

iFluor™ 610 Anti-human CD14 Antibody
HI221Catalog number: 101400D0, 101400D1
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
Isotype	Mouse IgM
Immunogen	CD14 (LPS-Receptor)
Clone	HI221
Conjugate	iFluor™ 610

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 610 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 610
Excitation Wavelength	610 nm
Emission Wavelength	628 nm

Applications

HI221 is an anti-human monoclonal antibody that is specific for the CD14 antigen. CD14 (sometimes referred to as myeloid cell-specific leucine-rich glycoprotein or LPS receptor) is a glycoprotein that is found on the surface of cells like granulocytes and macrophages. CD14 plays a role in vital cellular pathways, namely, the apoptotic signaling pathway, lipopolysaccharide-mediated signaling pathway and cell surface receptor

signaling pathway. In addition, in many organisms, it enhances interleukin-8 secretion, acts to positively regulate interferon-gamma production and positively regulates NIK/NF-kappaB signaling. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as Endotoxin and TLR4. CD14 is a very popular antibody target, with over 42000 publications in the last decade. CD14 has a variety of applications in neuroinflammation, neuroscience and cell biology 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™ 610 (ex/em = 610/628 nm). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).