

RBM39 antibody (pAb)

Catalog Nos: 61669, 61670

RRID: AB_2793727

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 68 kDa

Background: **RBM39** (RNA Binding Motif Protein 39) is a transcriptional coactivator for steroid nuclear receptors ESR1/ER-alpha and ESR2/ER-beta, and JUN/AP-1 (By similarity). May be involved in pre-mRNA splicing process.

Immunogen: This antibody was raised against a peptide within the N-terminal region of human RBM39.

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

Application Notes:

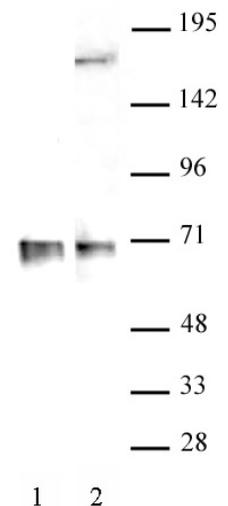
Applications Validated by Active Motif:

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

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

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.



RBM39 antibody (pAb) tested by Western blot.

Nuclear extract (20 μ g) of Raji cells (Lane 1) or nuclear extract (30 μ g) of HeLa cells (Lane 2) probed with RBM39 antibody at a 1:500 dilution.