

Instruction Manual

Catalog Number	PK-AB718-3915P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of avian influenza NS1
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 Avian Influenza Nonstructural Protein 1 peptide at -20°C, stable for one year.
Application	Avian influenza nonstructural protein 1 peptide is used for blocking the activity of avian influenza nonstructural protein 1 antibody.
References	Krug RM, Yuan W, Noah D, et al. Intracellular warfare between human influenza viruses and human cells: the role of the viral NS1 protein. <i>Virology</i> 2003; 309:181-9. Aivazian D, Serrano RL, and Pfeffer S. TIP47 is a key effector for Rab9 localization. <i>J. Cell Biol.</i> 2006; 173:917-26. Ganley IG, Carroll K, Bittova L, et al. Rab9 GTPase regulates late endosome size and requires effector interaction for its stability. <i>Mol. Biol. Cell</i> 2004; 15:5420-30.

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