

Slc22A17 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Slc22A17 (Solute carrier family 22 member 17)

Instruction Manual

Catalog Number	PK-AB718-4651
Synonyms	Slc22A17 Antibody: Solute carrier family 22 member 17, Brain-type organic cation transporter, BOCT, BOIT
Description	The Slc22 family of organic anion and cation transporters (OATs, OCTs, OCTNs) are transmembrane proteins expressed predominantly in kidney and liver. Each contain 12 predicted alpha-helical transmembrane domains (TMDs) and one large extracellular loop between TMDs 1 and 2. Transporters of the SLC22 family function in different ways such as uniporters that mediate facilitated diffusion in either direction (OCTs), as anion exchangers (OAT1, OAT3 and URAT1), and as Na(+)/l-carnitine cotransporter (OCTN2). Slc22 family members participate in the absorption and/or excretion of drugs, xenobiotics, and endogenous compounds in intestine, liver, and kidney, and perform homeostatic functions in brain and heart. Mutations in the Slc22 family may cause specific diseases such as primary systemic carnitine deficiency or idiopathic renal hypouricemia and may change drug absorption or excretion. Recent studies show the expression of Slc22A17 as receptor for Lipocalin 2 is relatively high in hematopoietic stem cells.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Slc22A17 antibody was raised in rabbits against a 14 amino acid peptide near the carboxy terminus of the human Slc22A17.
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, IF Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Slc22A17 antibody can be used for detection of Slc22A17 by Western blot at 1 µg/mL. Antibody can also be used for immunofluorescence starting at 20 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Koepsell H and Endou H. The SLC22 drug transporter family. <i>Pflugers Arch.</i> 2004; 447:666-76. Rizwan AN and Burckhardt G. Organic anion transporters of the SLC22 family: biopharmaceutical, physiological, and pathological roles. <i>Pharm. Res.</i> 2007; 24:450-70. Miharada K, Hiroyama T, Sudo K, et. al. Lipocalin 2-mediated growth suppression is evident in human erythroid and monocyte/macrophage lineage cells. <i>J. Cell Physiol.</i> 2008; 215:526-37. Devireddy LR, Gazin C, Zhu X, et al. A cell-surface receptor for lipocalin 24p3 selectively mediates apoptosis and iron uptake. <i>Cell</i> 2005; 123:1293-305.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4651P; Slc22A17 Peptide Cat. No. PK-AB718-1220; SK-N-SH Cell Lysate

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