

## Tbpl1 antibody (mAb)

---

**Catalog Nos:** 61331, 61332

**RRID:** AB\_2793594

**Clone:** 3TRF-2A1

**Isotype:** IgG1k

**Application(s):** IHC, WB

**Reactivity:** Human, Mouse

**Quantity:** 100 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 0.69 µg/µl

**Molecular Weight:** 22 kDa

---

**Background:** **Tbpl1** (TATA box-binding protein-like protein 1) does not bind to the TATA box and initiates transcription from TATA-less promoters. This gene plays a critical role in spermatogenesis, and single nucleotide polymorphisms in this gene may be associated with male infertility. Alternatively spliced transcript variants have been observed for this gene, and a pseudogene of this gene is located on the long arm of chromosome 3.

**Immunogen:** This antibody was raised against a recombinant protein corresponding to mouse Tbpl1.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 32 mM glycine, 0.07 mM EDTA, 30% glycerol, and 0.035% sodium azide.

**Application Notes:**

Published Applications:

IHC

WB

See references for more information. Individual optimization may be required.

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.