



## 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

## RABBIT ANTI MOUSE INTERLEUKIN-13

### With HRP-conjugated Secondary Antibody

**Catalog No.**  
DS-PB-01088

**Species**  
Mouse

**Isotype**  
Rabbit IgG

### **Preparation**

**Immunogen:** Recombinant murine IL-13

**Purification:** Antisera to mouse IL-13 were raised by repeated immunisations of rabbits with highly purified antigen. Purified IgG was prepared from whole serum by affinity chromatography.

### **Specificity**

DS-PB-01088 is specific for mouse Interleukin-13 (IL-13), a 12.3kDa immunoregulatory cytokine produced primarily by activated T helper type 2 (Th2) cells and also by mast cells and NK cells. The biological properties of IL-13 include stimulation of B cell proliferation and antibody production and studies suggest that IL-13 plays an important role in the induction of the pathophysiological features of allergic asthma, independently of Immunoglobulin E and eosinophils (1).

### **Formulation**

**Product Form:** Purified IgG - lyophilised

**Buffer Solution:** Phosphate buffered saline pH7.4

**Reconstitution:** Reconstitute with 0.1 ml distilled water

For long term storage, the addition of 0.09% sodium azide is recommended.

**Approx. Protein Concentrations:** IgG concentration 1.0 mg/ml

### **Storage**

12 months shelf life.

Prior to reconstitution store at +4°C.

Following reconstitution store at -20°C.

This product should be stored undiluted. Storage in frost-free freezers is not recommended. Should this product contain a precipitate we recommend microcentrifugation before use.

**Please avoid freeze-thaw cycles.**

**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 0.5ug/ml
Western Blotting	•			0.1-0.2ug/ml
Immunohistology - frozen			•	
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Functional Assays			•	
Immunofluorescence staining			•	
Neutralization			•	

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

### Secondary Antibody Applications

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Immunoassay (ELISA, Western blot)	•			1:5000-1:10000

### Reference

1. Wills-Karp, M. et al. (1998) Interleukin-13: central mediator of allergic asthma. Science. 282: 2258-2261.

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