

## TLR9 antibody (mAb)

**Catalog No:** 40956

**Clone:** 26C593

**Isotype:** IgG1

**Application(s):** WB

**Reactivity:** Human, Mouse, Rat

**Purification:** Affinity Purified

**Host:** Mouse

**Concentration:** 1 µg/µl

**Quantity:** 100 µg

**Molecular Weight:** 120 kDa

**Background:** TLR9 (Toll-Like receptor 9) 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. TLR9 forms a homodimer and recognizes unmethylated bacterial DNA. While primarily extracellular, TLR9 may be also localized intracellularly.

**Immunogen:** This TLR9 antibody was raised against a KLH-conjugated synthetic peptide corresponding to amino acid residues 268-284 of human TLR9.

**Buffer:** PBS containing 0.05% BSA and 0.05% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

Validated Applications:

WB: 1 - 2 µg/ml dilution

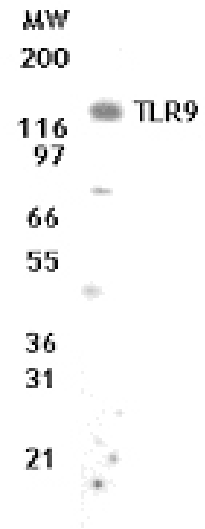
Published Applications: see references

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

**Storage and Guarantee:** Antibodies in solution can be stored at -20°C for 2 years.

Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

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



**TLR9 mAb tested by Western blot.**

TLR9 detection by Western blot. The analysis was performed using PBMC cell extract and TLR9 mAb at a 2 µg/ml dilution.