

PE/XFD610 Anti-human CD79b Antibody
***CB3-1, XFD610 Same Structure to Alexa**
Fluor™ 610*

Catalog number: 107911P0, 107911P1, 107911P2
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	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	CD79b (B29, IGB)
Clone	CB3-1
Conjugate	PE/AF610

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/AF610
Excitation Wavelength	567 nm
Emission Wavelength	627 nm

Applications

CB3-1 is an anti-human monoclonal antibody that recognizes the CD79b antigen. CD79b (sometimes called B29) is a 37 - 39 kD single-pass type I membrane protein that is located on the surface of cells such as B cells. CD79b has been closely linked to essential biological processes like immune response, particularly adaptive immune response. Additionally, it is a member of essential cellular pathways, for example, the B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as CD79a, CD19, CD22 and CD5. CD79b is a fairly uncommon antibody target, with a little more than 1700 publications in the last decade. Even

still, CD79b is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 610 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).