



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 Interleukin-2

Catalog No.
IP-01-102

Isotype
IgG2a

Synonyms

T-cell growth factor (TCGF), Aldesleukin, Lymphokine, IL-2.

Introduction

IL2 is a secreted cytokine that is important for the proliferation of T and B lymphocytes. The receptor of this cytokine is a heterotrimeric protein complex whose gamma chain is also shared by interleukin 4 (IL4) and interleukin 7 (IL7). The expression of this gene in mature thymocytes is monoallelic, which represents an unusual regulatory mode for controlling the precise expression of a single gene. The targeted disruption of a similar gene in mice leads to ulcerative colitis-like disease, which suggests an essential role of this gene in the immune response to antigenic stimuli.

Preparation

Product Type: Mouse Monoclonal Antibody
Immunogen: Recombinant Human IL-2.
Purification Method: Ion Exchange

Formulation

Shipping Condition: Antibody is shipped lyophilized at ambient temperature.
Protein Concentration: 1mg/ml in PBS (after reconstitution)
Reconstitution: Reconstitute with sterile H₂O. Mix gently, wash the sides of the vial and wait 30-60 seconds before use.

Titer

By direct ELISA: 1:10,000 dilution will yield 0.4 O.D using alkaline phosphatase conjugated rabbit anti-mouse Ig (Jackson Laboratories). 1:50 dilution will block 70-80% of CTLL proliferative response to 50 units of r.human IL-2.

Crossreactivity

A certain degree of functional cross reactivity with murine IL-2 was reported by some investigators.

**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

Storage

In lyophilized form, for long periods, store at 4°C in a dry environment. After reconstitution, if not intended for use within a month, aliquot and store at -20°C.

Stable for two years lyophilized, one month in solution at 4°C.

Applications

Table Summary of antibody applications and working conditions

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA(direct)	.			
Western Blotting	.			
Block T Cell Proliferation	.			
Immunoprecipitation	.			
Immunohistochemistry	.			

Note: 1. 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.**