

FITC Anti-human CD163 Antibody *GHI/61*Catalog number: 116301G0, 116301G1
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	CD163 (M130, GHI/61, RM3/1)
Clone	GHI/61
Conjugate	FITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

GHI/61 is an anti-human monoclonal antibody that forms an immune complex with the CD163 antigen. CD163 (alternatively called GHI/61, M130 or RM3/1) is a 134 kD glycoprotein that is found on the surface of cells such as macrophages. CD163 is associated with a variety of biologically interesting macromolecules/ligands, for instance, hemoglobin. CD163 is a fairly uncommon antibody target, with a little more than 9600 publications in the last decade. Even still, CD163 has a variety of applications in innate immunity, cell biology and immunology 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 FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 525/50 nm bandpass filter (for

example, as in the Miltenyi Biotec MACSQuant Analyzer 16).