

Instruction Manual

Catalog Number	PK-AB913-216
Quantity	50 µg
Description	Influenza A virus subtype H5N1 is a subtype of the Influenza A virus. Subtype H5N1 is commonly known as avian influenza or bird flu. H5N1 may cause more than one influenza pandemic as it is expected to continue mutating in birds. The dominant strain of HPAI A (H5N1) evolved creating the Z genotype. It has also been called Asian lineage HPAI/A/H5N1 which is divided into 2 antigenic clades. "Clade 1" includes human and bird isolates from Vietnam, Thailand, and Cambodia and bird isolates from Laos and Malaysia. "Clade 2" viruses were first identified in bird isolates from China, Indonesia, Japan, and South Korea before spreading westward to the Middle East, Europe, and Africa. Hybridoma clones have been derived from hybridization of Sp2/0 myeloma cells with spleen cells of Balb/c mice immunized with purified avian influenza virus type A H5N1.
Source / Host	Mouse
Immunogen	Influenza A hemagglutinin H5N1
Purification Method	Protein A affinity chromatography
Clone / IgG Subtype	clone IA-H5N1; Mouse IgG1
Specificity	Specifically recognizes Influenza A hemagglutinin H5N1.
Formulation	Sterile-filtered solution in PBS (pH 7.4) with 0.1 % NaN3
Reconstitution	Please Note: Always centrifuge product briefly before opening vial.
Storage & Stability	Shipped at 4°C. Stable at 4°C for up to 2 years.
Applications	Influenza A haemagglutinin H5N1 immunodetection in direct or indirect ELISA, and Western Blotting. Optimal dilution has to be determined by user. Antibody might also be suited for other applications not tested so far.

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