

## Recombinant Human GDF-2 (BMP-9)

Growth differentiation factor 2 (GDF-2, BMP-9) is a member of the bone morphogenetic protein (BMP) family that plays a role in cholinergic CNS neuron differentiation and response to the neurotransmitter acetylcholine, as well as the regulation of embryonic and adult tissue cell growth and differentiation. Potent circulatory inhibitor of angiogenesis, signaling through the type I activin receptor kinase ACVRL1.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1249</b>
<b>Quantity Supplied:</b>	1 vial(s), 10 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>	24100 g/mol
<b>Source:</b>	Recombinant
<b>Purity:</b>	≥ 95%
<b>Endotoxin Level:</b>	N/A
<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>	0.1 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)