

## Anti-STAT5

Rabbit Polyclonal Antibody

**Catalog # S56-63R**

Lot # H257-3

### Cited Applications

- Western blot (1:250)

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

### Specificity

Recognizes the STAT5 protein

### Cross Reactivity

- Western blot of human STAT5

*STAT5 from other species may also be detectable*

### Host

Rabbit

### Immunogen

Protein Code: P42229

### 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.

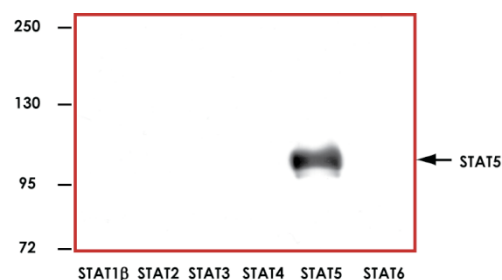
### Scientific Background

STAT5 (signal transducers and activators of transcription 5) is a member of the STAT family and this protein is phosphorylated by receptor associated protein kinases in response to cytokines and growth factors (1). The differentiation of T helper (Th) cells is regulated by members of the STAT family of signaling molecules. In BCR-ABL-positive cells, STAT5 is constitutively activated by tyrosine phosphorylation. STAT5 activation results in upregulation of Bcl-XL and increased resistance to induction of apoptosis. STAT5 is involved in expression and growth hormone-mediated sexually dimorphic regulation of cytochrome P450 3A10/lithocholic acid 6 beta-hydroxylase (2).

### References

1. Wang, D. et al: Naturally occurring dominant negative variants of Stat5. *Molec. Cell Biol.* 16: 6141-6148, 1996.
2. A Subramanian et al: STAT 5 and NF-Y are involved in expression and growth hormone-mediated sexually dimorphic regulation of cytochrome P450 3A10/lithocholic acid 6beta-hydroxylase. *Nucleic Acids Res.* 1998; 26(9): 2173-2178.

### Sample Data



Representative western blot with Anti-STAT5 (1:250) using 20 ng of GST tagged human recombinant STAT1β, STAT2, STAT3, STAT4, STAT5, and STAT6.

## Anti-STAT5

Rabbit Polyclonal Antibody

Catalog Number

S56-63R

Specific Lot Number

H257-3

Purification

Affinity chromatography

Concentration

0.25 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.**