

Recombinant Human PlGF

Recombinant human protein expressed in E. coli cells

Catalog # P829-40N

Lot # G3223-6

Product Description

Recombinant Human PlGF was expressed in E. coli cells. The protein accession number is [P49763](#)

Gene Aliases

PGF, PGLF

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 Citrate, phosphate and NaCl (pH 4.0)

Reconstitution Protocol

A quick spin of the vial followed by reconstitution in PBS to a concentration not less than 0.1 mg/mL. This solution can then be diluted into other buffers.

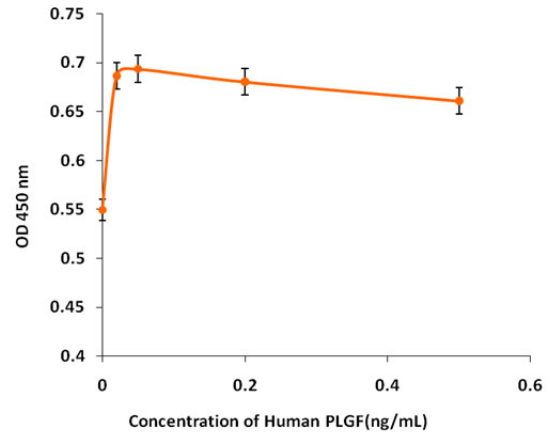
Storage and Stability

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.

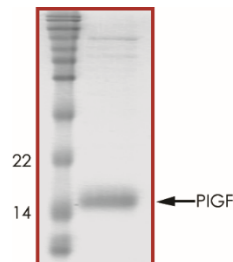
Scientific Background

Human PlGF, or Placental Growth Factor, is part of the PDGF/VEGF family and contains the 8 cysteine residues conserved in the family. It can be alternately spliced into three different mature forms. PGF is typically found as a disulfide linked dimer which is highly glycosylated. It is a key molecule involved in vasculogenesis and angiogenesis, especially during embryogenesis. It also induces the activation of monocytes, as well as their migration and production of inflammatory cytokines. It is primarily expressed by the placental trophoblast during pregnancy, but is also expressed in villous trophoblasts, decidual cells, erythroblasts, keratinocytes, and somewhat in endothelial cells. PGF signals by binding VEGF R1, and competes with VEGF for binding. Recombinant PlGF is a 15.0 kDa protein.

Specific Activity



Purity



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

Approx. MW **15.0 kDa**

Recombinant Human PlGF

Recombinant human protein expressed in E. coli cells

Catalog Number	P829-40N
Specific Lot Number	G3223-6
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.