



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 MOUSE INTERLEUKIN-1 ALPHA

With HRP-conjugated Secondary Antibody

Catalog No.
DS-PB-01060

Species
Mouse

Isotype
Rabbit IgG

Preparation

Immunogen: Recombinant mouse IL-1 alpha

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

Specificity

DS-PB-01060 is specific for mouse interleukin-1 alpha (IL-1 alpha).

Interleukin-1 alpha is a potent immuno-modulator, which mediates a wide range of immune and inflammatory responses.

Formulation

Product Form: Purified IgG - lyophilised

Buffer Solution: Phosphate buffered saline pH7.4

Reconstitution: Reconstitute with 0.1 ml distilled water.

For long term storage, the addition of 0.09% sodium azide is recommended.

For functional studies do not add azide.

Approx. Protein Concentrations: IgG concentration 1.0 mg/ml

Storage

12 months shelf life.

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. Should this product contain a precipitate we recommend microcentrifugation before use.

Please avoid freeze-thaw cycles.

**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 0.5ug/ml
Western Blotting	•			0.1ug/ml-0.2ug/ml
Immunohistology - frozen			•	
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Functional Assays	•			0.03ug/ml-0.05ug/ml
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.**