

# DDX3 (NT) antibody (pAb)

## Rabbit Anti-Human/Mouse Dead-box Polypeptide 3 X-linked (DBX)

### Instruction Manual

Catalog Number	PK-AB718-3757
Synonyms	DDX3 Antibody: Dead box polypeptide 3 X-linked, DBX, Helicase-like protein 2
Description	DDX3 contains all of the motifs of the DEAD-box family of RNA helicases, including the Asp-Glu-Ala-Asp sequence that gives the protein family its name and distinguishes it from other RNA helicases. DDX3 is localized to the X chromosome and has a highly conserved functional homolog (DBY) on the Y chromosome. DDX3 is thought to be involved in RNA splicing, RNA transport, and translation initiation. It has also been shown to be involved in cell growth control and is deregulated in hepatitis virus-associated hepatocellular carcinoma. Recent experiments suppressing DDX3 expression blocked HIV-1 RNA export from the nucleus, suggesting that DDX3 functions as a shuttling protein that transports the HIV-1 protein Rev and its cofactor CRM1 from the nucleus to the cytoplasm.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	DDX3 antibody was raised against a 16 amino acid peptide from near the amino terminus of human DDX3.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	DDX3 antibody will detect both DDX3 and DBY.
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, ICC, IF <span style="float: right;">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.</span> DDX3 antibody can be used for the detection of DDX3 by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Linder P, Lasko PF, Ashburner M, et al. Birth of the D-E-A-D box. Nature 1989; 337:121-2. Park SH, Lee SG, Kim Y, et al. Assignment of a human putative RNA helicase gene, DDX3, to human X chromosome bands p11.3-->p11.23. Cytogenet. Cell Genet. 1998; 81:178-9. Lahn BT and Page DC. Functional coherence of the human Y chromosome. Science 1997; 278:675-80. Abdelhaleem M. RNA helicases: regulators of differentiation. Clin. Biochem. 2005; 38:499-503.
Images	Available upon request.
Related Products	DDX3 Peptide, Cat. No. PK-AB718-3757P HepG2 Cell Lysate, Cat. No. PK-AB718-1211

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