

PRODUCT INFORMATION



PDGF-B (human, recombinant)

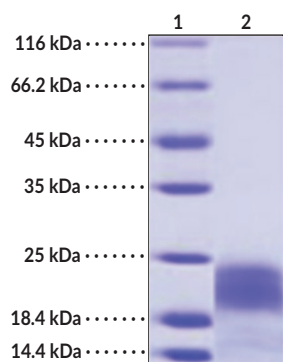
Item No. 44191

Overview and Properties

Synonyms:	Becaplermin, IBGC5, PDGF-2, PDGFB, Platelet-Derived Growth Factor B Chain, Platelet-Derived Growth Factor β Polypeptide, Proto-oncogene c-Sis, SIS, SSV
Source:	Active recombinant human N-terminal His-tagged PDGF-B expressed in yeast
Amino Acids:	82-190
Uniprot No.:	P01127
Molecular Weight:	14.3 kDa
Storage:	-80°C (as supplied)
Stability:	≥ 1 year
Purity:	$\geq 90\%$ estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile 30% acetonitrile and 1/1,000 trifluoroacetic acid
Endotoxin Testing:	< 1.0 EU/ μ g, determined by the LAL endotoxin assay
Concentration:	<i>batch specific</i> mg/ml
Activity:	<i>batch specific</i> U/ml
Specific Activity:	<i>batch specific</i> U/mg
Bioactivity:	See figures for details

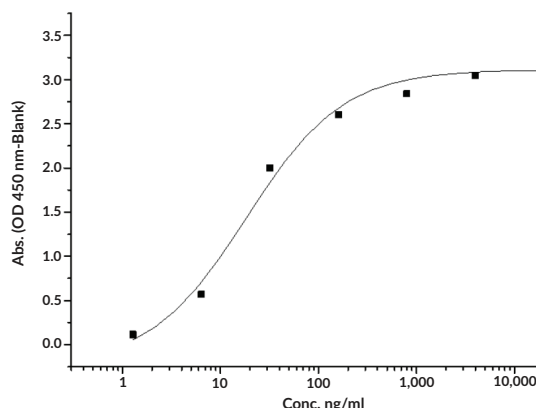
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: PDGF-B

SDS-PAGE Analysis of PDGF-B. This protein has a calculated molecular weight of 14.3 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 20 kDa due to glycosylation.



Immobilized recombinant human PDGFR β at 2 μ g/ml (100 μ l/well) binds to PDGF-B with an EC₅₀ value of 10-30 ng/ml.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/09/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

PDGF-B is a secreted growth factor and member of the PDGF family.¹ It is composed of a signal peptide, a pro-peptide sequence, a growth factor domain containing a cystine knot, and a C-terminal tail. PDGF-B is expressed at high levels in the heart, placenta, and fetal kidney and in the brain in the substantia nigra.² It exists as either a homodimer, PDGF-BB, which can activate PDGFR α or PDGFR β , or a heterodimer with PDGF-A, PDGF-AB, which can only activate PDGFR α , and induces signaling through multiple pathways, including ERK/MAPK, PI3K/Akt/mTOR, PLC/PKC, and JAK/STAT.^{1,3,4} PDGF-B is a potent mitogen for cells of mesenchymal origin and is necessary for vascular development and vascular smooth muscle cell proliferation in the CNS, lungs, heart, placenta, and skin.^{4,5} A fusion between the genes encoding PDGF-B and collagen 1A1, *PDGFB* and *COL1A1*, leads to overproduction of PDGF-B and is associated with various benign and malignant tumors, including dermatofibrosarcoma protuberans.⁴ Formulations containing PDGF-B have been used in the treatment of diabetic neuropathic foot ulcers. Cayman's PDGF-B (human, recombinant) protein can be used for binding assays. This protein consists of 124 amino acids and has a calculated molecular weight of 14.3 kDa.

References

1. Chen, P.-H., Chen, X., and He, X. Platelet-derived growth factors and their receptors: Structural and functional perspectives. *Biochim. Biophys. Acta* **1834**(10), 2176-2186 (2013).
2. LaRochelle, W.J., Jeffers, M., McDonald, W.F., *et al.* PDGF-D, a new protease-activated growth factor. *Nat. Cell Biol.* **3**(5), 517-521 (2001).
3. Heldin, C.-H. and Lennartsson, J. Structural and functional properties of platelet-derived growth factor and stem cell factor receptors. *Cold Spring Harb. Perspect. Biol.* **5**(8), a009100 (2013).
4. Roskoski, R., Jr. The role of small molecule platelet-derived growth factor receptor (PDGFR) inhibitors in the treatment of neoplastic disorders. *Pharmacol. Res.* **129**, 65-83 (2018).
5. Gaengel, K., Genové, G., Armulik, A., *et al.* Endothelial-mural cell signaling in vascular development and angiogenesis. *Arterioscler. Thromb. Vasc. Biol.* **29**(5), 630-638 (2009).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM