

**mFluor™ Blue 570 Anti-human CD114
Antibody *LMM741***Catalog number: 111400T0, 111400T1
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 IgG1 kappa
Immunogen	CD114 (CSF3R, GCSFR)
Clone	LMM741
Conjugate	mFluor™ Blue 570

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Blue 570 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Blue 570
Excitation Wavelength	553 nm
Emission Wavelength	565 nm

Applications

The LMM741 monoclonal antibody binds to human CD114, a 130 kD single-pass type I membrane protein typically found on the surface of endothelial cells, platelets, myeloid progenitor cells, neutrophils and granulocytes. CD114 acts in key cellular pathways, in particular, the cytokine-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with key

macromolecules/ligands such as jak2, jak1 and G-CSF. CD114 is a relatively rare antibody target, with fewer than 100 publications in the last decade. Even still, CD114 is essential for immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to mFluor™ Blue 570 (ex/em = 553/565 nm). It is compatible with the 561 nm laser and 577/15 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).