

TAZ / WWTR1 antibody (pAb)

Catalog Nos: 61265, 61266

RRID: AB_2614991

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Concentration: 1.15 μ g/ μ l

Molecular Weight: 55 kDa

Background: The TAZ / WWTR1 (Transcriptional coactivator with PDZ-binding motif, WW domain-containing transcription regulator protein 1) gene encodes a protein that is a transcriptional coactivator which acts as a downstream regulatory target in the Hippo signaling pathway that plays a pivotal role in organ size control and tumor suppression by restricting proliferation and promoting apoptosis. The core of this pathway is composed of a kinase cascade wherein MST1/MST2, in complex with its regulatory protein SAV1, phosphorylates and activates LATS1/2 in complex with its regulatory protein MOB1, which in turn phosphorylates and inactivates YAP1 oncoprotein and WWTR1/TAZ. WWTR1 enhances PAX8 and NKX2-1/TTF1-dependent gene activation. Regulates the nuclear accumulation of SMADS and has a key role in coupling them to the transcriptional machinery such as the mediator complex. Regulates embryonic stem-cell self-renewal, promotes cell proliferation and epithelial-mesenchymal transition.

Immunogen: TAZ / WWTR1 antibody was raised against a peptide corresponding to the C-terminal region of human TAZ / WWTR1.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

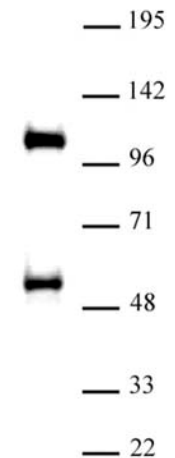
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 dilution

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.



TAZ / WWTR1 antibody (pAb) tested by Western blot

Nuclear extract of A-431 cells (30 μ g) probed with TAZ / WWTR1 antibody (1:500).