



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

Rat Anti Mouse B220 (CD 45R)

Catalog No.	Species	Isotype
IP-11-139	Mouse	Rat IgM

Synonyms

LCA, PTPRC, LY5, B220, CD45R, T200, GP180, PROTEIN-TYROSINE PHOSPHATASE, RECEPTOR-TYPE C

Type

Rat Anti Mouse Monoclonal

Introduction

B220 (CD45R) is a member of the protein tyrosine phosphatase (PTP) family. PTPs are known to be signaling molecules that regulate a variety of cellular processes including cell growth, differentiation, mitotic cycle, and oncogenic transformation. This PTP contains an extracellular domain, a single transmembrane segment and two tandem intracytoplasmic catalytic domains, and thus belongs to receptor type PTP. This gene is specifically expressed in hematopoietic cells. This PTP has been shown to be an essential regulator of T- and B-cell antigen receptor signaling. It functions through either direct interaction with components of the antigen receptor complexes, or by activating various Src family kinases required for the antigen receptor signaling. This PTP also suppresses JAK kinases, and thus functions as a regulator of cytokine receptor signaling. Four alternatively spliced transcripts variants of this gene, which encode distinct isoforms, have been reported.

B220 (CD45R) is required for t-cell activation through the antigen receptor. The first ptpase domain has enzymatic activity, while the second one seems to affect the substrate specificity of the first one

Immunogen

Purified mouse LN B cells

Purification Method

Ion exchange column

**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

Protein concentration

1mg/ml in PBS (after reconstitution)

Storage Procedures

Lyophilized: store at 4°C. After reconstitution, if not intended for use within a month, aliquot and store at -20°C.

Reconstitution

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

Stability / Shelf Life

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

Antibody Properties

Staining antibody. For staining, use 10 µl/10⁶ cells.

Available Conjugates

This antibody is only available as purified antibody.

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