



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

RABBIT ANTI HUMAN IP-10

With HRP-conjugated Secondary Antibody

Catalog No.
DS-PB-01174

Species
Human

Isotype
Rabbit IgG

Preparation

Immunogen: Recombinant human IP-10

Purification: Antisera to human IP-10 were raised by repeated immunisations of rabbits with highly purified antigen. Purified IgG prepared by affinity chromatography.

Specificity

DS-PB-01174 is specific for human IP-10 (10-kDa interferon-inducible protein), otherwise known as CXCL10, a 77 amino acid CXC chemokine which, like monokine induced by IFN-gamma (Mig) and IFN-inducible T-cell alpha chemoattractant (I-TAC), signals through the CXCR3 receptor.

IP-10 acts as a chemoattractant for Th1 lymphocytes and natural killer (NK) cells and is secreted by a variety of cells, including neutrophils, monocytes, keratinocytes and epithelial cells, in response to interferon gamma (IFNgamma) and lipopolysaccharide (LPS). IP-10 is thus associated with several clinical conditions linked with Th1 type immune responses and raised IFNgamma expression, such as psoriasis, sarcoidosis and viral meningitis. In addition, studies have demonstrated the ability of IP-10 to suppress the synergistic action of stimulatory cytokines on hematopoietic progenitor cell proliferation

Formulation

Product Form: Purified IgG - lyophilised

Buffer Solution: Phosphate buffered saline pH7.4

Reconstitution: Reconstitute with 0.1ml distilled water. For long term storage the addition of 0.09% sodium azide is recommended. For functional studies do not add azide.

Approx. Protein Concentrations: IgG concentration 1.0 mg/ml

**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

12 months shelf life.

Prior to reconstitution store at +4°C.

After reconstitution store at -20°C.

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.

Applications

Table Summary of antibody applications and working conditions

Options Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA (1)	•			0.5ug/ml
Western Blotting	•			0.1-0.2ug/ml
Immunohistology - frozen			•	
Immunohistology - paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Functional Assays	•			2.0-3.0ug/ml
Immunofluorescence staining			•	
Neutralization			•	

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

(1) Direct

**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

Secondary Antibody Applications

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

Reference

1. Cox, M.A. et al. (2001) Human interferon-inducible 10-kDa protein and human interferon-inducible T cell alpha chemoattractant are allotypic ligands for human CXCR3: Differential binding to receptor states. *Mol. Pharmacol.* 59: 707-715.
2. Sauty, A. et al. (1999) The T cell-specific CXC chemokines IP-10, Mig and I-TAC are expressed by activated human bronchial epithelial cells. *J. Immunol.* 162: 3549-3558.

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