

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](#).

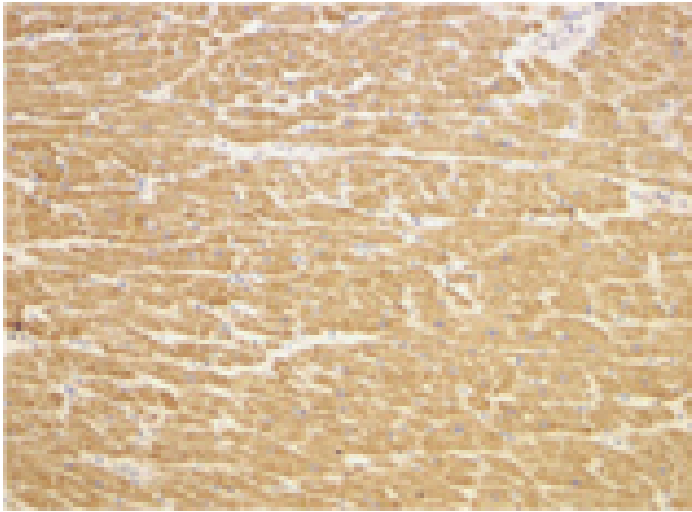
Ordering Information

[Order Online »](#)

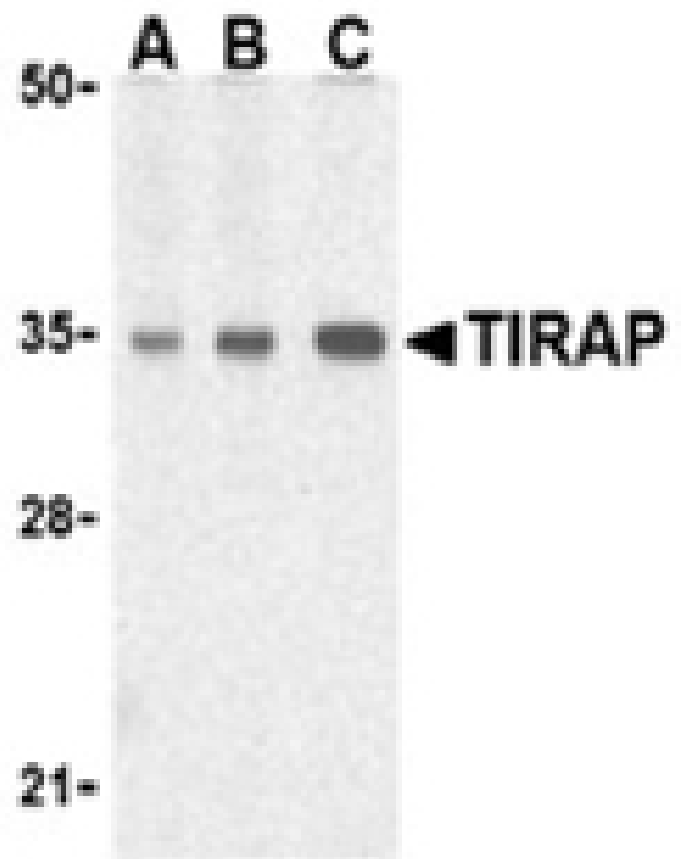
ADI-CSA-514-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



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



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



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: +1 940 948 9120
info@enzolifesciences.com
usa@enzolifesciences.com

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

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

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

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

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