



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

Anti-Human Interleukin-1 beta Antibody

Catalog No.

RB-IP-01-101

Species:

Human

Isotype:

Mouse Ig G

Description

IL-1 production is generally thought to be associated with inflammation, but it has been shown to be expressed during kidney development, thymocyte differentiation and cartilage degradation. IL-1 plays a critical role in the regulation of immune response and inflammation, acting as an activator of T and B lymphocytes and natural killer (NK) cells. In T cells, IL-1 stimulates the production of IL-2 and selectively inhibits IL-4 expression. IL-1 induces B cell proliferation and maturation, and immunoglobulin synthesis. NK cells require IL-1 β for production of the anti-pathogen IFN- γ . IL-1 has also been implicated in several pathological conditions including rheumatoid arthritis, inflammatory bowel disease and atherosclerosis.

Applications

Table Summary of antibody applications and working conditions

Options Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA	*			For Sandwich-based ELISA and membrane-based assay, detecting level is up to 1pg/ml when it is used for capture
Western Blotting	*			0.1-0.2 μ g/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.

**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

Preparation

Immunogen was recombinant protein derived from human IL-1 beta. This antibody was produced from a hybridoma resulting from the fusion of a mouse myeloma with B cells obtained from a mouse immunized with the immunogen. The IgG fraction of tissue culture supernatant was purified by Protein G affinity chromatography.

Specificity

This antibody was selected for its ability to detect human IL-1beta in Sandwich ELISA and Array. In ELISA, this antibody shows approximately 5% cross-reactivity with human IL-RI, and no cross-reactivity with other tested antibodies.

Reconstitution

Supplied as 100 µg of lyophilized purified antibody. It can be reconstituted with sterile PBS before use. **Please avoid freeze-thaw cycles.**

Storage

Store at -20⁰C, if not intended for use within a month. Keep it at 4⁰C and minimize freezing and thawing when use.

Reference

P.E., Webb, A.C., Rosenwasser, L.J., Mucci, S.F., Rich, A., Wolff, S.M. and Dinarello, C.A. 1985.
Nucleotide sequence of human monocyte interleukin-1 precursor cDNA. Proc. Natl. Acad. Sci.
USA 81: 7907-7911.

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