



6405 Mira Mesa Blvd, Ste 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

## PDGF-BB Recombinant

Catalog #: 90800  
Lot#: 230705

**Description:** Recombinant PDGF-BB is a disulfide-linked homodimeric protein consisting of a two 109 amino acid residue subunits, and migrates as an approximately 32 kDa protein under non-reducing and 16 kDa under reducing conditions.

**Source:** Optimized DNA sequence encoding human PDGF-BB mature chain was expressed in *Saccharomyces cerevisiae* cells.

**Formulation:** Lyophilized from a 0.2 µm filtered 0.02M acetic acid solution pH 6.5.

**Reconstitution:** A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/ml. This solution can then be diluted into other buffers.

**Storage:** The lyophilized protein is stable for at least 2 years from date of receipt when stored at -20°C. Upon reconstitution, store in working aliquots at 2 - 8° C for up to one month, or at -20°C for up to six months, in the presence of a carrier protein. Avoid repeated freeze/thaw cycles.

**Purity:** >98%, as determined by SDS-PAGE and HPLC.

**Endotoxin Level:** Endotoxin level was found to be < 0.1 ng/µg (1EU/µg), using the LAL gel clot method.

**Biological Activity:** Determined by the dose-dependent proliferation of BALB/c 3T3 cells to be 1.0-3.0 ng/ml, corresponding to a specific activity of  $\geq 1 \times 10^6$  IU/mg.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)