

LSH antibody (pAb)

Catalog Nos: 61595, 61596

RRID: AB_2793693

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 100 kDa

Background: LSH (lymphoid-specific helicase) plays an essential role in normal development and survival. Involved in regulation of the expansion or survival of lymphoid cells. Required for de novo or maintenance DNA methylation. May control silencing of the imprinted CDKN1C gene through DNA methylation. May play a role in formation and organization of heterochromatin, implying a functional role in the regulation of transcription and mitosis (By similarity).

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

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

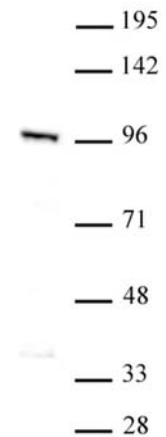
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:2,000 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.



LSH antibody (pAb) tested by Western blot.

Mouse testis nuclear cell extract (20 μ g) probed with LSH antibody at a dilution of 1:1,000.