

## APC/XFD750 Anti-human CD32 Antibody \*3D3\*

Catalog Number: 103211E0,  
103211E1, 103211E2  
Unit Size: 25 tests, 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	Lot specific (please consult certificate of analysis for given lot)
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	APC/AF750

### Biological Properties

---

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/AF750 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

### Spectral Properties

---

Conjugate	APC/AF750
-----------	-----------

Excitation Wavelength 651 nm

Emission Wavelength 785 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 APC/XFD750 (ex/em = 756/785 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).