

Anti-STAT3

Rabbit Polyclonal Antibody

Catalog # S54-63R

Lot # H257-2

Cited Applications

- Western blot (1:250)

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

Specificity

Recognizes the STAT3 protein

Cross Reactivity

- Western blot of human STAT3

STAT3 from other species may also be detectable

Host

Rabbit

Immunogen

Protein Code: P40763

Formulation

TBS, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) 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.

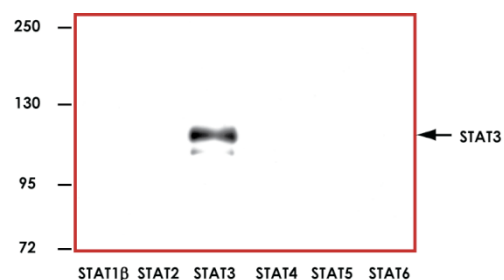
Scientific Background

STAT3 is a member of the signal transducers and activators of transcription (STAT) family of proteins that carry out a dual function: signal transduction and activation of transcription. STAT3 is widely expressed and becomes activated through phosphorylation on tyrosine as a DNA binding protein in response to a variety of stimuli such as EGF, IL-6, PDGF, IL-2 and G-CSF (1). This phosphoprotein forms homodimers as well as heterodimers with other members of the STAT family and translocate to the nucleus in order to modulate the transcription of various genes (2).

References

1. Zhong, Z. et al: Stat3: a STAT family member activated by tyrosine phosphorylation in response to epidermal growth factor and interleukin-6. *Science*. 1994 Apr 1;264(5155):95-8.
2. Tian, S. S. et al: Rapid activation of the STAT3 transcription factor by granulocyte colony-stimulating factor. *Blood*. 1994 Sep 15;84(6):1760

Sample Data



Representative western blot with Anti-STAT3 (1:250) using 20 ng of GST tagged human recombinant STAT1 β , STAT2, STAT3, STAT4, STAT5, and STAT6.

Anti-STAT3

Rabbit Polyclonal Antibody

Catalog Number

S54-63R

Specific Lot Number

H257-2

Purification

Affinity chromatography

Concentration

0.25 μ g/ μ L

Stability

1 yr 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.