

PDGF receptor β monoclonal antibody (42G12)

Mitogenic platelet derived growth factor (PDGF) signaling is mediated by the PDGF tyrosine kinase receptors PDGFR α and PDGFR β , which exist as homo- or heterodimer isoforms ($\alpha\alpha$, $\beta\beta$, $\alpha\beta$). Secretion of PDGFs and expression of PDGFRs in a broad distribution of cell types contributes to regulation of cellular processes including cell growth, actin reorganization, prevention of apoptosis, and chemotaxis. Structurally, the transmembrane PDGFRs are composed of an N-terminal extracellular domain with five immunoglobulin repeats, and an intracellular domain containing two spaced tyrosine kinase domains. PDGF ligands display preferential affinity to PDGFR α and PDGFR β , with PDGFR β activation restricted to binding by PDGF BB and PDGF DD. Growth factor binding initiates receptor dimerization and activates the cytosolic tyrosine kinase domains.

This antibody is covered by our [Worry-Free Guarantee](#).

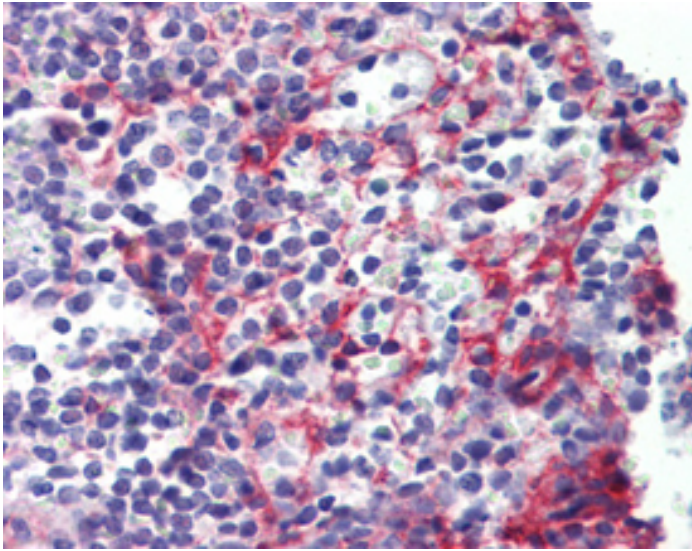
Ordering Information

[Order Online »](#)

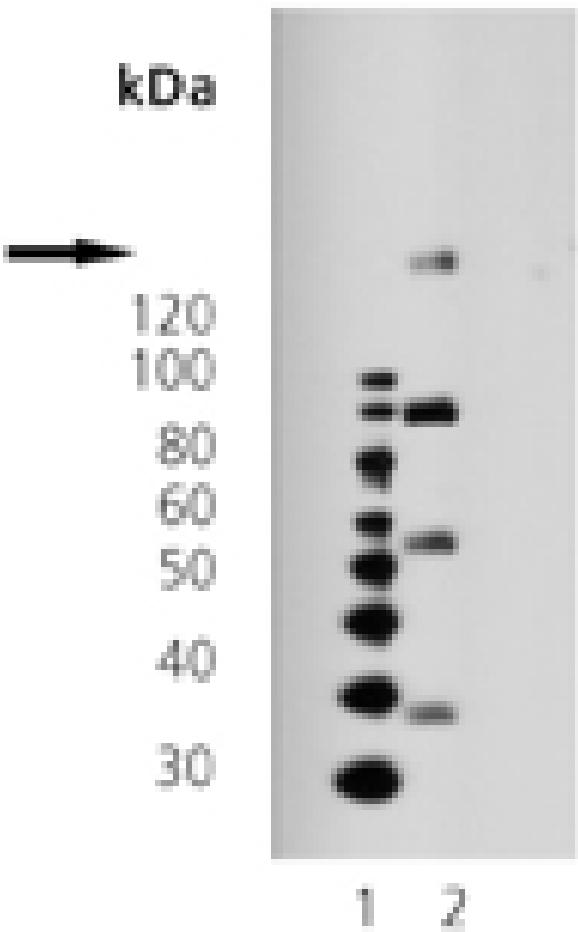
ADI-905-682-100	100 μ g
-----------------	-------------

Manuals, SDS & CofA

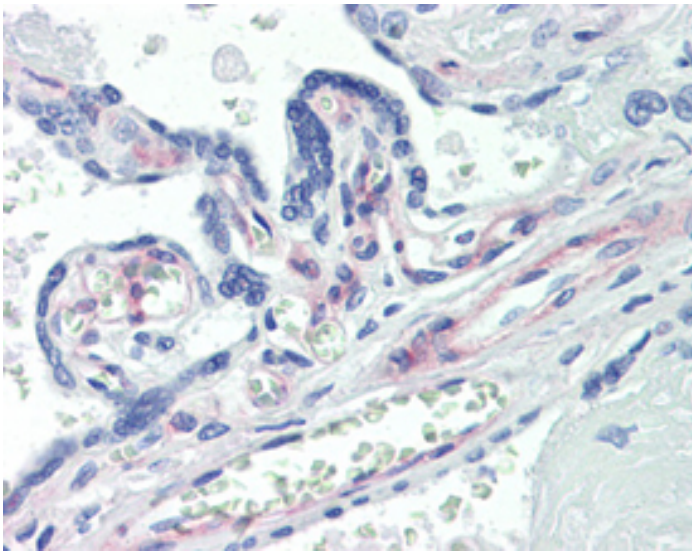
[View Online »](#)



IHC Analysis: Human spleen tissue stained with PDGF receptor β , mAb (42G12) at 20 μ g/ml.



Western blot analysis of MW marker (1) and mouse heart muscle lysate (2) probed with PDGF Receptor β mAb (42G12) (extra bands are secondary reactivity).



IHC Analysis: Human placenta tissue stained with PDGF receptor β , mAb (42G12) at 20 μ g/ml.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PDGFR, platelet-derived growth factor receptor
Application	IHC (PS), WB
Application Notes	Detects a band of ~125kDa by Western blot.
Clone	42G12
Crossreactivity	Weakly cross-reacts with human PDGF receptor β.
Formulation	Liquid. In PBS, pH 7.2, containing 0.09% sodium azide and 50% glycerol.
Host	Mouse
Immunogen	Recombinant human PDGF Receptor β.
Isotype	IgG1
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from mouse ascites.
Species Reactivity	Mouse, Rat
UniProt ID	P09619

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

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)