



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

Rat Anti-Mouse Interleukin-10 (rAmuIL-10)

Catalog No.
IP-01-113

Immunogen
r.Murine IL-10

Isotype:
Rat IgM

Accession No:
NM_000572

Description

Interleukin 10, or IL-10, is a 178 amino acid protein that is primarily secreted by TH2 clones. IL-10 has dual functions, the first of which is the suppression of cytokine production by TH1 clones responding to antigen presented by monocyte and macrophage antigen presenting cells (APCs). The second function consists of the inhibition of response of cytokine targeted cells, possibly by the downregulation of CD25 (the interleukin-2 receptor) on macrophages and B lymphocytes. Human and murine IL-10 exhibit 81% sequence identity at the amino acid level, and share 73% identity at the cDNA level. Both human and murine IL-10 are acid-labile and exist as non-covalently-linked homodimers in solution. IL-10 exerts its biological activity through the IL-10 receptor (IL-10R), a 560 amino acid, glycoprotein whose expression can be induced in cultured macrophages and fibroblasts by lipopolysaccharide (LPS) stimulation. IL-10 expression has been shown to be elevated in HIV-1 infected individuals and has been implicated in the progression of the disease.

Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Direct ELISA	•			1:5,000 dilution
Western Blotting	•			
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

Purification Method:

Boric Acid precipitation

Protein Concentration:

1 mg/ml in PBS (after reconstitution)

Storage Procedure:

Lyophilized: store at 4 C. After reconstitution, if not used within a month, aliquot and store at 20 C.

Reconstitution:

Reconstitute with 0.5 ml of H₂O. Mix gently, wash the sides of the vial and wait 30-60 seconds before use.

Stability:

Two years lyophilized, one month in solution at 4 C.

Activity:

This is a neutralizing antibody both in vitro and in vivo and will block the activity of IL-10 in a variety of assays

Reference:

Feng, L., Tang, W.W., Chang, J.C. and Wilson, C.B. 1993. Molecular cloning of rat cytokine synthesis inhibitory factor (IL-10) cDNA and expression in spleen and macrophages. *Biochem. Biophys. Res. Commun.* 192: 452-458.

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