

## Anti- SH2D1A (SAP)

Rabbit Polyclonal Antibody

**Catalog # S03-63R**

Lot # M132-9

### Cited Applications

- Western blot (1:1000)

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

### Specificity

Recognizes the SH2D1A (SAP) protein

### Cross Reactivity

- Western blot human SH2D1A (SAP)

SH2D1A (SAP) from other species may also be detectable

### Host

Rabbit

### Immunogen

Synthetic peptide corresponding to amino acids 113-128 of human SH2D1A (SAP)

### Formulation

TBS, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. 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.

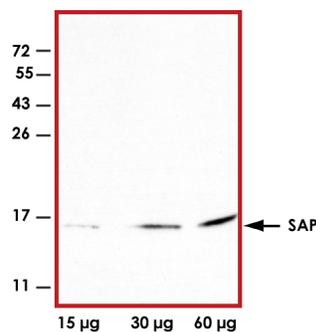
### Scientific Background

SH2D1A contains an SH2 domain and a short tail, and expresses at a high level in thymus and lung. It acts as an inhibitor of the SLAM (signaling lymphocyte-activation molecule) proteins by blocking the recruitment of the SH2-domain-containing signal-transduction molecule SHP-2 to its docking site. It modifies signal transduction pathways in activated T, B and NK cells. SH2D1A mediates interaction between FYN and SLAMF1 (1). Mutations in SH2D1A are a cause of lymphoproliferative syndrome X-linked type 1 (XLP1) (2).

### References

1. Chan B., et al: SAP couples Fyn to SLAM immune receptors. *Nat. Cell Biol.* 5:155-160(2003)
2. Dutz, J. P., et al: Lymphocytic vasculitis in X-linked lymphoproliferative disease. *Blood* 97: 95-100, 2001.

### Sample Data



Representative Western blot with Anti-SH2D1A (SAP) (1:1000) using 15 µg, 30 µg and 60 µg of Jurkat cell lysate.

## Anti-SH2D1A (SAP)

Rabbit Polyclonal Antibody

Catalog Number

S03-63R

Specific Lot Number

M132-9

Purification

Affinity chromatography

Concentration

1.0 µg/µL

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.**