

iFluor™ 430 Anti-human CD32 Antibody
IV.3Catalog number: 10320030, 10320031
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 igg2b
Immunogen	CD32 (FcγRII, Fc gamma RII)
Clone	IV.3
Conjugate	iFluor™ 430

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 430 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 430
Excitation Wavelength	433 nm
Emission Wavelength	498 nm

Applications

IV.3 is an anti-human monoclonal antibody that forms an immune complex with the CD32 antigen. CD32 (sometimes referred to as FCGR2A or FcγRII) is a 40 kD single-pass type I membrane protein that is expressed on the surface of cells such as . CD32 plays a role in 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 important macromolecules/ligands like . CD32 is a fairly uncommon antibody target, with a little more than 7000 publications in the last decade. Even still, CD32 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology and innate immunity. This antibody was purified through affinity chromatography and conjugated to iFluor™ 430 (ex/em = 433/498 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ III).