

## TCF7L1 / TCF3 antibody (pAb)

**Catalog Nos:** 61125, 61126**RRID:** AB\_2793517**Application(s):** ChIP, ChIP-Seq, WB**Reactivity:** Human**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 80 kDa

**Background:** TCF3 / TCF7L1 (T-cell factor 3, TCF7L1) is a member of the TCF/LEF family and a component of the Wnt signaling pathway and a dominant downstream effector in embryonic stem cells (ESCs). TCF3 binds to DNA and serves as both a repressor as well as an activator of transcription. TCF3 brings developmental signals directly to the core regulatory circuitry of ES cells to influence the balance between pluripotency and differentiation. TCF3 transcriptionally represses many genes important for maintaining pluripotency and self-renewal, as well as those involved in lineage commitment and stem cell differentiation. This effect is in part mediated by the corepressors transducin-like enhancer of split 2 (TLE2) and C-terminal Binding Protein (CtBP).

**Immunogen:** This TCF7L1 / TCF3 antibody was raised against a peptide from the N-terminus of human TCF7L1 / TCF3.

**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:

ChIP: 10 µl per ChIP

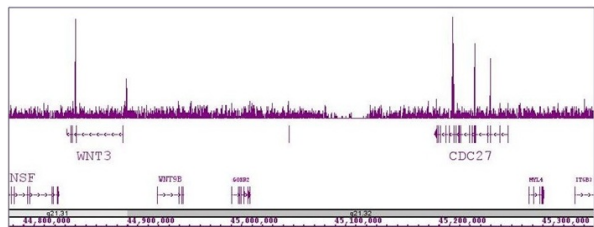
ChIP-Seq: 10 µl each

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

ChIP-Seq validation was performed by Active Motif's Epigenetics Services; the complete data set is available in the UCSC Genome Browser by clicking [here](#).

**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.



#### TCF7L1 / TCF3 antibody (pAb) tested by ChIP-Seq.

ChIP was performed using the ChIP-IT® High Sensitivity Kit (Cat. No. 53040) with 30 ug of chromatin from undifferentiated hESC cells and 7 µl of antibody. ChIP DNA was sequenced on the Illumina HiSeq and 26 million sequence tags were mapped to identify TCF7L1 / TCF3 binding sites. The image shows binding across a region of chromosome 17. You can view the complete data set in the UCSC Genome Browser, starting at this specific location, [here](#).



#### Western blot of TCF7L1 / TCF3 antibody.

Nuclear extract of PANC-1 cells (20 µg) probed with TCF7L1 / TCF3 antibody (1:500).