

Recombinant Human PDGFA

Recombinant human protein expressed in E. coli cells

Catalog # P841-40BN

Lot # G3223-3

Product Description

Recombinant Human PDGFA was expressed in E. coli cells. The protein accession number is [P04085](#)

Gene Aliases

PDGF subunit A, PDGF-1, Platelet-derived growth factor A chain, Platelet-derived growth factor alpha polypeptide, PDGFA, PDGF-AA, PDGF1, Recombinant Human Platelet-Derived Growth Factor Subunit A (PDGFA)

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

Formulation

Lyophilized from a 0.2 µm filtered solution in Tris and NaCl

Reconstitution Protocol

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.

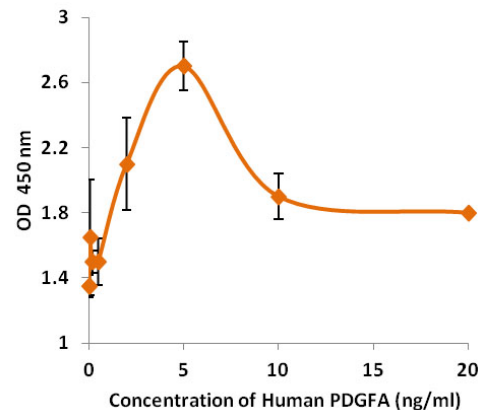
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

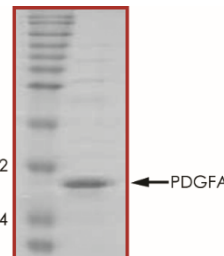
The PDGF family is comprised of five different disulphide-linked dimers of four different polypeptide chains: A, B, C and D (PDGF-AA, PDGF-BB, PDGF-AB, PDGF-CC and PDGF-DD). Synthesized mainly by megakaryocytes, PDGFs are stored in the alpha granules of platelets from which they are released following platelet activation. PDGF receptors are expressed in fibroblasts, osteoblasts, chondroblasts, smooth muscle cells, glial cells and endothelial cells. Functioning as an autocrine and paracrine growth factor, PDGFs are involved in a number of biological processes that include but not limited to hyperplasia, chemotaxis, embryonic neuron development, wound healing and respiratory tubule epithelial cell development. Aberrant expression of PDGFs is observed with vascular proliferative diseases such as atherosclerosis. PDGFs regulate the synthesis of their own receptor and also influence the expression of membrane receptors for IL1, EGF, 5-Hydroxytryptamine, LDL and transferrin. Recombinant Human PDGFA is a disulfide-linked homodimer of two A chains.

Specific Activity



The ED50 as determined by the dose-dependent proliferation of mouse embryonic fibroblast cells was found to be ≤ 5.0 ng/mL

Purity



Recombinant Human PDGFA resolved on a 17% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **14.5 kDa**

Recombinant Human PDGFA

Recombinant human protein expressed in E. coli cells

Catalog Number	P841-40BN
Specific Lot Number	G3223-3
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com

www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.