

## Recombinant Human BDNF

BDNF is widely expressed in the central nervous system and acts in an autocrine and paracrine manner on several classes of neurons. BDNF has also been shown to promote neuronal survival and differentiation, and to play a critical role in memory formation and synaptic regulation of neural tissues.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1024</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>	27 kg/mol
<b>Source:</b>	E. coli
<b>Purity:</b>	≥ 98% by SDS-PAGE and HPLC
<b>Endotoxin Level:</b>	< 0.1 ng/µg (<1EU/µg).
<b>Formulation:</b>	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
<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>	0.1 mg/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)