

TORC2 antibody (pAb)

Catalog Nos: 39779, 39780

RRID: AB_2615051

Isotype: Serum

Application(s): IP, WB

Reactivity: Human, Mouse

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 85 kDa

Background: TORC2 (Transducer of regulated cAMP response element-binding protein 2) (CRTC2 or CREB-regulated transcription coactivator 2) is a coactivator of transcription through the CREB1 signaling pathway, independent of phosphorylation at serine 133. Involved in metabolic signaling pathways, it is responsive to glucagon signaling and regulates gluconeogenesis. It is a potent activator of mitochondrial biogenesis in muscle cells. TORC2 is inactivated by phosphorylation at serine 171 (in response to both insulin signaling and the activation of AMPK), which facilitates recognition by 14-3-3 protein and subsequent translocation to the cytoplasm.

Immunogen: This TORC2 antibody was raised against a peptide at the C-terminus of human TORC2.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

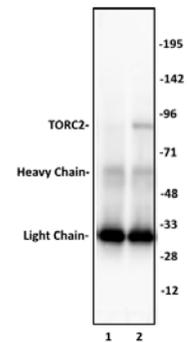
Applications Validated by Active Motif:

IP: 5 - 10 µl per IP

WB: 1:500 - 1:2,000 dilution

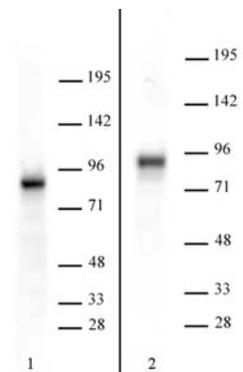
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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



TORC2 antibody (pAb) tested by Immunoprecipitation.

10 µl of TORC2 antibody was used to immunoprecipitate TORC2 from 400 µg of HEK293 whole cell extract (lane 2). 10 µl of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the TORC2 antibody at a dilution of 1:2,000.



TORC2 antibody (pAb) tested by Western blot.

Detection of TORC2 by Western blot analysis. Whole cell extract (20 µg) of NIH/3T3 (lane 1) and HEK293 (lane 2) blotted with TORC2 antibody at a dilution of 1:2,000.