

TNPO3 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat TNPO3 (Transportin-3, Importin-12)

Instruction Manual

Catalog Number	PK-AB718-4597
Synonyms	TNPO3 Antibody; Transportin-3, Importin-12, IPO12, TRNSR, MTR10A, TRN-SR
Description	TNPO3 is a member of the importin beta/transportin family of nuclear import receptors and is specific for the essential splicing factors termed SR proteins, binding specifically and directly to the arginine/serine (RS)-rich domains of these proteins. Further experiments demonstrated that this interaction occurs only when these RS domains are phosphorylated. More recently, TNPO3 was identified as an HIV dependency factor (HDF) and that depletion of TNPO3 resulted in HIV inhibition after reverse transcription but before integration, suggesting that TNPO3 may be an important drug target in HIV treatment. At least three isoforms of TNPO3 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TNPO3 antibody was raised in rabbits against an 18 amino acid peptide from near the amino terminus of human TNPO3.
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. TNPO3 antibody can be used for detection of TNPO3 by Western blot at 1 - 2 µg/mL.
Images	Available upon request.
References	Kataoka N, Bachorik JL and Dreyfuss G. Transportin-SR, a nuclear import receptor for SR proteins. <i>J. Cell Biol.</i> 1999; 145:1145-52. Lai M-C, Lin R-I, Huang S-Y, et al. A human importin-b family protein, transportin-SR2, interacts with the phosphorylated RS domain of SR proteins. <i>J. Biol. Chem.</i> 2000; 275:7950-7. Brass AL, Dykxhoorn DM, Benita Y, et al. Identification of host proteins required for HIV infection through a functional genomic screen. <i>Science</i> 2008; 319:921-6.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4597P; TNPO3 Peptide Cat. No. PK-AB718-1207; Raji Cell Lysate

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