

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4333P
<b>Quantity</b>	50 µg
<b>Source</b>	15 amino acids near the amino terminus of H5N1 influenza A M2
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	Store Avian Influenza A M2 peptide at -20°C, stable for one year.
<b>Application</b>	Avian Influenza A M2 peptide is used for blocking the activity of M2 antibody.
<b>References</b>	Lamb RA, Holsinger LJ, and Pinto LH. The influenza A virus M2 ion channel protein and its role in the influenza virus life cycle. In Cellular receptors of animal viruses. Edited by Wimmer E. Cold Spring Harbor, Cold Spring Harbor Laboratory; 1994; 303-321.

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