

SOX6 antibody (pAb)

Catalog Nos: 61749, 61750

RRID: AB_2793756

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 100 kDa

Background: **SOX6** (SRY (Sex Determining Region Y)-Box 6) is a transcriptional activator. Binds specifically to the DNA sequence 5-AACAAT-3. Plays a key role in several developmental processes, including neurogenesis and skeleton formation.

Immunogen: This antibody was raised against a recombinant protein corresponding amino acids 1-20 of human SOX6.

Buffer: Purified IgG in 70 mM Tris, 105 mM NaCl, 0.07 mM EDTA, 31 mM glycine, 30% glycerol, and 0.035% sodium azide. Sodium azide is highly toxic.

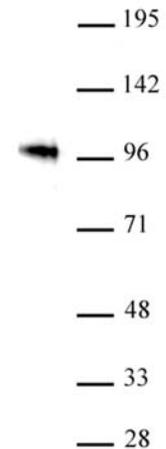
Application Notes:

Applications Validated by Active Motif:

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

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.



SOX6 antibody (pAb) tested by Western blot.
SOX6 antibody detection by Western blot. The analysis was performed using 20 μ g of mouse testes nuclear extract and SOX6 at a 1:500 dilution.