

Tbp antibody (mAb)

Catalog No: 61329

RRID: AB_2793593

Clone: 3TF1-3G3

Isotype: IgG1k

Application(s): ICC, IF, IHC, IP, WB

Reactivity: Human, Mouse

Quantity: 100 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 0.73 µg/µl

Molecular Weight: 40 kDa

Background: Tbp (TATA-box-binding protein) is a general transcription factor that functions at the core of the DNA-binding multiprotein factor TFIID. Binding of TFIID to the TATA box is the initial transcriptional step of the pre-initiation complex (PIC), playing a role in the activation of eukaryotic genes transcribed by RNA polymerase II. Component of the transcription factor SL1/TIF-IB complex, which is involved in the assembly of the PIC (preinitiation complex) during RNA polymerase I-dependent transcription.

Immunogen: This antibody was raised against a recombinant protein corresponding to human Tbp.

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:

Validated Applications:

ICC/IF: 0.5 µg/ml dilution

Published Applications:

IP

ICC/IF

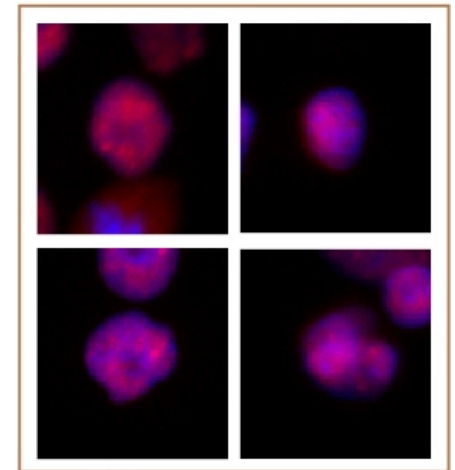
IHC

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.



Tbp antibody (mAb) tested by immunofluorescence.

Formaldehyde fixed HeLa cells stained with Tbp antibody at a 0.5 µg/ml dilution.