

IRAK-4 polyclonal antibody

IRAK (IL-1 associated serine/threonine kinase) is a critical regulator of IL-1 induced activation of the NF- κ B pathway. IL-1 stimulates the association of IRAK with the heterodimer of IL-IRI and IL-IRAcP via the adapter protein MyD88. In addition to its role in IL-1 signaling, IRAK is a component of the signal transduction pathway through which TNF-R1 activates NF- κ B dependent gene expression.

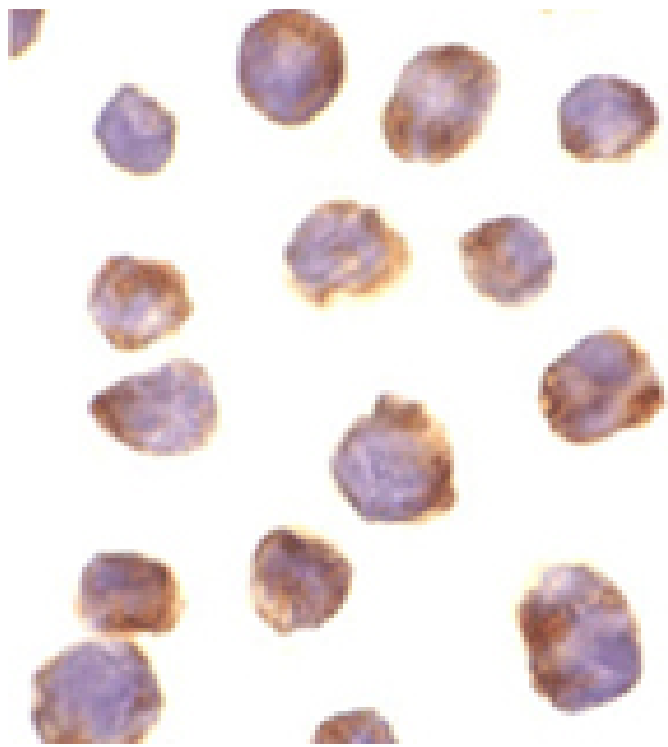
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 3

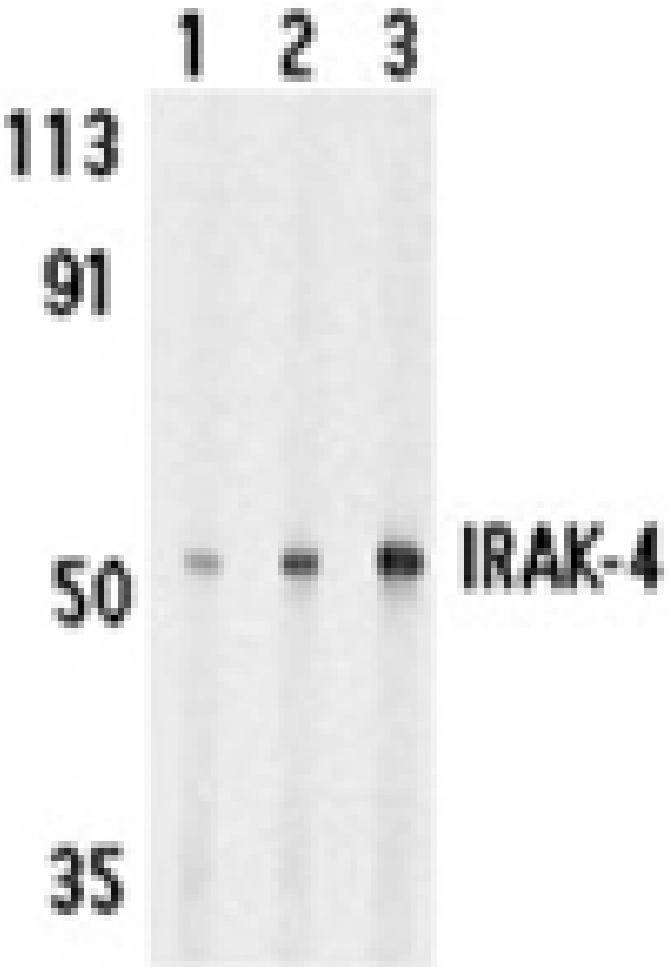
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Immunocytochemistry analysis of IRAK-4 in K562 cells with IRAK-4 antibody at 10µg/ml.



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

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Interleukin-1 receptor-associated kinase 4
Application	ICC, WB
Application Notes	Detects a band of ~50kDa by Western blot.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human IRAK-4.
Purity Detail	Peptide affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (10µg/ml)Western Blot (1-4µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q9NWZ3
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com