

Unc93b antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Unc93b

Instruction Manual

Catalog Number	PK-AB718-4553
Synonyms	Unc93b Antibody; Unc93b1, homolog of <i>C. elegans</i> Unc93
Description	The endoplasmic reticulum (ER) protein Unc93b, a human homolog of the <i>C. elegans</i> Unc93 gene, was initially identified by a forward genetic screen using N-ethyl-N-nitrosourea where a histidine-to-arginine substitution in Unc93b caused defects in Toll-like receptor (TLR) 3, 7 and 9 signaling. Unlike Unc93a, another homolog of the <i>C. elegans</i> Unc93 gene whose function is unknown, Unc93b specifically interacts with TLR3, 7 and 9; the histidine-to-arginine point mutation used to identify Unc93b abolishes this interaction. Mice carrying this point mutation are highly susceptible to infection with a number of viruses, indicating that Unc93b plays an important role in innate immunity. Multiple isoforms of Unc93a are known to exist. This antibody will not cross-react with Unc93a.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Unc93b antibody was raised in rabbits against a 19 amino acid peptide from near the amino terminus of human Unc93b.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	Multiple isoforms of Unc93a are known to exist. This antibody will not cross-react with Unc93a.
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. Unc93b antibody can be used for the detection of Unc93b by Western blot at 0.5 - 1 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	Tabeta K, Hoebe K, Janssen EM, et al. The Unc93b1 mutation 3d disrupts exogenous antigen presentation and signaling via Toll-like receptors 3, 7, and 9. <i>Nat. Immunol.</i> 2006; 7:156-64. 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. Brinkmann MM, Spooner E, Hoebe K, et al. The interaction between the ER membrane protein UNC93B and TLR3, 7, and 9 is crucial for TLR signaling. <i>J. Cell Biol.</i> 2007; 177:265-75.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4553P; Unc93b Peptide Cat.No. PK-AB718-1301; Human Heart Tissue Lysate

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