

PAX5 antibody (pAb)

Catalog Nos: 61565, 61566

RRID: AB_2614963

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 50 kDa

Background: PAX5 (paired box 5) is a member of the paired box family of transcription factors. PAX5 plays an important role in B-cell differentiation as well as neural development and spermatogenesis. Involved in the regulation of the CD19 gene, a B-lymphoid-specific target gene.

Immunogen: This antibody was raised against a peptide within the N-terminal region of human PAX5.

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

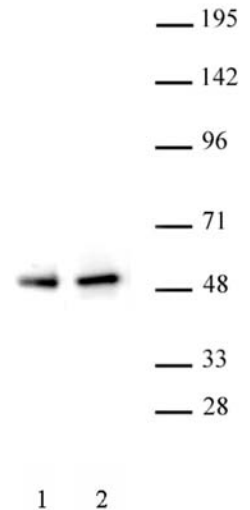
Application Notes:

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.



PAX5 antibody (pAb) tested by Western blot.
 Detection of PAX5 by Western blot. Lane 1: Raji nuclear cell extract (20 µg). Lane 2: Ramos nuclear cell extract (10 µg). Both probed with PAX5 antibody (pAb) at a 1:500 dilution.