

MyD88 antibody (pAb)

Catalog No: 40914

RRID: AB_2793435

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Concentration: 0.5 µg/µl

Molecular Weight: 35 kDa

Background: MyD88 is an adapter protein involved in the Toll-like receptor and IL-1 receptor signaling pathway in the innate immune response. MyD88 acts via IRAK1, IRAK2 and TRAF6, leading to NFκB (NFκB p50 & NFκB p65) activation, cytokine secretion and the inflammatory response. MyD88 increases IL-8 transcription, and may be involved in myeloid differentiation.

Immunogen: This MyD88 antibody was raised against a synthetic peptide corresponding to amino acid residues 233-248 of human MyD88.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

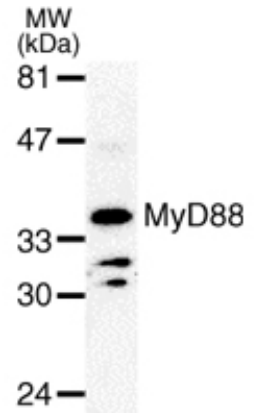
Applications Validated by Active Motif:

WB: 2 - 4 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



MyD88 pAb tested by Western blot.

MyD88 detection by Western blot. The analysis was performed using Jurkat whole-cell extract and MyD88 pAb at a 2 µg/ml dilution.