

Sox2 antibody (pAb)

Catalog Nos: 39823, 39824

RRID: AB_2793356

Isotype: Serum

Application(s): ICC, IF, IHC, WB

Reactivity: Mouse

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 35 kDa

Background: Sox2 (SRY related HMG BOX gene 2) is a DNA binding transcription factor and a member of the SOX family of proteins. SOX proteins have an HMG box that binds DNA. Sox2 forms a complex with Oct-4 and controls the expression of a number of genes involved in embryonic. Sox2 is critical for early embryogenesis and for embryonic stem cell pluripotency and thus can serve as a stem cell marker. Overexpression of Sox2 (along with Oct-4, KLF4 and c-Myc) can transform mouse fibroblasts into a state resembling embryonic stem cells (ES cells), referred to as Induced Pluripotency. Defects in Sox2 are the cause of microphthalmia syndromic type 3.

Immunogen: This Sox2 antibody was raised against a peptide corresponding to amino acids 111-140 mouse Sox2.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

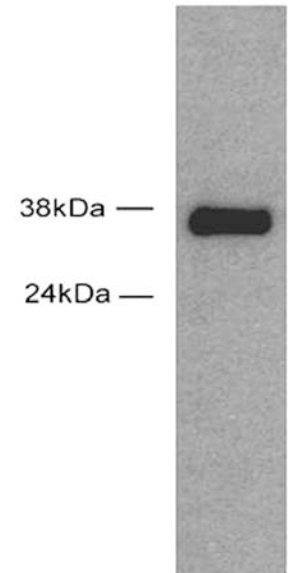
Application Notes:

Applications Validated by Active Motif:

WB: 1:2,000 - 1:10,000 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.



Sox2 antibody (pAb) tested by Western blot. HeLa cell extract expressing mouse Sox2 (50 µg per lane) probed with Sox2 antibody at a dilution of 1:5,000.