

# MyD88 polyclonal antibody

MyD88 (Myeloid differentiation primary response gene 88) is a general adapter protein for the Toll/IL-1R family of receptors that have been shown to induce growth arrest, terminal differentiation, and apoptosis. MyD88 plays an important role in the inflammatory response induced by cytokines, IL-1 and IL-18, and endotoxin. It has been shown that signaling via MyD88 is essential for an LPS response. MyD88 is a cytoplasmic protein, which is expressed in many tissues.

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

Citations: 1

[View Online »](#)

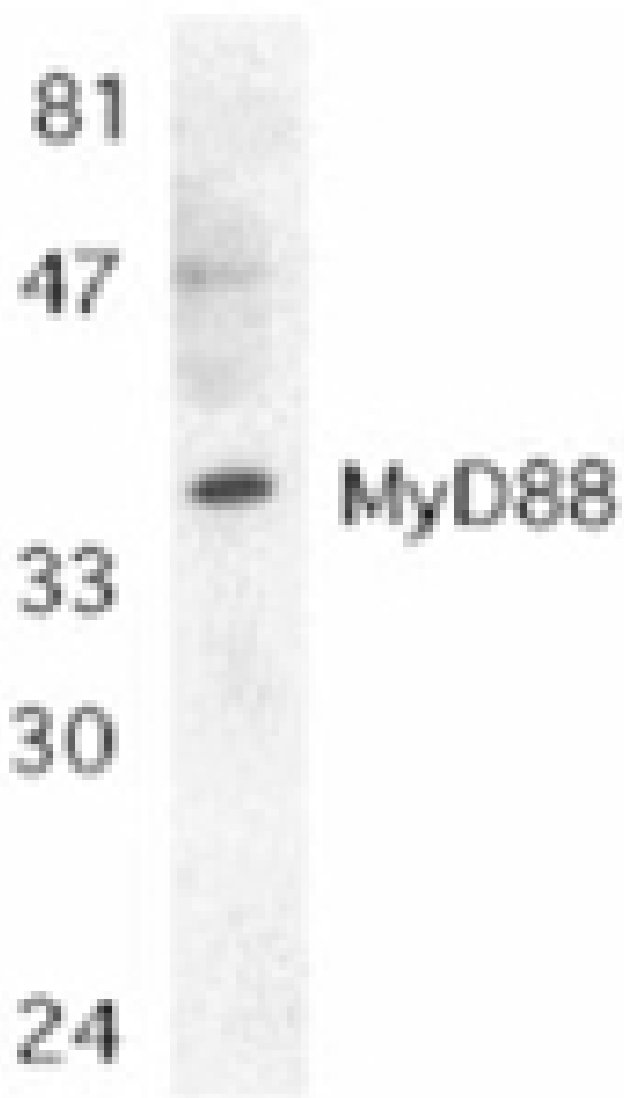
## Ordering Information

[Order Online »](#)

ADI-CSA-510-E	100µg
---------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of MyD88 in Jurkat whole cell lysate with MyD88 antibody at 1:500 dilution.

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

**Alternative Name**      Myeloid differentiation primary response gene

**Application**      WB

**Application Notes**      Detects a band of ~35kDa by Western blot.

**Formulation**      Liquid. In PBS containing 0.02% sodium azide.

**GenBank ID**      NM\_002468.1 (RefSeq)

**Host**      Rabbit

**Immunogen**      Synthetic peptide corresponding to a portion of human MyD88. The sequence is completely conserved in mouse.

**Purity Detail**      Peptide affinity purified.

**Recommendation  
Dilutions/Conditions**      Western Blot (1-2µg/ml, ECL)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**      Q99836

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



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-  
fr@enzolifesciences.com](mailto:info-fr@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)