



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Certificate of Analysis and Data Sheet

Influenza A Strain A/ Beijing/262/95 (H1N1)

Catalog No.
MD-05-0110P

Source: Allantoic fluid of 10 day old
embryonated egg

Description

H1N1 is a subtype specie 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 about 100nm 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.

Preparation

Allantoic fluid of 10 days old embryonated eggs, inoculated with influenza A virus, strain A/Beijing/262/95. The Influenza Virus was purified by Ultracentrifugation with 10-40 % sucrose gradient.

Formulation

Sterile Filtered colorless solution

The H1N1 A/Beijing/262/95 solution contains STE, 0.1 % sodium azide (NaN_3) and 0.005 % thimerosal. Greater than 90.0% purity as determined by SDS-PAGE.

Protein Concentration: 0.8mg/ml (BCA method)

Inactivation

Thimerosal and beta propiolactone treatment

This product has been treated in a manner consistent with methods of inactivation. Generally accepted good laboratory practices appropriate to microbiological/viral safe handling practices and techniques are required when handling this product.

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Storage

H1N1 A/Beijing/262/95 although stable 14°C for 4 weeks, should be stored desiccated below -18°C.
Please prevent freeze-thaw cycles.

Application

Serological studies of influenza A virus, immunogen for antibody production, has been tested with monoclonal antibodies in ELISA. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not necessarily be excluded.

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**