



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-2

With HRP-conjugated secondary antibody

Catalog No.

DS-PB-01101

Target Species

Mouse

Isotype

IgG

Preparation

Immunogen: Recombinant mouse IL-2

Purification: Antisera to mouse IL-2 were raised by repeated immunizations of rabbits with highly purified antigen. Purified IgG was prepared from whole serum by affinity chromatography.

Specificity

DS-PB-01101 is specific for mouse Interleukin-2 (IL-2), a potent lymphoid cell growth factor which primarily acts on T lymphocytes, in response to antigenic or mitogenic stimulation.

IL-2 stimulates the growth/differentiation of monocytes, macrophages, B lymphocytes, NK cells, and lymphokine activated killer cells which are used in experimental cancer therapy.

DS-PB-01101 may be used in an indirect ELISA or as a capture antibody in a sandwich ELISA with DS-PB-01102 as the detection reagent.

Formulation

Product Form: Purified IgG – lyophilized

Reconstitution: Reconstitute with 0.1 ml distilled water. Care should be taken during reconstitution as the protein may appear as a film at the bottom of the vial. RayBiotech recommends that the vial is gently mixed after reconstitution. For long term storage the addition of 0.09% sodium azide is recommended.

*For functional studies do not add sodium azide

Product Type: Polyclonal Antibody

Buffer Solution: Phosphate buffered saline pH7.4

Preservative Stabilizers: None Present

Approx. Protein Concentrations: IgG concentration 1.0 mg/ml after reconstitution

**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

Storage

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.

Shelf Life: 12 months from date of dispatch.

References

1. Smith, K.A. (1998) Interleukin-2: inception, impact, and implications. *Science*. 240: 1169-1176.

Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA	.			0.5 - 2.0ug/ml
Western Blotting	.			0.1 - 0.2ug/ml
Immunohistology - frozen			.	
Immunohistology – paraffin			.	
Immunohistology - resin			.	
Functional Assays	.			0.04 - 0.05ug/ml

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)	.			1:1000-1:2000

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