

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](#).

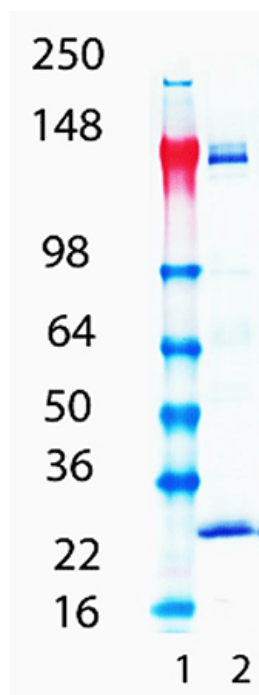
Ordering Information

[Order Online »](#)

ENZ-ABS688-0100	100 μ g
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of PDGF receptor β mAb (42G12) (Prod. No. ENZ-ABS688), Lane 1: MW marker, Lane 2: Rat Heart Cell Lysate.

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 Western Blot (1 μ g/ml, ECL) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Mouse, Rat

UniProt ID

P09619



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-fr@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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.