

## BRD3 antibody (pAb)

---

**Catalog Nos:** 61489, 61490

**RRID:** AB\_2737028

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

**Reactivity:** Human, Mouse

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Isotype:** IgG

**Molecular Weight:** 100 kDa

---

**Background:** BRD3 (Bromodomain Containing 3) is a member of the mammalian BET (Bromodomain Extra Terminal) family of proteins, binds hyperacetylated chromatin and has been reported to interact with acetylated Histone H4K5, Histone H4K12 and Histone H3K14. It plays a role in the regulation of transcription, probably by chromatin remodeling and interaction with transcription factors. Regulates transcription by promoting the binding of the transcription factor GATA1 to its targets (By similarity). Regulates transcription of the CCND1 gene.

**Immunogen:** This antibody was raised against a peptide within the C-terminal region of mouse BRD3.

**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: 10 µl per ChIP

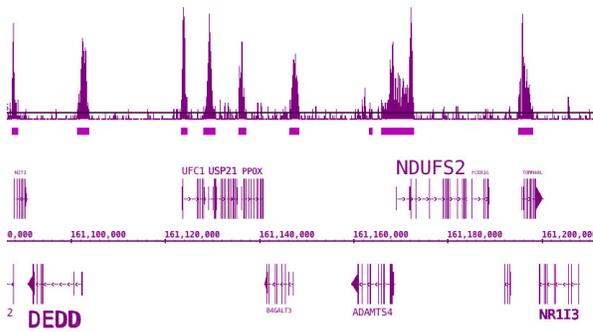
ChIP-Seq & ChIP-chip: 10 µl each

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

CUT&RUN: 1 µl per 50 µl reaction

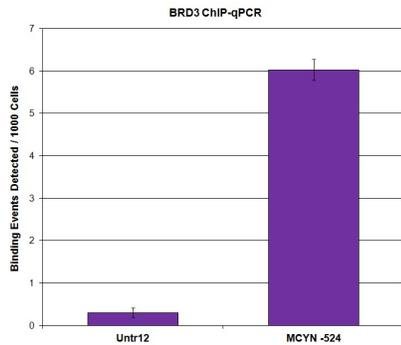
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



### BRD3 antibody (pAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30 µg of chromatin from human H526 cells and 10 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 16 million sequence tags were mapped to show the expected promoter localization.

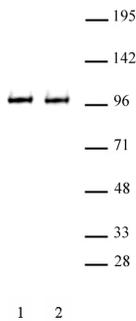


### BRD3 antibody (pAb) tested by ChIP.

Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT<sup>®</sup> High Sensitivity Kit (Cat. No. 53040) with 30 µg of chromatin from human H526 cells and 10 µl BRD3 antibody. ChIP DNA was used in qPCR with the control primer or gene-specific primer as indicated. Data are presented as Binding Events Detected per 1000 Cells using Active Motif's Epigenetic Services normalization scheme which accounts for primer efficiency and the amount of chromatin used in the ChIP reaction.

### BRD3 antibody (pAb) tested by Western blot.

Detection of BRD3 by Western blot analysis. Lane 1: Whole cell extract of HeLa cells (20 µg). Lane 2: Whole cell extract of NIH-3T3 cells. Both probed with BRD3 antibody (pAb) at a 1:500 dilution.



### BRD3 antibody (pAb) tested by CUT&RUN

CUT&RUN was performed using 500,000 K562 cells and sequenced using 38 base-pair, paired-end reads on the Illumina NovaSeq. Data was collected from 23 million reads, and BRD3 data is shown for Chromosome 19.

