

**APC/iFluor™ 750 Anti-human CD32  
Antibody \*IV.3\***Catalog number: 103201G0, 103201G1, 103201G2  
Unit size: 25 tests, 100 tests, 500 tests**Product Details**

|                    |                                                                             |
|--------------------|-----------------------------------------------------------------------------|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.                         |
| Expiration Date    | 12 months upon receiving                                                    |
| Concentration      | 0.1 mg/mL                                                                   |
| Formulation        | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA |

**Antibody Properties**

|                    |                             |
|--------------------|-----------------------------|
| Species Reactivity | Human                       |
| Class              | Primary                     |
| Clonality          | Monoclonal                  |
| Host               | Mouse                       |
| Isotype            | Mouse igg2b                 |
| Immunogen          | CD32 (FcγRII, Fc gamma RII) |
| Clone              | IV.3                        |
| Conjugate          | APC/iFluor™ 750             |

**Biological Properties**

|             |                                                                                                                |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 750 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                          |

**Spectral Properties**

|                       |                 |
|-----------------------|-----------------|
| Conjugate             | APC/iFluor™ 750 |
| Excitation Wavelength | 754 nm          |
| Emission Wavelength   | 793 nm          |

**Applications**

IV.3 is an anti-human monoclonal antibody that forms an immune complex with the CD32 antigen. CD32 (sometimes referred to as FCGR2A or FcγRII) is a 40 kD single-pass type I membrane protein that is expressed on the surface of cells such as . CD32 plays a role in essential cellular pathways, namely, the Fc-gamma receptor signaling pathway involved in phagocytosis. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like . CD32 is a fairly uncommon antibody target, with a little more than 7000 publications in the last decade. Even still, CD32 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of

cell types, particularly in the study of immunology and innate immunity. This antibody was purified through affinity chromatography and conjugated to APC/iFluor™ 750 (ex/em = 754/793 nm).