

## NKX2.5 antibody (pAb)

**Catalog Nos:** 61267, 61268

**RRID:** AB\_2793576

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 35 kDa

**Background:** The NKX2.5 (NK2 homeobox 5, CSX, Homeobox protein CSX, NKX2E, Homeobox protein NK-2 homolog E, NK2 transcription factor related, locus 5, CHNG5, cardiac-specific homeobox 1) gene encodes a homeobox-containing transcription factor that functions in heart formation and development. Alternative splicing results in multiple transcript variants. Implicated in commitment to and/or differentiation of the myocardial lineage. Acts as a transcriptional activator of ANF in cooperation with GATA4. Also interacts with JARID2 which represses its ability to activate transcription of ANF.

**Immunogen:** This NKX2.5 antibody was raised against a peptide within the N-terminal region of human NKX2.5.

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

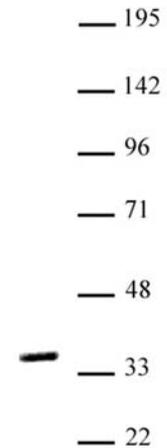
### Application Notes:

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.



### NKX2.5 antibody (pAb) tested by Western blot

Nuclear extract of human heart cells (20  $\mu$ g) probed with NKX2.5 antibody (1:1,000).