



## 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

## Parainfluenza Virus Type-1

**Catalog No.**  
IP-05-028P

### ***Introduction***

Parainfluenza virus, PIV, contains a group of four distinct serotypes of single-stranded RNA viruses belonging to the paramyxovirus family and two subtypes (4a and 4b). Parainfluenza virus is a negative-sense, single-stranded RNA virus that possess fusion and hemagglutinin-neuraminidase glycoprotein "spikes" on their surface. The virion varies in size (average diameter between 150 and 300 nm) and shape.

### ***Description***

Allantoic fluid of 10 days old embryonated eggs, inoculated with parainfluenza-1 virus, strain Sendai. The Parainfluenza-1 Virus was purified by Ultracentrifugation with 20-60 % sucrose gradient.

### ***Physical Appearance***

Sterile Filtered colorless solution.

### ***Formulation***

The Parainfluenza type-1 solution contains 0.1M NaCl, 10mM Tris Hcl, 1mM EDTA pH-8, and 0.1 % sodium azide (NaN<sub>3</sub>).

### ***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 at work.

### ***Immunological Activity***

Serological studies of Parainfluenza virus, immunogen for antibody production.

### ***Purity***

Greater than 90.0% as determined by:  
Analysis by SDS-PAGE.

### ***Stability***

Parainfluenza type-1 although stable 4°C for 4 weeks, should be stored desiccated below -18°C.  
**Please prevent freeze-thaw cycles.**

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