

Ets-1 antibody (pAb)

Catalog Nos: 39580, 39581

RRID: AB_2793266

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

Reactivity: Human, Mouse

Volumes: 200 µl, 10 µl

Purification: None

Host: Rabbit

Isotype: Serum

Molecular Weight: 54 kDa

Background: Ets-1 transcription factor is a member of the Ets family of proteins that share a unique binding domain. Ets1 (E26 transformation-specific sequence-1) is a proto-oncogene whose DNA binding activity is controlled by kinases and other transcription factors. Ets 1 has been associated with regulation of cell growth, differentiation and apoptosis.

Immunogen: This Ets-1 antibody was raised against a peptide corresponding to the C-terminal half of mouse Ets-1.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

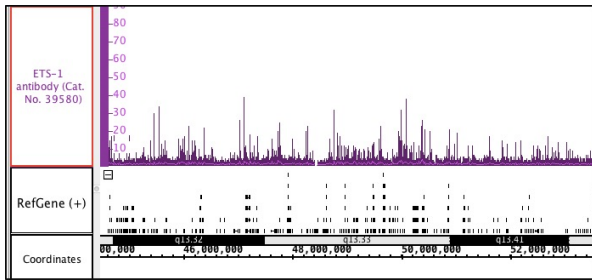
ChIP-Seq: 4 µg (7 µl) per ChIP

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

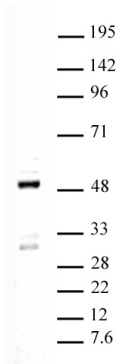
CUT&RUN: 1 µl per 50 µl reaction

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.

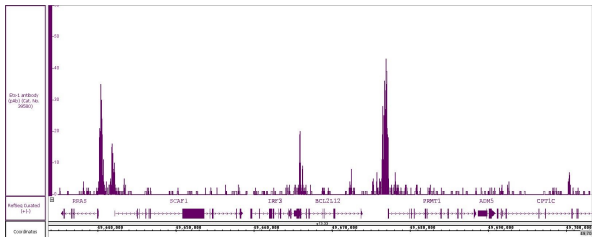


ETS-1 antibody (pAb) tested by ChIP-Seq.Chromatin immunoprecipitation (ChIP) was performed using the ChIP-IT® High Sensitivity Kit (Cat. No. 53040) with 30 µg of chromatin from human CD4+ T cell chromatin and 4 µg ETS-1 antibody. ChIP DNA was sequenced on the Illumina HiSeq and 7.9 million sequence tags were mapped to identify ETS-1 binding sites in a region of human chromosome 19.



Ets-1 pAb tested by Western blot.

Detection of Ets-1 by Western blot. The analysis was performed using CCRF-CEM nuclear extract and Ets-1 pAb at a 1:500 dilution.



Ets-1 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 49 million reads, and Ets-1 data is shown for Chromosome 19.