

TROY/TAJ antibody (pAb)

Catalog No: 40937**RRID:** AB_2793451**Application(s):** WB**Reactivity:** Human, Mouse**Quantity:** 100 µg**Purification:** Affinity Purified**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 46 kDa

Background: TROY / TAJ – Tumor Necrosis Factor Receptor Superfamily Member 19 is a member of the TNF-receptor superfamily. It is highly expressed during embryonic development and has been shown to interact with TRAF (TRAF2 and TRAF5) family members, and to activate the JNK signaling pathway when overexpressed in cells. This receptor can induce apoptosis by a caspase-independent mechanism, and it is thought to play an essential role in embryonic development.

Immunogen: This TROY/TAJ antibody was raised against a mixture of synthetic peptides corresponding to amino acid residues 29-44, 204-219, and 409-421 of human TROY/TAJ.

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

Application Notes:

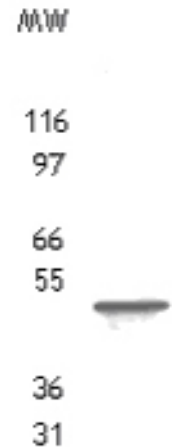
Applications Validated by Active Motif:

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. 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.



TROY/TAJ pAb tested by Western blot. TAJ/TROY detection by Western blot. The analysis of TAJ/TROY was performed using TROY/TAJ pAb in T98G, a human glioblastoma cell line extract. A protein band of approximately 46 kDa was detected.