



## 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](http://www.raybiotech.com)  
Email: [info@raybiotech.com](mailto:info@raybiotech.com)

---

# Certificate of Analysis and Data Sheet

## Influenza A Strain A/Wisconsin/67/05 (H3N2)

**Catalog No.**  
MD-14-0305P

### ***Description***

Influenza A Virus, Strain A/Wisconsin/67/05 (H3N2)

### ***Preparation***

**Purification:** >90% pure (SDS-PAGE). Ultracentrifugation with 10-40% sucrose gradient

**Source:** Allantoic fluid of 10 day old embryonated eggs, inoculated with influenza A virus  
**Strain:** A/Wisconsin/67/05

### ***Formulation***

**Format:** Purified, Liquid

**Buffer:** 0.05M Tris-HCl, pH 8.0 containing 0.1M Sodium chloride, 5mM EDTA

**Preservatives:** 0.1% Sodium azide, 0.005% Thimerosal

**Approx Protein Concentration:** Lot Specific

### ***Application***

Serological studies of Influenza A virus, immunogen for antibody production. 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.

### ***Inactivation***

Thimerosal and beta propiolactone treatment

**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](http://www.raybiotech.com)  
Email: [info@raybiotech.com](mailto:info@raybiotech.com)

---

### **Storage**

Store at -20°C. Avoid multiple freeze/thaw cycles.

### **Warnings**

This product contains sodium azide, which has been classified as Xn (Harmful) in European Directive 67/548/EEC in the concentration range of 0.1 – 1.0 %. When disposing of this reagent through lead or copper plumbing, flush with copious volumes of water to prevent azide build-up in drains.

No test guarantees a product to be non-infectious. All materials should be handled as if potentially infectious. Generally accepted laboratory practices appropriate for infectious materials should be employed when handling this product.

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