

## Recombinant Human GDNF

Glial-derived neurotrophic factor (GDNF) is a disulfide-linked, homodimeric neurotrophic factor protein that promotes the survival and differentiation of dopaminergic neurons in culture and able to prevent apoptosis of motor neurons induced by axotomy. GDNF specifically promotes dopamine uptake and survival, and morphological differentiation of midbrain neurons.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1299</b>
<b>Quantity Supplied:</b>	1 vial(s), 2 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>	30400 g/mol
<b>Source:</b>	Recombinant
<b>Purity:</b>	≥ 98%
<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>	0.020 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)