

PE/Cy5 Anti-human CD14 Antibody *61D3*Catalog number: 101411M0, 101411M1, 101411M2
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
Immunogen	CD14 (LPS-Receptor)
Clone	61D3
Conjugate	PE/Cy5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Cy5 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Cy5
Excitation Wavelength	565 nm
Emission Wavelength	666 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 PE/Cy5 (ex/em = 565/666 nm). It is compatible with the 561 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).