



## 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

## Rat Anti Mouse IFN-gamma

**Catalog No.**  
IP-01-166

**Immunogen**  
r.mouse IFN-gamma

**Isotype:**  
rat IgG1

**Accession No:**  
NM\_000619

### Description

Interferon (IFN)- $\gamma$  is an antiviral and antiparasitic agent produced by CD4<sup>+</sup>/CD8<sup>+</sup> lymphocytes and natural killer cells that undergo activation by antigens, mitogens or alloantigens. IFN- $\gamma$  production modulates T cell growth and differentiation and inhibits the growth of B cells. Synthesis of IFN- $\gamma$  is inducible by IL-2, FGF and EGF. The active form of IFN- $\gamma$  is a homodimer with each subunit containing six helices. The dimeric structure of human IFN- $\gamma$  is stabilized by non-covalent interactions through the interface of the helices. IFN- $\gamma$  translated precursor is 166 amino acids, including the 23 amino acid secretory sequence. Multiple forms exist due to variable glycosylation and under non-denaturing conditions due to dimers and tetramers.

### Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Direct ELISA	•			1:10,000
Western Blotting	•			
Immunohistology - frozen			•	
Immunohistology -paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation	•			
Flow Cytometry			•	
Immunofluorence staining			•	
Neutralization			•	
Intracellular Staining	•			

Note: 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.**



## 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)

---

### **Purification Method:**

Ion exchange column

### **Protein Concentration:**

1 mg/ml in PBS (after reconstitution).

### **Storage Procedure:**

Lyophilized: store at 4°C. After reconstitution, if not used within a month, aliquot and store at 20°C.

### **Reconstitution:**

Reconstitute with of H<sub>2</sub>O. Mix gently, wash the sides of the vial and wait 30-60 seconds before use.

### **Stability:**

Two years lyophilized, one month in solution at 4°C.

### **Reference:**

Young, H.A., et al. 1995. Role of IFN- $\gamma$  in immune cell regulation. J. Leukoc. Biol. 58: 373-381.

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