

ETO / RUNX1T1 antibody (pAb)

Catalog Nos: 61109, 61110

RRID: AB_2793509

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 68 kDa

Background: ETO / RUNX1T1 (AML1T1, CBFA2T1, MTG8, ZMYND2) belongs to a family of evolutionarily conserved nuclear factors that act as a transcriptional corepressor. ETO / RUNX1T1 does not interact directly with DNA but is recruited by transcription factors such as PLZF, BCL6, TAL1/SCL, Gfi1 and Heb recruit nuclear co-repressors such as NCoR, SIN3, and SMRT / NCoR2, which in turn interact with histone deacetylase (HDAC) to mediate transcriptional repression. ETO / RUNX1T1 contains four evolutionarily conserved functional domains called nervy homology regions (NHRs) involved in protein-protein interaction. ETO / RUNX1T1 is frequently found as a fusion partner of AML1 in acute myeloid leukemia with the t(8;21) translocation which gives rise to the AML-ETO fusion protein.

Immunogen: This ETO / RUNX1T1 antibody was raised against a peptide within the internal region of human ETO / RUNX1T1.

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:

IP: 10 µl per IP

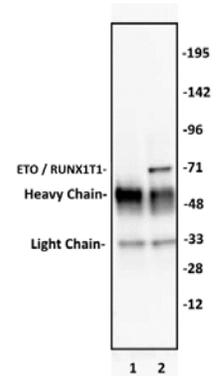
WB*: 1:500 - 1:2,000 dilution

For optimal results, primary antibody incubations should be performed at 4°C. Individual optimization may be required.

*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.

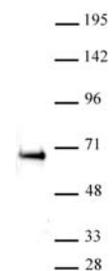
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



ETO / RUNX1T1 antibody (pAb) tested by Immunoprecipitation.

10 µl of ETO / RUNX1T1 antibody was used to immunoprecipitate ETO / RUNX1T1 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 ETO / RUNX1T1 antibody at a dilution of 1:500.



ETO / RUNX1T1 antibody (pAb) tested by Western Blot:

K-562 nuclear extract (20 µg per lane) probed with the ETO / RUNX1T1 antibody (pAb) at a dilution of 1:500.