



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-8

Catalog No.
RB-IP-01-111

Immunogen
Human IL-8

Isotype:
Mouse IgG1

Accession No:
NM_000584

Description

Interleukin-8, or IL-8, the prototypic member of the C-X-C, or α , family of chemokines, is a chemoattractant cytokine involved in the chemotaxis and activation of neutrophils. IL-8 expression has been correlated to a large number of chronic inflammatory diseases, including inflammatory bowel disease (IBD) and atherosclerosis. IL-8 is cleaved from a 99 amino acid precursor to a 72 amino acid, nonglycosylated, and biologically active protein. Whether IL-8 functions in vivo as a monomer or a hydrogen-bonded dimer has not yet been resolved. Two IL-8 receptors, designated IL-8RA and IL-8RB, have been described and share 77% sequence identity. Both are seven-transmembrane domain proteins (7TMD), similar to the G protein-coupled receptors, and, in addition to IL-8, serve as receptors for other members of the α and β chemokine families.

Applications

Table Summary of antibody applications and working conditions

Options Functions	YES	NO	Not determined	Recommended Work dilution or concentration
ELISA	•			For Sandwich-based ELISA and membrane-based assay, detecting level is up to 1pg/ml when it is used for capture
Western Blotting	•			0.1-0.2 μ g/ml
Immunohistology - frozen			•	
Immunohistology -paraffin			•	
Immunohistology - resin			•	
Immunoprecipitation			•	
Flow Cytometry			•	
Immunofluorence staining			•	
Neutralization			•	

Note: 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

Preparation

Immunogen was recombinant protein derived from human IL-8. This antibody was produced from a hybridoma resulting from the fusion of a mouse myeloma with B cells obtained from a mouse immunized with the immunogen. The IgG fraction of tissue culture supernatant was purified by Protein G/A affinity chromatography.

Specificity

This antibody was selected for its ability to detect human IL-8 in Sandwich ELISA and Array. In ELISA, this antibody shows approximately 5% cross-reactivity with human IL-1b, and no cross-reactivity with murine IL-8 and other tested proteins.

Reconstitution

Supplied as 100 µg of lyophilized purified antibody. It can be reconstituted with sterile PBS before use. **Please avoid freeze-thaw cycles.**

Storage Procedure

Lyophilized: store at 4°C. After reconstitution, if not used within a month, aliquot and store at 20°C.

Stability

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

Reference

Knall, C., Young, S., Nick, J.A., Buhl, A.M., Worthen, G.S., and Johnson, G.L. 1996. Interleukin-8 regulation of the Ras/Raf/mitogen-activated protein kinase pathway in human neutrophils. *J. Biol. Chem.* 271: 2832-2838.

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