

Anti-Human CD32
 Monoclonal Antibody
 Clone IV.3

Catalog #01470 0.1 mg



SCIENTIFIC BACKGROUND:

CD32 (Fc γ RII) is a 40,000 D receptor for the Fc region of Immunoglobulin G (IgG). CD32 is encoded by three genes, A, B, and C, and is expressed as 6 protein isoforms, IIa1, IIa2, IIb1, IIb2, IIb3 and IIc. All isoforms are expressed on monocytes/macrophages, the IIb isoform is also expressed on B cells, IIa on platelets, granulocytes and, weakly, on B cells. Isoform IIc is expressed on NK cells. CD32 binds weakly to the Fc portion of monomeric IgG, but efficiently to IgG aggregates and immune complexes. The Fc/FcR interactions may result in nonspecific staining in antibody-based detection and cell separation experiments.

SPECIFICITY:

The mouse monoclonal antibody binds to CD32 (Fc γ RII) on human cells.

CLONE: IV.3

ISOTYPE: IgG_{2b}

PREPARATION:

The antibody was generated in a CAF1 mouse immunized with the K562 cell line.

FORMAT:

1 mg/mL purified antibody in phosphate buffered saline.

STABILITY AND STORAGE:

Store at 2-8°C in undiluted aliquots. Stable for at least 12 months when kept sterile. Repeated freezing and thawing is not recommended.

APPLICATIONS AND DIRECTIONS FOR USE:

Centrifuge tube briefly before use to ensure recovery of entire contents.

Flow cytometry:

This antibody blocks the binding of the Fc region of IgG antibodies to CD32 (Fc γ RII). In flow cytometry experiments, it can be used to stain CD32⁺ cells or to block undesired CD32 / IgG Fc interactions. Recommended amount per 1x10⁶ cells in a volume of 100 μ L is <1 μ g (1 μ L). Larger volumes or numbers of cells may require more antibody for blocking or for staining.

Cell separation:

Blocking of nonspecific antibody binding to monocytes and other Fc γ RII⁺ cells in positive selection or depletion of Fc γ RII⁻ cells.

REFERENCES:

1. Van Sorge NM, van der Pol WL, van de Winkel JG: Fc γ RII polymorphisms: Implications for function, disease susceptibility and immunotherapy. *Tissue Antigens* 61:189, 2003
2. Looney RJ, Abraham GN, Anderson CL: Human monocytes and U937 cells bear two distinct Fc receptors for IgG. *J Immunol.* 136:1641, 1986
3. Van den Herik-Oudijk IE, Westerdal NAC, de Haas M, Kemper L, Capel PJA, van dem Borne AEGKr, Van de Winkel JGJ: Binding heterogeneity within the CD32 panel of mAb, in Schlossman SF, Boumsell L, Gilks W, Harlan JM, Kishimoto T, Morimoto C, Ritz J, Shaw S, Silverstein R, Springer T, Tedder TF, Todd RF (eds): *Leucocyte Typing V*, New York, Oxford University Press, 1995, p 832
4. Tomiyama Y, Kunicki TJ, Zipf TF, Ford SB, Aster RH: Response of human platelets to activating monoclonal antibodies: importance of Fc gamma RII (CD32) phenotype and level of expression. *Blood* 80:2261, 1992

**THIS REAGENT IS FOR RESEARCH ONLY.
 IT IS NOT TO BE ADMINISTERED TO HUMANS.**

StemCell Technologies

In North America

Tel: 1.604.877.0713
 Fax: 1.604.877.0704
 Toll Free Tel: 1.800.667.0322
 Toll Free Fax: 1.800.567.2899
 e-mail: info@stemcell.com
 www.stemcell.com

In the United Kingdom

Tel: +44.(0).20.7537.7565
 Fax: +44.(0).20.7515.5408
 Toll Free within United Kingdom:
 Tel: 0800.731.27.14
 Fax: 0800.731.27.13
 e-mail: info@stemcellgb.com
 www.stemcellgb.com

In Europe

Tel: +33.(0).4.76.04.75.30
 Fax: +33.(0).4.76.18.99.63
 e-mail: info@stemcellfrance.com

Version 1.0.0
 Revised:
 September 2005

Printed on recycled paper.