



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-4 (rAmuIL-4)

Catalog No.
IP-01-108

Immunogen
r.Murine IL-4

Isotype:
Rat IgG1

Accession No:
NM_172348

Description

Interleukin-4 (IL-4), also designated B cell stimulatory factor-1, is a glycosylated cytokine secreted by activated T lymphocytes, basophils and mast cells. The secreted IL-4 protein promotes the growth and differentiation of cells that participate in immune defense by favoring such events as the expansion of the Th2 lineage relative to Th1 cells. These T helper cell subsets are defined by their pattern of cytokine secretion: Th1 cells secrete IL-2, TNF β and IFN- γ , while Th2 cells secrete IL-4, IL-5 and IL-10. Another key immunological function of IL-4 is to induce immunoglobulin class switching. IL-4 has been shown to induce the production of IgE and enhance IgG4 secretion by B cells, but suppress the production of IgM, IgA, IgG1, IgG2 and IgG3. It has been determined that Stat6 is rapidly tyrosine phosphorylated following stimulation of IL-3 or IL-4, but is not detectably phosphorylated following stimulation with IL-2, IL-12 or erythropoietin.

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			•	
Immunofluorence 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:

Ion exchange column

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.

Reference:

Kamogawa, Y., et al. 1993. The relationship of IL-4- and IFN γ -producing T cells studied by lineage ablation of IL-4-producing cells. Cell 75: 985-995.

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