

## Recombinant Human PDGF-BB

Platelet-derived growth factor BB (PDGF-BB) is a homodimeric protein that is released upon platelet activation that induces proliferation, migration, and survival of endothelial, epithelial, and smooth muscle cells, as well as fibroblasts. Potent mitogen for mesenchymal cells. Able to bind all three of the PDGF receptors, and plays an important part in healing, regeneration, and angiogenesis and formation of smooth muscle during embryonic development. May facilitate tumor metastasis and vascularization.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1110</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>	24300 g/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>	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)