

# H5N1 NS2 / NEP antibody (pAb)

## Rabbit Anti-Nonstructural Protein 2 (Nuclear Export Protein)

### Instruction Manual

Catalog Number	PK-AB718-3917
Synonyms	Avian Influenza Nonstructural Protein 2 Antibody; Avian Influenza Nonstructural Protein 2, H5N1 NS2, Nonstructural protein 2, nuclear export protein, NEP
Description	Influenza A virus is a major public health threat, killing more than 30,000 people per year in the USA. Novel influenza virus strains caused by genetic drift and viral recombination emerge periodically to which humans have little or no immunity, resulting in devastating pandemics. Influenza A can exist in a variety of animals; however, it is in birds that all subtypes, including the so-called "avian flu" or H5N1, can be found. These subtypes are classified based on the combination of the virus coat glycoproteins hemagglutinin (HA) and neuraminidase (NA) subtypes. One of the less studied proteins encoded by the influenza virus is the nonstructural protein (NS) 2. NS2 binds to the influenza matrix protein M1 that is bound to the ribonucleoprotein (RNP) complex and mediates the contact between the M1/RNP complex and the cellular exportin CRM, but does not interact with nucleosomes.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	NS2 antibody was raised against a synthetic peptide corresponding to 14 amino acids near the middle of the NS2 protein. Efforts were made to use relatively conserved regions of the viral sequence as the antigen.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Virus
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 Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Avian Influenza Nonstructural Protein 2 antibody can be used for the detection of the Avian Influenza Nonstructural Protein 2 protein from influenza A in ELISA.
Images	NA
References	Thompson WW, Shay DK, Weintraub, et al. Mortality associated with influenza and respiratory syncytial virus in the United States. JAMA 2003; 289:179-186. Alexander DJ. A review of avian influenza. Proceedings of the European Society for Veterinary Virology (ESVV) Symposium on Influenza Viruses of Wild and Domestic Animals. Vet. Microbiol. 2000; 74:3-13. Neumann G, Hughes MT and Kawaoka Y. Influenza A virus NS2 protein mediates vRNP nuclear export through NES-independent interaction with hCRM1. EMBO J. 2000; 19:6751-8. Garcia-Robles I, Akarsu H, Muller CW, et al. Interaction of influenza virus proteins with nucleosomes. Virology 2005; 332:329-36.
Images	NA
Related Products	H5N1 NS1 Antibody, Cat. No. PK-AB718-3915; Hemagglutinin Antibody, Cat. No. PK-AB718-3425; Hemagglutinin Antibody, Cat. No. PK-AB718-3427; Neuraminidase Antibody, Cat. No. PK-AB718-3423; Neuraminidase Antibody, Cat. No. PK-AB718-3421

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