

Anti-phospho-STAT5A (Ser780)

Rabbit Polyclonal Antibody

Catalog # S56-365R

Lot # J1178-22

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 MEK1/2 protein phosphorylated at serine 780

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser780

Formulation

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

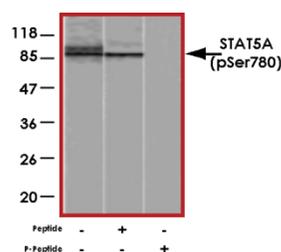
Scientific Background

Signal transducers and activators of transcription (STAT)5A is phosphorylated by receptor associated protein kinases in response to cytokines and growth factors (1). STAT proteins are responsible for the differentiation of helper T cells, thus playing key roles in regulating immune functions. In BCR-ABL-positive cells, STAT5 is constitutively activated by tyrosine phosphorylation. STAT5 activation results in upregulation of Bcl-XL and results in increased resistance to apoptosis. STAT5 is involved in expression and growth hormone-mediated sexually dimorphic regulation of cytochrome P450 3A10/lithocholic acid 6 beta-hydroxylase (2). STAT5A is phosphorylated at Ser780 in response to IL-2 stimulation (3).

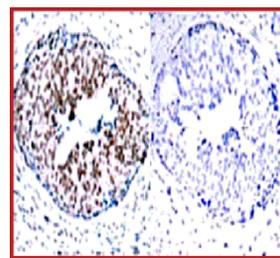
References

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- κ B 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.
3. Xue H H. et al: Serine phosphorylation of Stat5 proteins in lymphocytes stimulated with IL-2. *Int Immunol.* 2002 Nov;14(11):1263-71.

Sample Data



Western blot analysis of extracts from HeLa cells using anti-phospho-STAT5A (Ser780) antibody.



P-Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using anti-phospho-STAT5A (Ser780) antibody.

Anti-phospho-STAT5A (Ser780)

Rabbit Polyclonal Antibody

Catalog Number

S56-365R

Specific Lot Number

J1178-22

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

www.signalchem.com

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