

TLR8 antibody (mAb)

Catalog No: 40955

RRID: AB_2793456

Clone: 44C143

Isotype: IgG1

Application(s): WB

Reactivity: Human, Mouse

Quantity: 100 µg

Purification: Affinity Purified

Host: Mouse

Molecular Weight: 120 kDa

Background: TLR8 – Toll-Like receptor 8 is a member of the Toll-like receptor (TLR) family that plays a fundamental role in pathogen recognition and activation of innate immunity. TLRs are highly conserved from *Drosophila* to humans and share structural and functional similarities. They recognize specific molecular features associated with pathogens, and mediate the production of cytokines necessary for the development of effective immunity. Many of the TLRs heterodimerize with each other to modulate receptor function. TLR8 expression cells is elevated after exposure to both Gram-positive and Gram-negative bacteria. While primarily extracellular, TLR8 may be also localized intracellularly.

Immunogen: This TLR8 antibody was raised against a peptide corresponding to amino acid residues 750-850 of human TLR8.

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

Application Notes:

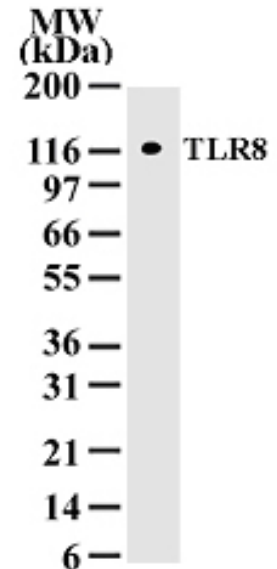
Applications Validated by Active Motif:

WB: 1 - 2 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. 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.



TLR8 mAb tested by Western blot.

TLR8 detection by Western blot. The analysis was performed using TLR8 mAb and nuclear extract from 293 cells that had been transfected with human TLR8 cDNA.