

Instruction Manual

Catalog Number	PK-AB718-3925P
Quantity	50 µg
Source	14 amino acids near the middle of avian influenza hemagglutinin 1
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 Hemagglutinin 1 peptide at -20°C, stable for one year.
Application	Avian influenza hemagglutinin 1 peptide is used for blocking activity of avian influenza hemagglutinin 1 antibody. It usually blocks antibody activity completely in ELISA assays by incubating peptide with equal volume of antibody for 30 min at 37°C.
References	Shortridge KF, Zhou NN, Guan Y, et al. Characterization of avian H5N1 influenza viruses from poultry in Hong Kong. <i>Virology</i> . 1998; 252:331-342.

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