

## TLR6 antibody (mAb)

**Catalog No:** 40954

**RRID:** AB\_2793455

**Clone:** 86B1153

**Application(s):** WB

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Affinity Purified

**Host:** Mouse

**Isotype:** IgG1

**Concentration:** 0.5 µg/µl

**Molecular Weight:** 92 kDa

**Background:** TLR6 – Toll-Like receptor 6 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. TLR6 forms a heterodimer with TLR2. TLR6 is thought to specify the sensitivity of TLR2 and contribute to its signaling capabilities through heterodimerization.

**Immunogen:** This TLR6 antibody was raised against a synthetic peptide corresponding to amino acid residues 408-424 of human TLR6.

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

### Application Notes:

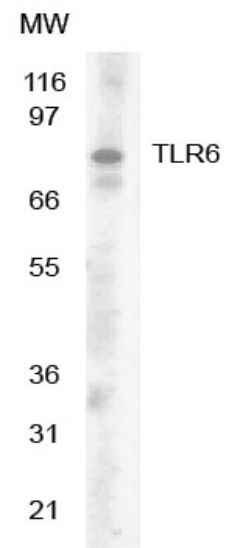
Applications Validated by Active Motif:

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.



### TLR6 mAb tested by Western blot.

TLR6 detection by Western blot. The analysis was performed using Daudi whole-cell extract and TLR6 mAb at a 2 µg/ml dilution.