

Sox11 antibody (pAb)

Catalog Nos: 61181, 61182

RRID: AB_2793541

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 65 kDa

Background: Sox11 (SRY-box containing protein 11) is a member of a family of transcription factors containing an HMG DNA-binding domain. Sox11 is important for the development of nervous system. Sox11 is upregulated in a variety of tumor types, including medulloblastomas, gliomas and epithelial ovarian tumors. In B cell lymphomas, information regarding the expression of Sox11 aids in diagnosis and prognosis. Sox11 also functions as a tumor suppressor in glioma-initiating cells.

Immunogen: This Sox11 antibody was raised against a peptide from the N-terminus of human Sox11.

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

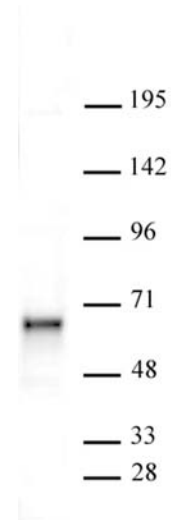
Application Notes:

Applications Validated by Active Motif:

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

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.



Sox11 antibody (pAb) tested by Western blot. Nuclear extract of Y-79 cells (20 µg) was probed with Sox11 pAb (1:500 dilution).