

**mFluor™ Green 620 Anti-human CD16
Antibody *3G8***Catalog number: 101610U0, 101610U1
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	CD16a (FCRIIIA)
Clone	3G8
Conjugate	mFluor™ Green 620

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Green 620
Excitation Wavelength	525 nm
Emission Wavelength	623 nm

Applications

The 3G8 monoclonal antibody recognizes human CD16a, a 50 - 65 kD transmembrane protein typically found on the surface of macrophages, natural killer cells, dendritic cells, monocytes and neutrophils. CD16 is a member of key cellular pathways, namely, the Fc-gamma receptor signaling pathway involved in phagocytosis. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like IgG Fc. CD16 is a very popular antibody target, with over 25000 publications in the last decade. CD16a is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This

antibody was purified through affinity chromatography and conjugated to mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 610/20 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).