# PDGF-AB (human), (recombinant)

Platelet-derived growth factor (PDGF) is an important regulator of cell growth, proliferation, and angiogenesis. PDGF synthesis is induced by IL-1, IL-6, TNF-a, TGF-b and EGF signaling. PDGF functions as a mitogenic growth hormone on cells of mesenchymal lineage, such as smooth muscle and glial cells. PDGF is also stored in the alpha-granules of platelets and is released upon adherence to traumatized tissues. PDGF is a dimeric glycoprotein formed by two A chains (AA), two B chains (BB), or as a heterodimer with an A and a B chain (AB). The PDGF dimer binds the cell surface receptor tyrosine kinases PDGFR-a and PDGFR-b.

## **Ordering Information**

**Order Online** »

**ENZ-PRT258-0100** 100µg

Manuals, SDS & CofA

**View Online** »

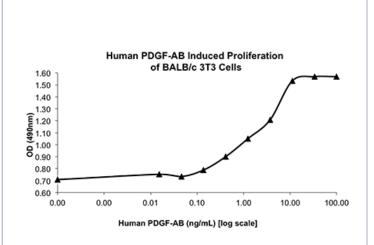


Figure 1: Serial dilutions of Human PDGF-AB, starting at 100 ng/mL, were added to 3T3 cells. Cell proliferation was measured after 46 hours and the linear portion of the curve can be used to calculate the ED50.

### **Handling & Storage**

**Handling** Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For

long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA

solution.

Long Term Storage -80°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Platelet-derived growth factor-AB, GDGF, ODGF

Endotoxin Content ≤1 EU/μg protein (LAL test)

Formulation Lyophilized from a sterile (0.2 micron) filtered aqueous

solution containing 0.1% Trifluoroacetic Acid (TFA).

**MW** ~Alpha14.4kDa, Beta 12.4/26.8kDa

Purity ≥95% (SDS-PAGE)

**Reconstitution** Reconstitute in sterile water at 0.1 mg/mL.

**Source** Produced in *E. coli*. Contains 126, 110/236 amino acids.

**UniProt ID** P04085/P01127

Last modified: May 29, 2024

