



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

Mouse Anti-Human Mouse Anti Human Interleukin-7 (mAHuIL-7)

Catalog No.
IP-01-110

Immunogen
r.Human IL-7

Isotype:
Mouse IgG1

Accession No:
NM_000880

Description

Interleukin-7 (IL-7) was originally described as a factor capable of inducing in vitro proliferation of pre-B cells from marrow cultures. The IL-7 gene encodes a protein 177 amino acids in length, with a molecular weight of 17.5 kDa. IL-7 exerts its biological function through the IL-7 receptor, which is expressed on pre-B cells, thymocytes and bone marrow-derived macrophages. The IL-7 receptor is composed of an IL-7 receptor-specific chain and the IL-2 receptor γ chain common to the IL-2, IL-4, IL-7, IL-9 and IL-15 receptors. IL-7 stimulation leads to the activation of Janus tyrosine kinase family members JAK1 and JAK3. Other studies have shown that in T cells, the IL-7 receptor-specific chain associates with the Src kinases family Lck and Fyn. IL-7 induces phosphorylation of insulin receptor substrate-1 (IRS-1) and insulin receptor substrate- 2 (IRS-2), originally called 4PS.

Applications

Table Summary of antibody applications and working conditions

Options Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Direct ELISA	•			1:10,000 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

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:

Goodwin, R.G., Lupton, S., Schmierer, A., Hjerrild, K.J., Jerzy, R., Clevenger, W., Gillis, S., Cosman, D. and Namen, A.E. 1989. Human interleukin-7: molecular cloning and growth factor activity on human and murine B-lineage cells. Proc. Natl. Acad. Sci. USA 86: 302-306.

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