

## TAC1 antibody (pAb)

**Catalog No:** 40936**RRID:** AB\_2793450**Isotype:** IgG**Application(s):** WB**Reactivity:** Human, Mouse**Quantity:** 100 µg**Purification:** Affinity Purified**Host:** Rabbit**Concentration:** 0.5 µg/µl**Molecular Weight:** 32 kDa

**Background:** TAC1 – Tachykinin 1 is a member of a family of peptide hormones that serve as neurotransmitters. The tachykinins excite neurons, evoke behavioral responses, are potent vasodilators and secretagogues, and stimulate contraction of smooth muscle cells.

**Immunogen:** This TAC1 antibody was raised against a synthetic peptide corresponding to amino acid residues 116-132 of human TAC1.

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

**Application Notes:**

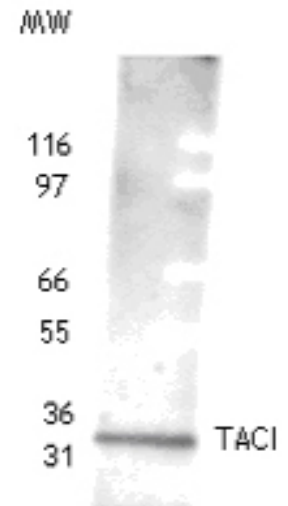
Applications Validated by Active Motif:

WB: 0.5 - 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.

**TAC1 pAb tested by Western blot.**

TAC1 detection by Western blot. The analysis of TAC1 was performed using 10 µg of human spleen nuclear extract and TAC1 pAb at 1 µg/ml. A protein band of approximately 32 kDa was detected.