

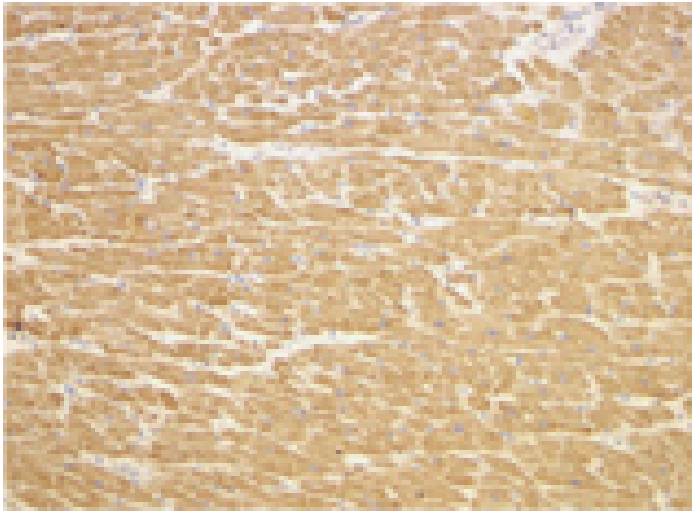
TIRAP polyclonal antibody

Toll-like receptors (TLRs) are signaling molecules that recognize different microbial products during infection and serve as an important link between the innate and adaptive immune responses. These proteins act through adaptor molecules such as TIRAP and MyD88 to activate various kinases and transcription factors.

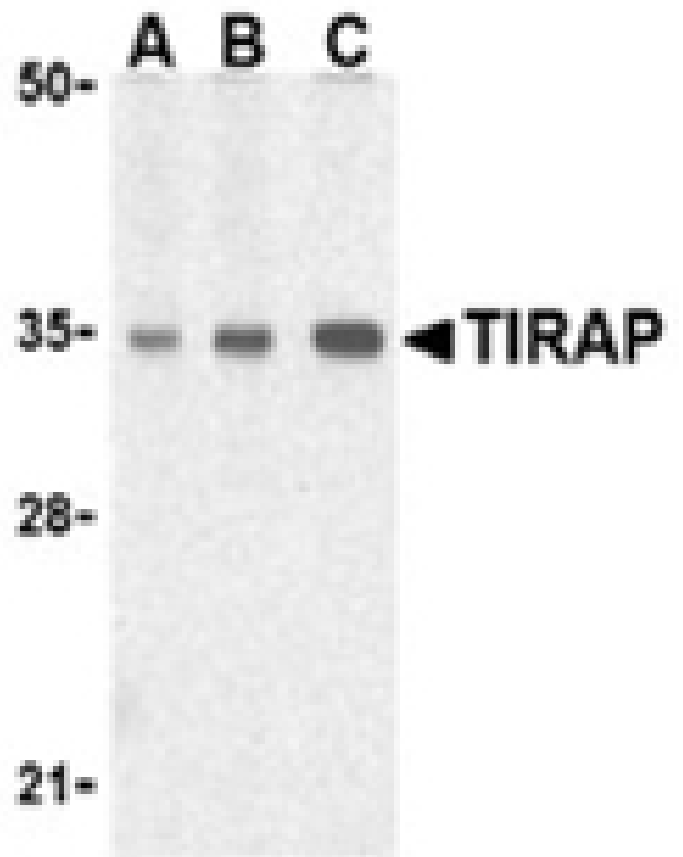
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human heart cells using TIRAP antibody at 2 μ g/ml.



Western blot analysis of TIRAP in human heart cell lysates with TIRAP antibody at (A) 0.5, (B) 1, and (C) 2 μ g/ml.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Toll-interleukin 1 receptor domain-containing adapter protein, MyD88 adapter-like protein, Mal
Application	IHC, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of mouse TIRAP.
Purity Detail	Peptide affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (2µg/ml)Western Blot (0.5-2.0µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q99JY1
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee