

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-7:Biotin

With HRP-conjugated Secondary Antibody

Catalog No.
DS-PB-01137

Target Species
Mouse

Isotype
Polyclonal IgG

Preparation

Antiserum Preparation: Antisera to mouse IL-7 were raised by repeated immunizations of rabbits with highly purified antigen. Purified IgG was prepared from whole serum by affinity chromatography.

Immunogen: Recombinant mouse IL-7

Specificity

DS-PB-01137 is specific for mouse interleukin-7 (IL-7).

The biological activity of IL-7 includes stimulating the proliferation of pre-B cells, pro-B cells and activated T cells and inducing the synthesis of other inflammatory mediators, such as IL-1, IL-6 and MIP.

Formulation

Product Type: Polyclonal Antibody

Product Form: Purified IgG conjugated to Biotin – lyophilized

Buffer Solution: Phosphate buffered saline

Preservative Stabilizers: None present

Approx. Protein Concentrations: IgG concentration 50 ug/ml

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

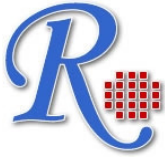
Storage

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 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

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			•	
Immunofluorence staining			•	
Functional Assays			•	

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.**