



## 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 BOVINE INTERLEUKIN-1 BETA:Azide Free With HRP-conjugated Secondary Antibody

Catalog No.	Species	Isotype
DS-PB-01074	Bovine	IgG

### Preparation

**Preparation:** Purified IgG prepared by affinity chromatography on Protein G

**Immunogen:** Recombinant bovine IL-1 beta.

### Specificity

DS-PB-01074 is specific for bovine interleukin-1 beta (IL-1 beta).

IL-1 beta is principally expressed by macrophages, and plays an important immunomodulatory role in the bovine immune response.

DS-PB-01074 recognizes both native and recombinant bovine IL-1 beta. No detectable cross-reactivity was observed with human, mouse or rat IL-1 beta.

### Formulation

**Product Type:** Polyclonal Antibody

**Product Form:** Purified IgG conjugated to Biotin - liquid

**Preservative Stabilizers:** 0.09% Sodium Azide

**Buffer Solution:** Phosphate buffered saline

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

### Storage

18 months from date of dispatch.

Store at 4°C or -20°C if preferred.

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.**

**Avoid repeated freezing and thawing as this may denature the antibody.**

**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)

### Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA	•			1ug/ml - 10ug/ml
Western Blotting			•	
Immunohistology - frozen			•	
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
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

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