



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 IP-10

With HRP-conjugated Secondary Antibody

Catalog No.
DS-MB-01718

Species
Human

Isotype
Mouse IgG2a

Preparation

Immunogen: Recombinant human IP-10

Purification: Purified IgG prepared by affinity chromatography on Protein G from tissue culture supernatant

Fusion Partners: Spleen cells from immunised Balb/c mice were fused with cells of the mouse P3X63-AgP-653 myeloma cell line

Specificity

DS-MB-01718 recognizes both natural and recombinant human IP-10, also known as CXCL10.

Formulation

Product Form: Purified IgG - liquid

Buffer Solution: Phosphate buffered saline pH7.4

Preservative & Stabilizers: 0.09% Sodium Azide

Approx. Protein Concentrations: IgG concentration 1.0 mg/ml

Storage

18 months shelf life. Store at +4°C or at -20°C if preferred.

This product should be stored undiluted. Storage in frost-free freezers is not recommended. Should this product contain a precipitate we recommend microcentrifugation before use.

Please avoid freeze-thaw cycles.

**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	•			1ug/ml-2ug/ml
Western Blotting	•			1/1000
Immunohistology - frozen			•	
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Immunofluorescence staining			•	
Neutralization			•	

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

Secondary Antibody Applications

Options / Functions	YES	NO	Not determined	Recommended Work dilution or concentration
Immunoassay (ELISA, Western blot)	•			1:5000-1:10000

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