

Boris / CTCFL antibody (pAb)

Catalog Nos: 39851, 39852

RRID: AB_2793368

Isotype: IgG

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

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 96 kDa

Background: BORIS (Brother of Regulator of Imprinted Sites) / CTCFL (CTCF-like) is a zinc-finger DNA binding protein that acts as a chromatin insulator and a paralog of the CTCF protein. Normally, BORIS / CTCFL is expressed only in primary spermatocytes and is thought to play a role in resetting DNA methylation marks during the differentiation of male germ cells. BORIS / CTCFL is also expressed in many cancers and may serve a pro-proliferative function.

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

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: 5 µl per ChIP

ChIP-Seq: 5 µl each

WB: 1:500 - 1:2,000 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](#).

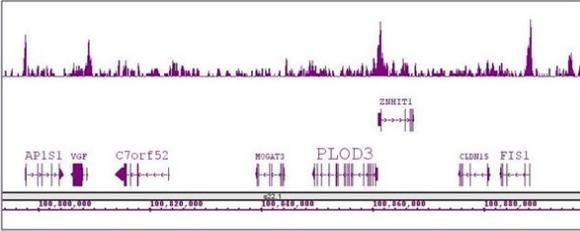
For CTCF, we also offer AbFlex® CTCF Recombinant Antibody (rAb). For details, see Catalog No. 91285.

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.

Boris / CTCFL 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 human testis and 2 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 20 million sequence tags were mapped to identify Boris / CTCFL binding sites. The image shows binding across a region of chromosome 7. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, here.



BORIS / CTCFL antibody (pAb) tested by Western blot.

20 µg of LNCaP nuclear extract was probed with BORIS / CTCFL antibody at 1:1,000.

