

iFluor™ 750 Anti-human CD14 Antibody
61D3Catalog number: 101410L0, 101410L1
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™ 750

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 750
Excitation Wavelength	757 nm
Emission Wavelength	779 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™ 750 (ex/em = 757/779 nm).