

Anti-phospho-JAK1 (Tyr1022)

Rabbit Polyclonal Antibody

Catalog # J01-65R

Lot # J1178-12

Cited Applications

IHC

Suggested Dilutions:

IHC: 1:50-1:100

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

Specificity

Recognizes the JAK1 protein phosphorylated at tyrosine 1022

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr1022

Formulation

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

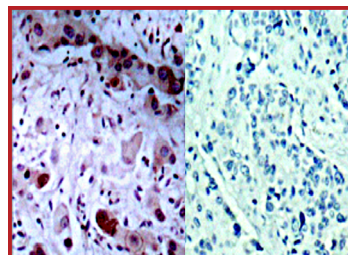
Scientific Background

JAK1 is a member of the Janus family of tyrosine kinases which are involved in cytokine receptor-mediated intracellular signal transduction (1). Upon stimulation by cytokines, JAK1 undergoes autophosphorylation. In turn, activated JAK1 is responsible for phosphorylating several signaling proteins, including STATs. Phosphorylated STATs dimerize and translocate to the nucleus where they regulate transcription of genes involved in growth, survival and differentiation. Phosphorylation of JAK1 at Tyr1022 and Tyr1023 is associated with cells that are undergoing proliferation (2).

References

1. Leonard, W J. et al : Jaks and STATs: biological implications. Annu Rev Immunol. 1998;16:293-322.
2. Trenary, M K. et al : JAK/STAT signaling and human in vitro myogenesis. BMC Physiol. 2011 Mar 9;11:6.

Sample Data



P-Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using anti-phospho-JAK1 (Tyr1022) antibody.

Anti-phospho-JAK1 (Tyr1022)

Rabbit Polyclonal Antibody

Catalog Number

J01-65R

Specific Lot Number

J1178-12

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.