



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

Mouse Anti Influenza-A Hemagglutinin H1N1

Catalog No.
IP-14-335

Isotype
IgG1

Introduction

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 are 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.

Description

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.

Preparation

Product Type: Mouse Monoclonal Antibody

Immunogen: Influenza A hemagglutinin H1N1.

Purification Method: Protein-A column.

Formulation

Shipping Condition: Antibody is shipped in liquid form with ice packs.

Protein Concentration: (7.3mg/ml) In PBS, pH 7.4 and 0.1 % NaN3.

Storage

For long periods, store at 4°C. Stable for two years at 4°C.

**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

Applications

Table Summary of antibody applications and working conditions

Options Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA(direct/indirect)	.			
Western Blotting	.			
Flow Cytometry			.	
Immunoprecipitation			.	
Immunohistochemistry			.	

Note: 1. Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.

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