

## FOXP1 antibody (mAb)

---

**Catalog Nos:** 61309, 61310

**RRID:** AB\_2615068

**Clone:** 1G1

**Isotype:** IgG1

**Application(s):** ChIP, EMSA, WB

**Reactivity:** Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein A Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 95, 70 & 50 kDa

---

**Background:** **FOXP1** (forkhead box P1) is a transcriptional repressor and belongs to subfamily P of the forkhead box (FOX) transcription factor family. It plays an important role in the specification and differentiation of lung epithelium. Can act with CTBP1 to synergistically repress transcription but CTPBP1 is not essential. Essential transcriptional regulator of B-cell development (By similarity).

**Immunogen:** This FOXP1 antibody was raised against a recombinant protein corresponding to amino acids 495-578 of mouse FOXP1. This antibody detects three isoforms.

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

**Application Notes:**

**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.