

**XFD350 Anti-human CD14 Antibody *61D3,
XFD350 Same Structure to Alexa Fluor™
350***Catalog number: 10141140, 10141141
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 | AF350 |

Biological Properties

| | |
|-------------|--|
| Appearance | Yellow liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with AF350 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|--------|
| Conjugate | AF350 |
| Excitation Wavelength | 343 nm |
| Emission Wavelength | 441 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 XFD350 (ex/em = 343/441 nm). XFD350 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 350 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).