

Anti-TYK2 (JTK1)

Rabbit Polyclonal Antibody

Catalog # T21-63R

Lot # O2121-52

Cited Applications

WB, ELISA

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

Specificity

Recognizes the TYK2 (JTK1) protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Rabbit polyclonal TYK2 antibody was raised against a 17 amino acid peptide near the amino terminus of human TYK2

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

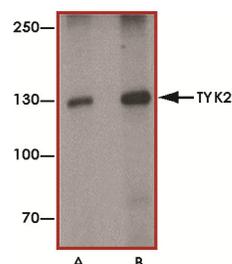
Scientific Background

TYK2 is a member of the tyrosine kinase and, more specifically, the Janus kinases (JAKs) protein families which associates with the cytoplasmic domain of type I and type II cytokine receptors and promulgate cytokine signals by phosphorylating receptor subunits. It is also component of both the type I and type III interferon signaling pathways. TYK2 play a role in anti-viral immunity and TYK2 plays an indispensable role in controlling responses to multiple cytokines in humans (1). Tyk2 contributes selectively to signals triggered by various biologic stimuli and cytokine receptors and TYK2 is an important regulator of lymphoid tumor surveillance (2).

References

1. Minegishi, Y. et.al: Human tyrosine kinase 2 deficiency reveals its requisite roles in multiple cytokine signals involved in innate and acquired immunity. *Immunity* 25: 745-755, 2006.
2. Stoiber, D. et.al: TYK2 is a key regulator of the surveillance of B lymphoid tumors. *J. Clin. Invest.* 114: 1650-1658, 2004.

Sample Data



Western blot analysis of TYK2 in human small intestine tissue lysate with TYK2 antibody at (A) 1 and (B) 2 µg/ml.

Anti-TYK2 (JTK1)

Rabbit Polyclonal Antibody

Catalog Number

T21-63R

Specific Lot Number

O2121-52

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.