

SFMBT1 antibody (pAb)

Catalog Nos: 61759, 61760

RRID: AB_2793761

Isotype: IgG

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

Reactivity: Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 110 kDa

Background: **SFMBT1** (Scm-Like With Four Mbt Domains 1) is a histone-binding protein, which is part of various corepressor complexes. Mediates the recruitment of corepressor complexes to target genes, followed by chromatin compaction and repression of transcription. Plays a role during myogenesis: required for the maintenance of undifferentiated states of myogenic progenitor cells via interaction with MYOD1. Interaction with MYOD1 leads to the recruitment of associated corepressors and silencing of MYOD1 target genes. Part of the SLC complex in germ cells, where it may play a role during spermatogenesis.

Immunogen: This antibody was raised against a peptide within the C-terminal region of mouse SFMBT1.

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

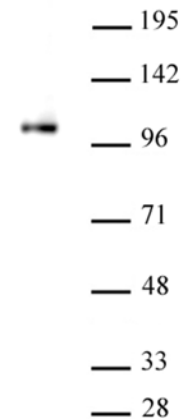
Application Notes:

Applications Validated by Active Motif:

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

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.



SFMBT1 antibody (pAb) tested by Western blot.

Detection of SFMBT1 antibody by Western blot. The analysis was performed using mouse testes nuclear extract (30 μ g) and SFMBT1 antibody at a 1:1,000 dilution.