

**iFluor™ 633 Anti-human CD14 Antibody**  
**\*61D3\***Catalog number: 101410E0, 101410E1  
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          | CD14 (LPS-Receptor) |
| Clone              | 61D3                |
| Conjugate          | iFluor™ 633         |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Black liquid                                                                                               |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 633 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 633 |
| Excitation Wavelength | 640 nm      |
| Emission Wavelength   | 654 nm      |

**Applications**

61D3 is an anti-human monoclonal antibody that targets the CD14 antigen. CD14 (sometimes referred to as myeloid cell-specific leucine-rich glycoprotein or LPS receptor) is a transmembrane protein that is located on the surface of cells such as macrophages. CD14 acts in critical cellular pathways, for example, the toll-like receptor signaling pathway, cell surface receptor signaling pathway and lipopolysaccharide-mediated signaling pathway. Moreover, in certain organisms, it promotes interleukin-8 secretion, is a positive regulator of tumor necrosis factor production and is a promoter of type I interferon production. From a research standpoint, it is of biological interest due to its association with

essential macromolecules/ligands like LY96. CD14 is a very popular antibody target, with over 42000 publications in the last decade. CD14 is essential for immunology, cell biology and neuroscience research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 633 (ex/em = 640/654 nm). It is compatible with the 640 nm laser and 660/20 nm bandpass filter (for example, as in the BD FACSJazz™).