

---

## SOX2 Polyclonal Antibody, Biotin Conjugated

(Catalog #A54443)

---

### 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, Biotin Conjugated. Biotin. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### 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; Recommended dilution: ELISA 1:500-1:1000

---