



## 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 CHICKEN INTERLEUKIN-1 BETA:Azide Free With HRP-Conjugated Second Antibody

Catalog No.	Species	Isotype
DS-PB-01075	Chicken	Polyclonal IgG

### Preparation

**Preparation:** Antisera to chicken Interleukin 1 Beta were raised by repeated immunizations of rabbits with highly purified antigen. Purified IgG prepared by affinity chromatography on protein G

**Immunogen:** Recombinant chicken IL-1

### Specificity

This product recognizes chicken Interleukin 1 beta (chIL-1beta), which shares homology with mammalian IL-1beta.

IL-1 is a major regulator in both innate and adaptive immune responses, plays an important role in the regulation of acute inflammation and has also been implicated in many other processes, including both fever and fibrinogen induction. The previous names by which IL-1 has been known includes lymphocyte activating factor, endogenous pyrogen, catabolin, hemopoietin-1, melanoma growth inhibition factor, and osteoclast activating factor, reflecting its widespread cell expression.

Studies using chicken models have indicated physiological similarities between chIL-1beta and mammalian IL-1beta in relation to its neural distribution and several factors have been shown to induce chIL-1beta secretion, including protein kinase C (PKC), Calmodulin-dependent protein kinase, lipopolysaccharide, phorbol esters, calcium and silica.

It neutralizes chicken IL-1 beta (1-2).

### Formulation

**Product Type:** Polyclonal Antibody

**Product Form:** Purified IgG - Liquid

**Buffer Solution:** Phosphate buffered saline

**Preservative Stabilizers:** None present

**Approx. Protein Concentrations:** IgG concentration 0.5 mg/ml

**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

18 months from date of dispatch.

Store at -20°C only.

Storage in frost-free freezers is not recommended. This product should be stored undiluted.

Avoid repeated freezing and thawing as this may denature the antibody. Should this product contain a precipitate we recommend microcentrifugation before use.

### Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA	•			
Western Blotting(*)	•			
Immunohistology - frozen			•	
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Immunofluorescence			•	
Functional Assays	•			

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.  
(\*) detects a band of approximately 29kDa in HD-11 cell lysates transfected with chIL-1.

### Secondary Antibody Applications

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

**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)

---

### **Reference**

1. Staeheli, P. et al. (2001) Cytokines of Birds: Conserved functions – A largely different look. *J. Interferon and Cytokine Research* 21: 993-1010.
2. Gyorfy, Z. et al. (2003) Truncated chicken interleukin-1beta with increased biologic activity. *J. Interferon and Cytokine Research* 23: 223-228.
3. Weining, K.C. et al. (1998) A chicken homolog of mammalian interleukin-1 beta: cDNA cloning and purification of active recombinant protein. *Eur. J. Biochem.* 258: 994-1000.
4. Bombara, C.J. et al. (1991) Signal transduction events in chicken interleukin-1 production. *Poult. Sci.* 70: 1372-1380.

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