

ANTI-MOUSE CD45R/B220

**Conjugated Rat Monoclonal Antibodies Against Mouse CD45R/B220
Clone RA3-6B2**

FITC-conjugate:	Catalog # 10711	100 mg
	Catalog # 10712	500 mg
PE-conjugate:	Catalog # 10811	100 mg
	Catalog # 10812	200 mg

SPECIFICITY:

Antibody RA3-6B2 recognizes an epitope on the extracellular domain of CD45 which is expressed on B lymphoid cells, at all stages from pro-B to mature B-cells, with lower expression on plasma cells and memory B cells. CD45R/B220 is also expressed on NK cells.

CLONE: RA3/6B2

ISOTYPE: IgG_{2a},κ (rat)

FORMAT:

FITC-conjugate: Catalog # 10711: 100 µg in 0.2 mL (0.5 mg/mL) diluent
Catalog # 10712: 500 µg in 1.0 mL (0.5 mg/mL) diluent
PE-conjugate: Catalog # 10811: 100 µg in 0.5 mL (0.2 mg/mL) diluent
Catalog # 10812: 200 µg in 1.0 mL (0.2 mg/mL) diluent
Diluent: Aqueous buffered solution, containing 0.09% (w/v) sodium azide.

STABILITY AND STORAGE:

Store at 4°C. Do not freeze. Product is stable for at least 6 months.

APPLICATIONS AND DIRECTIONS FOR USE:

Flow cytometry:

Recommended amount per 1x10⁶ cells in a volume of 100 µL:

- FITC-conjugate (Cat No. 10711,10712): ≤ 1 µg (2 µL)
- PE-conjugate (Cat No. 10811,10812): ≤ 1 µg (5 µL)

Appropriate conditions should be established for each application.

Cell separation:

Positive selection of murine CD45R⁺ cells with StemCell's reagents for immunomagnetic cell separation. Please contact us for more information.

**NOT FOR CLINICAL USE; INCLUDING IN VITRO DIAGNOSTIC USE, AND EX VIVO OR IN VIVO THERAPEUTIC USE IN CLINICAL TRIALS OR IN CLINICAL PRACTISE.
THIS REAGENT IS FOR RESEARCH ONLY**

Hazardous Ingredient: Sodium Azide. *Avoid exposure to skin and eyes, ingestion and contact with heat, acids and metals. Wash exposed skin with soap and water. Flush eyes with water. Dilute with running water before discharging into plumbing.*

REFERENCES:

1. Coffman RL, Weissman IL, B220: a B cell-specific member of the T200 glycoprotein family, 1981, Nature 289: 681-683.
2. Coffman RL, Surface antigen expression and immunoglobulin gene rearrangement during mouse pre-B cell development, 1982, Immunol Rev 69:5-23.
3. Johnson P, Greenbaum L, Bottomly K, Trowbridge IS. Identification of the alternatively spliced exons of murine CD45 (T200) required for reactivity with B220 and other T200-restricted antibodies, 1989, J Exp Med 169: 1179-1184.
4. Watanabe Y, Akaike T. Activation signal induces the expression of B cell-specific CD45R epitope (6B2) on murine T cells, 1994, Scand J Immunol 39:419-25.
5. Rolink A, ten Boekel E, Melchers F, Fearon DT, Krop I, Andersson J. A subpopulation of B220+ cells in murine bone marrow does not express CD19 and contains natural killer cell progenitors. J Exp Med. 1996, 183: 187-194.