

**iFluor™ 820 Anti-human CD163 Antibody**  
**\*GHI/61\***Catalog number: 116300P0, 116300P1  
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	iFluor™ 820

**Biological Properties**

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 820 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 820
Excitation Wavelength	822 nm
Emission Wavelength	850 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 iFluor™ 820 (ex/em = 822/850 nm).