

BRD8 / SMAP2 antibody (pAb)

Catalog Nos: 61007, 61008

RRID: AB_2615053

Isotype: IgG

Application(s): ChIP, ChIP-Seq, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 145 kDa

Background: BRD8 (Bromodomain-containing protein 8, SMAP2, TRCP120) is a member of the Bromodomain-containing family of chromatin associated proteins. BRD8 is a member of the NuA4 histone acetyltransferase complex, involved in the transcriptional activation of specific genes through the acetylation of histones H2A and H4. BRD8 contains two Bromodomains, a motif that has been demonstrated to facilitate interaction with acetylated histones.

Immunogen: This BRD8 / SMAP2 antibody was raised against a peptide corresponding to amino acids within the N-terminal region of human BRD8.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 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: 10 µl each

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

ChIP-Seq validation was performed by Active Motif's Epigenetics Services; the complete data set is available in the UCSC Genome Browser by clicking [here](#).

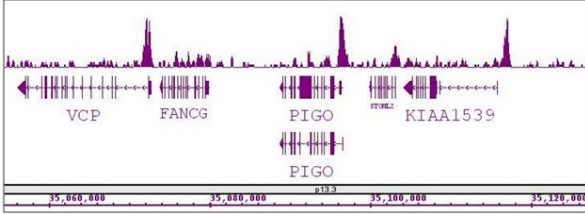
*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

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.

BRD8 / SMAP2 antibody (pAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) with 30 ug of chromatin from an adenocarcinoma cell line and 10 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify BRD8 / SMAP2 binding sites. The image shows binding across a region of chromosome 9. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, here.



Western blot of BRD8 / SMAP2 antibody.

Nuclear extract of HeLa cells (30 µg) probed with BRD8 / SMAP2 antibody (1:500).

