron survival and maintenance. Mutations have been associated with hereditary sensory and autonomic neuropathy. Extracellular ligand for the NTRK1 and NGFR receptors.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1283</b>
<b>Quantity Supplied:</b>	1 vial(s), 100 µg 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>	13200 g/mol
<b>Source:</b>	Recombinant
<b>Purity:</b>	≥ 97%
<b>Endotoxin Level:</b>	<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>	1 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)