## 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** 

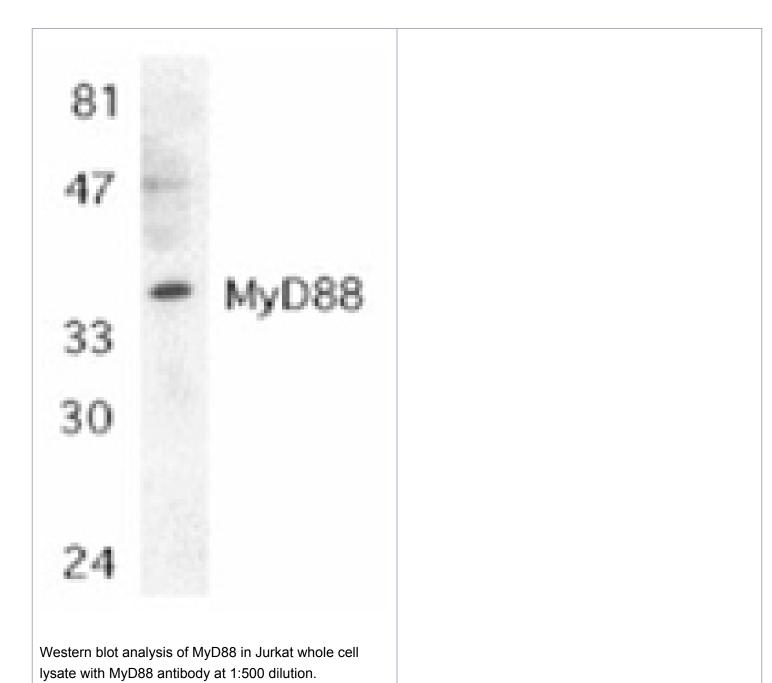
Order Online »

ADI-CSA-510-E

100µg

Manuals, SDS & CofA

**View Online** »



## **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

