

iFluor™ 450 Anti-human CD79b Antibody
CB3-1Catalog number: 10791040, 10791041
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	CD79b (B29, IGB)
Clone	CB3-1
Conjugate	iFluor™ 450

Biological Properties

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

Spectral Properties

Conjugate	iFluor™ 450
Excitation Wavelength	451 nm
Emission Wavelength	502 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 iFluor™ 450 (ex/em = 451/502 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).