

TCF7 antibody (pAb)

Catalog Nos: 61579, 61580

RRID: AB_2793685

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 52 & 55 kDa

Background: TCF7 (Transcription Factor 7) is a transcriptional activator involved in T-cell lymphocyte differentiation. Necessary for the survival of CD4(+) CD8(+) immature thymocytes. Isoforms lacking the N-terminal CTNNB1 binding domain cannot fulfill this role. Binds to the T-lymphocyte-specific enhancer element (5'-WWCAAAG-3') found in the promoter of the CD3E gene. May also act as feedback transcriptional repressor of CTNNB1 and TCF7L2 target genes. TLE1, TLE2, TLE3 and TLE4 repress transactivation mediated by TCF7 and CTNNB1.

Immunogen: This antibody was raised against a peptide within the N-terminal region of human TCF7.

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

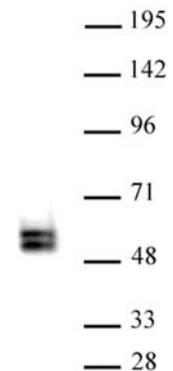
Application Notes:

Applications Validated by Active Motif:

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



TCF7 antibody (pAb) tested by Western blot.
TCF7 antibody detection by Western blot. The analysis was performed using 25 µg Molt-4 nuclear extract and TCF7 antibody at a 1:500 dilution.