

**PacBlue Anti-human CD16 Antibody \*3G8\***Catalog number: 101611J0, 101611J1  
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 (FCRIIIA) |
| Clone              | 3G8             |
| Conjugate          | PacBlue         |

**Biological Properties**

|             |                                                                                                        |
|-------------|--------------------------------------------------------------------------------------------------------|
| Appearance  | Light yellow liquid                                                                                    |
| Preparation | Antibody purified by affinity chromatography and then conjugated with PacBlue under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                            |

**Spectral Properties**

|                       |         |
|-----------------------|---------|
| Conjugate             | PacBlue |
| Excitation Wavelength | 404 nm  |
| Emission Wavelength   | 455 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 PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser

and 460/22 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).