

## Recombinant Human b-NGF

$\beta$ -NGF is a potent neurotrophic factor that signals through its receptor  $\beta$ -NGFR, and plays a crucial role in the development and preservation of the sensory and sympathetic nervous systems.  $\beta$ -NGF also acts as a growth and differentiation factor for B lymphocytes and enhances C cell survival in the body.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1023</b>
<b>Quantity Supplied:</b>	1 vial(s), 20 $\mu$ 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>	27 kg/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>	0.1 mg/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)