

Anti-phospho-STAT3 (Ser727)

Rabbit Polyclonal Antibody

Catalog # S54-365BR

Lot # J1178-34

Cited Applications

WB, IHC

Suggested Dilutions:

WB: 1:500-1:1000 IHC: 1:50-1:100

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

Specificity

Recognizes the GS3 beta protein phosphorylated at serine 727

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser727

Formulation

PBS (pH 7.4) 150mM NaCl, 0.02% sodium azide and 50% glycerol.

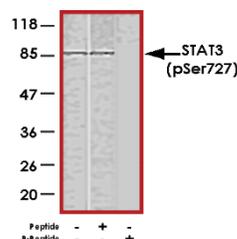
Scientific Background

STAT3 is a member of the signal transducers and activators of transcription (STAT) family of proteins that carry out dual functions. STAT3 is a DNA binding protein that is widely expressed and becomes activated upon phosphorylation on Tyr708, which leads to dimerization, intracellular translocation and DNA-binding. Phosphorylation at Ser727 regulates transcriptional activity. STAT3 is activated in response to a variety of stimuli, which includes EGF, IL-6, PDGF, IL-2 and G-CSF (1). STAT3 forms homo- and heterodimers with other STAT proteins to modulate the transcription of various genes (2).

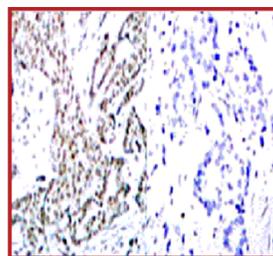
References

- 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.
- 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



Western blots from HeLa cells using anti-phospho-STAT3 (Ser727) antibody.



P-Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using anti-phospho-STAT3 (Ser727) antibody.

Anti-phospho-STAT3 (Ser727)

Rabbit Polyclonal Antibody

Catalog Number	S54-365BR
Specific Lot Number	J1178-34
Purification	Affinity chromatography
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.