

Instruction Manual

Catalog Number	PK-AB913-217
Quantity	50 µg
Description	Influenza-B virus is a genus in the virus family Orthomyxoviridae. The only species in this genus is called "Influenza-B virus". Influenza-B virus only infects humans and seals. This limited host range is apparently in contrast with those caused by the similar Influenza-A virus as both mutate by both genetic drift and reassortment. Influenza-B virus evolves slower than A viruses and faster than C viruses. Influenza-B virus mutates at a rate 2-3 times lower than type A. However, Influenza-B mutates enough that lasting immunity is not possible. The Influenza-B virus capsid is enveloped while its virion consists of a matrix protein + envelope + nucleoprotein complex + nucleocapsid, and a polymerase complex. Influenza-B is sometimes spherical and sometimes filamentous. Its 500 or so surface projections are made of hemagglutinin and neuraminidase. The Influenza B virus is 14648 nucleotides long and consists of eight segments of linear negative-sense, single-stranded RNA. The multipartite genome is encapsidated, each segment in a separate nucleocapsid, and the nucleocapsids are surrounded by one envelope. Hybridoma has been derived from hybridization of Sp2/O myeloma cells with spleen cells of Balb/c mice immunized with purified influenza virus type B strain B/Tokio/53/99.
Source / Host	Mouse
Immunogen	Influenza-B virus
Purification Method	Protein A affinity chromatography
Clone / IgG Subtype	clone PIB-633-HY; Mouse IgG1
Specificity	Specifically recognizes Influenza-B.
Formulation	Sterile-filtered solution (4.6 mg/ml) in PBS (pH 7.4) with 0.1% NaN ₃
Reconstitution	Please Note: Always centrifuge product briefly before opening vial.
Storage & Stability	Shipped at 4°C. Store at 4°C.
Applications	Western blotting, ELISA; can also be used in indirect immunofluorescence. 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.