



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

Certificate of Analysis and Data Sheet

RABBIT ANTI BOVINE INTERLEUKIN-1 BETA: Biotin With HRP-conjugated Secondary Antibody

Catalog No.	Species	Isotype
DS-PB-01073	Bovine	IgG

Preparation

Preparation: Preparation Purified IgG prepared by affinity chromatography on Protein G
Immunogen: Recombinant bovine IL-1 beta.

Specificity

This product 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.
It 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
 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	•			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.**