

SOX2 Polyclonal Antibody

(Catalog #A54440)

Background

Transcription factor that forms a trimeric complex with OCT4 on DNA and controls the expression of a number of genes involved in embryonic development such as YES1, FGF4, UTF1 and ZFP206 by similarity. Critical for early embryogenesis and for embryonic stem cell pluripotency. May function as a switch in neuronal development. Downstream SRRT target that mediates the promotion of neural stem cell self-renewal by similarity. Keeps neural cells undifferentiated by counteracting the activity of proneural proteins and suppresses neuronal differentiation

Description

SOX2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P48431

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transcription factor SOX-2 protein (1-317AA)

Storage

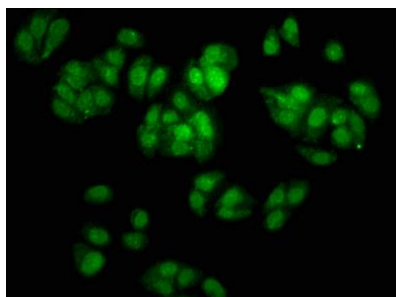
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

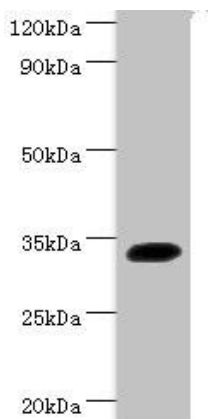
SOX2

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Immunofluorescence staining of HepG2 cells with SOX2 Polyclonal Antibody at 1:166, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).



Western blot

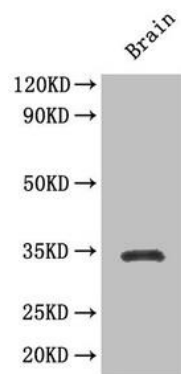
All lanes: Transcription factor SOX2 Polyclonal Antibody at 2µg/ml + Mouse kidney tissue

Secondary

Goat polyclonal to rabbit IgG at 1/15000 dilution

Predicted band size: 34 kDa

Observed band size: 34 kDa



Western Blot

Positive WB detected in: Mouse brain tissue

All lanes: SOX2 Polyclonal Antibody at 3.5µg/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 34 kDa

Observed band size: 34 kDa



Western Blot

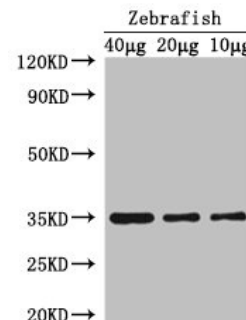
All lanes: Transcription factor SOX2 Polyclonal Antibody at 6µg/ml + 293T whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/15000 dilution

Predicted band size: 34 kDa

Observed band size: 34 kDa



Western Blot

Positive WB detected in: Zebrafish 40µg, 20µg, 10µg

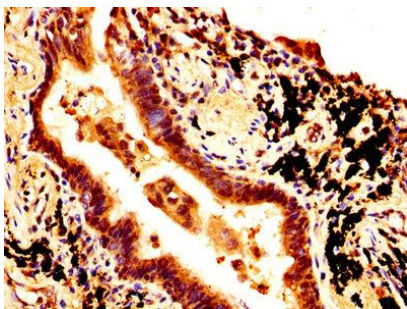
All lanes: SOX2 Polyclonal Antibody at 3µg/ml

Secondary

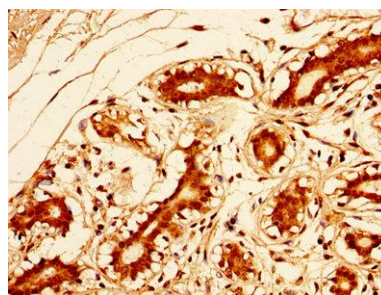
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 35 kDa

Observed band size: 35 kDa



IHC image of SOX2 Polyclonal Antibody diluted at 1:500 and staining in paraffin-embedded human lung cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



IHC image of SOX2 Polyclonal Antibody diluted at 1:500 and staining in paraffin-embedded human breast cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.