

# PDGF receptor $\beta$ monoclonal antibody (42G12)

Mitogenic platelet derived growth factor (PDGF) signaling is mediated by the PDGF tyrosine kinase receptors PDGFR $\alpha$  and PDGFR $\beta$ , 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 $\alpha$  and PDGFR $\beta$ , with PDGFR $\beta$  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](#).

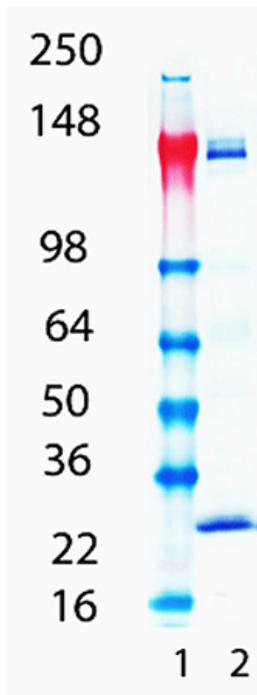
## Ordering Information

[Order Online »](#)

|                 |             |
|-----------------|-------------|
| ENZ-ABS688-0100 | 100 $\mu$ g |
|-----------------|-------------|

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of PDGF receptor  $\beta^2$  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  $\beta$ .

**Formulation** Liquid. In PBS, pH 7.2, containing 0.09% sodium azide and 50% glycerol.

**Host** Mouse

**Immunogen** Recombinant human PDGF Receptor  $\beta$ .

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (1 $\mu$ 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

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)