

**mFluor™ Violet 450 Anti-human CD32
Antibody *3D3***Catalog number: 103210Z0, 103210Z1
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	CD32 (FcγRII, Fc gamma RII)
Clone	3D3
Conjugate	mFluor™ Violet 450

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 450
Excitation Wavelength	406 nm
Emission Wavelength	445 nm

Applications

The 3D3 monoclonal antibody binds to human CD32, a 40 kD member of the Ig superfamily often located on the surface of platelets, monocytes, B cells, granulocytes and dendritic cells. CD32 is involved with essential 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 essential

macromolecules/ligands. CD32 is a fairly uncommon antibody target, with a little more than 7000 publications in the last decade. Even still, CD32 has been widely used in innate immunity 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 mFluor™ Violet 450 (ex/em = 406/445 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD FACSymphony™ A5).