

ARID3A antibody (pAb)

Catalog Nos: 61615, 61616

RRID: AB_2793703

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 80 kDa

Background: ARID3A (AT Rich Interactive Domain 3A (BRIGHT-Like)) is a transcription factor which may be involved in the control of cell cycle progression by the RB1/E2F1 pathway and in B-cell differentiation.

Immunogen: This antibody was raised against a peptide within the C-terminal region of human ARID3A.

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

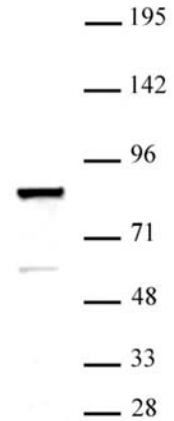
Application Notes:

Applications Validated by Active Motif:

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



ARID3A antibody (pAb) tested by Western blot.

Detection of ARID3A by Western blot. The analysis was performed using 30 μ g K562 whole-cell extract and ARID3A (pAb) at a 1:500 dilution.