

## Anti-IRAK1

Rabbit Polyclonal Antibody

**Catalog # 109-63R**

Lot # O2109-1

### Cited Applications

WB, ELISA, ICC, IF, IP

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

### Specificity

Recognizes the IRAK1 protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

IRAK1 antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human IRAK1

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

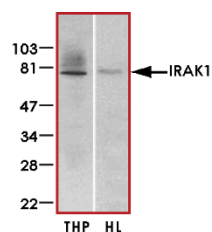
### Scientific Background

IRAK1 encodes the interleukin-1 receptor-associated kinase 1, one of two putative serine/threonine kinases that become associated with the interleukin-1 receptor (IL1R) upon stimulation. The exposure of HeLa cells or human embryonic kidney cells overexpressing IL1R to IL1 caused rapid association of IRAK with the IL1R complex and phosphorylation of IRAK (1). IRAK1 is partially responsible for IL1-induced upregulation of the transcription factor NF-kappa B. IRAK1 as a risk gene with critical role in the pathogenesis of systemic lupus erythematosus (2).

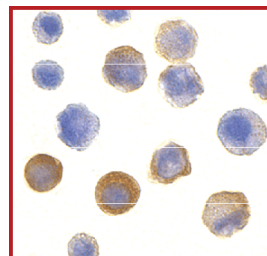
### References

1. Cao, Z. et al: IRAK: a kinase associated with the interleukin-1 receptor. *Science* 271: 1128-1131, 1996.
2. Jacob, C. O. et al: Identification of IRAK1 as a risk gene with critical role in the pathogenesis of systemic lupus erythematosus. *Proc. Nat. Acad. Sci.* 106: 6256-6261, 2009.

### Sample Data



Western blot analysis of IRAK1 in THP-1 (THP) and HeLa (HL) whole cell lysates with IRAK antibody at 1:2000 dilution.



Immunocytochemistry of IRAK1 in HeLa cells with IRAK1 antibody at 10 µg/ml.

## Anti-IRAK1

Rabbit Polyclonal Antibody

Catalog Number

**109-63R**

Specific Lot Number

**O2109-1**

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](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.**