

## Anti-PTPN11 (SHP2)

Rabbit Polyclonal Antibody

**Catalog # P38-63CR**

Lot # O2109-58

### Cited Applications

WB, ELISA, IF

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

### Specificity

Recognizes the PTPN11 protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

PTPN11 antibody was raised against a 14 amino acid synthetic peptide from near the carboxy terminus of human PTPN11.

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

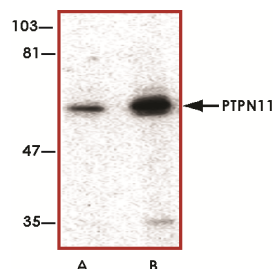
### Scientific Background

Mammalian PTPases can be subdivided into 1 of 2 broad categories: transmembrane receptor PTPases and intracellular PTPases. PTPN11 is one of the 2 closely related mammalian intracellular PTPases whose sequences encode 2 tandem SRC homology 2 (SH2) domains that are located at the amino-terminal side of a single PTPase catalytic domain (1). This PTP is widely expressed in most tissues and plays a regulatory role in various cell signaling events that are important for a diversity of cell functions, such as mitogenic activation, metabolic control, transcription regulation, and cell migration (2)

### References

1. Dechert, U. et al: Protein-tyrosine phosphatase SH-PTP2 (PTPN11) is localized to 12q24.1-24.3. Hum. Genet. 96: 609-615, 1995.
2. Ahmad, S. et al: A widely expressed human protein-tyrosine phosphatase containing src homology 2 domains. Proc. Nat. Acad. Sci. 90: 2197-2201, 1993.

### Sample Data



Western blot analysis of PTPN11 in mouse skeletal muscle tissue lysate with PTPN11 antibody at (A) 0.5 and (B) 1 ug/ml.

## Anti-PTPN11 (SHP2)

Rabbit Polyclonal Antibody

Catalog Number

P38-63CR

Specific Lot Number

O2109-58

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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**