

# 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- $\alpha$ , TGF- $\beta$  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.

**Manuals, SDS & CofA**

[View Online »](#)

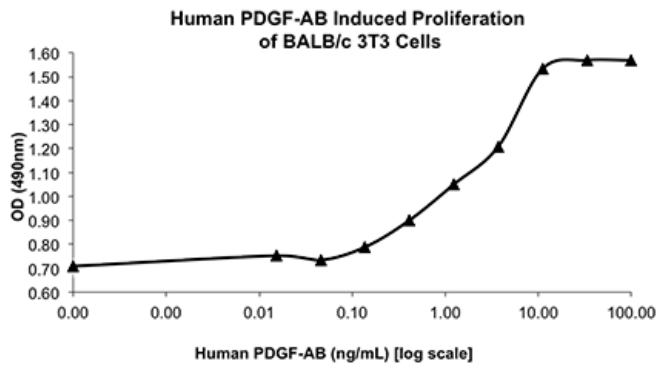


Figure 1: Serial dilutions of Human PDGF-AB, starting at 100 ng/mL, were added to 3T3 cells. Cell proliferation was measured after 46 hours and the linear portion of the curve can be used to calculate the ED50.

## Handling & Storage

<b>Handling</b>	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.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Platelet-derived growth factor-AB, GDGF, ODGF
<b>Endotoxin Content</b>	≤1 EU/μg protein (LAL test)
<b>Formulation</b>	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 0.1% Trifluoroacetic Acid (TFA).
<b>MW</b>	~Alpha14.4kDa, Beta 12.4/26.8kDa
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute in sterile water at 0.1 mg/mL.
<b>Source</b>	Produced in <i>E. coli</i> . Contains 126, 110/236 amino acids.
<b>UniProt ID</b>	P04085/P01127

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)