

LIMP2 antibody (pAb)

Rabbit Anti-Human/Mouse LIMP2 (Lysosomal integral membrane protein 2)

Instruction Manual

Catalog Number	PK-AB718-4655
Synonyms	LIMP2 Antibody: Lysosomal integral membrane protein 2, CD36L2, Scavenger receptor class B member 2, SR-B2, SCARB2
Description	The lysosomal integral membrane protein 2 (LIMP2) is a heavily glycosylated type III transmembrane protein, the majority of which exists in the lumen of the lysosome and a cytoplasmic domain of approximately 20 amino acids. A deficiency of LIMP2 in mice causes uretic pelvic junction obstruction, deafness, and peripheral neuropathy associated with impaired vesicular trafficking and distribution of apically expressed proteins. More recently, LIMP2 was shown to act as a receptor to bind b-glucocerebrosidase, the enzyme defective in Gaucher disease, a lysosomal storage disorder. LIMP2-deficient mice showed missorted as well as secreted b-glucocerebrosidase, suggesting that LIMP2 also functions as the mannose-6-phosphate-independent trafficking receptor. Despite its predicted molecular weight, LIMP2 runs at approximately 80 – 85 kDa in SDS-PAGE.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	LIMP2 antibody was raised in rabbits against an 18 amino acid peptide from near the carboxy terminus of human LIMP2.
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, IHC Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. LIMP2 antibody can be used for detection of LIMP2 by Western blot at 1 and 2 µg/mL. Despite its predicted molecular weight, LIMP2 runs at approximately 80 - 85 kDa in SDS-PAGE. Antibody can also be used for immunohistochemistry starting at 10 µg/mL.
Images	Available upon request.
References	Fujita H, Saeki M, Yasunaga K, et al. Isolation and sequencing of a cDNA clone encoding 85kDa sialoglycoprotein in rat liver lysosomal membranes. Biochem. Biophys. Res. Commun. 1991; 178:444-52. Gamp A, Tanaka Y, Lullmann-Rauch R, et al. LIMP-2/LGP85 deficiency causes uretic pelvic junction obstruction, deafness and peripheral neuropathy in mice. Hum. Mol. Genet. 2003; 12:631-46. Knipper M, Claussen C, Ruttiger L, et al. Deafness in LIMP2-deficient mice due to early loss of the potassium channel KCNQ1/KCNE1 in marginal cells of the stria vascularis. J. Physiol. 2006; 576:73-86. Reczek D, Schwake M, Schroder J, et al. LIMP-2 is a receptor for lysosomal mannose-6-phosphate-independent targeting of b-glucocerebrosidase. Cell 2007; 131:770-83.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4655P; LIMP2 Peptide

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