

Instruction Manual

Catalog Number	PK-AB718-2345
Synonyms	TP1 Antibody: TP1, Telomerase-associated protein 1, telomerase protein 1, TEP1
Description	Telomerase is an RNA-dependent DNA polymerase that uses an RNA component to add telomeric repeat sequences at the ends of chromosomes. Besides the RNA component which serves as the template that specifies the telomeric repeat, the telomerase complex contains a reverse transcriptase protein (TRT) and various accessory proteins including the telomerase-associated protein 1 (TP1). Telomerase activity is low in most somatic cells, causing the gradual shortening of telomeres which can ultimately lead to telomere fusion and cell death. High levels of telomerase activity are widely seen in cancerous cells and while recent experiments have suggested that telomerase may be a viable target in cancer therapy, expression levels of TP1 do not correlate with malignancy. At least two isoforms of TP1 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal TP1 antibody was raised against a 20 amino acid peptide from near the amino terminus of human TP1 (GenBank accession no. AAC51107).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. TP1 antibody can be used for detection of TP1 by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Greider CW. Telomere length regulation. Annu. Rev. Biochem. 1996; 274:92-7 Cong YS, Wright WE, and Shay JW. Human telomerase and its regulation. Microbiol. Mol. Biol. Rev. 2002; 66:407-25. Harrington L, McPhail T, Mar V, et al. A mammalian telomerase-associated protein. Science 1997; 275:973-7. Wong JMY and Collins K. Telomere maintenance and disease. Lancet 2003; 362:983-8.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2345P Human Kidney Tissue Lysate, Cat. No. PK-AB718-1305 TRT Antibody, Cat. No. PK-AB718-2341

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.