

BRD9 antibody (pAb)

Catalog Nos: 65705, 65905, 65706

Application(s): IP, WB

Reactivity: Human

Volumes: 100 µl, 50 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Molecular Weight: 85 kDa

Background: BRD9 (bromodomain containing 9) belongs to the BET subclass of proteins, which are characterized by two N-terminal bromodomains and one ET (Extra Terminal) domain. BRDs associate with chromatin through their bromodomains that recognize acetylated histone lysine residues. Bromodomains function as 'readers' of these epigenetic histone marks and regulate chromatin structure and gene expression by linking associated proteins to the acetylated nucleosomal targets. The ET domain functions as a protein binding motif and exerts atypical serine-kinase activity. The BET family consists of at least four members in mouse and human, BRD2 (also referred to as FSRG1, RING3), BRD3 (FSRG2, ORFX), BRD4 (FSRG4, MCAP/HUNK1), and BRDT (FSRG3, BRD6) that function in the regulation of transcriptional activation and chromatin remodeling. There are five isoforms of BRD9 that are produced by alternative splicing.

Immunogen: This antibody was raised against a peptide within the C-terminal region of human BRD9 (bromodomain containing 9).

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:

IP: 5 µl per IP

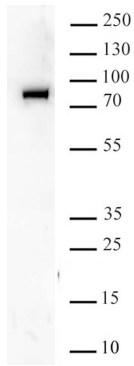
WB*: 1:250 - 1:1,000 dilution

*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.

For optimal IP results, the addition of 0.05% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

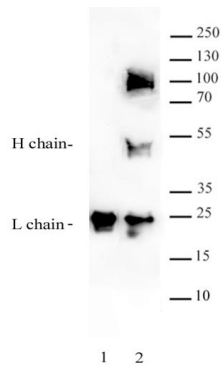
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



BRD9 antibody (pAb) tested by Western blot.

Nuclear extract (20 µg) of K562 probed with BRD9 antibody at a 2 µg/mL dilution.



BRD9 antibody (pAb) tested by Immunoprecipitation.

5 µl of BRD9 antibody was used to immunoprecipitate BRD9 from 500 µg of Jurkat WCE (lane 2). 5 µl of mouse IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the BRD9 antibody at a dilution of 1:1000.