



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 RAT INTERLEUKIN-10

With HRP-conjugated Secondary Antibody

Catalog No.	Species	Isotype
DS-PB-01078	Rat	Rabbit IgG

Preparation

Immunogen: Recombinant rat IL-10

Antisera to rat IL-10 were raised by repeated immunizations of rabbits with highly purified antigen. Purified IgG was prepared from whole serum by affinity chromatography.

Specificity

DS-PB-01078 recognizes rat interleukin-10 (IL-10). Rat IL-10 is a 18.7kD cytokine which exists as a homodimer.

Formulation

Product Form: Purified IgG - lyophilized

Buffer Solution: Phosphate buffered saline pH7.4

Approximate IgG Concentration: IgG concentration 1.0 mg/ml

Preservative / Stabilizer: none present

Reconstitution: Reconstitute with 0.1 ml distilled water

Storage

Prior to reconstitution, store at 4°C. After 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	•			0.5 µg/ml
Western Blotting	•			0.1 - 0.2 µg/ml
Immunohistology - frozen			•	
Immunohistology – paraffin*			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Immunofluorescence staining			•	
Functional Assays*	•			0.5 - 0.9µg/ml

Note: Other applications are not tested yet. Optimal dilutions should be determined by each laboratory for each application.

*For functional studies do not add sodium azide.

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.**