

Anti-IRAK2

Rabbit Polyclonal Antibody

Catalog # I10-63R

Lot # O2109-12

Cited Applications

WB, ELISA, ICC, IF

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

Specificity

Recognizes the IRAK2 protein and has no cross response to IRAK1

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

IRAK2 antibody was raised against a peptide corresponding to amino acids near the carboxy terminus of human IRAK2

Formulation

PBS + 0.02% sodium azide

Stability

1yr at -20°C from date of shipment

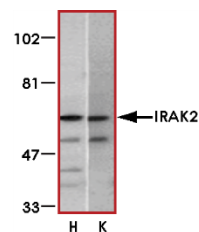
Scientific Background

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 (IRAK-1, IRAK-2, IRAK-4, and IRAK-M) (2). It is believed that IRAK-2 could be an additional therapeutic target for inhibiting IL-1-induced inflammation.

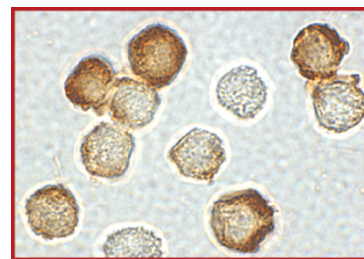
References

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 HeLa (H) and K562 (K) whole cell lysate with IRAK2 antibody (C2) at 2 µg/ml.



Immunocytochemical staining of HeLa cells using IRAK2 antibody at 2 µg/ml.

Anti-IRAK2

Rabbit Polyclonal Antibody

Catalog Number

I10-63R

Specific Lot Number

O2109-12

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.