

TRBP1 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse TRBP1

Instruction Manual

Catalog Number	PK-AB718-4011
Synonyms	TRBP1 Antibody: TAR RNA binding protein 1, TARBP1, TRP-185, TRP-1
Description	The human trans-activation response (TAR) RNA binding protein 1 (TRBP1) was initially identified as a protein that binds to the HIV-1 TAR RNA and activates the long terminal repeat (LTR) expression in the absence and presence of the viral trans-activator Tat. This binding is enhanced by the presence of co-factors such as elongation factor 1alpha (EF-1alpha), polypyrimidine tract-binding protein (PTB), and the chaperonin-like protein stimulator of TAR RNA-binding proteins (SRB). TRBP1 may act to disengage RNA polymerase II from TAR during transcriptional elongation if the RNA polymerase stalls at the TAR during transcriptional elongation.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TRBP1 antibody was raised in rabbits against a 17 amino acid peptide from near the center of human TRBP1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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 Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. TRBP1 antibody can be used for detection of TRBP1 by Western blot at 1 - 2 µg/mL.
Images	Available upon request.
References	Gatignol A, Buckler-White A, Berkhout B, et al. Characterization of a human TAR RNA-binding protein that activates the HIV-1 LTR. <i>Science</i> 1991; 251:1597-1600. Wu-Baer F, Lane WS, and Gaynor RB. Identification of a group of cellular cofactors that stimulate the binding of RNA polymerase II and TRP-185 to human immunodeficiency virus 1 TAR RNA. <i>J. Biol. Chem.</i> 1996; 271:4201-8. Wu-Baer F, Lane WS, and Gaynor RB. The cellular factor TRP-185 regulates RNA polymerase II binding to HIV-1 TAR RNA. <i>EMBO J.</i> 1995; 14:5995-6009. (07-01D)
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4011P; TRBP1 Peptide Cat.No. PK-AB718-1212; 3T3 Cell Lysate

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