

Anti-JAK2

Rabbit Polyclonal Antibody

Catalog # J02-63R

Lot # H283-4

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the JAK2 protein

Cross Reactivity

- Western blot JAK2 from human and mouse cells

JAK2 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 1005-1009 of human JAK2

Formulation

PBS pH 7.4, 0.02% sodium azide, 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.

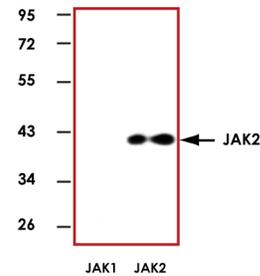
Scientific Background

JAK2 is a member of intracellular non-receptor tyrosine kinases that transduce cytokine-mediated signals via the JAK-STAT pathway (1). JAK2 has two near-identical phosphate-transferring domains. One domain exhibits the kinase activity while the other stabilizes the JAK conformational structure. JAK2 is the predominant JAK kinase activated in response to several growth factors and cytokines such as IL-3, GM-CSF and erythropoietin. JAK2 has been found to be constitutively associated with the prolactin receptor and is required for responses to gamma interferon (2).

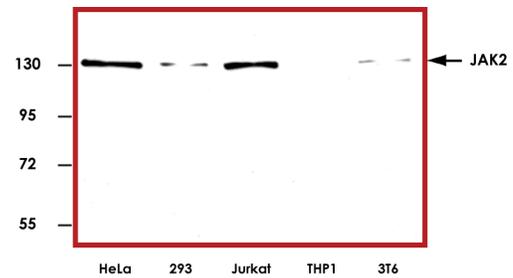
References

1. Lacronique, V. Et al : A TEL-JAK2 fusion protein with constitutive kinase activity in human leukemia. *Science*, 1997; 278: 1309-1312.
2. Ho, J.M. et al: TEL-JAK2 constitutively activates the extracellular signal-regulated kinase (ERK), stress-activated protein/Jun kinase (SAPK/JNK), and p38 signaling pathways. *Blood*, 2002; 100: 1438-1448.

Sample Data



Representative western blot with Anti-JAK2 (1:1000) using 20 ng of human recombinant JAK1 and JAK2.



Representative western blot with Anti-JAK2 (1:1000) using 30 µg of HeLa, HEK-293T, Jurkat, THP1, and 3T6 cell lysate.

Anti-JAK2

Rabbit Polyclonal Antibody

Catalog Number

J02-63R

Specific Lot Number

H283-4

Purification

Affinity chromatography

Concentration

1.0 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
www.signalchem.com