

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- α , TGF- β 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- α and PDGFR- β .

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------------|
| ENZ-PRT258-0100 | 100 μ g |
|-----------------|-------------|

Manuals, SDS & CofA

[View Online »](#)

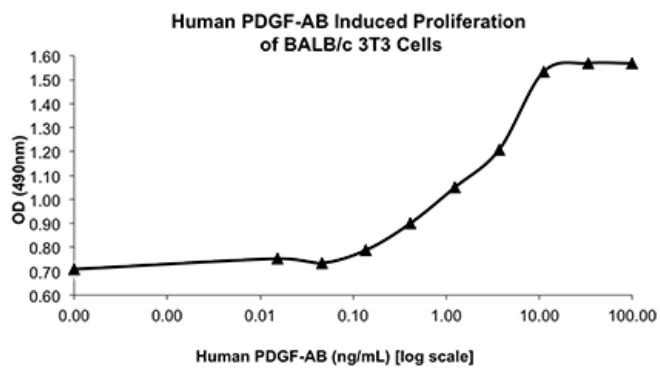


Figure 1: Serial dilutions of Human PDGF-AB, starting at 100 ng/mL, were added to 3T3 cells. Cell proliferation was measured after 46 hours and the linear portion of the curve can be used to calculate the ED50.

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 <i>E. coli</i> . Contains 126, 110/236 amino acids. |
| UniProt ID | P04085/P01127 |



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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