

Instruction Manual

Catalog Number	PK-AB718-4651P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human Slc22A17
Formulation	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 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	Store Slc22A17 peptide at -20°C, stable for one year.
Application	Slc22A17 peptide is used for blocking the activity of Slc22A17 antibody.
References	Koepsell H and Endou H. The SLC22 drug transporter family. Pflugers Arch. 2004; 447:666-76.

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