

## Anti-IRAK4

Rabbit Polyclonal Antibody

**Catalog # I12-63R**

Lot # O2109-23

### Cited Applications

WB, ELISA, ICC, IF

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

### Specificity

Recognizes the IRAK4 protein and has no cross response to other IRAKs

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

IRAK4 antibody was raised against a synthetic peptide corresponding to amino acids at the carboxy terminus of human IRAK4

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

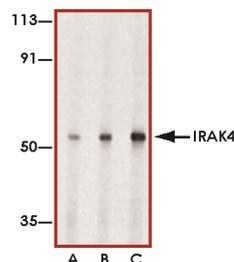
### Scientific Background

Interleukin-1 receptor-associated kinase 4 (IRAK4) is an important mediator in the signal transduction of Toll-like receptor (TLR) and IL1R family members (1). IRAK4 is involved in the Toll-like receptor signaling pathway leading to Apoptosis. IRAK4 has molecular functions like protein binding, ATP binding, kinase activity and magnesium ion binding. Toll/IL-1 receptor family members like IRAK4 are central components of host defense mechanisms in a variety of species (2). One well conserved element in their signal transduction is the Ser/Thr kinase activity which couple early signaling events in a receptor complex at the plasma membrane to larger signalosomes in the cytosol.

### References

1. Li, S. et al: IRAK-4: a novel member of the IRAK family with the properties of an IRAK-kinase. Proc. Natl. Acad. Sci. USA. 2002; 99:5567-72.
2. Suzuki, N. et al: IRAK-4 as the central TIR signaling mediator in innate immunity. Trends Immunol. 2002; 23:503-6.

### Sample Data



Western blot analysis of IRAK4 in HeLa cell lysate with IRAK4 antibody at 1 (lane 1), 2 (lane 2), and 4 (lane 3) µg/ml, respectively.



Immunocytochemistry of IRAK4 in K562 cells with IRAK4 antibody at 10 µg/ml.

## Anti-IRAK4

Rabbit Polyclonal Antibody

Catalog Number

I12-63R

Specific Lot Number

O2109-23

Purification  
Stability  
Storage & Shipping

Affinity chromatography  
1yr at -20°C from date of shipment  
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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