

# IRAK-1 polyclonal antibody

IRAK (IL-1 associated serine/threonine kinase) is a critical regulator of IL-1 induced activation of the NF- $\kappa$ 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- $\kappa$ B dependent gene expression.

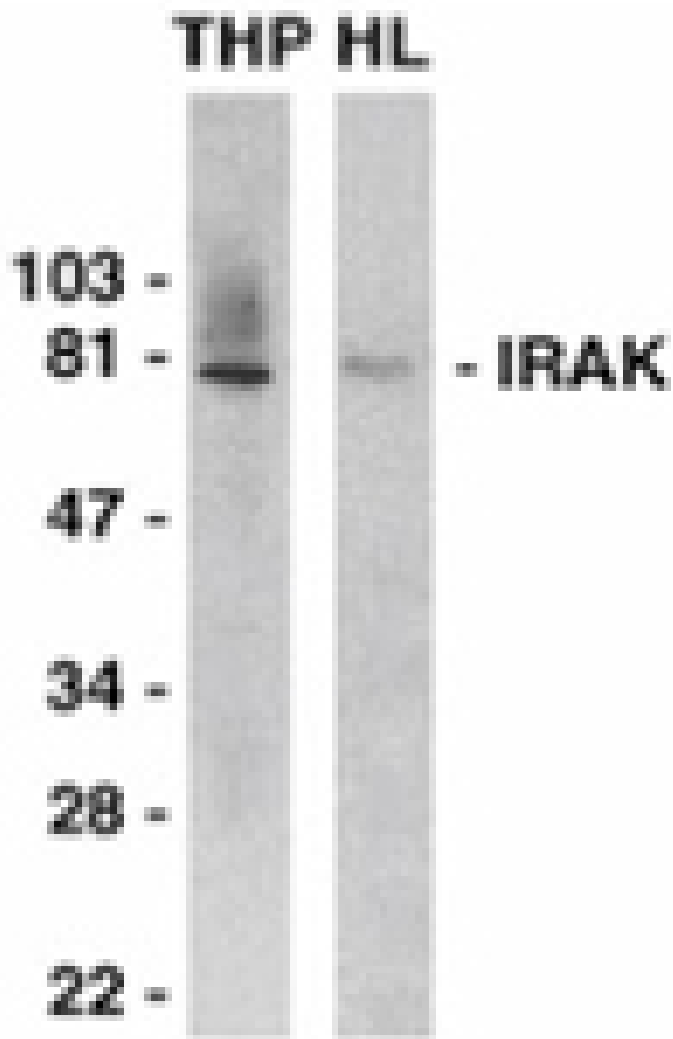
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 3

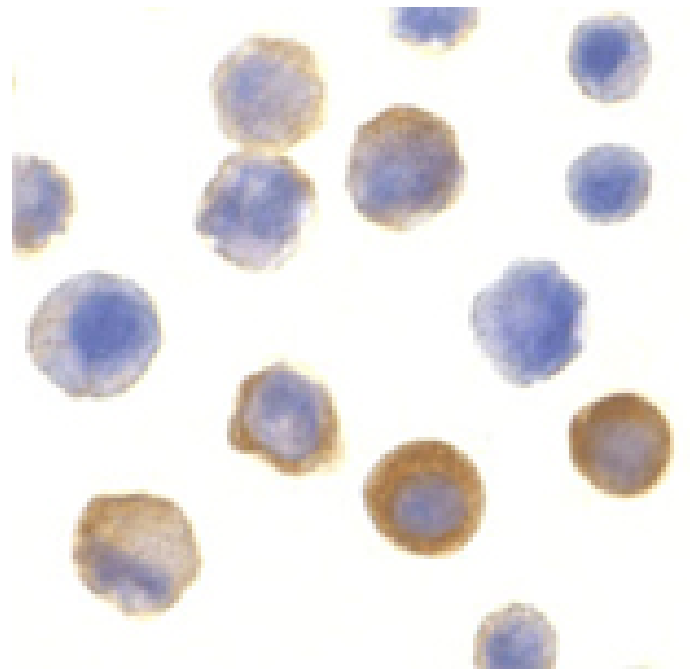
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



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



Immunocytochemistry analysis of IRAK in HeLa cells with IRAK antibody at 10 $\mu$ g/ml.

## 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 1
Application	ICC, IP, WB
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-1.
Purity Detail	DEAE purified.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P51617
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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