

Smyd3 antibody (pAb)

Catalog Nos: 61407, 61408

RRID: AB_2793623

Isotype: IgG

Application(s): IHC, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 50 kDa

Background: Smyd3 (SET and MYND domain containing 3) is a histone methyltransferase. Specifically methylates 'Lys-4' of histone H3, inducing di- and tri-methylation, but not monomethylation. Plays an important role in transcriptional activation as a member of an RNA polymerase complex. Binds DNA containing 5'-CCCTCC-3' or 5'-GAGGGG-3' sequences. Recently, Smyd3 was described as the methyltransferase responsible for a novel site of histone modification, H4K5 methylation.

Immunogen: This antibody was raised against a peptide within human Smyd3.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

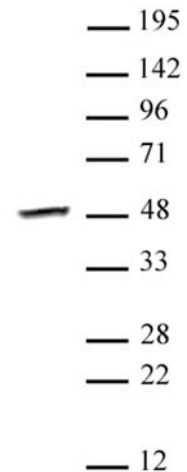
Applications Validated by Active Motif:

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

*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.



Smyd3 antibody (pAb) tested by Western blot.

K-562 nuclear extract (30 μ g) probed with Smyd3 antibody at a dilution of 1:500.