

JunD antibody (pAb)

Catalog Nos: 61403, 61404

RRID: AB_2793621

Application(s): ChIP-Seq, CUT&RUN, IP, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Molecular Weight: 40 & 45 kDa

Background: JunD (Jun D proto-oncogene) is a member of the JUN family and forms a component of the AP-1 transcription factor complex. JunD binds DNA as a dimer. Other AP-1 family members consist of c-Fos, Fra-1, Fra-2, FosB, c-Jun, and JunD.

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

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:

ChIP-Seq: 2 - 5 µl per ChIP

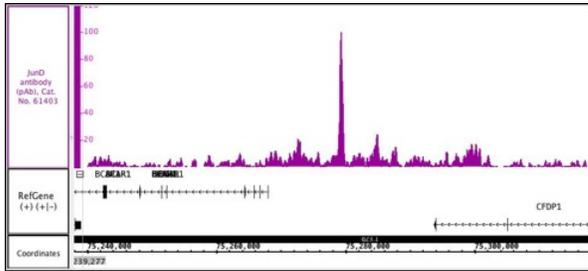
IP: 10 µl per IP

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

CUT&RUN: 1 µl per 50 µl reaction

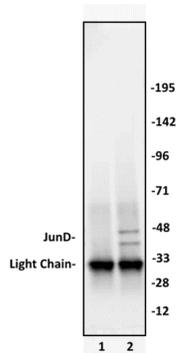
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



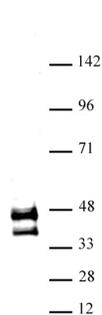
JunD antibody (pAb) tested by ChIP-Seq

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT® High Sensitivity Kit (Cat. No. 53040) with 30 µg of chromatin from human endometrial stromal cell chromatin and 4 µg of JunD antibody. ChIP DNA was sequenced on the Illumina HiSeq and 8.6 million sequence tags were mapped to identify JunD binding sites on chromosome 16. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, here.



JunD antibody (pAb) tested by Immunoprecipitation.

10 µl of JunD antibody was used to immunoprecipitate JunD from 250 µg of K562 nuclear 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 JunD antibody at a dilution of 1:1,000.



JunD antibody (pAb) tested by Western blot.

K-562 nuclear extract (20 µg) probed with JunD antibody at a dilution of 1:1,000.

JunD antibody (pAb) tested by CUT&RUN

CUT&RUN was performed using 500,000 K562 nuclei and sequenced using 38 base-pair, paired-end reads on the Illumina NovaSeq. Data was collected from 46 million reads, and JunD data is shown for Chromosome 19.

