



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 SHEEP INTERLEUKIN-1 BETA

With HRP-conjugated Secondary Antibody

Catalog No.	Species	Isotype
DS-PB-01071	Sheep	IgG

Preparation

Preparation: Antisera to ovine IL-1 beta were raised by repeated immunisation of rabbits with highly purified antigen.

Immunogen: Recombinant IL-1 beta

Specificity

This product recognizes ovine interleukin-1 beta (IL-1 beta). No cross-reactivity is seen with ovine IL-6, IL-8, MCP-1 or TNF alpha.

It neutralises the bioactivity of ovine IL-1 beta. Removal of sodium azide is recommended prior to use in functional assays.

Formulation

Product Type: Polyclonal Antibody

Product Form: Serum - liquid

Preservative Stabilizers: 0.09% Sodium Azide

Storage

18 months from date of despatch.

Store at +4°C or at -20°C if preferred.

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	.			1/500
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

Reference

Rothel, J.S. et al. (1997) Analysis of ovine IL-1 beta production in vivo and in vitro by EIA and immunohistochemistry. *Vet. Immunol. Immunopathol.* 57: 267-278.

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