

## Anti-IRAK2

Rabbit Polyclonal Antibody

**Catalog # I10-63BR**

Lot # O2109-43

### Cited Applications

WB, ELISA, ICC, IF

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

### Specificity

Recognizes the IRAK2 protein

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

IRAK2 antibody was raised against a peptide corresponding to 17 amino acids near the carboxy terminus of mouse IRAK2a which is common to all four isoforms

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

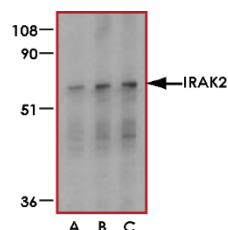
### Scientific Background

Interleukin-1 receptor-associated kinase 2 (IRAK2) is important downstream signaling components of Toll-like receptors (TLRs) (1). IRAKs were first described as signal transducers for IL-1 and later have been implicated in signal transduction of other members of the Toll/IL-1 receptor family. The interleukin-1 receptor (IL-1R) signaling pathway leads to NFκB activation in mammals. To date, four mammalian IRAKs have been identified (IRAK1, IRAK2, IRAK4, and IRAKM) (2). It is believed that IRAK-2 could be an additional therapeutic target for inhibiting IL-1-induced inflammation.

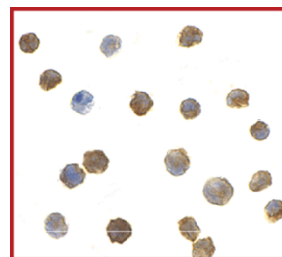
### References

1. Muzio, M. et al: IRAK (Pelle) family member IRAK-2 and MyD88 as proximal mediators of IL-1 signaling. *Science*. 1997;278:1612-5.
2. Wesche, H. et al: IRAK-M is a novel member of the Pelle/interleukin-1 receptor-associated kinase (IRAK) family. *J. Biol. Chem.* 1999; 274:19403-10.

### Sample Data



Western blot analysis of IRAK2 in RAW264.7 whole cell lysate with mIRAK2 antibody at (A) 0.5, (B) 1, and (C) 2 µg/ml.



Immunocytochemistry of IRAK2 in A-20 cells with IRAK2 antibody at 1 µg/ml.

## Anti-IRAK2

Rabbit Polyclonal Antibody

Catalog Number

**I10-63BR**

Specific Lot Number

**O2109-43**

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.**