

Anti-IL-21

Rabbit Polyclonal Antibody

Catalog # 1881-463R

Lot # Z2014-13

Cited Applications

WB, ELISA, ICC

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

Specificity

Recognizes the IL-21 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Anti-IL-21 antibody was raised against a synthetic peptide corresponding to 15 amino acids near the center of human IL-21 precursor.

Formulation

PBS containing 0.02% sodium azide.

Stability

1yr at -20°C from date of shipment

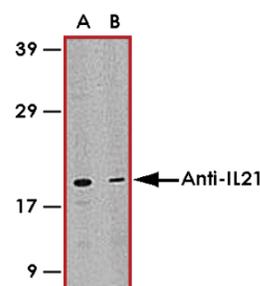
Scientific Background

A novel cytokine was recently identified in human and mouse and designated IL-21, which has significant homology to IL-2, IL-4, and IL-15. The receptor for IL-21 (IL-21R, also termed NILR for novel Interleukin receptor) is a member of the class I cytokine receptor family. IL-21R forms a complex with the common cytokine receptor gamma chain, gamma c, and mediates IL-21 signaling. IL-21 and its receptor activate JAK-STAT signaling pathway. IL-21 is expressed in activated T cells, and HL-60 and THP-1 cell lines. IL-21 plays a role in the proliferation and maturation of NK, B and T cell populations.

References

1. Parrish-Novak J, Dillon SR, Nelson A, et al. Interleukin 21 and its receptor are involved in NK cell expansion and regulation of lymphocyte function. *Nature*. 2000; 408(6808):57-63.
2. Asao H, Okuyama C, Kumaki S, et al. the common γ -chain is an indispensable subunit of the IL-21 receptor complex. *J Immunol*. 2001;167(1):1-5.

Sample Data



Western blot analysis of IL-21 expression in HL-60 cell lysate in the absence (lane A) or presence of blocking peptide (lane B) with Anti-IL-21 antibody at 1 µg/ml.



Immunocytochemistry of IL-21 in K562 cells with Anti-IL-21 antibody at 2 µg/mL.

Anti-IL-21

Rabbit Polyclonal Antibody

Catalog # **1881-463R**

Lot # **Z2014-13**

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.