

Instruction Manual

Catalog Number	PK-AB913-214
Quantity	50 µg
Description	H1N1 is a subtype species of Influenza A virus. H1N1 Influenza Virus has mutated into various strains such as the Spanish Flu strain, mild human flu strains, endemic pig strains, and various strains found in birds. The Influenza A Virus is a globular particle of about 100 nm in diameter, sheathed in a lipid bilayer derived from the plasma membrane of its host. Studded in the lipid bilayer are two integral membrane proteins: some 500 molecules of hemagglutinin ("H") and some 100 molecules of neuraminidase ("N"). Within the lipid bilayer are 3000 molecules of matrix protein and 8 pieces of RNA. Each of the 8 RNA molecules is associated with many copies of a nucleoprotein, several molecules of the three subunits of its RNA polymerase some "non-structural" protein molecules of uncertain function. Hybridoma clones have been derived from hybridization of Sp2/0 myeloma cells with spleen cells of Balb/c mice immunized with Influenza A/NewCaledonia/20/99 H1N1 derived from allantoic fluid of 10 days old embryonated eggs.
Source / Host	Mouse
Immunogen	Influenza A hemagglutinin H1N1
Purification Method	Protein A affinity chromatography
Clone / IgG Subtype	clone IA-H1N1; Mouse IgG1
Specificity	Specifically recognizes Influenza A hemagglutinin H1N1.
Formulation	Sterile-filtered solution (3.8 mg/ml) in PBS (pH7.4) with 0.1 % NaN3
Reconstitution	Please Note: Always centrifuge product briefly before opening vial.
Storage & Stability	Shipped at 4°C. Store at 4°C. Do not freeze.
Applications	Influenza A haemagglutinin 1 H1N1 immunodetection in direct or indirect ELISA, sandwich immunoassay, Western Blotting. Recommended pairs for Influenza A H1N1 in sandwich immunoassay (coating – conjugate): All pairs detect H1N1 of Influenza A as well as recombinant H1N1 with high specificity and do not interact with haemagglutinin H3N2. Antibody might also be suited for other applications not tested so far.

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