

**PE Anti-human CD64 Antibody \*10.1\***Catalog number: 106401L0, 106401L1, 106401L2  
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 IgG1   |
| Immunogen          | CD64 (FcR I) |
| Clone              | 10.1         |
| Conjugate          | PE           |

**Biological Properties**

|             |                                                                                                   |
|-------------|---------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                             |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | PE     |
| Excitation Wavelength | 566 nm |
| Emission Wavelength   | 574 nm |

**Applications**

The 10.1 monoclonal antibody binds with human CD64, a 72 kD single-pass type I membrane protein typically expressed on the surface of granulocytes, monocytes and dendritic cells. In some organisms, CD64 enhances protein tyrosine kinase activity. Moreover, it is a member of vital cellular pathways, for example, the interferon-gamma-mediated signaling pathway and 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 such as . CD64 is a fairly uncommon antibody target, with a little more than 4000 publications in the last decade. Even still, CD64 is often 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 PE (ex/em = 566/574 nm). It is compatible with the 561 nm laser

and 586/15 nm bandpass filter (for example, as in the BD FACSCelesta™).