

# LIMP2 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat LIMP2 (Lysosomal integral membrane protein 2, CD36L2)

## Instruction Manual

|                     |  |
|---------------------|--|
| Catalog Number      | PK-AB718-4621  |
| 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 approx. 80–85 kDa in SDS-PAGE. |
| Quantity            | 100 µg   |
| Source / Host       | Rabbit   |
| Immunogen           | LIMP2 antibody was raised in rabbits against a 16 amino acid peptide from near the center of human LIMP2.  |
| 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, IHC, IF<br>INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.<br><br>LIMP2 antibody can be used for detection of LIMP2 by Western blot at 1 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 2.5 µg/mL. For immunofluorescence start at 20 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.<br>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.<br>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.<br>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-AB718-4621P; LIMP2 Peptide; Cat. No. PK-AB718-1375; Human Skeletal Muscle Tissue Lysate   |

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