



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 RAT INTERLEUKIN-1 BETA:Biotin With HRP-conjugated Secondary Antibody

Catalog No.	Species	Isotype
DS-PB-01069	Rat	Rabbit IgG

Preparation

Antisera to rat IL-1 beta were raised by repeated immunisations of rabbits with highly purified antigen. Purified IgG was prepared from whole serum by affinity chromatography.

Immunogen

Recombinant rat IL-1 beta.

Specificity

AAR15B recognizes rat IL-1 beta.

Formulation

Product Type Polyclonal Antibody

Product Form Purified IgG conjugated to Biotin - lyophilized

Preservative Stabilisers None Present

Buffer Solution Phosphate buffered saline

Approx. Protein Concentrations IgG concentration 50 ug/ml

Reconstitution Reconstitute with 1.0 ml sterile PBS containing 0.1% Bovine Serum Albumin

Storage

12 months from date of reconstitution.

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.

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

**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)	.			0.15ug/ml - 0.30ug/ml
Western Blotting	.			0.1ug/ml - 0.2ug/ml
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.**