

Instruction Manual

Catalog Number	PK-AB718-4551
Synonyms	Unc93a Antibody; Unc93a, homolog of <i>C. elegans</i> Unc93
Description	Unc93a, a human homolog of the <i>C. elegans</i> Unc93 gene, was initially identified by the sequencing of chromosome 6q27 at D6S193 as loss at this region correlated with sporadic ovarian cancer. Further analysis however, showed no tumor suppressor role for Unc93a. Unlike Unc93b, another homolog of the <i>C. elegans</i> Unc93 gene that is crucial for Toll-like receptor (TLR) signaling, the function of Unc93a is unknown. It is expressed in testis, small intestine, spleen, prostate and ovary. Two isoforms of Unc93a are known to exist. This antibody will not cross-react with Unc93b.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Unc93a antibody was raised in rabbits against an 18 amino acid peptide from near the carboxy terminus of human Unc93a.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	Two isoforms of Unc93a are known to exist. This antibody will not cross-react with Unc93b.
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 I Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Unc93a antibody can be used for the detection of Unc93a by Western blot at 0.5 - 1 µg/mL.
Images	Available upon request.
References	Liu Y, Dodds P, Emilion G, et al. The human homolog of unc-93 maps to chromosome 6q27 – characterisation and analysis in sporadic epithelial ovarian cancer. <i>BioMed Central Genetics</i> 2002; 3:20. Massey DC and Parkes M. Genome-wide association scanning highlights two autophagy genes, ATG16L1 and IRGM, as being significantly associated with Crohn's disease. <i>Autophagy</i> 2007; 3:649-51.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4551P; Unc93a Peptide Cat.No. PK-AB718-1201; HeLa Cell Lysate

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