

## SOX2 antibody (rAb), 100 µg

**Catalog No:** 91729

**Clone:** 230225G1

**Application(s):** IF

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Protein A Chromatography

**Host:** Rabbit

**Isotype:** IgG

**Molecular Weight:** 34 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 antibody was raised against the SOX2 fusion protein.

**Buffer:** Purified IgG in PBS, pH 7.3, with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

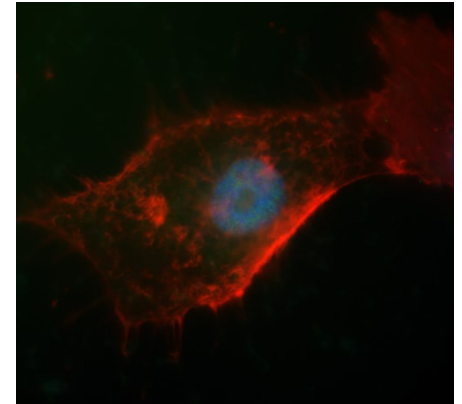
**Application Notes:**

Applications Validated by Active Motif:

IF: 1:150-1:600 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 (rAb) tested by immunofluorescence.** SOX2 was detected in U-251 cells using SOX2 antibody (rAb) at 1:300, followed by Multi-rAb CoraLite® Plus 488-Goat Anti-Rabbit Recombinant Secondary Antibody (H+L).