

**iFluor™ 750 Anti-human CD16 Antibody**  
**\*3G8\***Catalog number: 101610L0, 101610L1  
Unit size: 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          |
| Immunogen          | CD16a (FCRIIA) |
| Clone              | 3G8            |
| Conjugate          | iFluor™ 750    |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Dark blue liquid                                                                                           |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 750 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 750 |
| Excitation Wavelength | 757 nm      |
| Emission Wavelength   | 779 nm      |

**Applications**

The 3G8 monoclonal antibody recognizes human CD16a, a 50 - 65 kD transmembrane protein typically found on the surface of macrophages, natural killer cells, dendritic cells, monocytes and neutrophils. CD16 is a member of key 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 IgG Fc. CD16 is a very popular antibody target, with over 25000 publications in the last decade. CD16a is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This

antibody was purified through affinity chromatography and conjugated to iFluor™ 750 (ex/em = 757/779 nm).