



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-gamma (IFN-g Antibody)

Catalog No.
IP-01-123

Species
Human

Isotype
Mouse IgG2a

Synonyms

Immune Interferon, type II interferon, T cell interferon, MAF, IFNG, IFG, IFI, IFN-gamma.

Introduction

IFN-gamma produced by lymphocytes activated by specific antigens or mitogens. IFN-gamma, in addition to having antiviral activity, has important immunoregulatory functions, it is a potent activator of macrophages, and has antiproliferative effects on transformed cells and it can potentiate the antiviral and antitumor effects of the type I interferons.

Preparation

Immunogen: r.Human IFN-g

Purification: Ion exchange

Applications

Table Summary of antibody applications and working conditions

Options	YES	NO	Not determined	Recommended Work dilution or concentration
Functions				
ELISA (Direct)	.			
Western Blotting	.			
Immunoprecipitation	.			
Immunohistochemistry	.			
Intracellular staining.	.			
Flow Cytometry			.	

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.**



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

Formulation

Product Form: Mouse Anti Human Monoclonal

Shipping Conditions: Antibody is shipped lyophilized at ambient temperature.

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

Protein Concentrations: 1mg/ml in PBS (after reconstitution)

Titer

This antibody loses its antigen binding capacity when immobilized on plastic. Biological activity is determined by virus neutralization assay and MHC class II induction.

Storage

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

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.

Avoid multiple freeze/thaw cycles.

Centrifuge before opening to ensure complete recovery of vial contents.

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