



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 Interferon Alpha, Western blot

Catalog No.
IP-14-308

Isotype
IgG1

Introduction

IFN-alpha is produced by macrophages, IFN-alpha has antiviral activities and stimulates the production of two enzymes: a protein kinase and an oligoadenylate synthetase.

Preparation

Product Type: Mouse Monoclonal Antibody
Immunogen: Recombinant Human IFN-Alpha.
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 0.5 O.D using alkaline phosphatase conjugated rabbit anti-mouse Ig (Jackson Laboratories).

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.

**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(direct) (1)	.			
Western Blotting		.		
Block T Cell Proliferation	.			
Immunoprecipitation	.			
Immunohistochemistry	.			

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

(1) The antibody can be used for Capture ELISA together with RayBiotech's Mouse Anti Human IFN-alpha Neutralizing antibody.

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