

Anti-PDGFB

Rabbit Polyclonal Antibody

Catalog # P842-63R

Lot # O2121-55

Cited Applications

WB, ELISA

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the PDGFB protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Rabbit polyclonal PDGFB antibody was raised against a 19 amino acid peptide near the carboxy terminus of human PDGFB.

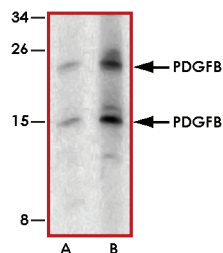
Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

Sample Data



Western blot analysis of PDGFB in rat liver tissue lysate with PDGFB antibody at (A) 1 and (B) 2 µg/ml.

Anti-PDGFB

Rabbit Polyclonal Antibody

Catalog Number

P842-63R

Specific Lot Number

O2121-55

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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.