

TBK1 antibody (mAb)

Catalog No: 40909**RRID:** AB_2793432**Clone:** 108A429**Application(s):** WB**Reactivity:** Human, Mouse, Rat**Quantity:** 100 µg**Purification:** Affinity Purified**Host:** Mouse**Isotype:** IgG1**Concentration:** 1 µg/µl**Molecular Weight:** 80 kDa

Background: TBK1 – TANK-binding kinase 1 is a regulator of the NFκB (NFκB p50 & NFκB p65) signaling pathway. NFκB signaling is controlled to a large extent by the sequestration of the NFκB complex in the cytoplasm by its association with one of the IκB family of proteins. IκB is phosphorylated by the IκB Kinase (IKK) complex, resulting in the degradation of IκB and the nuclear translocation of NFκB. TANK is an IκB family member that is specifically recognized and phosphorylated by TBK1, leading to the destruction of TANK and the activation of the NFκB pathway.

Immunogen: This TBK1 antibody was raised against a peptide corresponding to amino acid residues 563-577 of human TBK1.

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

Application Notes:

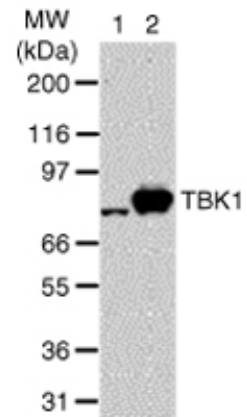
Applications Validated by Active Motif:

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. 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.

**TBK1 mAb tested by Western blot.**

TBK1 detection by Western blot. The analysis of TBK1 was performed using TBK1 mAb at a 2 µg/ml dilution and nuclear extract from 293 cells (lane 1) and nuclear extract from 293 cells that had been transfected with human TBK1 cDNA (lane 2).